



GEM5000 Gas Analyser

Operating Manual



Geotechnical Instruments (UK) Ltd

Sovereign House

Queensway

Leamington Spa

Warwickshire

CV31 3JR

England

Tel: +44 (0)1926 338111

Fax: +44 (0)1926 338110

Email: sales@geotech.co.uk

Website: www.geotechuk.com

Table of contents

1.0	MANUAL GUIDELINES	5
1.1	Hazard warnings and safety symbols	5
1.2	Notes.....	5
2.0	INTRODUCTION	6
2.1	Safety instructions.....	6
2.2	Instructions for safe use.....	7
2.3	MCERTS	9
2.4	CIRIA.....	9
2.5	EU Declaration of conformity.....	10
3.0	THE GEM5000 GAS ANALYSER	12
3.1	The GEM5000	12
3.2	GEM5000 standard product.....	13
4.0	GEM5000 OPTIONAL PRODUCTS AND ACCESSORIES.....	14
4.1	Optional products	14
4.1.1	Pitot tube (optional)	14
4.1.2	Orifice plate (optional).....	14
4.1.3	Temperature probe (optional).....	14
4.1.4	Anemometer (optional).....	14
4.1.5	H ₂ S filter (optional).....	15
4.1.6	Gas analyser manager – GAM (optional)	15
4.1.7	GPS (optional).....	15
4.1.8	Bluetooth.....	15
4.2	Instrument accessory products	16
5.0	GEM5000 INSTRUMENT FEATURES.....	17
5.1	Physical characteristics of the instrument panel.....	17
5.2	Analyser features and keys	18
5.3	Instrument connection points.....	19
6.0	GENERAL OPERATIONAL INSTRUCTIONS	20
6.1	Switching the instrument on	20
6.1.1	Power on self-test	20
6.2	Switching the analyser off	20
6.3	Instrument status icons.....	21
6.4	Instrument LED power states	22
6.5	Changing between parameters	22
6.6	Entering data	22
6.7	Instrument main gas read screen	23
6.8	Storage.....	23
6.9	Battery/charging	23
6.10	Cleaning instructions.....	24
6.11	Memory.....	25
6.12	Warning and error codes	25
6.13	Remote updates	25
6.13.1	What you need	25
6.13.2	Registration	25
6.13.3	Installation of the 5000 updater software.....	26
6.13.4	Installation of XP drivers.....	26
6.13.5	Installation of Vista and Win 7 drivers	26
6.13.6	Updating the analyser	26
6.13.7	Update from file.....	27
7.0	OPERATOR SETTINGS.....	28
7.1	Menu key	28

7.2	Device settings	28
7.2.1	Date and time	28
7.2.2	Bluetooth.....	29
7.2.3	Device information.....	30
7.2.4	Diagnostics	30
7.2.5	Navigation (optional).....	31
7.3	User settings	32
7.3.1	Operating language.....	32
7.3.2	Units of measurement	33
7.3.3	ID selection.....	34
7.3.4	Routes.....	36
7.3.4	Adjust backlight.....	36
7.3.5	Adjust volume	37
7.3.6	User Prompts	37
7.4	Operation settings	38
7.4.1	Timers.....	38
7.4.2	Gas Check	39
7.4.3	View data	39
7.4.4	Set alarms	40
7.4.5	Adjust flow fail	41
7.4.6	Technician login.....	42
7.5	Exit menu.....	43
8.0	TAKING READINGS.....	44
8.1	Preliminary checks before taking readings (best practice)	44
8.1.1	Creating an ID	45
8.2	Change screen layout	46
8.3	Answering site questions	46
8.4	Answering ID questions	46
8.5	Special action	47
8.5.1	Configuration of the data logging option	48
8.5.2	Profiling option	48
8.6	GEM analyser in GA mode	49
8.7	The gas flow measurement screen.....	50
8.8	How to use an anemometer (optional)	50
8.9	How to use a pitot tube (optional)	53
8.10	How to use an H ₂ S filter (optional).....	54
8.8.1	Cross gas effects on chemical cells.....	54
8.9.2	Cross-gas effects on methane, carbon dioxide and oxygen.....	55
8.11	How to use a temperature probe (optional).....	57
8.12	How to identify a borehole using the GPS feature (optional).....	57
8.13	Taking gas and flow measurement.....	61
9.0	CALIBRATION	62
9.1	Calibration introduction	62
9.2	Frequency of calibration – best practice.....	62
9.3	Calibration gases.....	63
9.4	Calibration set-up.....	64
9.5	Calibration equipment.....	64
9.6	Gas analyser.....	65
9.7	Calibration processes – best practice	65
9.7.1	Gas Check in fresh air	66
9.7.2	Calibration – mixtures 1, 2, & 3.....	67
9.8	Restore to factory.....	68
9.9	Calibration history	68
9.10	Calibration summary.....	68
10.0	PROBLEM SOLVING	69
11.0	WARNING AND ERROR DISPLAY.....	69

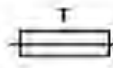
12.0	SERVICE	70
13.0	WARRANTY POLICY	71
14.0	IMPORTANT NOTICE TO ALL CUSTOMERS.....	72
15.0	APPENDICES – SAFETY INSTRUCTIONS.....	73
15.1	Instructions for safe use – Italian language.....	73
15.2	Instructions for safe use – German language	79
15.3	Instructions for safe use – French language	85
15.4	Instructions for safe use – Spanish language	91
15.5	Instructions for safe use – Portuguese language	97
15.6	Instructions for safe use – Chinese language	99
15.7	Instructions for safe use – Polish language	103
15.8	Instructions for safe use – Romanian language	106
17.0	GLOSSARY OF TERMS	109

1.0 Manual guidelines

1.1 Hazard warnings and safety symbols

 Warning	Information in this manual that may affect the safety of users and others is preceded by the warning symbol. Caution - Failure to follow the correct information may result in physical injury which in some cases could be fatal. If the equipment is used in a manner not specified by the manufacturer, the protection provided by the equipment may be impaired.
--	--

General product label symbols are listed as follows:

	CE conformity-The CE-marking is the manufacturer's statement to the EU authorities that the product complies with all relevant CE-marking Directives.		If the CSA mark appears with the indicator "US" or "NRTL" it means that the product is certified for the U.S. market, to the applicable U.S. standards.
	VDE mark is a symbol for electrical, mechanical, thermal, toxic, radiological and other hazards.		Separate collection, handling and disposal for waste electrical and electronic equipment and its components.
	Electric shock warning.		Refer to operators manual.
	Double insulated construction - does not require an Earth.		Specific marking of explosion protection (ATEX only).
II 2G	Equipment group and category. G = gases; the type of explosive atmosphere.		IECEx licenced mark (IECEx only).
	Fuse.		Equipment for indoor use only.

1.2 Notes

Important/useful information and instructions are shown clearly throughout the manual in a note format. For example:

 **Note:** For further information please contact Technical Support at Geotech on +44(0)1926 338111 or email technical@geotech.co.uk

2.0 Introduction

This manual explains how to use the GEM5000 landfill gas analyser. The GEM5000 has additional functionality to the GA5000 gas analyser. The GA5000 measures gas concentrations only, whereas the GEM5000 measures flow and gas concentrations. The GEM5000 measures flow and calculates the calorific values of the gas as well as being a useful tool for balancing the gas field. The GEM5000 gas analyser may be used to monitor, calculate, adjust and record the flow at each monitoring point.

This instrument may also be used in GA5000 mode of operation if required. The operator may change the analyser between a gas extraction monitor (a GEM5000 gas analyser) or a landfill gas analyser (a GA5000 gas analyser). The mode of operation can be changed from the 'Special Action' menu. Please refer to section '8.0 – Taking Readings' for further instruction.

The 5000 series of gas analysers complies with Part 15 of the FCC Rules. Operation is subject to the following two conditions:

- 1) This device may not cause harmful interference.
- 2) This device must accept any interference received, including interference that may cause undesired operation.

2.1 Safety instructions

 Warning	The 5000 series of gas analysers can be used for measuring gases from landfill sites and other sources as described in this manual. The operator may be exposed to harmful gases during the use of the instrument. Inhaling these gases may be harmful to health and in some cases may be fatal. It is the responsibility of the user to ensure that he/she is adequately trained in the safety aspects of the gases being used and appropriate procedures are followed. In particular, where hazardous gases are being used the gas exhausted from the analyser must be piped to an area where it is safe to discharge the gas. Hazardous gas can also be expelled from the instrument when purging with clean air. The instrument has been designed to be used in explosive atmospheres as defined by the classification. The instrument can be configured to measure low levels of several gases, but may not be certified for use in potentially explosive atmospheres of these gases. It is the responsibility of the operator to determine the protection concept and classification of equipment required for a particular application and whether these gases create a potentially explosive atmosphere.
--	---

 **Note:** Gas analysers are a sensitive piece of scientific equipment, and should be treated as such. If the equipment is used in a manner not specified by the manufacturer, the protection provided by the instrument may be impaired.

2.2 Instructions for safe use

For ATEX and IECEx the 5000 series of gas analysers are certified to Hazardous Area Classification

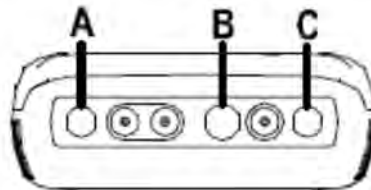
Ex II 2G Ex ib IIA T1 Gb (Ta = -10°C to +50°C)

It is vital instructions are followed closely. It is the responsibility of the operator to determine the protection concept and classification required for a particular application.

(Reference European ATEX Directive 94/9/EC, Annex II, 1.0.6)

The following instructions apply to equipment covered by certificate numbers SIRA 11ATEX2197X and IECEx SIR 11.0089X:

- The equipment may be used with flammable gases and vapours with apparatus group IIA and temperature class T1.
- The equipment can contain gas sensing heads for the detection of particular gases. The inclusion of a sensor does not infer that the equipment is suitable for the use of gases with a temperature class of less than T1.
- The equipment is only certified for use in ambient temperatures in the range -10°C to +50°C and should not be used outside this range.
- The equipment must not be used in an atmosphere of greater than 21% oxygen.
- Repair of this equipment shall be carried out in accordance with the applicable code of practice.
- When used in a hazardous area only use GF5.2 temperature probe (SIRA 11ATEX2197X and IECEx SIR11.0089X). For connector C, the GF5.4 anemometer (BVS 04ATEXE194) for use with ATEX only. The analyser should not be connected to any other devices in the hazardous area including the GF-USB lead (connector A) or GF3.9 battery charger (connector B) supplied with the analyser.



Do not charge, recharge or open in a potentially explosive atmosphere.
In hazardous area only use "Temperature Probe GF5.2" in Connector B.
Connector C (Uo=10V,Io=5mA,Po=50mW,Ci=0,Li=0,Co=100uF,Lo=1000mH),
Connector B (Uo=5V,Io=6mA,Po=7mW,Ci=0,Li=0,Co=100uF,Lo=1000mH)

MAXIMUM NON-HAZARDOUS SUPPLIES:

Connector A - Um=6V Connector B - Um=10.1V

- The safe area apparatus that is to be connected to the USB Port shall be a Safety Extra Low Voltage (SELV) or Protective Extra Low Voltage (PELV) circuit.
- Only a Geotechnical Instrument battery pack part number 20087 or 2011113 is permitted as a replacement. This battery pack shall only be changed in a safe area by Geotech personnel or authorised distributors.
- Only Battery Charger type GF3.9 shall be used to recharge the batteries via Connector 'B'.
- If the equipment is likely to come into contact with aggressive substances, e.g. acidic liquids or gases that may attack metals, or solvents that may affect polymeric materials, then it is the responsibility of the user to take suitable precautions, e.g. regular checks as part of routine inspections or establishing from the material's data sheet that it is resistant to specific chemicals that prevent it

from being adversely affected, thus ensuring that the type of protection is not compromised.

- The relative pressure range is +/-500 mbar. Note, however, that the input pressure should not exceed +/- 500 mbar relative to atmospheric pressure and the output pressure should not exceed +/- 100 mbar relative to atmospheric pressure.

For CSA (Canada) the 5000 series of gas analysers are certified to Hazardous Area Classification

CLASS 2258 03 - PROCESS CONTROL EQUIPMENT - Intrinsically Safe and Non-Incendive Systems - For Hazardous Locations



Ex ib IIA:

Model GA 5000, GEM 5000 and BIOGAS 5000 Methane Detectors; portable, battery powered with non-field-replaceable Battery Pack P/N 20087 or 2011113; intrinsically safe and providing intrinsically safe circuits ("ib" for Zone 1) to Model GF5.2 Temperature Probe (Connector B) and with entity output parameters as tabulated below; Temperature Code T1; $-10^{\circ}\text{C} \leq T_{\text{amb.}} \leq +50^{\circ}\text{C}$.

Connector	Entity Parameters						
	Uo (V)	Io (mA)	Po (mW)	Co (uF)	Lo (mH)	Ci (uF)	Li (mH)
B	5.0	6	7	100	1000	0	0
C	10.0	5	50	100	1000	0	0

Note: This device has been investigated for electrical safety features only.

For CSA (USA) the 5000 series of gas analysers are certified to Hazardous Area Classification

CLASS 2258 83 - PROCESS CONTROL EQUIPMENT - Intrinsically Safe and Non-Incendive Systems - For Hazardous Locations - CERTIFIED TO U.S. STANDARDS



AEx ib IIA:

Model GA 5000, GEM 5000 and BIOGAS 5000 Methane Detectors; portable, battery powered with non-field-replaceable Battery Pack P/N 20087 or 2011113; intrinsically safe and providing intrinsically safe circuits ("ib" for Zone 1) to Model GF5.2 Temperature Probe (Connector B) and with entity output parameters as tabulated below; Temperature Code T1; $-10^{\circ}\text{C} \leq T_{\text{amb.}} \leq +50^{\circ}\text{C}$.

Connector	Entity Parameters						
	Uo (V)	Io (mA)	Po (mW)	Co (uF)	Lo (mH)	Ci (uF)	Li (mH)
B	5.0	6	7	100	1000	0	0
C	10.0	5	50	100	1000	0	0

Note: This device has been investigated for electrical safety features only.

2.3 MCERTS

MCERTS is the UK Environment Agency's Monitoring Certification Scheme. The scheme provides a framework within which environmental measurements can be made in accordance with the Agency's quality requirements. The scheme covers a range of monitoring, sampling and inspection activities.

MCERTS promotes public confidence in monitoring data and provides industry with a proven framework for choosing monitoring systems and services that meet the Environment Agency's performance requirements.

The Environment Agency has established its Monitoring Certification Scheme (MCERTS) to deliver quality environmental measurements. The MCERTS product certification scheme provides for the certification of products according to Environment Agency performance standards, based on relevant CEN, ISO and national standards.

MCERTS certified instruments have been tested by an independent body to ensure that they meet certain performance requirements. In addition the manufacturer of an MCERTS product is regularly audited to ensure that the performance requirements of the certification are being continually met.

The 5000 series of gas analysers have been certified to Version 3.1 of the 'Performance Standards for Portable Emission Monitoring Systems'.

2.4 CIRIA

The CIRIA guideline 'Assessing the risks posed by hazardous ground gases to buildings' proposes that gas concentrations and flow rates should be monitored.

As an example methodology, they suggest using a gas analyser to first measure flow and pressure and then afterwards to measure gas concentration.

The logging profile option offers frequency of data to be collected within a timed period which, in return, identifies a gas profile of the sample point being monitored, information about whether the sample point is performing correctly, when the peaks occur and whether air is drawn in after a certain period. This logging option is available on firmware software version 1.6.5

Versions of the GA5000 analyser range with internal flow on firmware version 1.6.5 and above have the ability to take measurements according to the CIRIA guidelines, while still allowing other users to take the measurements as before.

2.5 EU Declaration of conformity



EU Declaration of Conformity

This Declaration of Conformity is issued under the sole responsibility of the manufacturer:

Geotechnical Instruments (UK) Ltd.
Sovereign House,
Queensway
Leamington Spa, Warwickshire
CV31 3JR
ENGLAND

Product: GA5000, GEM5000, BIOGAS 5000

Type of equipment:

GA5000 - Landfill Gas Analyser

GEM5000 - Landfill Gas Analyser and Extraction Monitor

BIOGAS 5000 – Anaerobic Digester Gas Analyser



The GA5000, GEM5000 & BIOGAS 5000 described above are in conformity with the relevant Union harmonisation legislation:

2014/34/EU: Equipment and protective systems intended for use in potentially explosive atmospheres (ATEX)

CSA Group (0518) performed assessment against:

- EN60079-0:2006
- EN60079-0:2009
- EN60079-11:2007

Issuing certificate number SIRA 11ATEX2197X

2014/30/EU: Electromagnetic capability (EMC)

- EN 301 489 Pt 1 (V1.9.1 – 2011-04)
- EN 301 489 Pt 17 (V2.1.1 – 2009-05)
- EN 301 489 Pt 19 (V1.2.1 – 2002-11)
- BS EN 61000-3-2:2006 + A2:2009
- BS EN 61000-3-3:2008

In addition, the following International requirements are met:



International Electrotechnical Commission system for certification to standards relating to equipment for use in explosive atmospheres (IECEx System)

CSA Group (0518) performed assessment against:

- IEC60079-0:2004 Ed4
- IEC60079-0:2007 Ed5
- IEC60079-11:2006 Ed5[list standards applied]

Issuing certificate number SIR 11.0089X



The CSA mark shows that our products have been certified by an accredited third party laboratory and have met applicable standards as required by North American law (cCSAus)

CSA International performed assessment against:

- C22.2 No. 0-10
- CAN/CSA-C22.2 No. 60079-0:07
- CAN/CSA-C22.2 No. 60079-1:07
- CAN/CSA-E60079-11:02
- ANSI/UL 60079-0:09
- ANSI/UL 60079-1:09
- ANSI/UL 60079-11:09

Issuing certificate number CSA 11 2445306

Signed for and on behalf of:

Name: Mr. Dean Kavanagh

Position: Programme Director

Done at: Geotechnical Instruments (UK) Ltd

On: 20th April 2016

3.0 The GEM5000 gas analyser

3.1 The GEM5000



GEM5000 gas analyser

The GEM5000 gas analyser is designed to monitor landfill gas extraction systems.

Benefits:

- Allows balancing of gas extraction site.
- Maximise power generation from site.
- Field proven.
- Standardises monitoring routines.
- Easy transfer of data.
- GPS for compliance.

Features:

- ATEX certified.
- MCERTS certified.
- H₂ compensated CO.
- Calculates flow (m³/hr) and calorific value (kW or BTU).
- Technician log-in.
- Event log.
- Two instruments in one (GA and GEM mode).
- Measures static and differential pressure.
- Simultaneous display of gases.
- Storage of changes in set-up of gas field.
- Data logging.

Applications:

- Gas extraction fields.
- Flare monitoring.
- Landfill sites.

3.2 GEM5000 standard product

GEM5000

Reference:

- A Hard carry case
- B In-line water trap tubing & filter
- C Gas analyser instrument
- D H2S filter (optional – if the compensated CO cell is fitted)
- E 5000 series safety manual
- F Mains battery charger and adaptors:
 - UK
 - Europe
 - US
 - Australia

4.0 GEM5000 optional products and accessories

4.1 Optional products

The GEM5000 gas analyser has a number of optional products for purchase which enhance the usability and enable further analysis of data and reading information.

✍ Note: For more information on the features listed in this section please contact Sales at Geotech on +44(0)1926 338111 or email Sales@geotech.co.uk.

4.1.1 Pitot tube (optional)

The GEM5000 gas analyser enables the use of a pitot tube to aid accurate flow measurement. The pitot tube is used for gas extraction systems and the pressure readings are taken in mbar. High gas flow is calculated in the analyser in m³/hr (metres cubed per hour).

4.1.2 Orifice plate (optional)

Geotech recommends the use of orifice plates as good practice when using the GEM5000 gas analyser, enabling repeatability in flow measurement via a fixed method. Orifice plates are a specialised product and configuration and Geotech does not currently manufacture or sell spare parts for this product. Please refer to a specialist supplier if the use of Orifice plates is required.

4.1.3 Temperature probe (optional)

The GEM5000 gas analyser has the facility to automatically display and record the borehole temperature via an optional temperature probe.

When a temperature probe is fitted the temperature reading will be displayed on the 'Main Gas Read Screen' and recorded with all other data.

✍ Note: Temperature probes with an Ex label are part of the GEM5000 Ex certification SIRA 11ATEX2197X and IECEx SIR11.0089X, and therefore certified for use under the same conditions as the analyser.

4.1.4 Anemometer (optional)

The GEM5000 gas analyser has the facility to automatically display and record high flow via an optional anemometer probe. It is designed to plug into the instrument and instantly provide a flow indication. An anemometer probe adds flow measurements to the professional reporting ability of the GEM5000 range along with gas concentrations, pressure and temperature.

The anemometer has a simple connection, a narrow diameter measurement head (11mm), a wide temperature operating range (up to 80°C) and indicates flows up to 40 m/sec.

When an anemometer probe is fitted to the analyser the flow will be displayed in the 'Main Gas Read Screen' and recorded with all other data.

Flow can be measured in either m/s (gas velocity) or m³/hr (volume flow rate). In order to calculate the volume flow rate the pipe diameter will need to be entered into the instrument, either manually or via the Gas Analyser Manager software.

✍ Note: The anemometer probe is ATEX certified for use in a potentially explosive atmosphere under Ex certificate BVS 04ATEXE194.

4.1.5 H₂S filter (optional)

The GEM5000 gas analyser has the capability to use an H₂S filter and is required as standard if the compensated CO cell is fitted and configured at the time the instrument is manufactured. H₂S gas can have a cross-gas effect on the CO reading. By using a filter, the H₂S is removed from the gas sample, therefore providing a more accurate CO reading.

The filter only needs to be used when you are trying to get rid of any possible cross gas effects H₂S might have on other gases. Do not use the filter on all boreholes.

4.1.6 Gas analyser manager – GAM (optional)

Gas Analyser Manager (GAM) enables the operator to maximise the operation of the gas analyser. It enables direct communication with the unit, features a simple upload and download facility and is fully compatible with the latest Microsoft™ operating systems.

Features:

- Organisation and transfer of borehole IDs and readings to and from the gas analyser.
- Configuration of the gas analyser.
- Flexible grouping of the IDs.
- Structured organisation of transferred data.
- Automatic detection of instrument type and available options.
- Secure data mode to prevent tampering.
- First time set-up wizard.
- Enable flow measurements for GEM5000 gas analysers.

4.1.7 GPS (optional)

An optional GPS feature is available for the GEM5000 gas analyser. It enables the site engineer to automatically locate borehole IDs using GPS satellite signal from predefined borehole IDs uploaded from Gas Analyser Manager or set on the analyser when out in the field prior to taking a reading. The GPS reading data is stored for each measurement reading providing an audit trail confirming that a reading was taken.

4.1.8 Bluetooth

The analysers are fitted with a Bluetooth receiver which enables the operator to download readings and upload IDs without the need to connect the analyser to a PC via a USB lead.

4.2 Instrument accessory products

Optional accessory and replacement parts must **ONLY** be purchased for the GEM5000 gas analyser direct from Geotech or your agent. Please contact sales@geotech.co.uk for further details on pricing and how to order.



Ref	Description	Order Code
A	Hard carry case	GF2.5
B	Soft carry case	GF5.1
C	Battery charger and adaptors	GF3.9
D	Anemometer (ATEX certified)	GF5.4
E	H ₂ S/Hydrocarbon filters H ₂ S/Hydrocarbon filters (pack of 5)	GA6.4 GA6.5
F	Temperature probe (ATEX certified)	GF5.2
G	Pitot tube kit	PITOT
H	In-line water trap filter & tubing	GF1.8
I	Water trap with barbed filters (pack of 10) Water trap with barbed filters (pack of 30)	GA4.9 GA4.9(30)
J	Gas ports connectors (pack of 10)	MC10
K	Sampling tube 5m length flexible PVC tubing (3/16" id)	GA2.3
L	Spare window for the soft carry case	GF5.3
M	Gas Analyser Manager (GAM) USB lead	GAM GFUSB
N	Check gas regulator and tubing for calibration gas. Calibration gas canister. Please contact sales@geotech.co.uk for gas canister concentrations	GA6.8

5.0 GEM5000 instrument features

5.1 Physical characteristics of the instrument panel

Front view:



Reference:

- A Main Gas Read Screen
- B Soft-keys
- C Backlight Key
- D Menu Key
- E Pump key
- F LED Light
- G On/Off Key
- H Assistance key
- I Enter Key
- J Scroll up key
- K Scroll down key
- L Key 0 – Space key

Back view:



Reference:

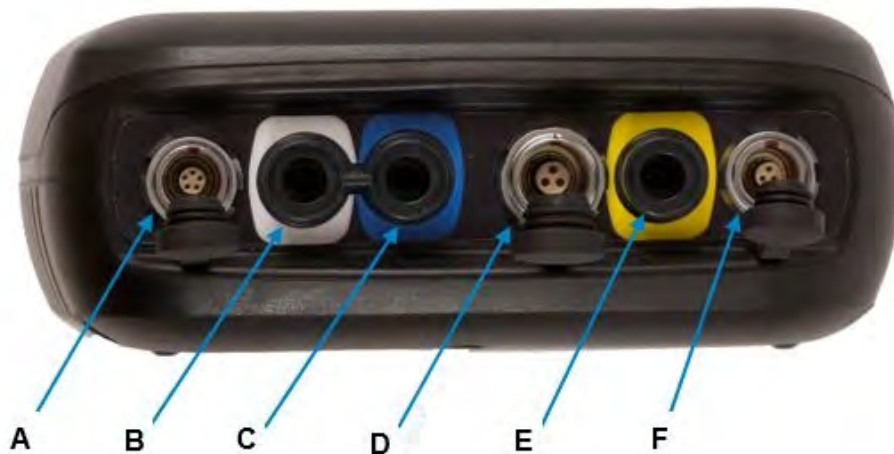
- M Model Number
- N Serial number
- O Part number
- P Certificate number
- Q Recalibrated date

5.2 Analyser features and keys

Ref	Feature	Function
A	Main Gas Read Screen	Start and end screen when using the instrument.
B	Soft-keys	The function of the three 'soft-keys' on the front of the instrument panel is determined by menu options taken. Functions vary from screen to screen.
C	Backlight key	Enables the operator to turn the backlight on/off on the analyser display panel.
D	Menu key	Press the 'Menu' key to view and maintain User, Device and Operation settings.
E	Pump key	Press the 'Pump' key to start or stop the pump.
F	LED light	LED power light is visible on the front of the analyser when the instrument is powered on.
G	On/Off key	Press the 'On/Off' key for 2 seconds to switch the instrument on and off.
H	Assistance key	Press for on-screen assistance and help.
I	Enter key	Use to accept changes, options, user inputted answers etc.
J	Key 2 – Page Up	Also 'Key 2'. Press scroll up to view further information on the instrument screen.
K	Key 8 – Page down	Also 'Key 8'. Enables the operator to scroll down to display more information.
L	Key 0 – Space key	Also 'Key 0'. Press to enter a space when entering text on the instrument screen.
M	Model Number	Instrument model type identification.
N	Serial Number	Unique identification for the instrument. Verification of the serial number will be required if Technical Support assistance is needed.
O	Part Number	Manufacturer's part number.
P	Certification Number	Displays instrument certification information.
Q	Recalibrated Date	The date displayed is the date the instrument is due to be calibrated.

5.3 Instrument connection points

Top view:



Ref:	Connection Point:	Function:
A	Connector A	Attach the USB lead for PC-to-analyser connectivity.
B	Inlet Port & Static Pressure Port (White port)	Attach the sample tube to take a gas sample. Also used to measure the static pressure.
C	Differential Pressure Port (Blue port)	Attach the sample tube to measure differential pressure.
D	Connector B	Attach the temperature probe and also used to attach the mains charger to the analyser for charging.
E	Gas Outlet Port (Yellow port)	The gas outlet port is the point at which the sample gas is expelled. Tubing may also be attached to the port.
F	Connector C	Attach the anemometer.

6.0 General operational instructions

6.1 Switching the instrument on

- 1) To switch on the analyser, press and hold the 'On/Off' key. The Geotech logo will display followed by the instrument warm up.
- 2) Following the instrument warm up, the 'Date and Time' screen is displayed prompting the technician to set the date and time and required format.
- 3) When complete select the soft-key to 'Exit' and the 'Power On Self-test' screen is displayed followed by instrument status. Instrument status displays the instruments service due date, serial number, options, service scheme and software version. Text will also display stating 'Self-test complete'.
- 4) Select the soft-key 'Next' to move onto the next screen and the 'Technician Login' screen is displayed.
- 5) Use the cursor keys to move through the list of ID's. Select either the required 'Technician ID' from the list followed by the soft-key 'Accept', or select 'Default' followed by the soft-key 'Accept' to continue to the 'Main Gas Read Screen'.

✍ Note: The selected technician ID is displayed at the top left corner of the Main Gas Read Screen.

6.1.1 Power on self-test

When switched on, the read-out will perform a pre-determined self-test sequence. During this time many of the analyser's functions are tested, including:

- General operation
- Gas flow measurement
- Calibration
- Battery charge level

During the self-test the following information is also displayed:

- Manufacturer's service due date
- The last gas check date
- Software version programmed
- Date format
- Serial number
- Operating language
- The currently enabled sales option

✍ Note: The self-test should only be done with the analyser sampling fresh air.

6.2 Switching the analyser off

- 1) To switch off the analyser, press and hold the 'On/Off' key, at which point a clean air purge will be carried out and the instrument will then switch off.
- 2) If for any reason the analyser 'locks up' and will not switch off in this manner, press and hold the 'On/Off' key for 15 seconds; this will force the instrument to switch off.

6.3 Instrument status icons

The following icons may be displayed on the instrument screen:

Icon	Description	Icon	Description
	Battery charge state Gives the operator an estimation of the battery charge state. For example 100% gives about 8 hours use in the field and 50% would mean that there is approximately 4 hours battery life remaining.		Battery charge state Indicates less than 2 hours of charge remaining.
	Pump status This icon is displayed along with a counter showing the pump run-time. This counts down where the operator has specified the pump run-time; if not it counts up; the icon turns red when stalled.		Pump stalled This icon is displayed when the pump stalls. The instrument's gas inlet (or outlet) may be blocked. This warning is most commonly caused by a water-logged or dirty sample filter. Change the sample filter and check for obvious blockages in the sample tubes. Alternatively, a small amount of adjustment can be made to the low flow detection point to compensate for minor changes in the performance of the pump fitted to the instrument.
	GPS signal strength This icon shows the signal strength the analyser's GPS module is able to provide. Full, okay and fair strength respectively.		GPS failure The GPS was unable to get a line of sight lock on enough satellites. Or, it may be that it hasn't had time to get a lock.
	This indicates when Bluetooth has been enabled. The colour changes from grey to blue when connected.		Language This icon indicates the currently selected operating language. This can be changed via the main menu.
	Data logging This icon indicates that the data logging feature is in operation.		Service overdue This icon indicates that the analyser is overdue for its service
	Legacy mode This icon indicates that the analyser is in legacy mode and hence is ready to connect to a PC.		USB disabled This icon indicates that the analyser has reached a battery critical state, and hence has turned off its USB connectivity.

6.4 Instrument LED power states

When the instrument is powered on a LED power light is visible on the front of the analyser, located above the 'On/Off' key. The following LED power light states are as follows:

Steady yellow	Unit turning on. This will extinguish when software has loaded correctly.
Flashing (rapid)	Unit is powering off.
Flashing (slow)	Power off is being delayed for purge/shutdown handling.
Flashing yellow	Unit is turning off due to power button being pressed.
Flashing red	Unit is turning off due to critically low battery.

✍ Note: Pressing and holding the power button for ~20s resets the analyser.

6.5 Changing between parameters

By default, the instrument displays the 'Main Gas Read Screen' (for gas measurement). The instrument will return to this screen after power on or when returning from the menus. The 'Scroll' keys can be used to switch to another measurement screen.

6.6 Entering data

During normal operation the operator may be prompted to enter data or information via the keypad, i.e. entering an ID code or setting an alarm level.

When entering data into the instrument all fields are fixed format and are populated from the left.

Text:

Entering text uses similar multi-tap functionality as a mobile phone. Key the numeric/alpha key pad the required number of times to select the appropriate letter. To key numeric data continue to press the numeric/alpha key until the required number is displayed.

Numeric data:

To enter a new date 09/11/11 the operator would type in 091111 using the numeric keypad in the following sequence:-

```
* 0 _/_/_
* 09/_/_
* 09/1 _/_
* 09/11/_
* 09/11/1_
* 09/11/11
```

Press the 'Enter' key to confirm/accept data keyed.

Any mistakes can be corrected using the soft-key 'Delete' which will delete the last digit typed. Alternatively, the sequence can be retyped before the 'Enter' key is pressed and the existing numbers will be pushed off the screen.

✍ Note: The instrument will not allow invalid data to be entered; this should be deleted and re-entered.

6.7 Instrument main gas read screen

The 'Main Gas Read Screen' is considered to be the normal operating screen and all operations are carried out from this starting point.

The actual data shown on this display will depend on the version of the instrument and the options that have been selected. In general, all of the main readings will be shown.



6.8 Storage

The analyser should not be exposed to extreme temperature. For example, do not keep the analyser in a hot car. When not in use analysers should be kept in a clean, dry and warm environment, such as an office and protect the analyser with either the soft carry case or store in the hard carry case provided with the instrument.

The instrument should be discharged and fully charged at least once every four weeks, regardless of indicated charge state.

6.9 Battery/charging

The battery used in the 5000 series of gas analysers is nickel metal hydride and manufactured as a pack from six individual cells. This type of battery is not so susceptible to the top-up charging 'memory effects' as nickel cadmium batteries, although it is not recommended that the unit is given small top-up charges.

Note: To reach optimum charge, it is recommended that the instrument is switched off when being charged and remains switched off during the charging process.

A full charge will take approximately 4 hours from a fully discharged battery.



Warning

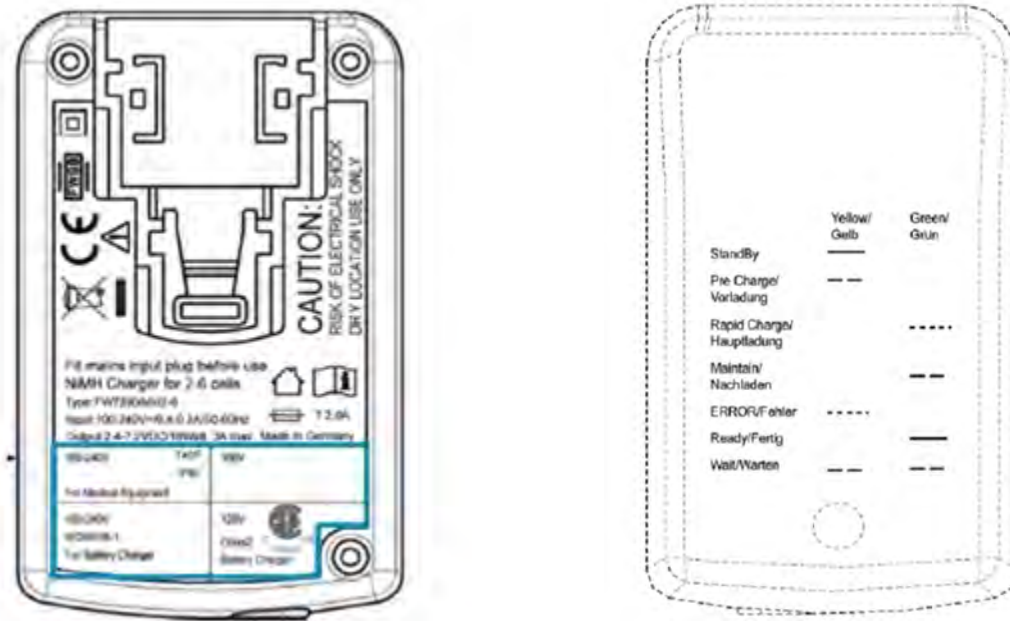
The battery charger is NOT covered by the Ex certification. The battery must be charged only in a safe area.

The battery charger is intelligent and will indicate when the unit is charging and charged.

The instrument must be charged **ONLY** using the battery charger supplied with the instrument. The battery charger supplied is intended for indoor use only. Please ensure adequate ventilation whilst charging. Typically, a fully charged battery will last 7-8 hours. A quick 30 minute charge can be used to give approximately one hours use in the field but this may shorten the battery life. Temperature can dramatically affect the battery life; please take this into account when estimating battery life.

✍ Note: Connect the charger to the mains attaching the appropriate adaptor.

Power supply front and back drawing:



Charger: Input voltage: 100-240V AC +/- 10%
 Input frequency: 50-60Hz +/- 10%
 Input current: 0.4A@100VAC .. 0.2A@240VAC

Output voltage: 10.1VDC max
 Output current: 1.5A max

✍ Note: This charger has been internally restricted to 1.5A

6.10 Cleaning instructions

Do NOT use any cleaning agents to clean the analyser or battery charger as they may have an adverse effect on the safe use of these devices.

6.11 Memory

The analyser's memory is stored in a readings and configuration database. The analyser will prompt when its memory is full, and you will not be able to store any further readings. Please download your readings via GAM or the Basic Download Software and then clear the memory.

✍ Note: The analyser should never be stored for prolonged periods with valuable data in its memory. It is advisable to download all readings to GAM at the end of each day's monitoring. To clear the memory, please refer to the Gas Analyser Manager (GAM) operating manual.

6.12 Warning and error codes

When switched on the instrument will perform a predetermined self-test sequence taking approximately ten seconds. During this time many of the instrument's working parameters and settings are checked. If any operational parameters are out of specification or if the pre-programmed recommended calibration/service date has passed, errors or warnings may be displayed.

✍ Note: For further information please refer to section '10.0 Problem Solving'.

6.13 Remote updates

The 5000 range of analysers (GA5000, GEM5000 and BIOGAS 5000) can be updated to the latest version of operating software using the free Geotech 5000 updater software. This allows customers to update their analyser's operating software at their convenience rather than waiting to return the analyser for service. The Geotech 5000 updater software needs to be installed on to a PC by a user with administrator rights, it is then ready to be used by the end-user. A GFUSB download cable is required along with a PC with access to a reasonably fast internet connection. This allows the users to check and download the latest data files from the web.

6.13.1 What you need

- Administrator rights (or access to IT person)
- Access to the internet and e-mail address
- Geotech's GFUSB download lead

6.13.2 Registration

- Go to <http://www.geotechuk.com/analyser-update-registration.aspx>
- Enter the registration information including the analyser's serial number (on the rear label)
- Wait for an e-mail with your access information and link to the files
- Select the embedded link which takes you to a new web page
- Enter the login details (email address and password) provided in the email you received from us when you registered (not your own email address)
- Select the "5000 series updater software.zip" file from available downloads.
- Select "Save as" and store it locally on your desktop or similar (this may take some time dependant on internet connection speed)

6.13.3 Installation of the 5000 updater software

This requires administrator rights.

- Unzip the file (extract contents), note its location (do NOT run from the compressed folder)
 - To do this, right click the zip folder and select 'extract all'
 - Save the new unzipped folder to a location on your computer
- Open this folder and double click the setup.exe (administration rights will be required)
- Follow the on-screen instructions, accepting where necessary. (Please note the progress bar doesn't move for about a minute)

6.13.4 Installation of XP drivers

This requires administrator rights.

- Locate the folder that contains the downloaded files (noted from before)
- Unzip the folder XPRNDIS_drivers.zip
- Turn the analyser on and wait for it to warm up
- Press the menu key
- Select device settings
- Select 3 for 'device information'
- Change the comms mode to "GA5K" by pressing the comms mode soft key
- Connect the analyser to the PC via the USB Lead
- Windows will start the "Found New Hardware Wizard"
- Select "No, not at this time" click next
- Select "Install from a list or specific location (advanced)" and then click next
- Unselect "search removable media (floppy, CD-ROM...)"
- Select "Include this location in the search:"
- Browse to the location of the XPRNDIS_drivers which were previously unzipped and click next
- Select continue or accept if any notifications appear.
- The drivers will now be installed

6.13.5 Installation of Vista and Win 7 drivers

This requires administrator rights.

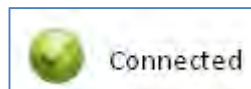
- Turn the analyser on and wait for it to warm up
- Press the menu key
- Select device settings
- Select 3 for 'device information'
- Change the comms mode to "GA5K" by pressing comms mode soft key
- Connect the analyser to the PC via the USB Lead
- Follow the onscreen instructions and accept all of the prompts
- The drivers will now be installed

6.13.6 Updating the analyser

Once all of the above operations are completed the software is fully installed and the instrument is ready to be updated. This can be done by a user who does not have administration rights on their PC.

- Ensure the analyser is fully charged - you do NOT want the analyser turning off during the update process
- Turn the analyser on and wait for it to start-up

- Navigate to the main read screen and press the menu key.
- Select device settings.
- Press 3 for 'device information'
- Change the comms mode to GA5K by pressing "Comms mode" soft key (**Please note:** if you have just installed the drivers the analyser may be in GA5K mode already)
- Connect the analyser to the PC via the Geotech GFUSB lead.
- Start the Geotech 5K updater software
- Press the "Check internet" button
- The software will now attempt the download the data files to the PC
- This may take several minutes depending on your internet speed
- Due to individual company's firewalls and I.T.' policies the update cannot always be done via this option. You will receive an error message at this point if unsuccessful. Refer to section 6.13.7 for an alternative method
- It may also take a minute at this point for the analyser to be detected as connected, please wait for the software to show the analyser as connected



- Once connected the software will be ready to update, press "update" which will copy the files to the analyser memory
- The analyser will then start to update, and may restart several times. Please wait for the analyser to show back on the update software as "connected". Do not turn off or disconnect the USB lead until this point
- The analyser is now fully updated. Please note that you may need to adjust the backlight on the analyser: (menu key > user settings > 4 Adjust Backlight > press 6 to make it lighter until you can see the screen. Remember to hit "save").
- Now set your analyser back to "legacy mode". (menu key > device settings > 3 Device information > press "Comms mode" soft key.) Your analyser will not work with GAMS unless it is in "legacy mode"

6.13.7 Update from file

If the update from "Check Internet" option does not work you can try the following steps. Each company may have certain IT policies in place that stop the download of certain files which may stop the successful download of the update file. You may need help from you I.T. department to download this.

- Click [here](http://www.geotechuk.com/G5000Files/G5000_UPDATE_FILE.5kc) to download the update file (http://www.geotechuk.com/G5000Files/G5000_UPDATE_FILE.5kc)
- When prompted select the save as function, and save the file to a location you can access later. Do not select open
- This may take several minutes depending on your internet speed
- Follow the instructions as per section 6.13.6 - Updating the analyser
- However, chose "Select file" instead of "Check internet"
- Navigate to where you saved the file (G5000_UPDATE_FILE.5kc)
- Click open to select the file
- The rest of the process is the same as 6.13.6 - Updating the analyser

7.0 Operator settings

7.1 Menu key



The 'Menu' key enables the operator to select options to set up specific parameters and perform operational tasks prior to sample readings being taken or to view data and information stored in the instrument.

- 1) Select the 'Menu' key on the front of the analyser and the following screen is displayed:



Device Settings

- 2) Press the relevant numeric key on the analyser keypad to select the required option.
- 3) To exit this menu, select the soft-key 'Exit' on the front of the analyser and the operator is returned to Main Gas Read Screen.

7.2 Device settings

7.2.1 Date and time



Date and Time

This option enables the operator to set the instrument date and time or to receive and update the settings automatically from satellite signal.

- 1) Select the 'Menu' key on the front of the analyser to display the 'Device Settings' menu followed by 'Key 1 – Date and Time' and the following screen is displayed:



- 2) Select 'Key 1 – Set Date' and key in the required date. Type the date using the numeric keypad. Press the soft-key 'Date Format' to toggle and select the required date format i.e. dd/mm/yy. Press the 'Enter' key to confirm and update the date setting.
- 3) Select 'Key 2 – Set Time' and key in the required time (hh:mm). Type the time using the numeric keypad and press the 'Enter' key to confirm the update.
- 4) The operator may also change the default time zone. Selecting the 'Key 4 Scroll-left' or 'Key 6 – Scroll right' to move through the different time zones. Press the 'Enter' key to confirm your default setting.
- 5) Select 'Key 3' to toggle between 'Manual Update' and 'Automatic Update' in order to choose how the date and time is set if updating from satellite signal.

Manual	Used to manually obtain and update the date and time from the satellite signal when requested. Select soft-key 'Set now' to set date and time from satellite when available.
--------	--

Automatic	Used to automatically update the date and time received from the satellite signal when available. This option is only available when the GPS option is fitted to the analyser at the time of manufacture.
-----------	---

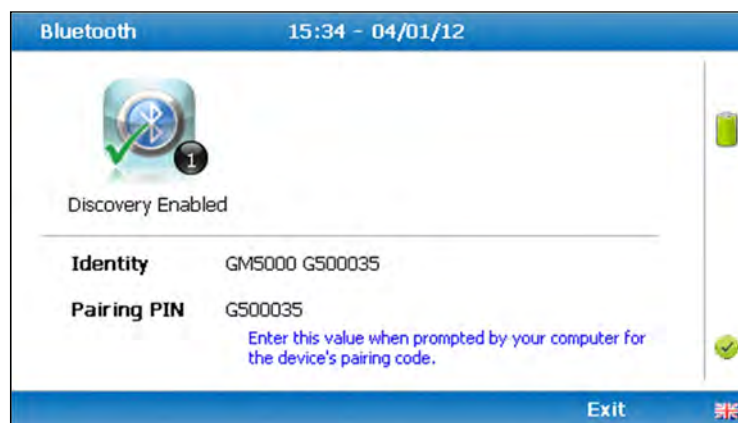
- 6) Select the soft-key 'Exit' to exit and return to the 'Device Settings' menu.

7.2.2 Bluetooth



This option enables the operator to utilise Bluetooth technology. This may be useful when downloading gas readings from the analyser to the PC instead of connecting the analyser to a PC via a USB lead.

- 1) Select the 'Menu' key on the front of the analyser to display the 'Device Settings' menu.
- 2) Select 'Key 2 – Bluetooth' and the following screen is displayed:



Bluetooth options

- 3) Enter the 'Pairing PIN' value when prompted by your computer for the device's pairing code.
- 4) Select soft-key 'Exit' to exit the screen and return to the 'Device Settings' menu.

7.2.3 Device information



This option displays default instrument information and settings such as serial number, service due date, last zero calibration date and last span calibration date.

- 1) Select the 'Menu' key on the front of the analyser to display the 'Device Settings' menu.
- 2) Select 'Key 3 – Device Information' and the following screen is displayed:

Device Information		12:23 – 17/06/14	
Serial Number	G500037	ID Count	7/2000
Version Number	1.11.9	Readings Count	142/4000
SB Version	1.19.400	Comms Mode	Serial
Last Check	---	SB Hardware	100
Last Cal	27/03/14	A5F Bias	Enabled
Manufacturer	--/--/--	A5F Status	Good
Agent Service	--/--/--		
This instrument is approved by the UK Environment Agency's Monitoring Certification Scheme, MCERTS.			
Comms Mode		Exit	

Device information

- 3) The information displayed on this screen is informational only and cannot be edited by the operator. The operator may be asked serial number, service due date and version number information when contacting Geotech.

✍ Note: The comms setting 'Legacy' mode is for use with GAM $\geq$ v1.5. GA5K mode is reserved for future applications – use with 5000 range updater tool

- 4) Select soft-key 'Exit' to exit the screen and return to the 'Device Settings' menu.

7.2.4 Diagnostics



This option enables the Geotech Technical Support Team to identify and resolve issues with the instrument and settings. If required, the operator may be asked to confirm the diagnostics displayed.

- 1) Select the 'Menu' key on the front of the analyser to display the 'Device Settings' menu.

- 2) Select 'Key 4 – Diagnostics' and the following screen is displayed:

Diagnostics		15:19 - 05/01/12				G500046	
Channel		ADC	Filt	Lin	Linz	Status	
Ref	(N/A)	10138	10134	+10134	+10134	✓	
CH ₄	(%)	-1	9944	0.2	0.2	✓	
CO ₂	(%)	5343	5343	2.6	2.6	✓	
O ₂	(%)	43511	43507	13.7	13.7	✓	
S4Cell	()					✗	
CO	(ppm)	32836	32836	0	0	✓	
H ₂	(ppm)	32833	32833	0	****	✓	
H ₂ S	(ppm)	32866	32859	3	3	✓	
PID	()					✗	
Next		Previous				Exit	

Diagnostics

- 3) Select soft-key 'Next' to display the next screen, 'Previous' to return to the previous screen, or select soft-key 'Exit' to exit this screen and return to the 'Device Settings' menu.

 Note: For further information please contact Technical Support at Geotech on +44(0)1926 338111 or email technical@geotech.co.uk.

7.2.5 Navigation (optional)



This option enables the operator to switch the 'GPS Navigation' functionality on or off. (This is optional and dependent upon purchasing the navigation option).

- 1) Select the 'Menu' key and the 'Device Settings' menu is displayed.
- 2) Select 'Key 5 – Navigation On' to switch on the GPS navigation functionality or 'Key 5 – Navigation Off' to switch the GPS navigation functionality off.



 Note: For further information please refer to section '8.0 – Taking Readings'.

7.3 User settings

To access the 'User settings' menu, select the 'Menu' key on the front of the analyser to display the 'Device Settings' menu followed by the soft-key to display 'User Settings' menu. The following menu is displayed:



User settings menu

To exit the user settings menu select the soft-key 'Exit'.

7.3.1 Operating language



This option enables the operator to specify the operating language displayed for the instrument.

- 1) Select 'Key 1 – Operating Language' and the following screen is displayed:



Set language

- 2) Set the required language for the gas analyser by selecting the appropriate function key. Choose from:

Key 1 English

Key 2	Spanish
Key 3	French
Key 4	German
Key 5	Italian
Key 6	Portuguese

Use the soft-keys to move to the next page for further language options, including simplified Chinese

- 3) To exit this option, select the soft-key 'Exit' and the operator is returned to the 'User Settings' menu.

7.3.2 Units of measurement



This option enables the operator to specify the default units of measurement for the instrument.

- 1) Select 'Key 2 – Units of Measurement' and the following screen is displayed:



Units of measurement menu

- 2) To set the required units of measurement toggle and choose from the following:

Key 1	Temperature	°C °F
Key 2	Flow	scfm m3/hr
Key 3	Measurement	Inches Millimetres
Key 4	Pressure	mb "H2O

Key 5 Balance Balance
Residual N2

Key 6 Energy BTU/h
kW

3) Select soft-key 'Exit' to exit this screen and return to the 'User Settings' menu.

7.3.3 ID selection

The ID selection screen allows the operator to scroll through all IDs, including those uploaded from GAM and added directly onto the instrument, and then make a selection. Detailed information regarding the currently selected ID, such as flow device type and pump runtime, are displayed below:-



By selecting 'Key 5' the operator can toggle between showing 36 IDs and showing 5 IDs with more detailed information relating to the chosen ID.



The technician can scroll between the IDs using the following keys on the instrument keypad:

Two (2) and eight (8) move the selection up/down
Four (4) and six (6) move the selection left/right on the list view
One (1) and three (3) move the selection left/right a page in the list view only
Seven (7) and nine (9) move the selection to first/last ID

Five (5) toggles between the 'ID with information' and 'ID list'
Return/enter key selects the desired ID and proceeds to the navigation or reading screen.

If there are no IDs present the technician can either add a new ID or press the enter key on the instrument keypad to return to the previous screen.

Soft keys:

Left - Select 'No ID' and go to the purge/reading screen.

Centre - Enabled when there is a list of IDs, allowing the technician to dynamically filter the IDs displayed in the list.

Right - Allows the technician to add a new ID to the instrument 'in the field'.

 **Note:** If your analyser has firmware version v1.12 or greater, used IDs will have a strikethrough.

Changing the sort order

By default the IDs are sorted in the order in which they were transferred to the instrument. To change the sort order between unsorted, sort by name or sorted by distance to travel press Key 0.



Sorted by original order (not sorted)



Sorted alphabetically



Sorted by distance to travel

 **Note:** Only available when GPS is enabled

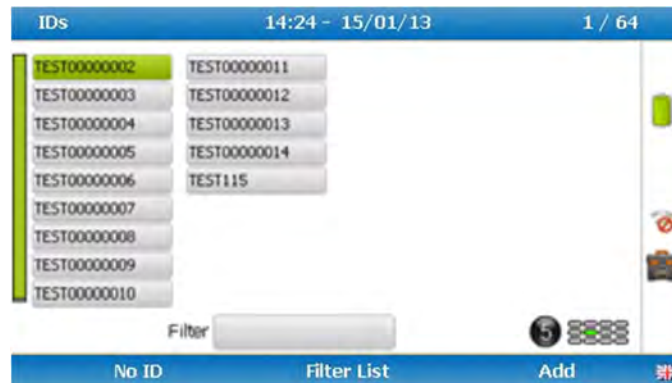
For analysers with firmware v1.12 and above:

- 1) Press the menu key
- 2) Press the middle soft key for 'User Settings'
- 3) Press key 3 – 'ID options'
 - a. Key 1 to change the sort order
 - b. Key 2 to change how the IDs are displayed
 - c. Key 3 to clear the line through on the current ID being used
 - d. Key 4 to remove the line through on all IDs

7.3.4 Routes

Using the optional Gas Analyser Manager Software (GAM) you can upload a route, this is a predefined list of IDs that you can go through in an order. This feature is useful if you have a certain order to complete your IDS in.

Furthermore if your GEM5000 has v1.12 or higher firmware it will strike through your IDs once you have used them.



 **Note:** If the analyser is set to "route mode" (v1.12 and above), the filter box will no longer be there, as you cannot filter in this mode. It will be replaced by "Route" and the route name.

7.3.5 Adjust backlight



This option enables the operator to adjust the backlight (brightness). Setting this to a darker setting will help preserve the battery power.

- 1) Select 'Key 4 – Adjust Backlight' and the following screen is displayed:



- 2) Manual adjustment of the instrument backlight is available via this option and can be carried out with use of 'Key 4 - Scroll left' (<) darker and 'Key 6 - Scroll right' (>) lighter.
- 3) Select the soft-key 'Save' to store the setting or select soft-key 'Exit' to exit the screen without saving the change.
- 4) The operator will return to the 'User settings' menu.

 **Note:** The manually set contrast setting is retained when the read-out is switched off and may require resetting when next switched on.

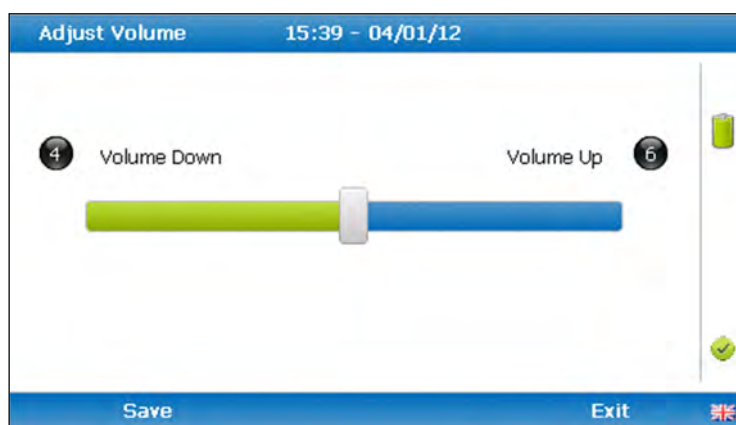
7.3.6 Adjust volume



Adjust Volume

This option enables the operator to adjust the volume for the internal speaker, for example the alarm tone. A lower setting will help preserve the battery power.

- 1) Select the 'Menu' button on the front of the analyser to display the 'Device Settings' menu. Press the soft-key to display 'User Settings'.
- 2) Select 'Key 5 – Adjust Volume' and the following screen is displayed:



Adjust volume

- 3) Manual adjustment of the volume is available via this option and can be carried out with use of 'Key 4 - Scroll left' (<) volume down and 'Key 6 - Scroll right' (>) volume up.
- 4) Select the soft-key 'Save' to store the setting or select soft-key 'Exit' to exit the screen without saving the change.
- 5) The operator will return to the 'User settings' menu.

7.3.7 User Prompts



User Prompts On

This option enables the operator to either turn on or off the context-sensitive user prompts which are displayed during the gas sample process. The analyser will have the user prompts on when it is first used, so if they are no required they can be switched off by selecting soft-key '6' and this will now be its default setting. Prompts can be switched back on at any time by returning to this menu and selecting soft-key '6'.

7.4 Operation settings

To access the 'Operation settings' menu, select the 'Menu' key on the front of the analyser to display the 'Device Settings' menu followed by the soft-key to display the 'Operation Settings' menu. The following menu is displayed:



Operation settings menu

7.4.1 Timers



Data Logging

The timers function enables the operator to set standard purge times and set auto-power off if the unit is untouched for the period of time specified.

- 1) Select 'Key 1 – Timers' and the following screen is displayed:



- 1) Select 'Key 1' to edit the purge time. Enter the 'Pump Running Time' in seconds; this is the length of time you wish to run the pump to draw the sample, e.g. key in 030 then press the 'Enter' key to accept.
- 2) Select 'Key 2' to edit the auto power off time. Enter the 'Auto power off' in minutes; the instrument will automatically power off to preserve the battery life after the specified time if no activity has occurred on the instrument. Press the 'Enter' key to accept.
- 3) Select the soft-key 'Exit' key to exit the screen and return to the 'Operation settings' menu.

-  **Note:** Setting the purge time and auto power off functions to zero, disables the option. It is not recommended to reduce the purge time to below 30 seconds.

7.4.2 Gas Check



This option displays the 'Gas Check' menu and enables the operator to zero and span the gas channels on the instrument. Historical/previous gas checks data can also be viewed and factory settings can be restored.

- 1) Select the 'Menu' button on the front of the analyser to display the 'Device Settings' menu. Press the soft key to display 'Operation Settings'.
- 2) Select 'Key 2 – Gas Check' and the following menu is displayed:



Gas Check

- 3) For more information about the Gas Check Menu please refer to section 9.0 – Calibration.
- 4) Select soft-key 'Exit' to exit operation settings and return to the main screen.

7.4.3 View data



This option enables the operator to view the readings collected and stored on the instrument. Readings may be downloaded to the optional Gas Analyser Manager (GAM) software if further analysis is required.

- 1) Select the 'Menu' button on the front of the analyser to display the 'Device Settings' menu. Press the soft-key to display 'Operation Settings'.

- 2) Select 'Key 3 – View Data' and the following screen is displayed:

View Data				10:49 - 09/01/12	3 / 3
ID: EEEE		06/01/12 11:36:37			
CH ₄	(%)	23.1	PEAKCH4	(%)	23.1
CO ₂	(%)	0.1	PEAKCO2	(%)	0.1
O ₂	(%)	17.0	MINO2	(%)	17.0
CO	(ppm)	0	SysP	(mb)	0.00
H ₂	(ppm)	----	Baro	(mb)	982
H ₂ S	(ppm)	2	Temp	(°C)	----
Bal	(%)	59.7	Anemo	(m/s)	----
			Flow	(m3/h)	----
Filter		Delete		Exit	

View data

- 3) Toggle through the reading by selecting 'Key 4 – Scroll left' and 'Key 6 – Scroll right' on the analyser. Select 'Key 2 – Page up' and 'Key 8 – Page down' to page through the auxiliary channels listed.
- 4) Select the soft-key 'Filter' to filter the data by sample point ID, or specify before or after date. Press the soft-key 'Exit' to exit the filter menu and return to the 'View Data' screen.

Filter				12:35 - 09/01/12
				
Sample Point ID	After Date	Before Date		
ANAW001	00/00/00	00/00/00		
Delete				ABC

Filter data

- 5) Select the soft-key 'Delete' followed by the appropriate soft-key to delete a single reading or all filtered readings. Press soft-key 'Cancel' to cancel the deletion request.
- 6) Select the soft-key 'Exit' to exit the view data screen.

7.4.4 Set alarms



This option enables the operator to define the conditions for which an alarm will be triggered. These conditions apply to the general operation of the instrument and are not ID specific. A summary of the alarm settings can be found in 'Key 3 – Summary'.

Types of alarms

Common Alarms – Are non-ID specific alarms which apply to all the readings taken with the analyser.

ID specific alarms – Are ID specific, i.e. they will only trigger when a certain Id is being used.

Tuning/targets – You can also set targets for your gas channels, these will highlight gas channels green as oppose to when they alarm (yellow). These can be common or ID specific.

Setting up alarms/targets

- 1) Select the 'Menu' button on the front of the analyser to display the 'Device Settings' menu. Press the soft key to display 'Operation Settings'.
- 2) Select 'Key 4 – Set Alarms' and the following menu is displayed:



Alarms Summary		12:57 - 04/05/16	
Common			
	Channel	Alarm	Low Limit High Limit
1	CH ₄ (%)	Disabled	-- --
2	CO ₂ (%)	Disabled	-- --
3	O ₂ (%)	Disabled	-- --
4	CO (ppm)	Disabled	-- --
5	H ₂ S (ppm)	Disabled	-- --
0	Disable All		
Initial		ID Alarms	Exit

Set primary alarms

- 3) Select the corresponding key to select the gas for which you wish to set an alarm/target trigger for, followed by 'Key 1' to change the trigger condition of an alarm.
- 4) To manually adjust the alarm/target set press (<) 'Key 4 – Scroll left' or 'Key 6 – Scroll right' (>) and enter the trigger value. Once you are happy, press the middle soft key for 'save'.
- 5) For pressure, temperature and flow alarms, press the left soft key for 'Secondary' and then select the corresponding key to select the channel for which you wish to set an alarm trigger for, followed by 'Key 1' to change the trigger condition of an alarm/target. Once you are happy, press the middle soft key for 'save'.
- 6) To disable all alarm settings select key 0 – 'Disable All'

 Note: ID specific alarms cannot be added/edited on the analyser, to add/edit ID specific alarms, please use the optional Gas Analyser Manager (GAM) Software.

7.4.5 Adjust flow fail



This option enables the operator to adjust the flow fail tolerance of the instrument, i.e. the operator can adjust the sensitivity for when the pump will stop operating on the presence of a blockage or low flow.

- 1) Select the 'Menu' button on the front of the analyser to display the 'Device Settings' menu. Press the soft-key to display 'Operation Settings'.
- 2) Select 'Key 5 – Adjust Flow Fail' and the following screen is displayed:



Adjust flow fail

- 3) Manual adjustment of the flow fail is available via this option and can be carried out with use of 'Key 4 – Scroll left' (<) less sensitive and 'Key 6 – Scroll right' (>) more sensitive.
- 4) Select the soft-key 'Save' to store the setting or select soft-key 'Exit' to exit the screen without saving the change.
- 5) The operator will return to the 'Operation settings' menu.

 **Note:** The default setting displays the bar in the centre. BEFORE altering this setting, please contact Technical Support at Geotech on +44(0)1926 338111 or email technical@geotech.co.uk

7.4.6 Technician login

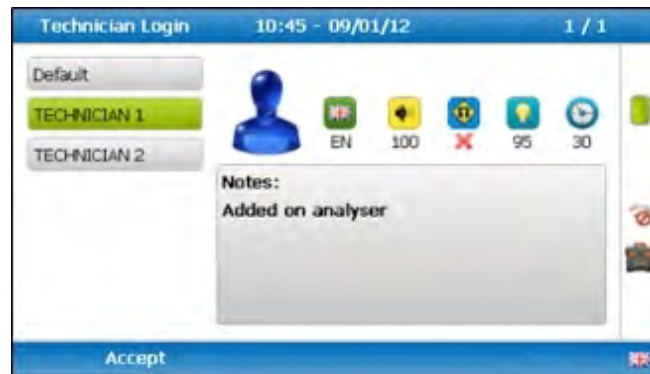


Technician Login

This option enables the operator to select or change a pre-defined technician login and all subsequent readings will be tagged with this Technician Login ID. The technician ID must already have been created using the Gas Analyser Manager (GAM) software and uploaded to the instrument.

- 1) Select the 'Menu' button on the front of the analyser to display the 'Device Settings' menu. Press the soft key to display 'User Settings'.

- 2) Select 'Key 6 – Technician login' and the following screen is displayed:



- 3) Use the cursor keys to move throughout the list of IDs displayed; select the 'Enter' key to select choice of ID, default if no IDs are listed or soft-key 'Skip' to skip the selection.
- 4) The operator will return to the 'User settings' menu.

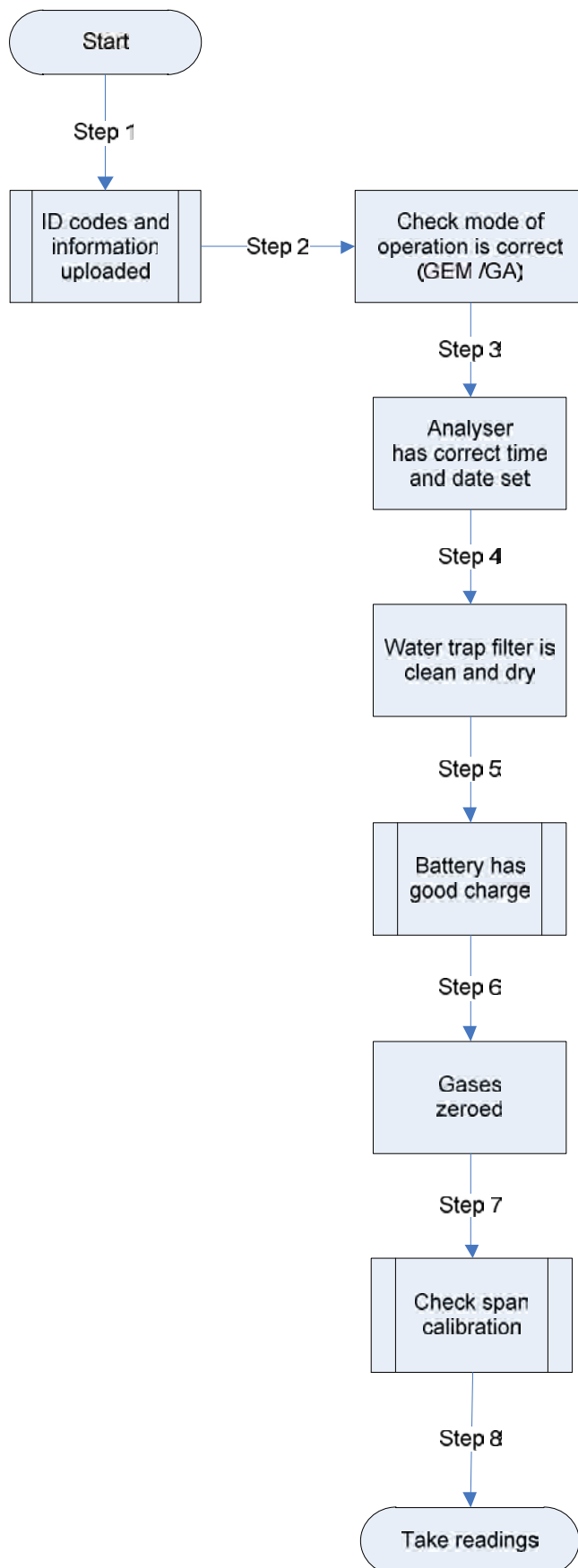
 Note: If no technicians are loaded via GAM this section is skipped during start up and the 'Technician ID' icon is removed from the menu.

7.5 Exit menu

- 1) Press the 'Menu' button on the front of the analyser to exit settings.

8.0 Taking readings

8.1 Preliminary checks before taking readings (best practice)



Prior to use, it is good practice to ensure that:

- Step 1** If using GAM - all necessary ID codes and information have been uploaded from GAM to the analyser. Please see section 8.1.1 for more information on this.
- Step 2** Check the 'Mode of Operation' is correct. Choose either GEM5000 for gas extraction monitoring analyser or GA5000 for landfill gas analyser. Change using 'Special Actions'.
- Step 3** The instrument has the correct time and date set.
- Step 4** The water trap filter is fitted and is clean and dry.
- Step 5** The battery has a good charge (minimum 25% charge, even if only a few readings are required).
- Step 6** The gas channels have been zeroed, without gas concentration present.
- Step 7** If necessary check the span calibration with a known concentration calibration gas.
- Step 8** Take readings.

**Warning**

Inhaling hydrogen sulphide gas (H₂S) or other harmful gases can cause death. It is the responsibility of the user to ensure that he/she is adequately trained in the safety aspects of using H₂S and other harmful gases. In particular, where hazardous gases are being used the gas exhausted from the analyser must be piped to an area where it is safe to discharge the gas. Hazardous gas can also be expelled from the instrument when purging with clean air.

Good practice

- Travel to site with the gas analyser in the vehicle's interior - not in the boot, where it may be subjected to extremes of temperature and possible shock damage. Do not place the gas analyser against anything hot (e.g. gas extraction pipe, car body or in an unattended car during the summer) as this will cause a temperature increase in the gas analyser and may cause erroneous readings.
- When moving around a site, protect the gas analyser from strong direct sunlight and heavy rain.
- Always use the water trap! If the water trap becomes flooded, change the filter and ensure all tubes are clear of moisture before re-use.

 **Note:** If the exhaust of a 5000 series gas analyser is connected to a pressurised system then this results in a flow of gas out of the inlet flow port.

8.1.1 Creating an ID

There are two different methods to creating an ID, either via Gas Analyser Manager or via the instrument.

If created on the analyser, you can only fill out basic information, such as ID code, description and ID type. Whereas on GAM, you can assign site and ID questions (please see below), assign flow devices, input GPS coordinates etc.

To create an ID on GAM please consult the GAM operating manual. To create an ID on the instrument:

- 1) Press the left blue arrow key for 'Next ID'
- 2) Press the right blue arrow key for 'Add'
- 3) Input an ID code using the keypad
- 4) Press enter
- 5) Using the corresponding number to input different properties of the ID
- 6) Once you are happy, press the middle key for 'add'

8.2 Change screen layout

Use key 2 to toggle between 2 different ways of viewing the gas reading screen:

***** 09:17 - 06/26/14 Baro: 997 mb				
Channel	Adjusted	Initial	Previous	Units
CH ₄		81.7	81.3	%
CO ₂		1.6	1.7	%
O ₂		16.7	17.0	%
S5Cell		0	----	ppm
Bal		0.0	0.0	%
Static.P			----	mb
Diff.P			----	mb
Temp.			----	°F
Flow			----	m3/h
Power			----	KW
Next ID		Special Action		Start

Default view

***** 15:18 - 06/17/14 Baro: 29.81 %				
Channel	Adjusted	Initial	Previous	Units
CH ₄	80.8	80.8	0.0	%
CO ₂	1.7	1.7	0.1	%
O ₂	17.5	17.5	20.6	%
Bal	0.0	0.0	79.3	%
Next ID		Special Action		Start

Alternative view

You can use key 4 and key 6 to switch between the gases page and the pressure & flow page. This feature is only available on analysers running on firmware version 1.11 or later.

8.3 Answering site questions

Prior to taking the readings at a particular site, the site questions should be populated, this is only necessary when using site questions, if not, please proceed with the reading as normal. This is accessed via the 'Special Action' menu. The answers to these questions are then stored and appended to each reading stored thereafter, until the site questions are updated for another site. You will need Gas Analyser Manager (GAM) software to create site questions and if you are using Gas Analyser Manager (GAM), this data will be uploaded to GAM along with the reading data.

8.4 Answering ID questions

Prior to, or after, you have taken the reading, the ID questions should be populated, this is only necessary when using ID questions, if not, please proceed with the reading as normal. When you have uploaded IDs with ID questions assigned to them (Gas Analyser Manager (GAM) software required), you will be prompted to answer these questions at the end of the reading, this means that the answer only equates to that reading.

8.5 Special action

This menu enables the operator to perform the additional following functions out of sequence if so desired.

- 1) From the 'Main Gas Read Screen' select the soft-key 'Special Action' and the following menu is displayed:



Special action menu

Note: The list of special action options displayed on the special action menu is dependent upon device type and sequence.

The following actions may be available:

Action	Function
Key 1 – Simple Gas	This action enables the operator to take a quick gas reading. The pump will start running automatically when this key is selected. The operator can stop the pump by pressing the pump key on the keypad at any time and the reading can then be stored by selecting soft-key 'Store'.
Key 2 – Site Questions	This action enables the operator to update site questions prior to taking a reading.
Key 3 – ID Questions	This action enables the operator to update ID questions specific to sample points prior to taking a reading.
Key 4 – Flow	This action enables the operator to measure internal flow first when taking a reading. Connect the blue hose to the sample point. The yellow hose can be vented a safe distance from the sample point or re-circulated back into the system. Select either soft-key 'Zero Flow', 'Flow Options' or 'Start' to commence internal flow. Select soft-key 'Store' to store and record the reading.
Key 5 – Enter Temperature	This action enables the operator to manually enter a temperature reading if not using a temperature probe prior to taking a gas measurement.
Key 6 – Start Logging	This action enables the operator to leave the analyser unattended to take samples at a predetermined time. The reading interval and pump run times may be edited prior to commencing the logging cycle.



8.5.1 Configuration of the data logging option



- 1) Connect the blue hose to the sample point. The yellow exhaust hose can be vented a safe distance from the sample point; do NOT re-circulate back into the system.
- 2) By selecting 'Next ID' the operator can select the ID which is being sampled at present.
- 3) Once the ID has been chosen the analyser will commence & complete its clean air purge cycle.
- 4) To gain access to the data logging option the operator will be required to select the 'Special Action' key to obtain the special user options. The data logging option can then be selected via 'Key 6' to configure the logging parameters.
- 5) Once the operator has confirmed the logging parameters, select soft-key 'Start Logging'.
- 6) Once the logging function has been activated the analyser will carry out a 30 second warm-up (displayed below the temperature read out at the right of the main gas read screen) and begin the first sample.
- 7) If for any reason during the logging cycle the inlet port becomes blocked, the analyser will sense this as a 'Flow Fail' and the pump will automatically retry until the reading can be obtained. As such care must be taken when positioning the sample tubing to ensure water/moisture ingress does not occur.
- 8) Select soft-key 'Stop Logging' to stop logging if required.

8.5.2 Profiling option

- 1) The 'Logging Mode' centre soft-key toggles between 'Logging Mode' and 'Profiling Mode' and pressing it will change the mode to the one the soft-key describes. For example, when on the profiling page the key will display as 'Logging Mode' and when on the logging page the key will display as 'Profiling Mode'.



- 2) To edit the parameter the operator will be required to select 'Key 3' to select the number of reading required. Once the number of readings has been updated press the return key to confirm parameter setting.
- 3) By selecting 'Key 2' the operator can edit the logging interval of their logging preferences and then confirm the amendments by pressing the return key.
- 4) Once the logging parameters are confirmed, commence the logging by selecting the 'Start Logging' key.
- 5) If for any reason during the logging cycle the inlet port becomes blocked, the analyser will sense this as a 'Flow Fail' and the pump will automatically retry until the reading can be obtained. As such care must be taken when positioning the sample tubing to ensure water/moisture ingress does not occur.
- 6) Select soft-key 'Stop Logging' to stop logging if required.

Select the soft-key 'Exit' to exit this menu and return to the 'Main Gas Read Screen'.

8.6 GEM analyser in GA mode

The operator may toggle between GEM (a gas extraction monitor) and GA (a landfill gas analyser) mode of operation if required. From the 'Special Action' menu, select the middle soft-key to toggle between GEM and GA Mode.



GEM analyser in GA mode - main gas read screen

Note: For operating instructions on how to use the GA5000 mode of operation, please refer to the GA5000 gas analyser operating manual.

8.7 The gas flow measurement screen

The GEM5000 gas analyser enables gas and flow measurements to be recorded by using:

- A flow device i.e. a PITOT tube, Orifice plate, Accu-Flo or anemometer.
- A temperature probe.

The GEM5000 instrument has the extra functionality to measure the calculated calorific value of the gas. The site engineer uses the information recorded by the analyser to make adjustments to the gas flow for specific boreholes. The initial and adjusted flow rates are recorded.

8.8 How to use an anemometer (optional)



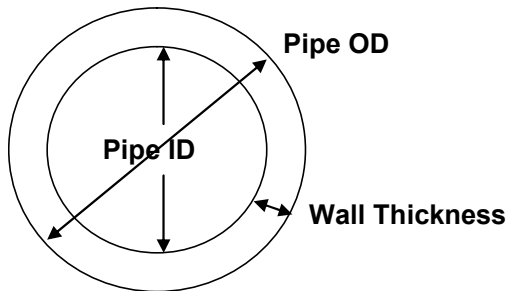
The GEM5000 gas analyser has the facility to attach an anemometer device enabling the site engineer to measure the flow of gas within an extraction system. The anemometer can be set to display two values; m/s (meters per second) and m³/hr (metres cubed per hour).

It is best practice to take the gas reading first before taking the flow reading with the anemometer attached.

If using a 'borehole ID' the internal pipe diameter can be predetermined in the optional Gas Analyser Manager (GAM) software. Once set, the site engineer cannot edit the pipe diameter setting.

If the site engineer is not using a borehole ID or the pipe diameter is not set in GAM the operator will be prompted to enter a pipe diameter with a new ID on the analyser. Select soft-key 'Next ID' from the Main Gas Read screen, followed by soft-key 'Add' and add a new borehole location.

In order to use the anemometer it is important to know the internal diameter (ID) of the pipe if you want to calculate the flow in m³/hr (metres cubed per hour). This must be the internal diameter not the outer diameter (OD) i.e. pipe outer diameter minus twice the pipe wall thickness.

For example:

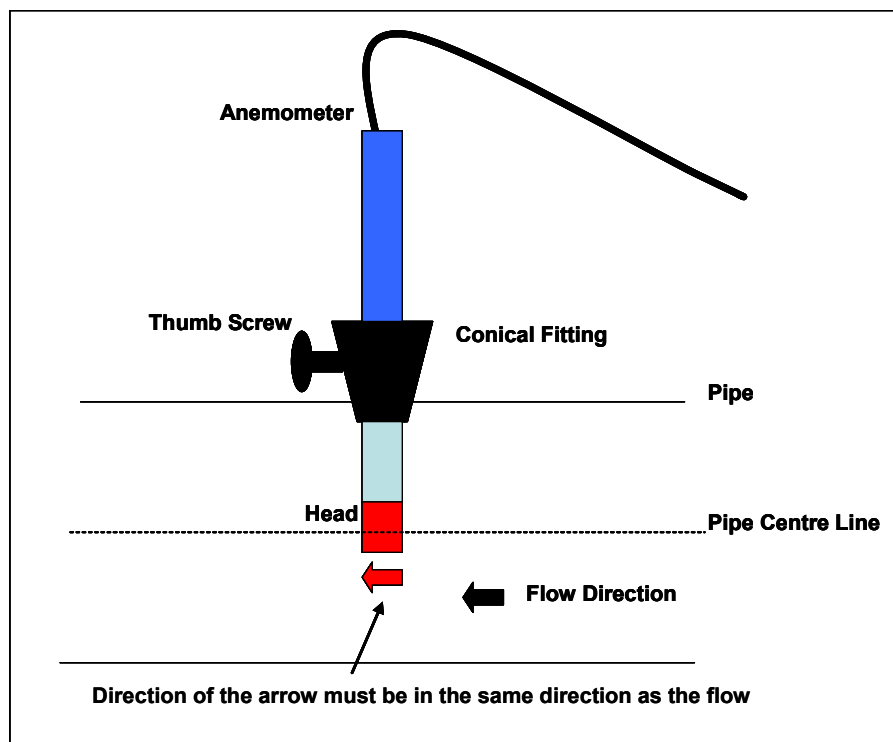
If you do not have any suitable monitoring points you will need to drill (tap of $\frac{3}{4}$ " BSP pipe thread) a hole in the piping of between 25mm and 30mm in diameter to seat the conical fitting on the anemometer (which is roughly between 20mm to 34mm). When not in use the hole can be re-sealed with a $\frac{3}{4}$ " BSP male bung.

 **Note:** When the anemometer is not in use the conical fitting should be placed over the probe to protect it.

The anemometer must fit centrally (the conical fitting must be set on the probe to half the pipe ID before insertion). The arrow on the tip of the probe must point in the direction of the gas flow.

 **Note:** Use the thumb screw to help align the direction of the probe into the gas stream.

Flow readings are most accurate when there is laminar flow (not turbulent). Turbulence can be caused by a change in pipe direction or restriction. Ideally, upstream you want at least 20 times the pipe ID along the length of the pipe without restriction or bend. Downstream, you want at least five times the pipe ID along the length of the pipe i.e. for a 100mm ID you need 2000mm of clear pipe upstream, 500mm downstream.

Example to show anemometer fitting into the sample point:

Instructions for use:

- 1) Attach the anemometer to 'Connector C' (refer to section 5.3 – Instrument connection points).
- 2) Place the anemometer into the pipe (sample point) ready to take the reading.
- 3) To take a flow measurement when using an anemometer, follow the instructions displayed on the analyser. When the reading has stabilised press the 'Enter' key to store the reading.

Anemometer cleaning instructions:**General handling tips:**

- Protect the probes against severe vibration.
- Do not kink the connector cable (risk of cable breakage).
- Never allow hard objects to contact rotating impellers.
- Always carry out probe cleaning according to the cleaning instructions.
- Never immerse probes in solvent.
- Never blow probes through with compressed air.
- Allow hot probes to cool slowly, never cool by plunging them in cold water etc.

Cleaning instructions:

- Instrument and probe must be switched off or disconnected prior to cleaning.

Vane probes:

- As the probes are highly sensitive measuring instruments, they must be cleaned with great care.
- Fibres or other foreign bodies can be carefully removed with fine tweezers. When doing so, take care not to bend or otherwise damage the vanes or the spindle.
- The adjustment of the bearing screws must never be changed. This can result in an erroneous measurement.
- Never allow hard objects to contact rotating impellers.
- Cleaning agents that extract the plasticizer from the plastic are never to be used for plastic probes (practically all solvents).

Cleaning the probes – best practice:

✍ Note: Use soapy water.

Cleaning example:

- 1) Carefully, swish the top part of the impeller back and forth in clean soapy water for approximately 10 minutes. Then swish the top part of the impeller back and forth in clean soapy water. If soapy water is used as a cleaning agent it is advisable to wash out the soap solution thoroughly with distilled water.
- 2) After cleaning the probe, rub it dry with a clean, dry cloth.

8.9 How to use a pitot tube (optional)

The GEM5000 gas analyser enables the site engineer to take gas measurements using a pitot tube. The pitot tube is used for gas extraction systems and the pressure readings are taken in mbar. Gas flow is calculated in the analyser in m³/hr (metres cubed per hour).



Example of a pitot tube

Fitting a pitot tube to the analyser:

- 1) It is important to seat the pitot tube into the monitoring point with the tip facing into the gas flow. The pitot tube should also be housed half way down the monitoring pipe. (Please refer to the anemometer instructions, which detail how to calculate the pipe diameter correctly).
- 2) Make sure that the sample tubing attached to the pitot tube fits correctly.
- 3) Attach the sample tube from the top connection of the pitot tube to the 'blue port' (the differential port) on the analyser making sure that the gas port connector secure into place.
- 4) Attach the sample tube from the side of the pitot tube to the 'white port' (inlet port/static pressure port) on the analyser making sure that the gas port connector secures into place.
- 5) House the pitot gland correctly onto the monitoring point.
- 6) When taking a gas reading and flow measurement follow the instruction as displayed on the front of the instrument.

 **Note:** Please refer to the following section for a diagram showing how to fit the tubing to the pitot tube.

8.10 How to use an H₂S filter (optional)**8.8.1 Cross gas effects on chemical cells**

Measurements of CO are important in landfill management. The GEM5000 analyser incorporates an improved CO measurement.

Measurements of CO can be affected by two other gases that can be found in landfill gas – hydrogen and hydrogen sulphide.

To reduce the effect of hydrogen, the GEM5000 analyser uses a technique that is hydrogen compensated. Hydrogen compensation is achievable up to a level of around 2000ppm. Above this level the CO reading will not be compensated for.

In order to assist the operator the GEM5000 instrument also indicates the level of hydrogen present as low, medium or high. If a high hydrogen reading is present then the CO reading may be affected.

The effect of hydrogen sulphide is eliminated by the use of a H₂S filter.

CO measurement

The CO measurement is sensitive to hydrogen sulphide. The presence of hydrogen sulphide can cause the CO reading to elevate (not to be the true value due to the interfering gas). If the presence of hydrogen sulphide is suspected to be causing false CO readings, then it is recommended that the external hydrogen sulphide filter is used whilst obtaining the CO measurement.

The H₂S filter only needs to be used when you are trying to remove of any possible cross gas effects H₂S might have on other gas channels. Do not use the filter on all boreholes.

When using the H₂S filter you will need to increase the gas sample and clean air purge run-time, as using an H₂S filter decreases the response time of the analyser.

 **Note:** There is an internal H₂S filter incorporated in the chemical cell that removes the H₂S; however this has a limited life span.

The electrochemical cells used to measure H₂S and CO do suffer from cross-gas effects. Such effects are not accurately specified. However, the following table may be useful as a guide; it represents how many ppm would be read by a cell if 100 ppm of the interfering gas were applied, with no other cross-contaminates being present in the sample.

GEM5000

		Interfering Gas				
		CO	H ₂ S	SO ₂	NO ₂	H ₂
Chemical Cell	CO (H ₂ compensated)	100	0 / ~300*	0	0	<1
	CO	100	<4 / ~300*	0	-20 to +5	<60
	H ₂ S	<4	100	20	<-25	<0.2

*after internal filter depleted.

 **Note:** Other gases could cause cross-gas effects. If you suspect a cross sensitivity problem please contact the Technical Support Team at Geotech on +44(0)1926 338111.

8.9.2 Cross-gas effects on methane, carbon dioxide and oxygen

Methane is measured using dual beam infrared absorption. Analysers are calibrated using certified methane mixtures and will give correct readings provided there are no other hydrocarbon gases present within the sample (e.g. ethane, propane, butane, etc.). If there are other hydrocarbons present, the methane reading will be elevated (never lower) than the actual methane concentration being monitored.

The extent to which the methane reading is affected depends upon the concentration of the methane in the sample and the concentration of the other hydrocarbons. The effect is totally non-linear and difficult to predict.

Note: The effect can be reduced by using an H₂S filter as it can reduce higher order hydrocarbons. When using the H₂S filter you will need to increase the gas sample and clean air purge run-time, as using an H₂S filter increases the response time of the analyser.

Carbon dioxide is measured by infrared absorption at a wavelength specific to carbon dioxide. Therefore, the carbon dioxide reading will not be affected by any other gases usually found on landfill sites.

The oxygen sensor is a galvanic cell type and suffers virtually no influence from CO₂, CO, H₂S, NO₂, SO₂ or H₂, unlike many other types of oxygen cell.

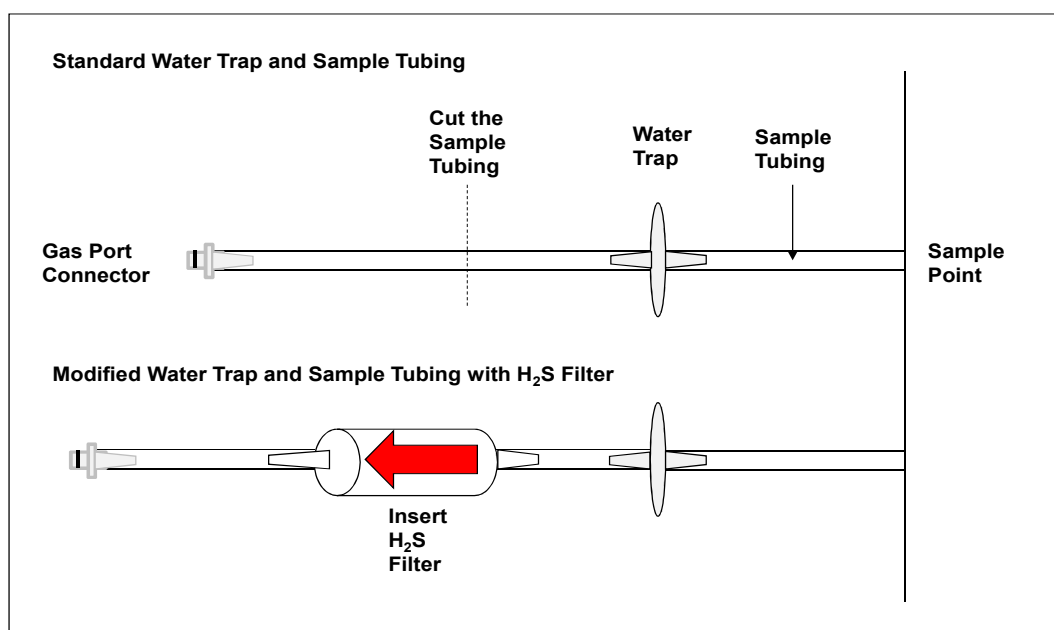
The infrared sensors will not be 'poisoned' by other hydrocarbons and will revert to normal operation as soon as the gas sample has been 'purged'.

H₂S filter:



Instructions for use:

The following diagram shows how to modify the standard water trap and sample tubing to fit the H₂S filter.

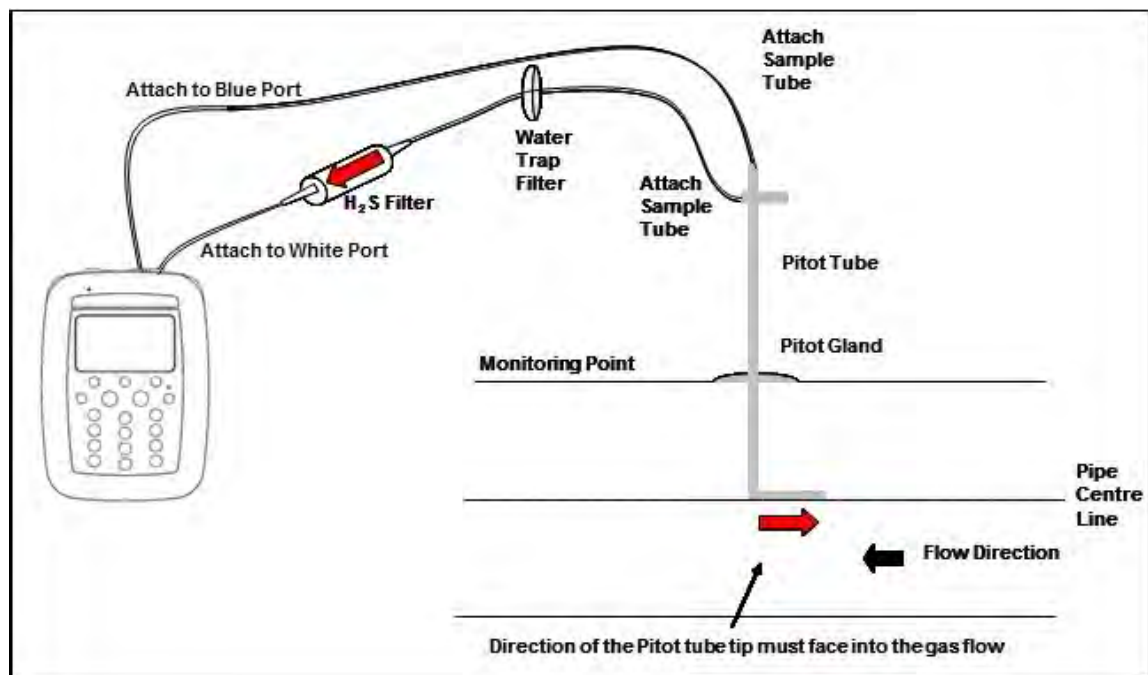


✍ Note: When onsite the site engineer must have an unmodified water trap assembly in addition to the modified water trap with a H₂S filter in order to take readings with and without a filter.

- 1) Once the H₂S filter is fitted follow the instructions as detailed on the front of the gas analyser displayed when taking readings using a H₂S filter.

✍ Note: This is only for analysers fitted with the CO compensated cell.

Fitting a pitot tube to a sample point when using an H₂S filter:



- 2) Make sure the water trap filter is fitted as close as possible to the pitot tube. This will help protect and trap moisture before it reaches the H₂S filter. Make sure that the water trap filter is clean and dry.
- 3) Attach the sample tubing into the analyser.
- 4) To take flow readings using a pitot tube, please follow the instructions displayed on the front of the instrument. The user will be prompted to complete each step and should follow the instructions on screen.

✍ Note: It is important to fit the pitot tube central and parallel with the pipe.

8.11 How to use a temperature probe (optional)

The temperature probe enables the site engineer to measure the temperature of the gas within a sample point. The GEM5000 gas analyser uses the temperature of the gas to give more accurate flow measurement readings as part of the instrument calculation.



GEM5000 & temperature probe

Instructions for use:

- 1) The temperature probe reading is taken along with the gas measurement reading.
- 2) The analyser must be at the 'Main Gas Read Screen'.
- 3) Attach the temperature probe to 'connector B' (refer to section 5.3 – Instrument connections points).
- 4) Insert the temperature probe into the sample point (borehole) at the same time as you attach the sample tube to the sample point (two sample points are required for the borehole).
- 5) Follow the instructions on the front of the instrument when taking your gas and measurement reading.
- 6) At the point in which the operator presses 'Enter' to store the gas reading the temperature is recorded.

 **Note:** Temperature probe readings can be analysed further when downloaded to Gas Analyser Manager (GAM).

8.12 How to identify a borehole using the GPS feature (optional)

An optional GPS feature is available for the GEM5000 gas analyser. It enables the site engineer to automatically locate borehole IDs using GPS satellite signal from predefined borehole IDs uploaded from Gas Analyser Manager. The GPS reading data is stored for each measurement reading.

 **Note:** Borehole IDs may be uploaded from GAM with or without location

information. If location details are not uploaded the location longitude and latitude coordinates can be stored when the borehole is located and downloaded to GAM with the reading measurements.

Screen navigation:

- 1) Switch on the analyser and wait for the self-test warm-up to complete and the analyser will display the 'Main Gas Read Screen'.
- 2) In order to use the navigation function if configured, you must switch 'Navigation - On' on the analyser. Select the 'Menu' key followed by 'Key 5' to toggle navigation to on. Select the 'Menu' key to exit and return to the 'Main Gas Read Screen'.
- 3) Select the soft-key 'Next ID', then select a borehole ID from the list displayed and press the 'Enter' key to continue.

Warning

Before entering the GPS Navigation Screen for the first time the following health and safety message will be displayed.

"Please be aware of the terrain when using this screen. You are responsible for your own safety whilst walking on-site!"

- 4) After reading the user warning message, select the soft-key 'Dismiss'. Use the tracking display to locate the borehole

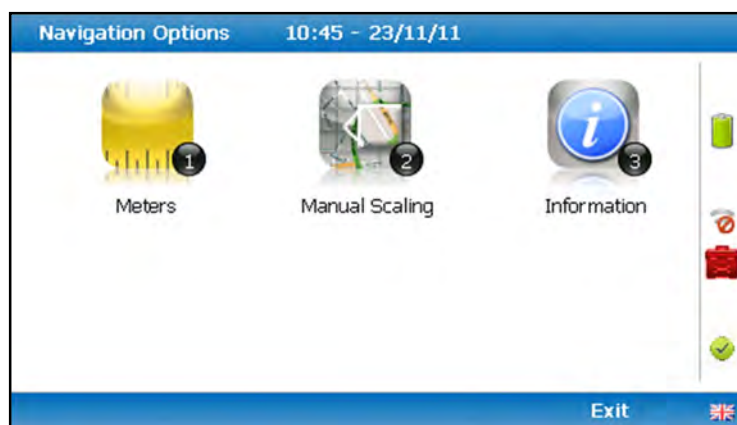
 **Note:** There is often a wait time frame of between 30 seconds to two minutes whilst getting a satellite signal. Be aware that heavy rain, trees overhead etc. will give a bad fix.

- 5) Once the operator selects a borehole ID the 'GPS Navigation' screen is displayed.



Navigation

- 6) If required select soft-key 'Options' to go to the 'Navigation Options' menu and the following screen is displayed:



Navigation options

Key 1	Key 1 to toggle between 'Meters', 'Feet' or 'Yards' to define the unit of measurement.
Key 2	Key 2 to toggle between 'Manual Scaling' and 'Automatic Scaling' to locate a sample point using a satellite signal.
Key 3	Key 3 to display sample point information:
Bearing:	Direction.
UTC time:	'Universal Time Code' received from the satellite and displays GMT.
Latitude:	Latitude displays as degrees, minutes, seconds and decimal seconds. The equator is 0.
Longitude:	E (East) or W (West) displays the longitude as degrees, minutes, seconds and decimal seconds. The Greenwich meridian defines the zero point.
Altitude:	Altitude displays in meters, feet or yards.
Satellite:	This is the number of satellites that the system can select. Four satellites will give reasonable position accuracy; eight or more satellites will increase accuracy.
Horizontal:	Estimate of horizontal error on the indicated position. The location is accurate to horizontal error 1.6m.
Vertical:	Estimate of vertical error on the indicated position.
HDOP:	'Horizontal Dilution of Precision' which measures the accuracy of the indicated position.

Changing the selected ID

It is possible to change the currently selected ID (identified by green icon ) by pressing the

2 and 8 and keys to move the cursor up and down the list of visible IDs. To change the selection press Enter. Notice that the selected ID moves to the top of the list and its icon turns green.

Changing the scale

You can change the display's scale by using the 4 and 6 keys to zoom in and out.

GPS signal strength:



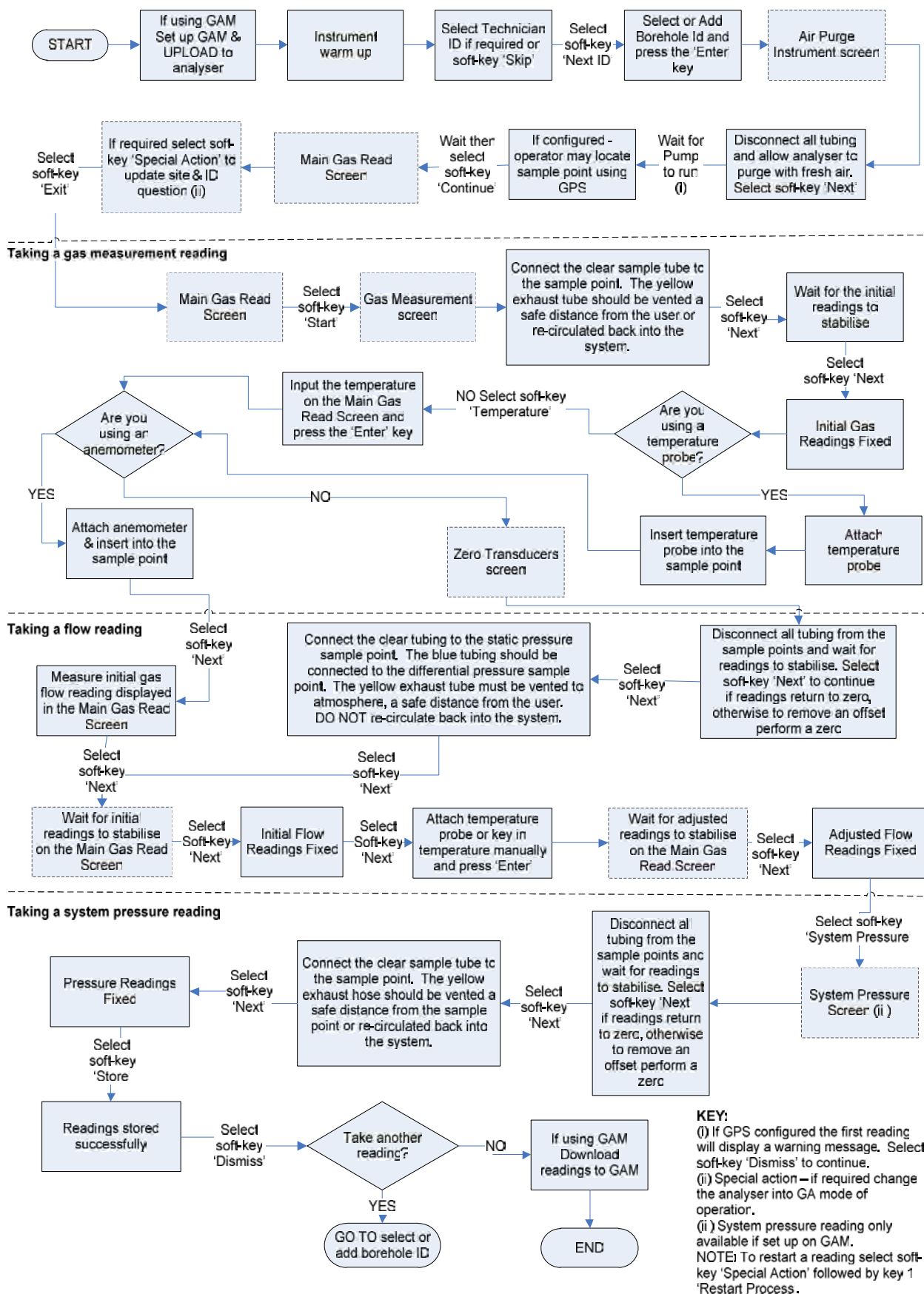
This icon shows the signal strength the analyser's GPS module is able to provide. Full, okay and fair strength respectively.



GPS failure - the GPS was unable to get a line of sight lock on enough satellites. Or, it may be that it hasn't had time to get a lock.

- 7) Select soft-key 'Continue' and the operator is returned to the 'Main Gas Read Screen'.

8.13 Taking gas and flow measurement



9.0 Calibration

9.1 Calibration introduction

The GEM5000 gas analyser is carefully calibrated at manufacture and when returned for service. However, it is sometimes desirable to be able to carry out a calibration process between services.

This section outlines the correct procedures to enable the site engineer to field calibrate the gas analyser.

 **Note:** This does not replace the factory service and calibration. If this calibration is completed incorrectly it may decrease the accuracy of the gas analyser.

CH₄, CO₂ and O₂ can be measured by GEM5000 gas analyser as standard; these channels can be user calibrated. The analysers have other gas channel options that are specified at manufacture; these too can be calibrated. This section will describe in detail how to calibrate the three standard gas channels plus the CO channel.

The GEM5000 instrument can have a H₂ compensated CO channel. This option requires that H₂ is used in the calibration process and is also set out within this section.

For the other gas channel options contact Geotech for advice.

Two important terms that are used within this section are 'Zero' and 'Span'.

Zero: The point at which the gas analyser is calibrated when there is none of the target gases present.

Span: The point at which the gas analyser is calibrated when a known quantity of the target gas is present.

9.2 Frequency of calibration – best practice

The GEM5000 gas analysers can be checked against a known concentration of gas, to give confidence that the analyser is operating as expected at the time and conditions in which it is being used.

It is recommended that the instrument is regularly serviced and calibrated by Geotech in accordance with the due date on the instrument.

When defining the frequency of user calibration, the following are factors to be considered:

- The frequency of use of the analyser. (daily?/monthly?)
- The level of confidence and accuracy required for readings to be taken.
- Historical user calibration data.
- Site specific requirements or conditions.
- Historical understanding of expected readings on site.

Zeroing of the gas analyser should be undertaken at the start of each day's monitoring.

Use historical data to drive your frequency of calibration.

If there is no historical data a good starting point for a daily monitoring round is performing a calibration once every week or every other week.

The results of the calibrations will need to be recorded to monitor over time whether the frequency of calibration needs to be increased or decreased relative to the confidence

required.

The confidence required will be driven by the site specific / user requirements.

When undertaking the monitoring with an understanding of the history of the gas levels of that site, a calibration check could be triggered if the readings measured are different to what is expected.

 **Note:** For assistance please contact Technical Support at Geotech on +44(0)1926 338111

9.3 Calibration gases

User calibration of a gas analyser will greatly improve the data accuracy in the range of the calibration gases used. This may cause less accurate readings of concentrations outside this calibrated range. Users should select the correct calibration gas for the expected gas levels on their particular application.

- To improve calibration at lower levels requires the use of gas mixtures 1 and 2.
- To improve higher levels use gas mixture 3.
- For standard CO only 100ppm CO gas is needed.
- For CO (H₂ compensated) both CO 100ppm and H₂ 1000ppm gases are needed.

The following table indicates the different gas mixture canisters used for calibration:

Calibration gas	CH ₄	CO ₂	O ₂
Mixture 1	5%	5%	6%
Mixture 2	5%	10%	0%
Mixture 3	60%	40%	0%

Calibration targets for gas cells are dependent on the gas/range and type of cell fitted. Contact Technical Support for assistance.

These are for general use but other gas concentrations can be used.

 **Note:** The above gases and most other gas concentrations can be supplied by Geotech. For further information please contact Sales on +44(0)1926 338111 or email sales@geotech.co.uk.

 Warning	Calibration gases can be dangerous. For each gas used the appropriate material safety data sheet must be read and understood before proceeding.
--	---

9.4 Calibration set-up**⚠ Warning**

Do NOT attach the gas supply to the gas analyser before putting the analyser into the 'Gas Check' screen. Select 'Check Spans' from the 'Operation Settings' menu.

The regulator supplied with the calibration kit has been configured to deliver a fixed flow.

As the regulator's flow is factory set, it only requires a few turns to open, no adjustment is necessary.

⚠ Warning**Exhaust port**

When the gas analyser is being calibrated, there are two possible exits for the gas, via the usual manner out of the exhaust (yellow) port of the analyser or in cases of over-pressurisation the 1/16" port on the red pressure relief valve located on the regulator.

It is recommended that both ports have exhaust tubing attached.

The exhaust tubing must emerge in a well-ventilated area. Ensure there are no leaks in the tubing and connections.

The calibration of the gas analyser should be carried out in a safe area with all necessary precautions taken when using potentially dangerous, explosive or toxic gases.

✍ Note: There is also potential for gas to expel from the internal flow (blue) port of the gas analyser (applies to the GA5000 only).

9.5 Calibration equipment

The diagram below displays the regulator and tubing equipment for user calibration:



- Certified calibration gas in 110 litre gas canisters are supplied with the Geotech calibration kit. Please refer to the Geotech website www.geotech.co.uk for further information.
- The regulator supplied with the calibration kit is pre-set for flow and pressure rates that are factory set.

9.6 Gas analyser

For the GEM5000 gas analyser the calibration options can be found by selecting the 'Menu' key followed by soft-key 'Operation Settings'. Select 'Key 2 – Gas Check' then follow the instructions on the analyser screen by selecting 'Key 2 – Gas Check'.

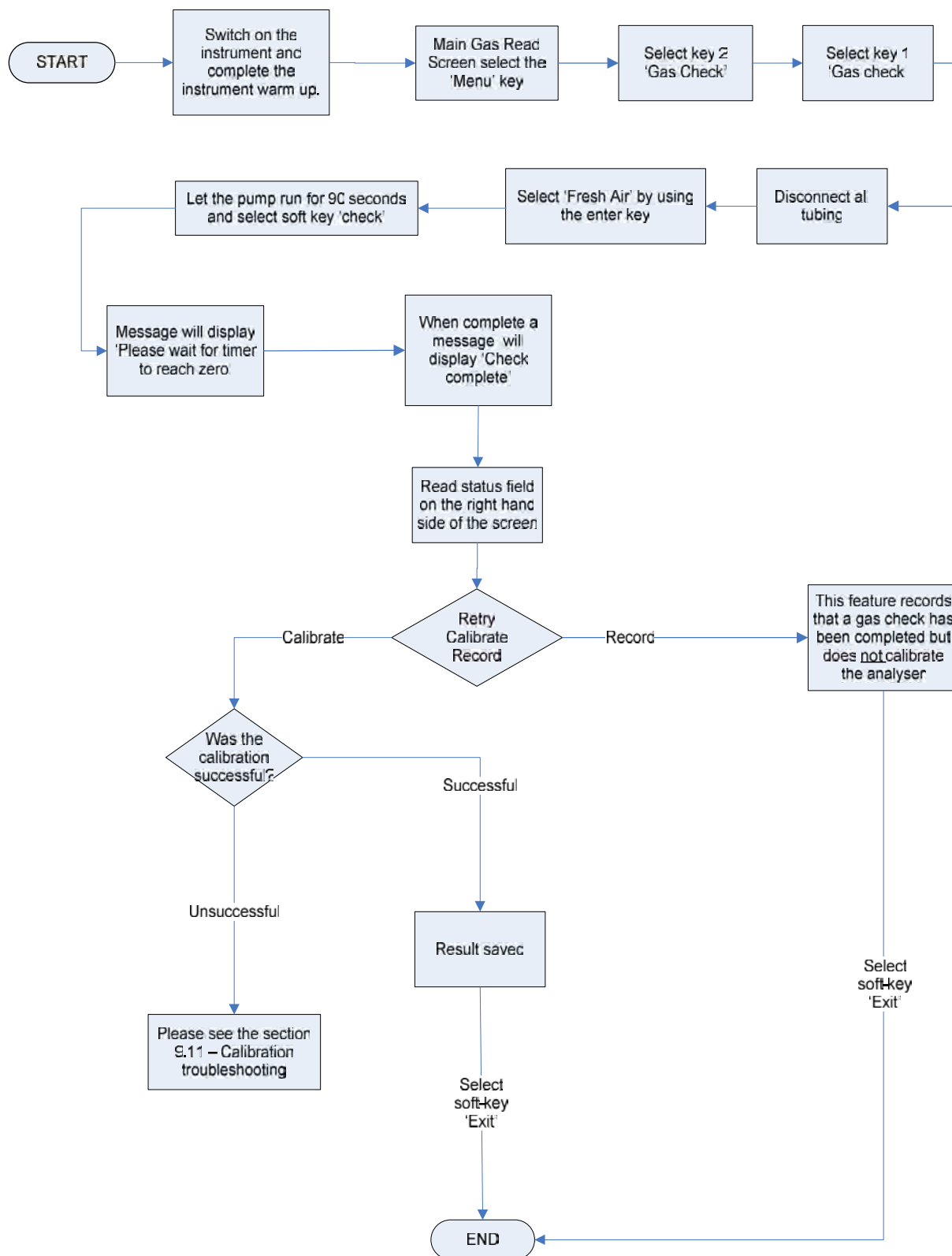


9.7 Calibration processes – best practice

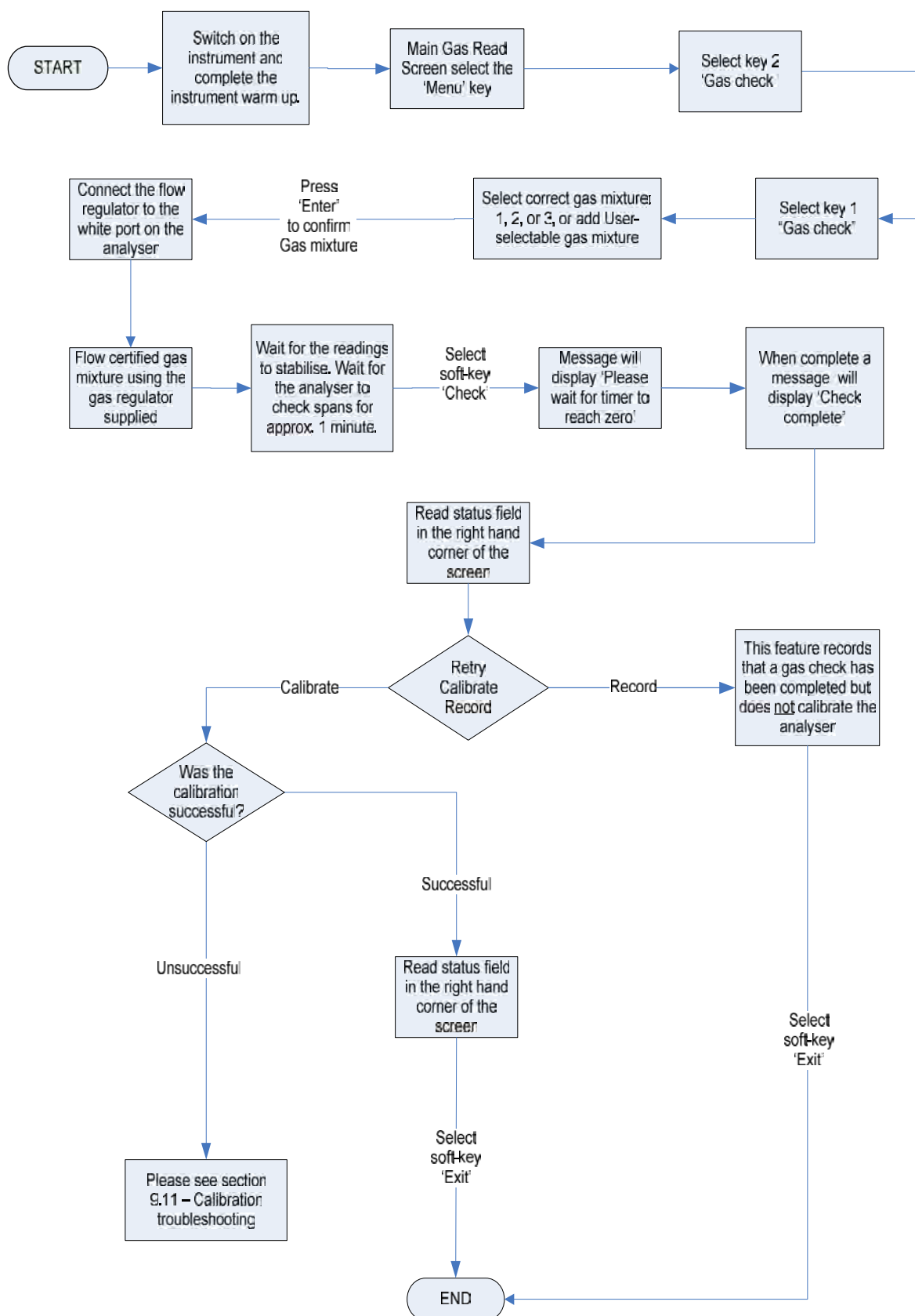
The following process diagrams outline the calibration steps.



9.7.1 Gas Check in fresh air



9.7.2 Calibration – mixtures 1, 2, & 3



9.8 Restore to factory



This option will reset the gas analyser to all of its factory programmed calibration settings and will clear ALL the user defined calibration points. It will not affect or remove ID's or readings from the analyser.

- 1) Select 'Key 2 - Restore to factory' followed by the soft-key 'Confirm' or 'Cancel'.
- 2) A validation message is displayed 'Reset user calibration?' Press the soft-key 'Confirm' to continue with the factory settings or soft-key 'Cancel' to cancel the operation and return to the Gas Check menu.

9.9 Calibration history

The GEM5000 gas analyser logs user calibrations in 'History' application. This can be used as an aid to ensuring that gas measurements are valid and accurate. Both good and failed calibration results are recorded for each channel calibrated.

- 1) Select 'Key 3 – History'.
- 2) The operator may view the calibration data stored. Use the soft-key 'Filter' to add a sort filter to the history enquiry.

9.10 Calibration summary

The GEM5000 gas analyser has the facility to log the history of user calibrations.

- 1) Select 'Key 4 – Summary'.
- 2) The operator may view the calibration data history stored by ID, technician, timestamp, type and calibration result. Use the soft-key 'Exit' to exit and return to the 'Gas Check' menu.

10.0 Problem solving

This section outlines various warning and error messages which the operator may receive during general operation of the instrument. For further assistance please contact Technical Support at Geotech on +44(0)1926 338111 or email technical@geotech.co.uk.

11.0 Warning and error display

When switched on the instrument will perform a pre-determined self-test sequence taking approximately 15 seconds. During this time many of the instrument's working parameters and settings are checked.

If any operational parameters are out of specification or the pre-programmed recommended calibration/service date has passed, errors or warnings may be displayed.

Use the 'Scroll up' and 'Scroll down' keys to move through the list if required.

Only three warnings/errors can be displayed at any time.

To ascertain if more errors have occurred use 'Key 8' – Scroll down' and 'Key 2' - Scroll up' through the list.

Warnings displayed:

All warnings displayed will be prefixed by the word WARNING followed by a relevant description.

There are two types of warning that may be displayed:

1. General warnings that may not affect the instrument's function and those where the self-test has detected a function that is outside the usual programmed operating criteria, e.g. battery charge low, memory nearly full.
2. Operational parameters that could affect the performance of the analyser: Cell out of calibration, CH₄ out of calibration, CO₂ out of calibration.

The most likely reason for the errors is either an incorrect user calibration or sensor failure. If an incorrect user calibration has caused the warning it should be correctable by way of returning the instrument to factory settings, zeroing or carrying out a user calibration as necessary for the relevant function.

12.0 Service

The GEM5000 gas analyser should be regularly serviced to ensure correct and accurate operation. Geotech recommends a service and recalibration every **6 months**.

The GEM5000 range is ATEX certified for use in potentially explosive areas. As such it should be serviced only by qualified engineers. Failure to observe this will result in the warranty becoming invalid and could invalidate the ATEX certification.

 Warning	If the GEM5000 is serviced by unqualified engineers the ATEX certification may be invalidated and the instrument may be unsafe for use in a potentially explosive atmosphere.
--	---

User serviceable parts:

There are no user serviceable parts inside the instrument.

The following parts can be user serviced:

In-line water filter	This should be regularly inspected for obstructions, moisture or damage and changed if needed. The instrument should never be operated without the in-line water filter as this may result in water entering the instrument.
Sample tubing	Always ensure that sample tubes are not contaminated or damaged.
Gas port connectors	Periodically check that the O-rings on the gas port connectors are not damaged. A damaged O-ring can let air into the sample gas and result in incorrect readings. If the O-ring is damaged the complete gas port connector should be replaced.
H ₂ S filter material	When the filter material changes colour to a <i>light grey</i> colour the filter should be replaced.

13.0 Warranty policy

This instrument is guaranteed, to the original end user purchaser, against defect in materials and workmanship for a period of **3 years** from the date of the shipment to the user.

During this period Geotech will repair or replace defective parts on an exchange basis.

The decision to repair or replace will be determined by Geotech.

To maintain this warranty, the purchaser must perform maintenance and calibration as prescribed in the operating manual.

Normal wear and tear, and parts damaged by abuse, misuse, negligence or accidents are specifically excluded from the warranty.

 Note: Please contact Technical Support at Geotech for further information.

14.0 Important notice to all customers**WEEE COMPLIANT**

The wheelie bin symbol now displayed on equipment supplied by Geotechnical Instruments (UK) Limited signifies that the apparatus must not be disposed of through the normal municipal waste stream but through a registered recycling scheme.

The Waste Electrical and Electronic Equipment directive (WEEE) makes producers responsible from July 1st 2007 in meeting their obligations, with the fundamental aim of reducing the environmental impact of electrical and electronic equipment at the end of its life.

Geotech is now registered with the Environmental Agency as a producer and has joined a recycling scheme provider who will manage and report on our electrical waste on our behalf.

Our Producer Registration Number is WEEE/NYN680

When your instrument is at the end of its life, please contact the Sales team at Geotech who will advise you on the next step in order to help us meet our obligations.

15.0 Appendices – safety instructions

15.1 Instructions for safe use – Italian language

Istruzioni per la sicurezza

 Avvertenza	Gli analizzatori di gas serie 5000 possono essere utilizzati per misurare i gas provenienti da scariche o da altre fonti, come descritto in questo manuale. L'operatore può essere esposto a gas nocivi durante l'utilizzo dello strumento. L'inalazione di questi gas può danneggiare la salute e in alcuni casi essere letale. Spetta all'utente controllare di essere sufficientemente informato sugli aspetti riguardanti la sicurezza dei gas utilizzati e di seguire le procedure appropriate. In particolare nel caso di gas pericolosi, quelli scaricati dall'analizzatore devono essere convogliati in un'area in cui lo scaricatore operazione possa essere effettuata in condizioni di sicurezza. È possibile che lo strumento scarichi gas pericolosi anche durante lo spurgo quando viene utilizzato per spurgare con aria pulita.
---	---

✎ **Nota:** gli analizzatori di gas sono apparecchi scientifici delicati e vanno trattati come tali. Utilizzando l'apparecchio in modo diverso da quanto specificato dalla casa produttrice, l'apparecchio stesso potrebbe non fornire più la protezione prevista.

Gli analizzatori di gas serie 5000 sono conformi alla Parte 15 delle norme FCC. L'utilizzo è soggetto alle seguenti due condizioni:

- 1) il dispositivo non deve causare interferenze dannose
- 2) il dispositivo deve accettare le interferenze che riceve, anche se possono causare effetti indesiderati per il suo funzionamento.

Gli analizzatori di gas della serie 5000 sono certificati con riferimento alla classificazione delle aree pericolose secondo ATEX e IECEx.



II 2G Ex ib IIA T1 Gb (Ta = da -10°C a +50°C)

È essenziale seguire scrupolosamente le istruzioni. Spetta all'operatore definire il concetto di protezione e la classificazione richiesta per una data applicazione.

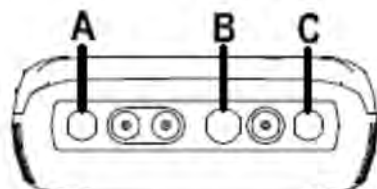
Istruzioni per l'uso in sicurezza - Italiano

(Riferimento alla Direttiva europea ATEX 94/9/CEC, Allegato II, 1.0.6.)

Le seguenti istruzioni si applicano agli apparecchi coperti dai certificati SIRA 11ATEX2197X e SIR 11.0089X della Direttiva IECEx:

- Gli apparecchi possono essere utilizzati con gas e vapori infiammabili di gruppo IIA e temperature di classe T1.
- Gli apparecchi sono certificati solo per l'uso a temperatura ambiente compresa tra -10°C e +50°C e non vanno utilizzati al di fuori di questo intervallo.

- Gli apparecchi non vanno utilizzati in un'atmosfera che contenga più del 21% di ossigeno.
- Le riparazioni di questi apparecchi vanno effettuate in conformità al codice professionale rilevante.
- In aree pericolose, utilizzare solo la sonda di temperatura GF5.2 (SIRA 11ATEX2197X e IECEx SIR11.0089X). Per il connettore C, l'anemometro GF5.4 (BVS 04ATEXE194) da usare esclusivamente con apparecchi ATEX. Nella zona pericolosa, l'analizzatore non va collegato ad altri dispositivi in dotazione, come il cavo GF-USB (per il connettore A) o il caricabatteria GF3.9 (per il connettore B).



**Non caricare, ricaricare o aprire in un'atmosfera potenzialmente esplosiva.
In aree pericolose, utilizzare solo la "sonda di temperatura GF5.2" con il connettore B.**

**Connettore C ($U_o=10V, I_o=5mA, P_o=50mW, C_i=0, L_i=0, C_o=100\mu F, L_o=1000mH$),
Connettore B ($U_o=5V, I_o=6mA, P_o=7mW, C_i=0, L_i=0, C_o=100\mu F, L_o=1000mH$)**

**ALIMENTAZIONI MASSIME NON PERICOLOSE:
Connettore A - $U_m=6V$ Connettore B - $U_m=10,1V$**

- Se esiste la possibilità che l'apparecchio potrebbe venire in contatto con sostanze aggressive, ad esempio liquidi acidi o gas che possono attaccare i metalli o solventi che possono agire su materiali polimerici, spetta all'utente adottare le precauzioni necessarie, ad es. controlli regolari come da programma o verifica della resistenza a sostanze chimiche specifiche consultando la scheda tecnica, per evitare effetti negativi e non compromettere il tipo di protezione di cui è dotato l'apparecchio.
- L'intervallo di pressione relativa è di ± 500 mbar. Si noti tuttavia che la pressione in entrata non deve superare ± 500 mbar relativamente alla pressione atmosferica e la pressione in uscita non deve superare ± 100 mbar relativamente alla pressione atmosferica.

Gli analizzatori di gas della serie 5000 sono certificati con riferimento alla classificazione delle aree pericolose secondo CSA (Canada)

CLASSE 2258 03 - APPARECCHIATURE DI CONTROLLO DEI PROCESSI - Sistemi intrinsecamente sicuri e ignifughi - Per luoghi pericolosi.



Ex ib IIA

Rilevatori di metano Modello GA 5000, GEM 5000 e BIOGAS 5000; portatili, a batteria, con pacco batterie non sostituibile sul campo, cod. parte 20087 o 201113; intrinsecamente sicuri, per circuiti intrinsecamente sicuri ("ib") per Zona 1) fino alla sonda di temperatura Modello GF5.2 (Connettore B) e con parametri di entità in uscita come indicato nella tabella sottostante; Codice temperatura T1; $-10\text{ }^{\circ}\text{C} \leq T_{amb.} \leq +50^{\circ}\text{C}$.

Connettore	Parametri entità
------------	------------------

	Uo (V)	Io (mA)	Po (mW)	Co (uF)	Lo (mH)	Ci (uF)	Li (mH)
B	5.0	6	7	100	1000	0	0
C	10,0	5	50	100	1000	0	0

✎ Nota: questo dispositivo è stato controllato solo per quanto riguarda le caratteristiche di sicurezza elettrica.

Gli analizzatori di gas della serie 5000 sono certificati con riferimento alla classificazione delle aree pericolose secondo CSA (USA)

CLASSE 2258 83 - APPARECCHIATURE DI CONTROLLO DEI PROCESSI - Sistemi intrinsecamente sicuri e ignifughi - Per luoghi pericolosi - CERTIFICATE PER USA NORME



AEx ib IIA

Rilevatori di metano Modello GA 5000, GEM 5000 e BIOGAS 5000; portatili, a batteria, con pacco batterie non sostituibile sul campo, cod. parte 20087 o 2011113; intrinsecamente sicuri, per circuiti intrinsecamente sicuri ("ib") per Zona 1) fino alla sonda di temperatura Modello GF5.2 (Connettore B) e con parametri di entità in uscita come indicato nella tabella sottostante; Codice temperatura T1; $-10^{\circ}\text{C} \leq T_{\text{amb.}} \leq +50^{\circ}\text{C}$.

Connettore	Parametri entità						
	Uo (V)	Io (mA)	Po (mW)	Co (uF)	Lo (mH)	Ci (uF)	Li (mH)
B	5,0	6	7	100	1000	0	0
C	10,0	5	50	100	1000	0	0

✎ Nota: questo dispositivo è stato controllato solo per quanto riguarda le caratteristiche di sicurezza elettrica.

MCERTS

MCERTS sta per Monitoring Certification Scheme, il programma di certificazione della Environment Agency, l'ente istituzione britannico per la protezione ambientale. Il programma offre uno schema in base al quale effettuare le misurazioni ambientali rispettando i requisiti qualitativi dell'Environment Agency e comprende varie attività di monitoraggio, campionatura e ispezione.

Lo strumento GEM5000 ha la certificazione MCERTS solo se:

- Dopo l'accensione iniziale, sullo schermo compare il logo MCERTS.

✎ Nota: MCERTS - Su questo strumento non si sono effettuati test di sensibilità incrociata con l'utilizzo di idrogeno solforato. Pertanto gli utenti devono accertare la presenza di H_2S , che potrebbe avere un effetto interferenziale.

MCERTS promuove fiducia nel monitoraggio dei dati e offre al settore una struttura sperimentata per scegliere sistemi e servizi di monitoraggio che soddisfino i requisiti dell'Environment Agency.

L'Environment Agency ha istituito il Monitoring Certification Scheme (MCERTS) per conseguire misurazioni ambientali qualitativamente valide. Il programma MCERTS offre la certificazione dei prodotti secondo gli standard qualitativi dell'Environment Agency, in base alle norme CEN, ISO e nazionali.

Gli strumenti con certificazione MCERTS sono stati testati da un organismo indipendente per verificare che soddisfino determinate caratteristiche di prestazione. Inoltre le aziende dei prodotti MCERTS vengono regolarmente controllate per accertare che le caratteristiche di prestazione da attestare nel certificato siano sempre conseguite.

Gli analizzatori di gas della serie 5000 sono stati certificati secondo la versione 3.1 delle 'Caratteristiche di prestazione dei sistemi portatili di monitoraggio delle emissioni'.

Batteria e ricarica

La batteria utilizzata negli analizzatori di gas della serie 5000 è al nichel-idruro metallico e viene prodotta come pacco con sei celle individuali. Anche se questo tipo di batteria non è soggetta all'effetto memoria della ricarica come quelle al nichel cadmio, non è consigliabile raccomandabile effettuare piccole ricariche parziali.

Il caricabatteria va scollegato solo quando indica la carica completa.



Avvertenza

Il caricabatteria NON è coperto dalla certificazione Ex. La batteria va caricata solo in un'area sicura.

Il caricabatteria è intelligente e indica lo stato di carica in corso o avvenuta.

Lo strumento va caricato utilizzando ESCLUSIVAMENTE il caricabatteria fornito in dotazione, che è inteso solo per l'uso in un locale interno. Effettuare la ricarica in un locale ventilato.

Caricabatteria:	Tensione d'ingresso:	100-240V CA +/- 10%
	Frequenza d'ingresso:	50-60Hz +/- 10%
	Corrente d'ingresso:	0,4A@100VCA .. 0,2A@240VCA
	Tensione d'uscita:	10,1VCC max
	Corrente d'uscita:	1,5A max

✍ Nota: collegare il caricabatteria all'elettricità di rete utilizzando l'adattatore necessario. Per ulteriori informazioni rivolgersi alla casa produttrice.

Istruzioni per la pulizia

NON utilizzare detergenti per pulire l'analizzatore o il caricabatteria, perché potrebbero avere un effetto negativo sulla sicurezza del loro uso.

Prassi ottimali per rilevare le letture



Avvertenza

L'inalazione del gas di solfuro d'idrogeno (H₂S) o di altri gas nocivi può causare la morte. Spetta all'utente accertarsi di essere sufficientemente addestrato negli aspetti della sicurezza relativi all'uso di H₂S e altri gas nocivi. In particolare nel caso di gas pericolosi, quelli scaricati dall'analizzatore devono essere convogliati in un'area in cui lo scaricotale operazione possa essere effettuata in condizioni di sicurezza. È possibile che lo strumento scarichi gas pericolosi anche durante lo spurgo quando viene utilizzato per spurgare con aria pulita.

Prassi ottimali

- Nel recarsi al luogo di utilizzo, collocare l'analizzatore di gas nell'abitacolo del veicolo, non nel bagagliaio, dove potrebbe essere soggetto a estremi di temperatura e danneggiarsi per eventuali urti. Non appoggiare l'analizzatore di gas contro superfici calde (ad es. tubo di aspirazione del gas, carrozzeria di un'autovettura o interno di un'autovettura incustodita in estate), perché ciò fa aumentare la temperatura dell'analizzatore di gas e può falsare le letture.
- Negli spostamenti nell'area di utilizzo dell'analizzatore di gas, proteggerlo dalla luce diretta del sole e dagli scrosci di pioggia.

Usare sempre il sifone per lo scarico della condensa!. Se il sifone si allaga, cambiare il filtro e controllare che tutti i tubi siano liberi dalla condensa prima di riutilizzarlo

✎ Nota: se lo scarico di un analizzatore di gas serie GA5000 è collegato ad un sistema pressurizzato, ciò fa uscire un flusso di gas dall'entrata.

Taratura

 Avvertenza	I gas di taratura possono essere pericolosi. Per ciascun gas da utilizzare, leggere preventivamente la relativa scheda tecnica di sicurezza, accertandosi di comprenderne il contenuto.
---	--

Il regolatore fornito col kit di taratura è stato configurato per permettere l'erogazione di un flusso fisso.

Dato che il flusso del regolatore è impostato in fabbrica, può essere aperto con una semplice rotazione, senza effettuare regolazioni.

 Avvertenza	Luce di scarico Durante la taratura dell'analizzatore, vi sono due possibili uscite per il gas: come di norma dalla luce di scarico (gialla) dell'analizzatore o, in caso di sovrappressione, dalla luce di 1/16" della valvola limitatrice di pressione rossa situata sul regolatore. Si raccomanda che entrambe le luci siano dotate di tubi di scarico, che devono sboccare in un'area ben ventilata. Controllare che i tubi e i raccordi non presentino perdite. La taratura dell'analizzatore di gas va eseguita in un'area idonea, adottando tutte le precauzioni necessarie per l'utilizzo di gas potenzialmente pericolosi, esplosivi o tossici.
---	---

✎ Nota: è possibile anche che il gas esca dall'uscita del flusso interno (blu) dell'analizzatore di gas (vale solo per GA5000).

Manutenzione

Gli analizzatori di gas della serie 5000 devono essere sottoposti a regolare controllo di manutenzione per accertare che il funzionamento sia corretto e preciso. Geotech raccomanda che la manutenzione e la taratura siano effettuate ogni 6 mesi.

Gli analizzatori di gas della serie 5000 hanno la certificazione ATEX per l'utilizzo in aree

potenzialmente esplosive. In quanto tali, la loro manutenzione deve essere affidata esclusivamente a tecnici qualificati. La mancata osservanza di questa regola invalida la garanzia ed eventualmente anche la certificazione ATEX.

 Avvertenza	Se la manutenzione dell'analizzatore di gas viene effettuata da tecnici non qualificati, la certificazione ATEX può venirne compromessa, pregiudicando così l'utilizzo in condizioni di sicurezza dello strumento in un'atmosfera potenzialmente esplosiva.
---	--

Parti riparabili dall'utente:

Non vi sono parti riparabili dall'utente all'interno dello strumento.
Le parti seguenti possono essere riparate:

Filtro acqua in linea	Va controllato regolarmente per evidenziare eventuali ostruzioni, condensa o danneggiamenti e sostituito se necessario. Lo strumento non va mai usato senza il filtro in linea, per evitare possibili infiltrazioni d'acqua.
Tubi di campionamento	Controllare sempre che i tubi di campionamento non siano contaminati o danneggiati.
Raccordi del gas	Controllare periodicamente che i gommini O-ring dei raccordi del gas non siano danneggiati. Se danneggiati, gli O-ring lasciano entrare aria nei tubi di campionamento, sfalsando le letture. Se l'O-ring risulta danneggiato, è necessario sostituire tutto il raccordo.
Materiale del filtro H ₂ S	Quando il materiale del filtro cambia colore e diventa <i>grigio chiaro</i> , il filtro va sostituito.

15.2 Instructions for safe use – German language

Sicherheitsvorschriften

 Warnhinweise	Die Gasanalysatoren der 5000er Serie können zum Messen der Gase von Deponien und anderen Quellen wie in diesem Handbuch beschrieben verwendet werden. Der Bediener kann bei Verwendung des Geräts schädlichen Gasen ausgesetzt werden. Das Einatmen dieser Gase kann gesundheitsschädlich und in manchen Fällen sogar tödlich sein. Es liegt in der Verantwortung des Benutzers sicherzustellen, dass er/sie angemessen über die Sicherheitsaspekte der eingesetzten Gase geschult ist und geeignete Verfahren befolgt werden. Vor allem beim Einsatz gefährlicher Gase muss das vom Analysator ausströmende Gas in einen Bereich geleitet werden, in dem das Gas sicher abgeführt werden kann. Gefährliches Gas kann ebenso vom Gerät ausgestoßen werden, wenn es mit sauberer Luft gereinigt wird.
---	--

 **Hinweis:** Gasanalysatoren sind empfindliche wissenschaftliche Geräte und sollten entsprechend behandelt werden. Wenn das Gerät anders als vom Hersteller spezifiziert verwendet wird, kann der vom Gerät gebotene Schutz beeinträchtigt werden.

Die Gasanalysatoren der 5000er Serie erfüllen Abschnitt 15 der FCC-Vorschriften. Der Betrieb unterliegt den folgenden zwei Bedingungen:

- 1) Dieses Gerät darf keine schädlichen Funkstörungen verursachen.
- 2) Dieses Gerät muss mögliche empfangene Funkstörungen und dadurch verursachte Funktionsstörungen dulden.

Für ATEX und IECEx sind die Gasanalysatoren der 5000er Serie für den Einsatz in explosionsgefährdeten Bereichen zertifiziert.



II 2G EX ib IIA T1 Gb (Ta = -10°C bis +50°C)

Die Vorschriften müssen unbedingt genau befolgt werden. Es liegt in der Verantwortung des Betreibers, das Schutzkonzept und die erforderliche Schutzklasse für eine bestimmte Anwendung festzulegen.

Vorschriften zur sicheren Verwendung - Deutsch

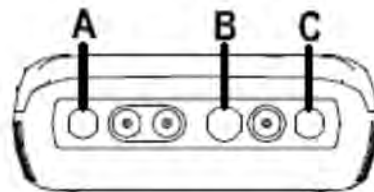
(Siehe Europäische ATEX-Richtlinie 94/9/EC, Anhang II, 1.0.6.)

Die folgenden Vorschriften gelten für Geräte, die in den Zertifikaten Nr. SIRA 11ATEX2197X und IECEx Richtlinie SIR 11.0089X behandelt werden:

- Die Geräte dürfen mit brennbaren Gasen und Dämpfen mit Apparategruppe IIA und Temperaturklasse T1 eingesetzt werden.
- Die Geräte sind nur für den Einsatz bei Umgebungstemperaturen im Bereich von -10 °C bis +50 °C zertifiziert und sollten nicht außerhalb dieses Bereichs eingesetzt werden.
- Die Geräte dürfen nicht in einer Atmosphäre mit mehr als 21 % Sauerstoffgehalt

eingesetzt werden.

- Die Reparatur dieser Geräte darf nur entsprechend der maßgeblichen Anleitung durchgeführt werden.
- Bei Einsatz in einem explosionsgefährdeten Bereich darf nur Temperaturfühler GF5.2 (SIRA 11ATEX2197X und IECExSIR11.0089X) verwendet werden. Anemometer GF5.4 (BVS 04ATEXE194), nur für den Einsatz mit ATEX, in Anschluss C. Der Analysator darf nicht an andere Geräte im explosionsgefährdeten Bereich angeschlossen werden, einschließlich des im Lieferumfang enthaltenen GF-USB-Kabels (Anschluss A) bzw. Ladegeräts GF3.9 (Anschluss B).



In einer explosionsgefährdeten Atmosphäre nicht laden, wieder aufladen oder öffnen.

In einem explosionsgefährdeten Bereich nur „Temperaturfühler GF5.2“ in Anschluss B verwenden.

**Anschluss C ($U_o=10\text{ V}$, $I_o=5\text{ mA}$, $P_o=50\text{ mW}$, $C_i=0$, $L_i=0$, $C_o=100\text{ uF}$, $L_o=1000\text{ mH}$),
Anschluss B ($U_o=5\text{ V}$, $I_o=6\text{ mA}$, $P_o=7\text{ mW}$, $C_i=0$, $L_i=0$, $C_o=100\text{ uF}$, $L_o=1000\text{ mH}$)**

MAXIMALE VERSORGUNG IN NICHT EXPLOSIONSGEFÄHRDETEN BEREICHEN:

Anschluss A - $U_m=6\text{ V}$ Anschluss B - $U_m=10,1\text{ V}$

- Falls die Möglichkeit besteht, dass die Geräte mit aggressiven Substanzen in Berührung kommen, z. B. mit sauren Flüssigkeiten oder Gasen, die Metalle angreifen können, oder mit Lösungsmitteln, die Polymerwerkstoffe schädigen können, liegt es in der Verantwortung des Benutzers, geeignete Sicherheitsvorkehrungen zu treffen, z. B. regelmäßige Kontrollen als Teil der Routineinspektionen oder die Prüfung des Materialdatenblatts darauf, ob das Gerät mit speziellen Chemikalien kompatibel ist, die es vor Schäden schützen, um zu gewährleisten, dass diese Art des Schutzes nicht beeinträchtigt wird.
- Der relative Druckbereich ist $\pm 500\text{ mbar}$. Es ist jedoch zu beachten, dass der Eingangsdruck $\pm 500\text{ mbar}$ relativ zum Atmosphärendruck nicht überschreiten darf und dass der Ausgangsdruck $\pm 100\text{ mbar}$ relativ zum Atmosphärendruck nicht überschreiten darf.

Für CSA (Kanada) sind die Gasanalysatoren der 5000er Serie für den Einsatz in explosionsgefährdeten Bereichen zertifiziert.

KLASSE 2258 03 - PROZESSKONTROLLGERÄT - Eigensichere und nicht zündgefährliche Systeme - Für explosionsgefährdete Standorte



Ex ib IIA:

Methandektoren Modell GA 5000, GEM 5000 und BIOGAS 5000; tragbar, batteriebetrieben mit nicht im Feld austauschbarem Akkupack Teilnr. 20087 oder 2011113; eigensicher, bietet eigensichere Kreise („[ib]“ für Zone 1) für Temperaturfühler Modell GFS.2 (Anschluss B), mit Entitätsausgabenparameter wie unten aufgeführt; Temperaturcode T1; $-10\text{ °C} \leq T_{amb.} \leq +50\text{ °C}$.

Anschluss	Entitätsparameter						
	Uo (V)	Io (mA)	Po (mW)	Co (uF)	Lo (mH)	Ci (uF)	Li (mH)
B	5,0	6	7	100	1000	0	0
C	10,0	5	50	100	1000	0	0

☞ Hinweis: Dieses Gerät wurde nur auf elektrische Sicherheitsfunktionen untersucht.

Für CSA (USA) sind die Gasanalysatoren der 5000er Serie für den Einsatz in explosionsgefährdeten Bereichen zertifiziert.

KLASSE 2258 83 - PROZESSKONTROLLGERÄT - Eigensichere und nicht zündgefährliche Systeme - Für explosionsgefährdete Standorte - NACH US-AMERIKANISCHEN NORMEN ZERTIFIZIERT



AEx ib IIA:

Methandektoren Modell GA 5000, GEM 5000 und BIOGAS 5000; tragbar, batteriebetrieben mit nicht im Feld austauschbarem Akkupack Teilnr. 20087 oder 2011113; eigensicher, bietet eigensichere Kreise („ib“) für Zone 1) für Temperaturfühler Modell GFS.2 (Anschluss B), mit Entitätsausgabenparameter wie unten aufgeführt; Temperaturcode T1; -10 °C ≤ Tamb. ≤ +50 °C.

Anschluss	Entitätsparameter						
	Uo (V)	Io (mA)	Po (mW)	Co (uF)	Lo (mH)	Ci (uF)	Li (mH)
B	5,0	6	7	100	1000	0	0
C	10,0	5	50	100	1000	0	0

☞ Hinweis: Dieses Gerät wurde nur auf elektrische Sicherheitsfunktionen untersucht.

MCERTS

MCERTS ist das Monitoring Certification Scheme (Zertifizierungsprogramm für Überwachungsgeräte) der britischen Umweltagentur. Das Programm bietet einen Rahmen, in dem Umgebungsmessungen gemäß den Qualitätsanforderungen der Agentur durchgeführt werden können. Es umfasst eine Reihe von Überwachungs-, Probenahme- und Prüfaufgaben.

Das Instrument GEM5000 ist nur MCERTS-zertifiziert, falls:

- Das MCERTS-Logo nach dem erstmaligen Einschalten auf dem Bildschirm erscheint.

☞ Hinweis: MCERTS - Bei diesem Gerät wurden keine Störempfindlichkeitsprüfungen mit Schwefelwasserstoff durchgeführt. Daher sollten Benutzer wissen, ob H₂S vor Ort vorhanden ist, da dies eine Störwirkung zur Folge haben könnte.

MCERTS fördert das öffentliche Vertrauen in Überwachungsdaten und liefert der Industrie einen erprobten Rahmen zur Auswahl von Überwachungssystemen und -dienstleistungen, die die Leistungsanforderungen der Umweltagentur erfüllen.

Die britische Umweltagentur hat MCERTS (Monitoring Certification Scheme) initiiert, um hochwertige Umweltmessungen bereitzustellen. Das MCERTS

Produktzertifizierungsprogramm ermöglicht die Zertifizierung von Produkten gemäß den Leistungsstandards der Umweltagentur basierend auf den entsprechenden CEN-, ISO- und nationalen Normen.

MCERTS-zertifizierte Geräte wurden durch eine unabhängige Stelle geprüft, um zu gewährleisten, dass bestimmte Leistungsanforderungen erfüllt werden. Darüber hinaus wird der Hersteller eines MCERTS-Produkts regelmäßig geprüft, um zu gewährleisten, dass die Leistungsanforderungen der Zertifizierung durchgehend erfüllt werden.

Die Gasanalysatoren der 5000er Serie wurden gemäß Version 3.1 der „Leistungsanforderungen für tragbare Emissionsüberwachungssysteme“ zertifiziert.

Akku und Aufladen

Bei dem Akku, der in den Gasanalysatoren der 5000er Serie verwendet wird, handelt es sich um einen Nickel-Metallhydrid-Akku, der als Akkupack aus sechs einzelnen Zellen hergestellt wird. Dieser Akkutyp ist weniger stark für den Memoryeffekt anfällig als Nickel-Kadmium-Akkus. Trotzdem wird davon abgeraten, das Gerät mit kleinen Ladungen nachzuladen.

Das Ladegerät sollte nur getrennt werden, wenn komplette Ladung angezeigt wird.



Warnhinweise

Das Akkuladegerät ist NICHT in der ATEX-Zertifizierung enthalten. Der Akku darf nur in einem sicheren Bereich aufgeladen werden.

Das Akkuladegerät ist intelligent und zeigt an, wenn die Einheit aufgeladen wird bzw. wenn sie vollständig aufgeladen ist.

Das Gerät darf NUR mit dem mitgelieferten Akkuladegerät aufgeladen werden. Das Akkuladegerät ist nur für den Gebrauch in Innenräumen vorgesehen. Bitte sorgen Sie für ausreichende Belüftung während des Aufladens.

Ladegerät:	Eingangsspannung:	100-240 V AC +/- 10 %
	Eingangsfrequenz:	50-60 Hz +/- 10 %
	Eingangsstromstärke:	0,4 A bei 100 VAC .. 0,2 A bei 240 VAC
	Ausgangsspannung:	Max. 10,1 VDC
	Ausgangsstromstärke:	Max. 1,5 A

✍ Hinweis: Schließen Sie das Ladegerät durch Verbinden des entsprechenden Adapters an das Stromnetz an. Wenden Sie sich für weitere Informationen an den Hersteller.

Reinigungsanweisungen

Verwenden Sie KEINE Reinigungsmittel zum Reinigen des Analysegeräts oder Akkuladegeräts, da sie die sichere Verwendung dieser Geräte beeinträchtigen können.

Good Practice beim Ablesen



Warnhinweise

Das Einatmen von Schwefelwasserstoffgas (H₂S) oder anderer schädlicher Gase kann tödlich sein. Es liegt in der Verantwortung des Benutzers sicherzustellen, dass er/sie angemessen über die Sicherheitsaspekte beim Einsatz von H₂S und anderer schädlicher Gase geschult ist. Vor allem beim Einsatz gefährlicher Gase muss das vom Analysator ausströmende Gas in einen Bereich geleitet werden, in dem das

	Gas sicher abgeführt werden kann. Gefährliches Gas kann ebenso vom Gerät ausgestossen werden, wenn es mit sauberer Luft ausgeblasen wird.
--	---

Good Practice

- Transportieren Sie den Gasanalysator im Inneren des Fahrzeugs an den Einsatzort - nicht auf der Ladefläche, wo er Temperaturschwankungen und möglichen Stößen ausgeliefert ist. Platzieren Sie den Gasanalysator nicht direkt an oder auf etwas Heißem (z. B. Gasleitung, Karosserie oder in einem unbeaufsichtigten Auto während des Sommers), da dies einen Temperaturanstieg im Gasanalysator verursacht und zu fehlerhaften Messwerten führen kann.
- Schützen Sie den Gasanalysator am Einsatzort vor starkem, direktem Sonnenlicht und starkem Regen.

Verwenden Sie stets den Wasserabscheider! Wenn der Wasserabscheider überschwemmt wird, tauschen Sie den Filter aus, und stellen Sie sicher, dass alle Schläuche frei von Feuchtigkeit sind, bevor Sie sie erneut verwenden.

 Hinweis: Wenn der Auslass eines Gasanalysators der Serie GA5000 an ein druckbeaufschlagtes System angeschlossen ist, dann führt dies zu einem Gasstrom aus dem Einlassstromanschluss.

Kalibrieren

 Warnhinweise	Eichgase können gefährlich sein. Vor dem Verfahren müssen die Materialsicherheitsdatenblätter aller verwendeten Gase gelesen und verstanden werden.
---	--

Der mit dem Kalibrierset gelieferte Regulator wurde so konfiguriert, dass er einen unveränderlichen Durchfluss liefert.

Da der Durchfluss des Regulators werkseitig eingestellt ist, lässt der Regulator sich mit nur wenigen Drehungen öffnen; eine Einstellung ist nicht erforderlich.

 Warnhinweise	Austrittsöffnung Wenn der Gasanalysator kalibriert wird, gibt es zwei mögliche Ausgänge für das Gas: auf die übliche Art über die Austrittsöffnung (gelb) des Analysators oder bei Überdruck über die 1/16" Öffnung des roten Druckentlastungsventils, das sich am Regler befindet. Für beide Öffnungen wird das Anbringen von Auslassrohrleitungen empfohlen. Die Auslassrohrleitungen müssen in einen gut belüfteten Bereich führen. Stellen Sie sicher, dass es in den Rohrleitungen und an den Verbindungen keine Undichtigkeiten gibt. Das Kalibrieren des Gasanalysators muss in einem sicheren Bereich unter Anwendung aller notwendigen Sicherheitsvorkehrungen durchgeführt werden, wenn möglicherweise gefährliche, explosive oder toxische Gase
---	---

	eingesetzt werden.
--	--------------------

- ☞ Hinweis: Es besteht außerdem die Möglichkeit, dass Gas aus dem internen Strömungsanschluss (blau) des Gasanalysators ausgestoßen wird (gilt nur für GA5000).

Wartung

Die Gasanalysatoren der 5000er Serie müssen regelmäßig gewartet werden, um den korrekten und genauen Betrieb zu gewährleisten. Geotech empfiehlt ein Wartungs- und Neukalibrierungsintervall von 6 Monaten.

Die Gasanalysatoren der 5000er Serie sind für den Einsatz in explosionsgefährdeten Bereichen ATEX-zertifiziert. Daher dürfen sie nur von qualifizierten Technikern gewartet werden. Im Falle der Nichtbefolgung erlischt die Garantie, und die ATEX-Zertifizierung kann ihre Gültigkeit verlieren.

 Warnhinweise	Falls der Gasanalysator von unqualifizierten Technikern gewartet wird, kann die ATEX-Zertifizierung ihre Gültigkeit verlieren, und das Gerät ist möglicherweise für den Einsatz in einer explosionsgefährdeten Atmosphäre nicht mehr sicher.
---	--

Teile, die durch den Benutzer gewartet werden können:

Im Gerät gibt es keine Teile, die durch den Benutzer gewartet werden können.

Die folgenden Teile können durch den Benutzer gewartet werden:

Wasserleitungsfilter	Er muss regelmäßig auf Verstopfungen, Feuchtigkeit bzw. Beschädigungen untersucht und, falls erforderlich, ausgetauscht werden. Das Gerät darf nie ohne den Wasserleitungsfilter betrieben werden, da dies dazu führen könnte, dass Wasser in das Gerät eindringt.
Probenrohrleitungen	Stellen Sie stets sicher, dass die Probenrohrleitungen weder verunreinigt noch beschädigt sind.
Gasausgangsanschlüsse	Prüfen Sie periodisch, ob die O-Ringe an den Gasausgangsanschlüssen beschädigt sind. Durch einen beschädigten O-Ring kann Luft in das Messgas eindringen und zu falschen Messwerten führen. Falls der O-Ring beschädigt ist, muss der gesamte Gasausgangsanschluss ausgetauscht werden.
H ₂ S Filtermaterial	Sobald die Farbe des Filtermaterials <i>leicht grau</i> wird, muss der Filter ausgetauscht werden.

15.3 Instructions for safe use – French language

Instructions concernant la sécurité

 Avertissement	Les analyseurs de gaz de la série 5000 sont conçus pour mesurer les gaz des sites d'enfouissement et d'autres sources, comme le décrit le présent manuel. L'opérateur risque d'être exposé à des gaz nocifs pendant l'utilisation de l'instrument. L'inhalation de ces gaz peut être nuisible à la santé et, dans certains cas, mortelle. Il incombe à l'utilisateur de s'assurer qu'il a reçu une formation adaptée aux aspects de la sécurité des gaz utilisés et de s'assurer du respect des procédures appropriées. En particulier, lors de l'utilisation de gaz dangereux, les gaz en sortie de l'analyseur doivent être évacués dans une zone où ils ne présentent aucun danger. Des gaz dangereux peuvent être également expulsés de l'instrument lors d'une purge à l'air propre.
--	---

☞ Remarque : Les analyseurs de gaz sont des instruments scientifiques sensibles qu'il convient de traiter en conséquence. Toute utilisation du matériel non conforme aux instructions du fabricant risque d'amoindrir la protection assurée par l'instrument.

Les analyseurs de gaz de la série 5000 sont conformes à l'article 15 de la réglementation FCC (Federal Communications Commission - Conseil supérieur de l'audiovisuel américain). Son utilisation est soumise aux deux conditions suivantes :

- 1) Cet appareil ne doit pas provoquer d'interférences nuisibles.
- 2) Cet appareil doit accepter toutes les interférences reçues, y compris celles qui pourraient provoquer un fonctionnement indésirable.

Dans le cadre des certifications ATEX et IECEx, l'analyseur de gaz de la série 5000 est certifié pour la catégorie zone dangereuse.

 **II 2G Ex ib IIA T1 Gb (Ta = -10°C à +50°C)**

Il est absolument indispensable de respecter les instructions contenues dans ce manuel. Il incombe à l'utilisateur de déterminer le type et la classification de protection requise pour une application spécifique.

Instructions pour une utilisation sûre– Français

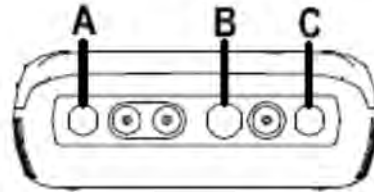
(Référence Directive européenne ATEX 94/9/CE, Annexe II, 1.0.6.)

Les instructions suivantes s'appliquent au matériel couvert par les numéros de certificat SIRA 11ATEX2197X et IECEx Directive SIR 11.0089X :

- Le matériel est utilisable avec des gaz et des vapeurs inflammables et des appareils de groupe IIA et de classe de température T1.
- Le matériel est certifié uniquement pour une utilisation à température ambiante entre -10°C et +50°C et ne doit pas être utilisé en dehors de cette plage.
- Le matériel ne doit pas être utilisé dans une atmosphère contenant plus de 21%

d'oxygène.

- Ce matériel devra être réparé conformément au code de pratique applicable.
- Lors d'une utilisation en zone dangereuse, utiliser exclusivement une sonde de température GF5.2 (SIRA 11ATEX2197X et IECEx SIR11.0089X). Pour le connecteur C, l'anémomètre GF5.4 (BVS 04ATEXE194) ne s'utilise qu'avec les dispositifs certifiés ATEX. L'analyseur ne doit être raccordé à aucun autre dispositif dans la zone dangereuse, ni au câble GF-USB (connecteur A) ni au chargeur de batterie GF3.9 (connecteur B) fourni avec l'analyseur.



**Ne pas charger, recharger ni ouvrir en atmosphère potentiellement explosive.
Dans une zone dangereuse, utiliser exclusivement la sonde de température GF5.2
branchée dans le connecteur B.**

Connecteur C ($U_o=10\text{ V}$, $I_o=5\text{ mA}$, $P_o=50\text{ mW}$, $C_i=0$, $L_i=0$, $C_o=100\text{ uF}$, $L_o=1\text{ 000 mH}$),

Connecteur B ($U_o=5\text{ V}$, $I_o=6\text{ mA}$, $P_o=7\text{ mW}$, $C_i=0$, $L_i=0$, $C_o=100\text{ uF}$, $L_o=1\text{ 000 mH}$)

ALIMENTATIONS MAXIMALES NON DANGEREUSES :

Connecteur A - $U_m=6\text{ V}$ Connecteur B - $U_m=10,1\text{ V}$

- Si le matériel est amené à être en contact avec des substances corrosives, par exemple des liquides ou des gaz acides susceptibles d'attaquer les métaux, ou des solvants pouvant affecter des polymères, il incombe alors à l'utilisateur de prendre des précautions appropriées, par exemple des contrôles réguliers dans le cadre d'inspections systématiques, ou des vérifications sur la fiche technique de la résistance du matériau à des produits chimiques spécifiques, ceci afin de préserver l'intégrité de la protection.
- La plage de pression relative est de $\pm 500\text{ mbar}$. Cependant, il convient de noter que la pression d'entrée ne doit pas dépasser $\pm 500\text{ mbar}$ par rapport à la pression atmosphérique et la pression de sortie ne doit pas dépasser $\pm 100\text{ mbar}$ par rapport à la pression atmosphérique.

Pour CSA (Canada), l'analyseur de gaz de la série 5000 est certifié pour la catégorie zone dangereuse

CLASSE 2258 03 - ÉQUIPEMENT DE CONTRÔLE DES PROCÉDÉS - Appareils à sécurité intrinsèque et appareils non incendiaires pour emplacements dangereux



Ex ib IIA :

Détecteurs de méthane Modèle GA 5000, GEM 5000 et BIOGAS 5000 ; appareils portatifs, avec batterie alimentée par bloc de batterie (numéro de pièce 20087 ou 2011113) non remplaçable sur place ; à sécurité intrinsèque et fournissant des circuits à sécurité intrinsèque (« [ib] » pour zone 1) pour sonde de température modèle GF5.2 (connecteur B) et avec paramètres de sortie comme indiqué au tableau ci-dessous ; code de température T1; $-10\text{ °C} \leq T_{amb.} \leq +50\text{ °C}$.

Connecteur	Paramètres						
	Uo (V)	Io (mA)	Po (mW)	Co (uF)	Lo (mH)	Ci (uF)	Li (mH)
B	5,0	6	7	100	1 000	0	0
C	10,0	5	50	100	1 000	0	0

✍ Remarque : Les tests effectués sur cet appareil concernaient uniquement la sécurité électrique.

Pour CSA (États-Unis), l'analyseur de gaz de la série 5000 est certifié pour la catégorie zone dangereuse

CLASSE 2258 83 - ÉQUIPEMENT DE CONTRÔLE DES PROCÉDÉS - Appareils à sécurité intrinsèque et appareils non incendiaires pour emplacements dangereux - CERTIFICATION AUX NORMES AMÉRICAINES (ÉTATS-UNIS)



AEx ib IIA :

Détecteurs de méthane Modèle GA 5000, GEM 5000 et BIOGAS 5000 ; appareils portatifs, avec batterie alimentée par bloc de batterie (numéro de pièce 20087 ou 2011113) non remplaçable sur place ; à sécurité intrinsèque et fournissant des circuits à sécurité intrinsèque (« [ib] » pour zone 1) pour sonde de température modèle GF5.2 (connecteur B) et avec paramètres de sortie comme indiqué au tableau ci-dessous ; code de température T1; $-10\text{ }^{\circ}\text{C} \leq T_{\text{amb.}} \leq +50\text{ }^{\circ}\text{C}$.

Connecteur	Paramètres						
	Uo (V)	Io (mA)	Po (mW)	Co (uF)	Lo (mH)	Ci (uF)	Li (mH)
B	5,0	6	7	100	1 000	0	0
C	10,0	5	50	100	1 000	0	0

✍ Remarque : Les tests effectués sur cet appareil concernaient uniquement la sécurité électrique.

MCERTS

MCERTS est le programme de certification de la surveillance, établi par l'agence britannique à l'Environnement. Ce programme forme le cadre dans lequel des mesures environnementales peuvent être effectuées conformément aux exigences de qualité de l'agence. Il couvre un ensemble d'activités de surveillance, d'échantillonnage et d'inspection.

L'instrument GEM5000 n'est certifié MCERTS que si :

- Le logo MCERTS s'affiche à l'écran après la mise sous tension de l'instrument.

✍ Remarque : MCERTS - Cet instrument n'a pas fait l'objet de tests de sensibilité croisée utilisant le sulfure d'hydrogène. Par conséquent, les utilisateurs doivent être conscients qu'en cas de présence de H_2S sur les sites, il peut y avoir un effet d'interférence.

Le programme MCERTS contribue à renforcer la confiance du public vis-à-vis des données de surveillance et donne à l'industrie des paramètres sûrs pour le choix de systèmes et de services de surveillance répondant aux exigences de performance de l'agence.

L'agence à l'Environnement a établi ce programme MCERTS (Monitoring Certification

Scheme) pour fournir des valeurs mesurées environnementales de qualité. La certification MCERTS concerne les produits aux normes de performance de l'agence à l'Environnement, sur la base des normes nationales, CEN et ISO pertinentes.

Les instruments certifiés MCERTS sont testés par un organisme indépendant pour assurer leur conformité à certaines exigences de performance. En outre, le fabricant de produits MCERTS fait l'objet d'audits réguliers pour s'assurer du respect continu des exigences de performance de ses produits aux fins de certification.

Les analyseurs de gaz de la série 5000 ont été certifiés conformément à la version 3.1 des « Normes de performance des systèmes portatifs de surveillance des émissions de substances dans l'atmosphère ».

Batterie et mise en charge

Les analyseurs de gaz de la série 5000 sont dotés d'une batterie au nickel-métal-hydrure, à six cellules individuelles intégrées. Ce type de batterie n'est pas aussi sensible à « l'effet de mémoire » qui affecte les capacités des piles au nickel cadmium. Il est cependant déconseillé de recharger la batterie par à-coups.

Débrancher le chargeur uniquement après indication de pleine charge.



Avertissement

Le chargeur de batterie n'est PAS couvert par la certification Ex. La batterie ne doit être chargée que dans un endroit sûr.

Le chargeur de batterie est intelligent et indique que la batterie est en cours de chargement ou chargée.

La batterie doit être chargée UNIQUEMENT à l'aide du chargeur fourni avec l'instrument. Le chargeur de batterie doit être utilisé uniquement à l'intérieur. Veiller à assurer une ventilation adéquate pendant le chargement.

Chargeur :	Tension d'entrée :	100-240 V CA +/- 10 %
	Fréquence d'entrée :	50-60 Hz +/- 10 %
	Courant d'entrée :	0,4 A@100 V CA.. 0,2 A@240 V CA
	Tension de sortie :	10,1 V CC max
	Courant de sortie :	1,5 A max

✍ Remarque : connecter le chargeur à l'alimentation secteur à l'aide de l'adaptateur approprié. Contacter le fabricant pour de plus amples informations.

Instructions de nettoyage

Ne PAS utiliser d'agents nettoyants pour nettoyer l'analyseur ou le chargeur de batterie car ces produits risquent d'avoir un effet préjudiciable sur l'utilisation sûre de ces appareils.

Bonnes pratiques lors des mesures**Avertissement**

L'inhalation de sulfure d'hydrogène (H₂S) ou d'autres gaz dangereux peut entraîner la mort. Il incombe à l'utilisateur de s'assurer qu'il a reçu la formation adaptée aux aspects sécurité de l'utilisation de H₂S et d'autres gaz dangereux. En particulier, lors de l'utilisation de gaz dangereux, les gaz en sortie de l'analyseur doivent être évacués dans une zone où ils ne présentent aucun danger. Des gaz dangereux peuvent aussi être expulsés de l'instrument lors d'une purge à l'air propre.

Bonnes pratiques

- Lors d'un déplacement vers un site d'utilisation, transporter l'instrument dans l'habitacle du véhicule, et non pas dans le coffre où il pourrait être soumis à des températures extrêmes, voire à des chocs risquant de l'endommager. Ne pas placer l'analyseur de gaz contre des surfaces chaudes (conduite d'extraction de gaz, carrosserie d'une voiture ou voiture laissée sans surveillance en été, par exemple) car les effets de la chaleur sur l'analyseur risquent d'entraîner des valeurs mesurées incorrectes.
- Lors du déplacement sur site, protéger l'analyseur de gaz de la lumière solaire directe et des fortes intempéries.

Toujours utiliser le piège à eau ! Si le piège à eau déborde, changer le filtre et vérifier que tous les tuyaux ne présentent aucune trace d'humidité avant toute utilisation.

✍ Remarque : si l'évacuation d'un analyseur de gaz de la série GA5000 est connectée à un système sous pression, un flux de gaz s'échappera du port de débit d'entrée.

Étalonnage**Avertissement**

Les gaz d'étalonnage peuvent être dangereux.
Pour chaque gaz utilisé, il convient de lire et de comprendre la fiche de données de sécurité correspondante avant de poursuivre.

Le régulateur fourni avec le kit d'étalonnage a été configuré pour assurer un débit fixe.

Le débit du régulateur étant réglé en usine, il ne faut que quelques tours pour l'ouvrir et aucun réglage n'est nécessaire.

**Avertissement****Orifice d'évacuation**

Lors de l'étalonnage de l'analyseur de gaz, le gaz peut être évacué par deux orifices : par l'orifice d'évacuation normal (jaune) de l'analyseur ou, dans les cas de surpression, par l'orifice de 1/16 de pouce de la soupape de surpression rouge située sur le régulateur.

Il est recommandé de raccorder des tuyaux d'évacuation à ces deux orifices.

Le tuyau d'évacuation doit laisser les gaz s'échapper dans un endroit bien ventilé. Vérifier que les tuyaux et les

	raccords ne présentent aucune fuite. L'étalonnage de l'analyseur de gaz doit s'effectuer dans un endroit sûr, en observant toutes les précautions nécessaires en présence de gaz potentiellement dangereux, explosifs ou toxiques.
--	--

✍ Remarque : le gaz peut être aussi expulsé au niveau du port de débit interne (bleu) de l'analyseur de gaz (applicable uniquement au modèle GA5000).

Entretien

Pour un fonctionnement correct et précis, l'analyseur de gaz de la série 5000 doit faire l'objet d'un entretien régulier. Geotech recommande un entretien et un réétalonnage tous les 6 mois.

Les analyseurs de gaz de la série 5000 sont certifiés ATEX pour l'utilisation en environnements potentiellement explosifs. En conséquence, leur entretien doit être effectué uniquement par des techniciens qualifiés. Le non-respect de cette exigence entraînera l'annulation de la garantie, voire de la certification ATEX.

 Avertissement	Si l'analyseur de gaz fait l'objet d'un entretien par des techniciens non qualifiés, la certification ATEX risque d'être annulée et l'appareil peut ne pas être sûr en cas d'utilisation dans un environnement potentiellement explosif.
--	---

Composants pouvant faire l'objet d'un entretien par l'utilisateur :

L'appareil ne contient aucun composant interne pouvant faire l'objet d'un entretien par l'utilisateur.

Les composants suivants peuvent faire l'objet d'un entretien par l'utilisateur :

Filtre à eau en ligne	Contrôler ce filtre régulièrement pour rechercher la présence d'obstructions, d'humidité ou de dommages ; le remplacer si besoin est. L'appareil ne doit jamais être utilisé sans le filtre à eau en ligne pour prévenir la pénétration d'eau dans l'appareil.
Tuyau d'échantillonnage	Toujours vérifier que les tuyaux d'échantillonnage ne sont ni contaminés ni endommagés.
Raccords d'orifices de gaz	Contrôler périodiquement les joints toriques des raccords d'orifices de gaz pour s'assurer qu'ils ne sont pas endommagés. Un joint torique endommagé peut laisser passer l'air dans le gaz d'échantillonnage et entraîner des valeurs mesurées incorrectes. Remplacer le raccord complet si le joint torique est endommagé.
Matériau du filtre à H ₂ S	Remplacer le filtre lorsque le matériau du filtre change de couleur et devient <i>gris clair</i> .

15.4 Instructions for safe use – Spanish language

Instrucciones de seguridad

 Advertencia	La serie 5000 de analizadores de gas puede usarse para medir gases de vertederos y otras fuentes de la forma descrita en este manual. El operario puede estar expuesto a gases perjudiciales durante el uso del instrumento. La inhalación de estos gases puede ser nociva para la salud y, en algunos casos, incluso mortal. El usuario es responsable de garantizar que está debidamente formado en los aspectos de seguridad de los gases utilizados y que se respetan los procedimientos adecuados, especialmente en los lugares en los que se usan gases peligrosos, en los cuales el gas emitido por el analizador debe conducirse por un tubo hasta una zona en la que pueda liberarse con seguridad. El instrumento también puede emitir gases peligrosos si se purga con aire limpio.
--	--

✎ **Nota:** los analizadores de gas son instrumentos especialmente delicados del equipamiento científico y deben tratarse con especial cuidado. Un uso del equipo no conforme a las especificaciones del fabricante podría afectar al sistema de protección.

La serie 5000 de los analizadores de gas cumple las estipulaciones establecidas en la Parte 15 de la normativa de la FCC. El funcionamiento depende de las dos condiciones siguientes:

- 1) El instrumento no debe causar interferencias perjudiciales.
- 2) El instrumento debe admitir cualquier interferencia que pueda recibir, incluidas aquellas que podrían causar un funcionamiento no deseado.

En cuanto a las directivas ATEX e IECEx, la serie 5000 de analizadores de gas ha recibido la certificación de clasificación de área peligrosa



II 2G Ex ib IIA T1 Gb (Ta = de -10 °C a +50 °C)

Es de vital importancia que se sigan rigurosamente las instrucciones. El operario es responsable de determinar la noción de protección y la clasificación necesaria para cada aplicación específica.

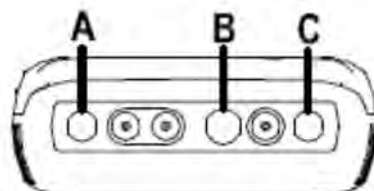
Instrucciones de seguridad (español)

(Directiva europea ATEX de referencia 94/9/CE, anexo II, 1.0.6.)

Las siguientes instrucciones se aplican a los instrumentos cubiertos por los números de certificado SIRA 11ATEX2197X y la norma IECEx SIR 11.0089X:

- El instrumento puede utilizarse con gases y vapores inflamables con el grupo de aparatos IIA y la clase de temperatura T1.
- El instrumento sólo está certificado para su uso a temperatura ambiente, comprendida entre -10 °C y +50 °C, y no debe utilizarse a diferentes temperaturas.
- El instrumento no debe usarse en lugares con una concentración de oxígeno superior al 21 %.

- La reparación del instrumento ha de realizarse de acuerdo con el código profesional aplicable.
- Si se utiliza en un área peligrosa, utilice solo la sonda de temperatura GF5.2 (SIRA 11ATEX2197X e IECEx SIR11.0089X). En cuanto al conector C, utilice el anemómetro GF5.4 (BVS 04ATEXE194) únicamente con ATEX. En las mencionadas zonas de riesgo, el analizador no debe conectarse a ningún otro aparato en el área peligrosa, incluidos el cable GF-USB (conector A) o el cargador de batería GF3.9 (conector B) suministrados con el propio analizador.



No cargar, recargar o abrir en lugares en los que exista riesgo de explosión. En áreas peligrosas, utilice solo la "sonda de temperatura GF5.2" en el conector B. Conector C ($U_o=10\text{ V}$, $I_o=5\text{ mA}$, $P_o=50\text{ mW}$, $C_i=0$, $L_i=0$, $C_o=100\text{ uF}$, $L_o=1000\text{ mH}$), Conector B ($U_o=5\text{ V}$, $I_o=6\text{ mA}$, $P_o=7\text{ mW}$, $C_i=0$, $L_i=0$, $C_o=100\text{ uF}$, $L_o=1000\text{ mH}$)

SUMINISTROS MÁXIMOS NO PELIGROSOS:
Conector A - $U_m=6\text{ V}$ Conector B - $U_m=10,1\text{ V}$

- En caso de riesgo de contacto con sustancias agresivas, p. ej. líquidos o gases ácidos que pueden atacar a los metales o disolventes que pueden afectar a los materiales poliméricos, el usuario es responsable de adoptar las medidas de precaución adecuadas, como las comprobaciones periódicas que sean necesarias como parte de las inspecciones rutinarias o determinar, a partir de la hoja de especificaciones del material, su resistencia a determinados productos químicos que puedan anular sus cualidades de protección, gracias a lo cual se asegura que el tipo de protección no queda comprometido.
- El intervalo de presión relativa es $\pm 500\text{ mbar}$. Tenga en cuenta que, sin embargo, la presión de entrada no debería exceder $\pm 500\text{ mbar}$ ni la presión de salida $\pm 100\text{ mbar}$ en relación a la presión atmosférica.

Con respecto a la CSA (Canadá), la serie 5000 de analizadores de gas cuenta con la certificación de clasificación de área peligrosa

CLASE 2258 03 - EQUIPO DE CONTROL DE PROCESOS - Sistemas intrínsecamente seguros y no inflamables - Para ubicaciones peligrosas



Ex ib IIA:

Detectores de metano modelos GA 5000, GEM 5000 y BIOGAS 5000; paquete de batería portátil, alimentado por batería no recambiable en el lugar de uso N/P 20087 o 2011113; intrínsecamente seguro y con circuitos intrínsecamente seguros ("ib" para zona 1) a la sonda de temperatura (conector B) modelo GF5.2 y con los parámetros de salida de entidad como se indica en la siguiente tabla; código de temperatura T1; temperatura ambiente de -10 °C a $+50\text{ °C}$.≤≤

Conector	Parámetros de entidad						
	U_o	I_o	P_o	C_o	L_o	C_i	L_i

	(V)	(mA)	(mW)	(uF)	(mH)	(uF)	(mH)
B	5,0	6	7	100	1000	0	0
C	10,0	5	50	100	1000	0	0

✍ Nota: Se ha investigado este instrumento únicamente en relación a las características de seguridad eléctrica.

Con respecto a la CSA (EE. UU.), la serie 5000 de analizadores de gas cuenta con la certificación de clasificación de área peligrosa

CLASE 2258 83 - EQUIPO DE CONTROL DE PROCESOS - Sistemas intrínsecamente seguros y no inflamables - Para ubicaciones peligrosas - CERTIFICADO PARA LA NORMATIVA DE EE. UU.



AEx ib IIA:

Detectores de metano modelos GA 5000, GEM 5000 y BIOGAS 5000; paquete de batería portátil, alimentado por batería no recambiable en el lugar de uso N/P 20087 o 2011113; intrínsecamente seguro y con circuitos intrínsecamente seguros ("ib" para zona 1) a la sonda de temperatura (conector B) modelo GF5.2 y con los parámetros de salida de entidad como se indica en la siguiente tabla; código de temperatura T1; temperatura ambiente de -10 °C a +50 °C.≤

Conector	Parámetros de entidad						
	Uo (V)	Io (mA)	Po (mW)	Co (uF)	Lo (mH)	Ci (uF)	Li (mH)
B	5,0	6	7	100	1000	0	0
C	10,0	5	50	100	1000	0	0

✍ Nota: Se ha investigado este instrumento únicamente en relación a las características de seguridad eléctrica.

MCERTS

MCERTS es el Esquema de certificación de control de la Agencia del Medio Ambiente del Reino Unido. Dicho esquema constituye el marco en el que realizar las mediciones medioambientales de acuerdo con los requisitos de calidad de la Agencia, y engloba toda una serie de actividades de control, recogida de muestras e inspección.

El instrumento GEM5000 cuenta con la certificación de MCERTS solo si:

- Aparece el logotipo de MCERTS en la pantalla al encenderlo.

✍ Nota: MCERTS - no se han llevado a cabo en este instrumento pruebas de sensibilidad cruzada con sulfuro de hidrógeno. Por lo tanto, el usuario debe tener en cuenta la posibilidad de interferencias en caso de presencia de H₂S.

El MCERTS proporciona fiabilidad y confianza a la supervisión de datos y ofrece a la industria un marco contrastado para la elección de los sistemas de supervisión y servicios conformes con las exigencias de rendimiento de la Agencia del Medio Ambiente.

El objetivo de la creación del Esquema de certificación de control (MCERTS) de la Agencia del Medio Ambiente es realizar controles medioambientales de calidad. El MCERTS ofrece certificaciones de productos de acuerdo con las normas de rendimiento de la Agencia del Medio Ambiente, tomando como base normas relevantes CEN, ISO y nacionales.

Los instrumentos con la certificación MCERTS han sido evaluados por un organismo independiente con el fin de garantizar el cumplimiento con determinadas exigencias de rendimiento. Además, el fabricante de un producto con certificación MCERTS está sometido a auditorías periódicas que garantizan el constante cumplimiento con las exigencias de rendimiento del certificado.

La serie 5000 de analizadores de gas cuenta con la certificación de la versión 3.1 de las Normas de rendimiento para sistemas portátiles de control de emisiones.

Batería y carga

La batería de la serie 5000 de analizadores de gas es un paquete de hidruro metálico de níquel compuesto por seis células individuales. Este tipo de batería es menos susceptible a los "efectos de memorización" en las cargas máximas que las baterías de níquel-cadmio, si bien no es recomendable cargar la unidad al máximo.

No desconecte el cargador hasta que se indique que el instrumento está totalmente cargado.

 Advertencia	NO se aplica la certificación Ex al cargador de la batería. Cargue siempre la batería en lugares seguros.
--	--

El cargador inteligente de la batería indica el estado de carga de la unidad y el fin de la carga.

El instrumento debe cargarse SÓLO con el cargador suministrado con el instrumento. El cargador de batería está concebido para su uso en interior. Asegúrese de que la ventilación es adecuada mientras carga la unidad.

Cargador: Voltaje de entrada: 100-240 V CA +/- 10%
Frecuencia de entrada: 50-60 Hz +/- 10%
Corriente de entrada: 0,4 A a 100 VCA .. 0,2 A a 240 VCA

Tensión de salida: 10,1 VCC máx.
Corriente de salida: 1,5 A máx.

 **Nota:** Conecte el cargador a la red eléctrica con el adaptador apropiado. Para más información, póngase en contacto con el fabricante.

Instrucciones de limpieza

NO utilice agentes limpiadores para limpiar el analizador o el cargador de la batería, ya que podrían tener efectos adversos en el uso seguro de los dispositivos.

Consejos de utilidad para efectuar lecturas

 Advertencia	La inhalación de sulfuro de hidrógeno (H ₂ S) o de otros gases nocivos puede ser mortal. El usuario es responsable de garantizar que está debidamente formado en aspectos de seguridad en cuanto al uso de H ₂ S y de otros gases nocivos especialmente en los lugares en los que se usan gases peligrosos, en los cuales el gas emitido por el analizador debe conducirse por un tubo hasta una zona en la que pueda liberarse con seguridad. El instrumento también puede emitir gases peligrosos si se purga con aire limpio.
--	--

Consejos de utilidad

- En los desplazamientos, el analizador de gas debe llevarse en el interior del vehículo

(nunca en el maletero), con objeto de evitar daños por temperaturas extremas o posibles impactos. No coloque el analizador de gas sobre superficies calientes (p. ej., tuberías de extracción de gas, carrocerías o vehículos aparcados a altas temperaturas); el aumento de la temperatura del analizador de gas afectaría a la precisión de las lecturas.

- Cuando se mueva por una zona, proteja el analizador de gas de la luz solar directa y la lluvia intensa.

Utilice siempre el colector de agua. Si rebosa, cambie el filtro y asegúrese de que ningún tubo esté húmedo antes de volver a usarlo.

✍ Nota: Si el escape de un analizador de gas de la serie GA5000 se conecta a un sistema presurizado, el gas se saldrá del orificio de entrada.

Calibrado

 Advertencia	Los gases de calibrado pueden ser peligrosos. Lea atentamente las especificaciones de seguridad de cada gas utilizado antes de proceder al calibrado.
--	---

El regulador suministrado con el juego de calibrado está configurado para proporcionar un caudal fijo.

Dado que el caudal del regulador se ajusta en fábrica, solo hay que girarlo unas vueltas para abrirlo; no es necesario ajustarlo.

 Advertencia	Orificios de escape Durante el calibrado del analizador de gas, existen dos salidas posibles para el gas: por la vía habitual, es decir, el orificio de escape del analizador (amarillo) o, en caso de sobrepresión, el orificio de 1,58 mm de la válvula ^{roja} de alivio de la presión situado en el regulador. Se recomienda conectar tubos de escape en ambos orificios. Los tubos de escape deben llegar a un área con ventilación suficiente. Asegúrese de que no haya fugas en los tubos ni las conexiones. El calibrado del analizador de gas debe realizarse en un área segura con todas las precauciones necesarias en el uso de gases potencialmente peligrosos, explosivos o tóxicos.
--	--

✍ Nota: También es posible que salga gas del orificio de caudal interno (azul) del analizador de gas (solo se aplica al GA5000).

Mantenimiento

Debe realizarse regularmente el mantenimiento de la serie 5000 de analizadores de gas con el fin de garantizar el funcionamiento correcto y preciso. Geotech recomienda que se lleve a cabo el mantenimiento y el recalibrado cada 6 meses.

La serie 5000 de analizadores de gas cuenta con la certificación ATEX para su uso en lugares con riesgo de explosión. Por lo tanto, únicamente los ingenieros cualificados pueden realizar el mantenimiento. En caso contrario, la garantía quedará anulada y podría

invalidar la certificación ATEX.

 Advertencia	Toda operación de mantenimiento del analizador de gas realizada por personal no cualificado puede dar lugar a la invalidación de la certificación ATEX, pudiendo igualmente afectar a la seguridad del uso del instrumento en lugares con riesgo de explosión.
--	---

Mantenimiento realizable por el usuario:

Este instrumento no contiene ninguna pieza o componente interior de cuyo mantenimiento pueda encargarse el usuario.

No obstante, el usuario puede encargarse del mantenimiento de los siguientes elementos:

Filtro de agua en línea	Comprobar periódicamente obstrucciones, humedad o daños y cambiar en caso necesario. No usar el instrumento sin el filtro de agua en línea; de lo contrario, podría penetrar agua en el interior.
Tubo de muestreo	Comprobar que no están contaminados ni dañados.
Conectores de los orificios de gas	Comprobar periódicamente que las juntas tóricas de los conectores de los orificios de gas no están dañadas. Una junta tórica dañada puede dejar pasar el aire en el gas de muestra y afectar a la precisión de las lecturas. En caso de daños en la junta tórica, sustituir todo el conector.
Material filtrante de H ₂ S	Si observa que el material filtrante cambia a un color <i>gris claro</i> , sustituir el filtro.

15.5 Instructions for safe use - Portuguese language

Instruções de Segurança

 Aviso	A série 5000 de analisadores de gás pode ser utilizada para a medição de gases de aterros e de outras fontes, tal como descrito no manual. O operador pode ser exposto a gases prejudiciais durante a utilização do instrumento. A inalação destes gases pode ser prejudicial para a saúde e, em alguns casos pode ser fatal. É de responsabilidade de o usuário assegurar que ele / ela seja devidamente treinado nos aspectos de segurança dos gases que estão sendo utilizados e os procedimentos apropriados sejam seguidos. Em particular, quando gases perigosos serão utilizados, os gases da exaustão do analisador devem ser canalizados para uma área onde é seguro para descarregar o gás. Gases perigosos também podem ser expelidos do instrumento quando purga com ar limpo.
--	---

Nota: Os Analisadores de gases são equipamentos científicos sensíveis, e devem ser tratados como tal. Se o equipamento for utilizado de uma maneira não especificada pelo fabricante, a proteção fornecida pelo instrumento pode ser prejudicada.

Instruções para o uso seguro

Para ATEX e IECEx a série 5000 de analisadores de gases são certificados para Classificação de área perigosa



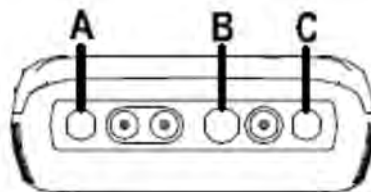
II 2G Ex ib IIA T1 Gb (Ta = +14°F para +122°F)

É vital que as instruções sejam seguidas cuidadosamente. É da responsabilidade do operador determinar o conceito de proteção e de classificação necessários para uma aplicação particular.

(Referência Europeia segundo a diretiva ATEX 94/9/CE, Anexo II, 1.0.6.).

As seguintes instruções se aplicam a equipamentos abrangidos pelos números de certificados SIRA 11ATEX2197X e Diretiva IECEx SIR 11.0089X:

- O equipamento pode ser usado com gases e vapores inflamáveis e com aparelhos grupo IIA e classe de temperatura T1.
- O equipamento está apenas certificado para uso em temperatura ambiente no intervalo (-10 °C a +50 °C) e não deve ser utilizado fora desta faixa.
- O equipamento não deve ser utilizado numa atmosfera de oxigénio superior a 21%.
- A reparação deste equipamento deve ser realizada em conformidade com o código de prática aplicável.
- Quando usado em uma área de risco somente utilizar sonda de temperatura TP-5000 (SIRA 11ATEX2197X e IECEx SIR11. 0089X).
- O analisador não deve ser ligado a outros dispositivos na área de risco, incluindo o cabo 5000-USB (conector A) ou carregador de bateria BC GEM5000 (conector B) fornecido com o analisador.



Não carregue, recarregue ou abra o equipamento em locais explosivos.

Em áreas perigosas somente utilize a "sonda de temperatura TP-5000" no conector B.

Conector A (Uo=10V, Io=5mA, Po=50mW, Ci=0, Li=0, Co=100uF, Lo=1000mH), **Conector B** (Uo=5V, Io=6mA, Po=7mW, Ci=0, Li=0, Co=100uF, Lo=1000mH) **FONTES MÁXIMAS NÃO PERIGOSAS** Conector A – Um=6V Conector B – UM=10.1V.

- Se o equipamento é suscetível a entrar em contato com substâncias nocivas, por exemplo, líquidos ou gases ácidos que podem atacar metais ou solventes e que possam afetar materiais poliméricos, então é de responsabilidade do usuário tomar as precauções adequadas, por exemplo: verificações regulares como parte das inspeções de rotina ou o estabelecimento de folha de dados do material resistente a produtos químicos específicos que impeçam que ele seja prejudicado, garantindo assim que a proteção não é comprometida.
- A faixa de pressão relativa é de + / -200 em H₂O (+ / -500 mbar). Note, no entanto, que a pressão de entrada não deve exceder + / -200, em H₂O (+ / - 500 mbar) em relação à pressão atmosférica e a pressão de saída não deve exceder os 40 em H₂O (+ / - 100 mbar) em relação à pressão atmosférica.

Para CSA (Canadá) a série 5000 de analisadores de gases são certificados para Classificação de área perigosa

CLASSE 2258 03 - EQUIPAMENTOS DE CONTROLE DE PROCESSO - sistemas intrinsecamente seguros e à Prova de Incêndio - Para Áreas Classificadas

Ex ib IIA:

 Modelo GA 5000, GEM5000 e Detector de metano BIOGAS 5000; bateria portátil, alimentado com Bateria não substituível em campo P / N 20087 ou 2011113; intrinsecamente seguras e fornecimento de circuitos de segurança intrínseca ("ib") para Zona 1) de modelo sonda de temperatura TP-5000 (conector B) e com parâmetros de saída como entidade tabulados abaixo; Código de Temperatura T1, 14 ° F ≤ Tamb. ≤ F 122 ° (-10 ° C ≤ Tamb. ≤ 50 ° C).

Conector	Uo (V)	Io (mA)	Po (mW)	Co (uF)	Lo (mH)	Ci (uF)	Li (mH)
B	5.0	6	7	100	1000	0	0
C	10.0	5	50	100	1000	0	0

Nota: Este dispositivo foi testado apenas para recursos de segurança elétrica.

Para CSA (USA) a série 5000 de analisadores de gases são certificados para a classificação de área de perigosa.

CLASSE 2258 83 - PROCESSO DE CONTROLE DE EQUIPAMENTO - Sistemas intrínsecos de segurança, e à Prova de Incêndio - Para Áreas Classificadas como perigosas - certificadas segundo as normas Norte-Americanas.



AEx ib IIA:

Modelo GA 5000, GEM5000 e Detectores de metano BIOGAS 5000; bateria portátil, alimentado com não substituível em campo Bateria P / N 20087 ou 2011113; intrinsecamente seguras e com fornecimento de circuitos de segurança intrínseca ("ib") para Zona 1) de modelo sonda de temperatura TP-5000 (conector B) e com parâmetros de saída como entidade tabulados abaixo; Código de Temperatura T1, 14 ° F < Tamb. > F122 ° F (-10 ° C < Tamb. > 50 ° C)

Conector	Uo (V)	Io (mA)	Po (mW)	Co (uF)	Lo (mH)	Ci (uF)	Li (mH)
B	5.0	6	7	100	1000	0	0
C	10.0	5	50	100	1000	0	0

Nota: Este dispositivo foi testado apenas para recursos de segurança elétrica.

15.6 Instructions for safe use - Chinese language

 Warning 警告	The 5000 series of gas analyzers' can be used for measuring gases from landfill sites and other sources as described in this manual. 5000系列气体分析仪可用来测量本手册中描述的填埋气及其它来源的气体成分。 The operator may be exposed to harmful gases during the use of the instrument. Inhaling these gases may be harmful to health and in some cases may be fatal. 操作员在使用仪器的过程中可能会接触到有毒有害气体。 吸入这些气体可能有害健康，在某些情况下这种损害可能是致命的。 It is the responsibility of the user to ensure that he/she is adequately trained in the safety aspects of the gases being used and appropriate procedures are followed. 操作员的责任是在使用仪器前，确保他 / 她在安全使用气体的安全性方面受过足够的培训，并且操作时按照正确和适当的程序。 In particular, where hazardous gases are being used the gas exhausted from the analyser must be piped to an area where it is safe to discharge the gas. 特别要注意的是，当测量有害气体浓度时，仪器排出的尾气必须通过气管排出到安全的区域。 Hazardous gas can also be expelled from the instrument when purging with clean air. 用泵抽入干净空气可以把仪器内残留的有害气体排走。 The instrument has been designed to be used in explosive atmospheres as defined by the classification. 本仪器设计可在危险区域定义中的易爆环境下使用。 The instrument can be configured to measure low levels of several gases, but may not be certified for use in potentially explosive atmospheres of these gases. 本仪器设计可用于测量低量程的气体浓度，但并没有认证证明可以用于测量潜在爆炸性环境中气体使用。 It is the responsibility of the operator to determine the protection concept and classification of equipment required for a particular application and whether these gases create a potentially explosive atmosphere. 操作员的责任是，确定一个对于特定的应用环境所需的设备保护模式和分类，以及这些气体是否会产生一个潜在的爆炸性气体环境。
--	---

 **Note:** Gas analysers are sensitive pieces of scientific equipment, and should be treated as such. If the equipment is used in a manner not specified by the manufacturer, the protection provided by the instrument may be impaired.

注意：气体分析仪是敏感性的科学器材，如果不按照制造商制定的方式使用仪器，可能会给仪器带来损害。

仪器安全使用指南

For ATEX and IECEx the 5000 series of gas analysers are certified to Hazardous Area Classification

5000系列气体分析仪已被ATEX及IECEx认证可用于危险性的爆炸区域。



II 2G Ex ib IIA T1 Gb (Ta = -10°C to +50°C)

It is vital instructions are followed closely. It is the responsibility of the operator to determine the protection concept and classification required for a particular application.

遵守操作要求非常重要。为特定的应用环境选定不同的保护模式及分类是操作员的责任。

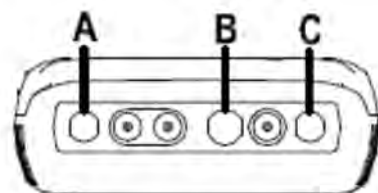
(Reference European ATEX Directive 94/9/EC, Annex II, 1.0.6.)

(参考欧洲ATEX指令94/9/EC, 附件II, 1.0.6)

The following instructions apply to equipment covered by certificate numbers SIRA 11ATEX2197X and IECEx Directive SIR 11.0089X:

以下的说明适用于已取得SIRA 11ATEX2197X 认证号码及IECEx Directive SIR 11.0089X认证号码的设备。

- The equipment may be used with flammable gases and vapours with apparatus group IIA and temperature class T1.
- 本设备可用于IIA级易燃气体和蒸气领域，以及T1级温度分类级别。
- The equipment can contain gas sensing heads for the detection of particular gases. The inclusion of a sensor does not infer that the equipment is suitable for the use of gases with a temperature class of less than T1.
- 本设备可装配检测特种气体的传感器探头。所装配的气体传感器探头并不意味着该设备适用于小于T1温度级别的气体检测。
- The equipment is only certified for use in ambient temperatures in the range -10°C to +50°C and should not be used outside this range.
- 本设备只适用于-10°C至+50°C的环境温度下使用，不要在这个温度范围之外环境下使用。
- The equipment must not be used in an atmosphere of greater than 21% oxygen.
- 本设备不适用于氧气浓度大于21%环境下使用。
- Repair of this equipment shall be carried out in accordance with the applicable code of practice.
- 本设备的维修维护需要按照适用的使用手册要求。
- When used in a hazardous area only use GF5.2 temperature probe (SIRA 11ATEX2197X and IECEx SIR11.0089X). For connector C, the GF5.4 anemometer (BVS 04ATEXE194) for use with ATEX only. The analyser should not be connected to any other devices in the hazardous area including the GF-USB lead (connector A) or GF3.9 battery charger (connector B) supplied with the analyser.
- 在危险区域能使用GF5.2型号的温度探头 (SIRA 11ATEX2197X 认证号码和 IECEx SIR11.0089X认证号码)。GF5.4型号风速仪(BVS 04ATEXE194认证号码)，只适用于ATEX认证区域使用。除上述配件外，本设备不应在危险区域下连接其它外部装置，包括制造商提供的USB数据连接线及充电器。



**Do not charge, recharge or open in a potentially explosive atmosphere.
In hazardous area only use "Temperature Probe GF5.2" in Connector B.
Connector C ($U_o=10V, I_o=5mA, P_o=50mW, C_i=0, L_i=0, C_o=100\mu F, L_o=1000mH$),
Connector B ($U_o=5V, I_o=6mA, P_o=7mW, C_i=0, L_i=0, C_o=100\mu F, L_o=1000mH$)**

**MAXIMUM NON-HAZARDOUS SUPPLIES:
Connector A - $U_m=6V$ Connector B - $U_m=10.1V$**

不要在易爆环境下充电或打开设备。

在危险区域只能使用连接在B接头的GF5.2温度探头。

C连接头($U_o=10V, I_o=5mA, P_o=50mW, C_i=0, L_i=0, C_o=100\mu F, L_o=1000mH$)

B连接头($U_o=5V, I_o=6mA, P_o=7mW, C_i=0, L_i=0, C_o=100\mu F, L_o=1000mH$)

最大安全性连接

A连接头 $U_m=6V$ B连接头 $U_m=10.1V$

- The safe area apparatus that is to be connected to the USB Port shall be a Safety Extra Low Voltage (SELV) or Protective Extra Low Voltage (PELV) circuit.
- 设备如需要连接USB线，需要在安全低电压或保护低电压的安全区下连接。
- Only a Geotechnical Instrument battery pack part number 20087 is permitted as a replacement. This battery pack shall only be changed in a safe area.
- 只有Geotechnical公司所配置电池组（编号20087）允许用于替换设备电池组。电池组的更换需要在安全区域下进行。
- Only Battery Charger type GF3.9 shall be used to recharge the batteries via Connector 'B'.
- 只允许使用GF3.9型号充电器通过B连接头给设备充电。
- If the equipment is likely to come into contact with aggressive substances, e.g. acidic liquids or gases that may attack metals, or solvents that may affect polymeric materials, then it is the responsibility of the user to take suitable precautions, e.g. regular checks as part of routine inspections or establishing from the material's data sheet that it is resistant to specific chemicals that prevent it from being adversely affected, thus ensuring that the type of protection is not compromised.
- 如果设备有可能接触到腐蚀性物质，例如酸性液体或气体，可能会对设备金属有腐蚀影响，或溶剂可能会影响到设备的高分子材料。用户有责任采取适当的预防措施，例如把定期检查作为常规检查的一部分，或者建立材料数据表，注明能对某些特殊化学物质有防腐功能，以防受到不利影响。
- The relative pressure range is +/-500 mbar. Note, however, that the input pressure should not exceed +/- 500 mbar relative to atmospheric pressure and the output pressure should not exceed +/- 100 mbar relative to atmospheric pressure.
- 压差范围为+/-500mbar。注意：输入压力不能超过大气压力的+/-500mbar范围，输出压力不能超过大气压力的+/-100mbar范围。

For CSA (Canada) the 5000 series of gas analysers are certified to Hazardous Area Classification

CLASS 2258 03 - PROCESS CONTROL EQUIPMENT - Intrinsically Safe and Non-Incendive Systems - For Hazardous Locations

CSA（加拿大）5000系列气体分析仪的危险区域等级认证：

CLASS225803 - 过程控制设备 - 本质安全和非易燃系统 - 用于危险场所



Ex ib IIA:

Model GA 5000, GEM 5000 and BIOGAS 5000 Methane Detectors; portable, battery powered with non-field-replaceable Battery Pack P/N 20087; intrinsically safe and providing intrinsically safe circuits ("ib" for Zone 1) to Model GF5.2 Temperature Probe (Connector B) and with entity output parameters as tabulated below; Temperature Code T1; $-10^{\circ}\text{C} \leq T_{\text{amb}} \leq +50^{\circ}\text{C}$.

型号GA5000, GEM5000和Biogas5000沼气分析仪；便携式，内置电池组（P/N 20087）；本质安全，并提供本安电路（"ib" 1区）GF5.2型号温度探头（B连接头），实际输出参数如下表；温度代码T1； $-10^{\circ}\text{C} \leq \text{温度} \leq +50^{\circ}\text{C}$ 。

Connector 连接头	Entity Parameters 实际参数						
	Uo (V)	Io (mA)	Po (mW)	Co (uF)	Lo (mH)	Ci (uF)	Li (mH)
B	5.0	6	7	100	1000	0	0
C	10.0	5	50	100	1000	0	0



Note: This device has been investigated for electrical safety features only.

注意：本设备已通过电气安全功能性检查。

or CSA (USA) the 5000 series of gas analysers are certified to Hazardous Area Classification

或CSA（USA）5000系列气体分析仪的危险区域等级认证。

CLASS 2258 83 - PROCESS CONTROL EQUIPMENT - Intrinsically Safe and Non-Incendive Systems - For hazardous Locations - CERTIFIED TO U.S. STANDARDS

CLASS 2258 83—过程控制设备 - 本质安全和非易燃系统 - 用于危险场所 - 美国标准认证

15.7 Instructions for safe use - Polish language

Instrukcje bezpieczeństwa

 Ostrzeżenie	Analizatory serii 5000 mogą być używane wyłącznie do pomiarów gazów na składowiskach oraz w innych miejscach, opisanych w instrukcji. Operator może mieć kontakt ze szkodliwymi gazami. Wdychanie tych gazów może nieść ze sobą poważne skutki dla zdrowia. Operator odpowiada za dostęp do analizatora wyłącznie dla osób posiadających odpowiednie kwalifikacji i wiedzę, szczególnie w przypadku kontaktu z gazami szkodliwymi. Miejsce wylotu gazów z analizatora powinno być odpowiednio wentylowane. Szkodliwe gazy mogą wydostać się z analizatora również podczas przepłukiwania czystym powietrzem. Urządzenie może być użytkowane w odpowiedniej strefie zagrożenia wybuchem. Urządzenie jest dostosowane do pomiarów niektórych gazów wybuchowych tylko na niskich stężeniach. Operator odpowiada za określenie i zapewnienie środków ochrony osobistej podczas pracy w strefie zagrożenia wybuchem.
---	--

✍ Uwaga: Analizatory gazu posiadają delikatne element, które należy odpowiednio traktować. Stosowanie urządzenie niezgodnie z przeznaczeniem powoduje powstawanie ryzyka innych niebezpieczeństw.

Instrukcje dot bezpieczeństwa

Analizatory posiadają poniższy certyfikat ATEX.



II 2G Ex ib IIA T1 Gb (Ta = -10°C to +50°C)

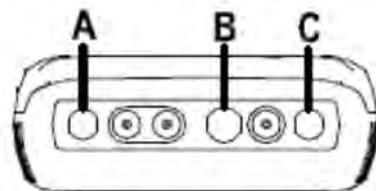
Poniższych instrukcje należy bezwzględnie przestrzegać. Do operatora należy określenie dodatkowych środków bezpieczeństwa w przypadku specyficznych zastosowań urządzenia.

(Europejska dyrektywa ATEX - 94/9/EC, Annex II, 1.0.6.)

Poniższe instrukcje dotyczą urządzeń z certyfikatami SIRA 11ATEX2197X oraz IECEx Directive SIR 11.0089X:

- Palne gazy oraz ich pary mogą mierzyć wyłącznie urządzenia z grupy IIA oraz klasy temperaturowej T1.
- Urządzeniem nie można mierzyć gazów o klasie temperaturowej gorszej niż T1.
- Bezpieczny zakres temperatury pracy urządzenia to -10°C do +50°C.
- Urządzenie nie może być użytkowane w atmosferze o większej zawartości tlenu niż 21%.
- Serwis urządzenia powinien być przeprowadzony zgodnie z wytycznymi producenta oraz najlepszą praktyką.

- Do pomiarów temperatury w strefie niebezpiecznej służy wyłącznie termopara o nr katalogowym GF5.2 (SIRA 11ATEX2197X and IECEx SIR11.0089X). Do pomiarów przepływu w strefie niebezpiecznej służy wyłącznie anemometr o nr katalogowym GF5.4 (BVS 04ATEXE194). Nie należy podłączać innych urządzeń w strefie niebezpiecznej, w tym używać portów A (port USB) oraz B do ładowania ładowarką GF3.9.



Nie ładować ani nie otwierać urządzenia w strefie wybuchowej.
W strefie wybuchowej używać wyłącznie portu B do pomiarów temperatury.
Port C (Uo=10V, Io=5mA, Po=50mW, Ci=0, Li=0, Co=100uF, Lo=1000mH),
Port B (Uo=5V, Io=6mA, Po=7mW, Ci=0, Li=0, Co=100uF, Lo=1000mH)

Maks. Napięcie dla portów:
Port A - Um=6V Port B - Um=10.1V

- Strefy w których można skorzystać z portu USB to Safety Extra Low Voltage (SELV) oraz Protective Extra Low Voltage (PELV).
- Należy używać wyłącznie oryginalnych baterii od producenta.
- Należy używać wyłącznie oryginalnej ładowarki od producenta, podłączając do portu B.
- Jeżeli analizator może mieć kontakt z substancjami agresywnymi, tj. Kwasy, ciecze korodujące, rozpuszczalniki polimerów, do użytkownika należy podjąć odpowiednie środki bezpieczeństwa.
- Ciśnienie względne w układzie wynosi +/- 500 mbar. Ciśnienie na porcie wylotowym nie powinno przekraczać +/- 100 mbar.

Kanada: Analizatory serii 5000 posiadają certyfikat dla poniższej klasy zagrożenia eksplozja.

Class 2258 03 – WYPOSAŻENIE KONTROLNO-POMIAROWE – samoistnie bezpieczny w strefie zagrożenia



Ex ib IIA:

Modele Biogas 5000, GA 5000, GEM 5000: przenośny, zasilany baterią niewymienialną w strefie zagrożenia (Battery Pack P/N 20087 lub 2011113), samoistnie bezpieczny i kompatybilny do termopary (Model GF 5.2). Klasa temperaturowa: T1; $-10^{\circ}\text{C} \leq T_{\text{amb}} \leq +50^{\circ}\text{C}$. Parametry dla konkretnych gniazd w tabeli poniżej.

Gniazdo	Nominalne parametry						
	Uo (V)	Io (mA)	Po (mW)	Co (uF)	Lo (mH)	Ci (uF)	Li (mH)
B	5.0	6	7	100	1000	0	0
C	10.0	5	50	100	1000	0	0

 Uwaga: Urządzenie zostało sprawdzone pod kątem bezpieczeństwa elektrycznego.

USA: Analizatory serii 5000 posiadają certyfikat dla poniższej klasy zagrożenia eksplozją.

CLASS 2258 83 – WYPOSAŻENIE KONTROLNO-POMIAROWE - samoistnie bezpieczny w strefie zagrożenia - certyfikowany wg amerykańskich norm.



AEx ib IIA:

Modele Biogas 5000, GA 5000, GEM 5000: przenośny, zasilany baterią niewymienialną w strefie zagrożenia (Battery Pack P/N 20087 lub 2011113), samoistnie bezpieczny i kompatybilny do termopary (Model GF 5.2). Klasa temperaturowa: T1; $-10\text{ °C} \leq T_{amb.} \leq +50\text{ °C}$. Parametry dla konkretnych gniazd w tabeli poniżej.

Gniazdo	Nominalne parametry						
	Uo (V)	Io (mA)	Po (mW)	Co (uF)	Lo (mH)	Ci (uF)	Li (mH)
B	5.0	6	7	100	1000	0	0
C	10.0	5	50	100	1000	0	0

 Uwaga: Urządzenie zostało sprawdzone pod kątem bezpieczeństwa elektrycznego.

15.8 Instructions for safe use - Romanian language**Instructiuni de protectie**

 Atentionare	Analizoarele de gaz din seria 5000 pot fi utilizate pentru masurarea gazelor din instalatiile din gropile de gunoi si alte surse precum cele descrise in acest manual. Operatorul poate fi expus gazelor nocive in timpul utilizarii instrumentului. Inhalarea acestor gaze poate fi periculoasa pentru sanatate si in unele cazuri poate fi fatala. Este responsabilitatea utilizatorului sa se asigure ca el/ea este instruit(a) corespunzator in privinta gazelor utilizate si ca procedurile adecvate sunt urmate. In cazuri particulare, cand sunt utilizate gaze periculoase, gazul evacuat din analizor trebuie eliminat intr-o zona unde este sigura eliberarea gazului. Gazul periculos poate fi de asemenea evacuat din instrument atunci cand se face purjarea cu aer curat. Instrumentul a fost proiectat pentru utilizarea in atmosfere explozive dupa cum este definit de clasificare. Instrumentul poate fi configurat sa masoare nivelurile joase ale catorva gaze, dar este posibil sa nu fie certificat pentru utilizarea in atmosferele explozive a acestor gaze. Este responsabilitatea utilizatorului sa determine conceptul de protectie si clasificarea echipamentului necesara pentru o anumita aplicatie si daca aceste gaze creeaza o potentiala atmosfera exploziva.
--	--

⚠ Nota: Analizoarele de gaz sunt componente sensibile ale echipamentului stiintific si trebuie tratate ca atare. Daca echipamentul este utilizat intr-o maniera nespecificata de producator, protectia oferita de instrument poate fi afectata.

Instructiuni pentru utilizarea in siguranta

Pentru ATEX si IECEx, analizoarele de gaz din seria 5000 sunt certificate Clasificarii de Zona Periculoasa



II 2G Ex ib IIA T1 Gb (Ta = -10°C to +50°C)

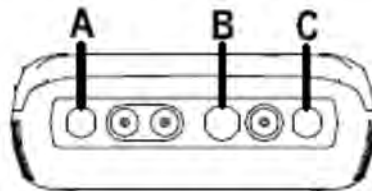
Este vital ca instructiunile sa fie urmate corespunzator. Este responsabilitatea operatorului sa determine conceptul de protectie si clasificarea necesara pentru o anumita aplicatie.

(Referinta Directiva Europeana ATEX 94/9/EC, Anexa II, 1.0.6.)

Urmatoarele instructiuni se aplica echipamentului cu numerele de certificat SIRA 11ATEX2197X si Directiva IECEx SIR 11.0089X:

- Echipamentul poate fi utilizat cu gaze explozive si vapori cu aparate grupa IIA si clasa de temperatura T1.
- Echipamentul poate contine senzori de gaz pentru detectarea anumitor gaze. Includerea unui senzor nu presupune ca echipamentul este adecvat pentru utilizarea gazelor cu o clasa de temperatura mai mica de T1.
- Echipamentul este certificat doar pentru utilizarea in temperaturi ambientale in intervalul de la -10°C la +50°C si nu ar trebui utilizat in afara acestui interval.
- Echipamentul nu trebuie utilizat in atmosfera cu mai mult de 21% oxygen.
- Repararea acestui echipament se realizeaza conform codului de practica in vigoare.

- Cand aparatul este folosit intr-o zona cu risc de explozive, utilizati doar sonda de temperatura GF5.2 (SIRA 11ATEX2197X si IECEx SIR11.0089X). Pentru conectorul C, se foloseste anemometrul GF5.4 (BVS 04ATEXE194) doar in zona ATEX. Analizorul nu trebuie conectat unui alt dispozitiv in zona exploziva inclusiv cablul GF-USB (conector A) sau incarcatorul GF3.9 (conector B) furnizat cu echipamentul.



Nu incarcati, reincarcati sau deschideti intr-o atmosfera cu potential exploziv.
In zona exploziva utilizati doar "Sonda de temperatura GF5.2" in Conector B.
Connector C (Uo=10V,Io=5mA,Po=50mW,Ci=0,Li=0,Co=100uF,Lo=1000mH),
Connector B (Uo=5V,Io=6mA,Po=7mW,Ci=0,Li=0,Co=100uF,Lo=1000mH)

ALIMENTARI MAXIME NEPERICULOASE:
Connector A - Um=6V Connector B - Um=10.1V

- Echipamentul pentru zone sigure care urmeaza a fi conectat la portul USB va fi cu circuit Safety Extra Low Voltage (SELV) sau Protective Extra Low Voltage (PELV).
- Se va utiliza doar bloc de acumulatori Geotechnical Instrument cu numar de serie 20087 sau 2011113 ca piese de schimb. Acest acumulator se va schimba doar in zona sigura.
- Doar incarcatorul tip GF3.9 se va utiliza pentru reincarcarea bateriilor prin Conectorul 'B'.
- Daca exista riscul ca echipamentul sa intre in contact cu substante agresive, ex. lichide acide sau gaze ce pot ataca metale, sau solventi ce pot afecta materiale polimerice, atunci este responsabilitatea utilizatorului sa ia masuri adecvate, cum a fi verificari regulate ca parte a inspectiei de rutina sau sa stabileasca din fisa de securitate a materialului ca este rezistent la chimicale specifice ce il impiedica sa fie afectat negativ, asigurandu-se astfel ca tipul de protectie nu este compromis.
- Intervalul presiunii relative este +/-500 mbar. Retineti, totusi, ca presiunea de intrare pentru GA5000 nu trebuie sa depaseasca +/- 500 mbar in raport cu presiunea atmosferica, si presiunea de intrare pentru GEM5000 si BIOGAS 5000 nu trebuie sa depaseasca +/- 250 mbar in raport cu presiunea atmosferica, presiunea de iesire nu trebuie sa depaseasca +/- 100 mbar in raport cu presiunea atmosferica.

Pentru CSA (Canada) analizoarele din seria 5000 sunt certificate Clasificarii de Zona Periculoasa

CLASS 2258 03 – ECHIPAMENTE DE CONTROL PROCES - Siguranta Intrinseca si Sisteme Antideflagrante – Pentru Locatii Explozive



Ex ib IIA:

Modele Detectoare Metan GA 5000, GEM 5000 si BIOGAS 5000; portabile, alimentate cu baterii cu bloc de acumulatori care nu pot fi inlocuiti in teren P/N 20087 sau 2011113; siguranta intrinseca si furnizare de circuite cu siguranta intrinseca ("ib") pentru Zona 1) pentru Modelul de Sonda de Temperatura GF5.2 (Conector B) si cu parametri de iesire in conformitate cu tabelul de mai jos; Cod Temperatura T1; -10 °C ≤ Tamb. ≤ +50°C.

Conector	Parametri						
	Uo (V)	Io (mA)	Po (mW)	Co (uF)	Lo (mH)	Ci (uF)	Li (mH)
B	5.0	6	7	100	1000	0	0
C	10.0	5	50	100	1000	0	0



Nota: Acest dispozitiv a fost verificat doar pentru caracteristici de siguranta electrica.

Pentru CSA (SUA) analizoarele din seria 5000 sunt certificate Clasificarii de Zona Periculoasa

CLASS 2258 83 - ECHIPAMENTE DE CONTROL PROCES - Siguranta Intrinseca si Sisteme Antideflagrante – Pentru Locatii Explozive – CERTIFICAT STANDARDELOR SUA



AEx ib IIA:

Modele Detectoare Metan GA 5000, GEM 5000 si BIOGAS 5000; portabile, alimentate cu baterii cu bloc de acumulatori care nu pot fi inlocuiti in teren P/N 20087 sau 2011113; siguranta intrinseca si furnizare de circuite cu siguranta intrinseca ("ib") pentru Zona 1) pentru Modelul de Sonda de Temperatura GF5.2 (Conector B) si cu parametri de iesire in conformitate cu tabelul de mai jos; Cod Temperatura T1; -10 °C Tamb. +50°C.

Conector	Parametrii Entitatii						
	Uo (V)	Io (mA)	Po (mW)	Co (uF)	Lo (mH)	Ci (uF)	Li (mH)
B	5.0	6	7	100	1000	0	0
C	10.0	5	50	100	1000	0	0



Nota: Acest dispozitiv a fost investigat doar pentru caracteristici de siguranta electrica.

17.0 Glossary of terms

5000 series	The 5000 series refers to the GA5000, GEM5000 and the Biogas5000 gas analysers.
Accu-Flo	A flow device used to aid accurate flow measurement.
Analyser error messages	Operational errors are prefixed on the analyser by the word ERROR followed by an error code. Refer to the list of standard error codes for more information.
Analyser warning	Analyser warnings are prefixed by the word WARNING followed by a relevant description. There are two types of warning messages displayed; general warnings that may not necessarily affect the instrument's function (for example, battery power low) and operational parameters that could affect the performance of the analyser (for example, CH₄ out of calibration).
Anemometer probe	Device for measuring velocity of gas in the pipe. The GEM5000 analyser can be set to convert into a flow. See also flow measurement.
ATEX certification	The GEM5000 is ATEX certified to zone 1 & 2 areas above ground not in mines.
Auxiliary channel	This refers to the channels where external devices will be connected or displayed.
Backlight	The analyser has a built-in backlight for low ambient light conditions. This can be toggled on/off using the backlight key.
Barometric pressure	The atmospheric pressure at the given location.
Borehole	Typical location from which a gas sample is obtained.
Calibration	The gas analyser is carefully calibrated against known standards.
Calibration record	The GEM5000 instrument has the facility to log user calibrations as a validation tool.
CH₄	Methane
Chemical cells	A method of gas detection that works on the basis of a chemical reaction with the target gas.
Clean air purge	Process used to clear out gas from the sample tube and analyser prior to taking a new

	reading.
CO	Carbon monoxide
CO₂	Carbon dioxide
Data logging	Data logging enables the operator to leave the analyser unattended to take samples at predetermined intervals for a set period of time
Download	Terminology used for the movement of data from the analyser to the GAM application on the PC.
Dual beam infrared absorption	Method of gas detection by measuring how much infrared is absorbed by the target gas.
Event log	Used as an aid to monitoring the use of the analyser. It can also be used as a diagnostic tool. The event log can be viewed via Gas Analyser Manager. It <u>cannot</u> be viewed on the analyser screen.
Exhaust port	The usual manner for the gas to exit the analyser is via the exhaust port located on the top side of the analyser. This port should have an exhaust tube attached.
Exhaust tube	Clear plastic tubing used to expel gases from the exhaust port.
Factory settings	Default settings preset at time of manufacture or service.
Firmware	Firmware is the term by which the internal analyser software is known and is not accessible by the client. This firmware is updated to the latest version when the analyser is returned in UK for servicing.
Flow measurement	Flow can be measured by either gas velocity m/s or volume flow rate m ³ /hr. This measurement of flow relates to the use of the anemometer and not the internal flow measurement technique.
Flow port	For the measurement of gas flow at the sample point.
Gas Analyser Manager	Also referred to as GAM. PC based software which enables the operator to upload and download information to/from the analyser. Gas Analyser Manager enables operators to maximise the operation of their gas analyser. It features a simple upload and download facility and is fully compatible with the latest

	Microsoft™ operating systems.
	This is optional.
Gas channels	The gases that are analysed by the instrument.
Gas velocity	The positional rate of change of the gas. Measured using the optional anemometer.
General warnings	Displayed throughout the documentation with a warning symbol. Warning information may affect the safety of operators.
H₂	Hydrogen
H₂S	Hydrogen sulphide
H₂S filter	Filter required for removal of H ₂ S. When the filter material changes colour to a light grey colour or if H ₂ S values are displayed, then the filter should be replaced.
Hydro-carbons	Organic compound consisting of only hydrogen and carbon.
In-line water filter	The component used to help protect the instrument from water ingress.
LCD display	Liquid Crystal Display
LEL	Lower Explosive Limit. Lower explosive limit of methane in air. 5% methane in air is the point at which it becomes explosive. 100% LEL equates to 5% methane.
m/s	Meters per second – measurement of gas velocity.
m³/hr	Meters cubed per hour – volumetric flow rate measurement.
Main Gas Read Screen	The main analyser screen for normal operations and all operations are carried out from this screen.
Material data sheet	Document from which information about a certain substance can be obtained.
MCERTS certification	MCERTS is the UK Environment Agency's Monitoring Certification Scheme. The scheme provides a framework within which environmental measurements can be made in accordance with the Agency's quality requirements. The scheme covers a range of monitoring, sampling and inspection activities.
Memory	Location where data and ID information is stored. The analyser memory should not be

	used as a permanent storage medium. Stored data should be regularly transferred using the GAM download software.
NO₂	Nitrogen dioxide
Operating language	The operator can choose the default operating language for the analyser. Choices are English, German, Spanish, French and Italian.
PPM	Parts per million
Pump	Used to draw the gas sample from the sample point to the analyser. Select the pump key  on the analyser to activate.
ID	The user definable identification tag allocated to a sample point.
Relative pressure	The pressure at the sample point 'relative' to atmospheric (barometric) pressure.
Relative pressure transducer	The internal component used to measure the relative pressure.
Residual N₂	The calculation for the residual N ₂ used on the latest version of the GEM5000 platform is as follows: Residual N ₂ = Balance – (O ₂ % x 3.76) Where, Balance = 100% - (CH ₄ % + CO ₂ % + O ₂ %) and 3.76 is the ratio of O ₂ to N ₂ in ambient air (79/21)
Sample tube	The tube used to obtain a sample of gas from the sample point to the analyser.
Span	The point at which the gas analyser is calibrated when a known quantity of the target gas is present.
Span multi gas	Term by which the span calibration of the three main gas channels is known. This option must only be used when the calibration gas being used is a combination of CH ₄ CO ₂ O ₂ .
Technician ID	An alpha-numeric code tagged to each gas reading. Facility only available via Gas Analyser Manager (GAM). This is an optional feature.
Temperature probe	External device used to measure the gas temperature at the sample point. This is optional.
Update site data	Enables the operator to answer pre-defined questions relating to the site, environment etc.

	These questions are defined via 5GAM software.
Upload	Terminology used for the movement of data from the PC via GAM software application to the analyser.
Volume flow rate	The volume of a gas that passes through a given surface per unit of time e.g. m ³ /hr
Warm-up self-test	Pre-determined self-test sequence to test the analyser functions which takes place after the analyser is switched on.
Warranty	The instrument is under guarantee against defect in materials and workmanship for a period of 3 years from the date of shipment to the operator and is subject to the recommended service and recalibration requirements.
Water trap	Device used to protect the instrument from water or moisture ingress.
Zero	The point at which the gas analyser is calibrated when there is none of the target gas present.
Zero transducers	This option allows the relative pressure transducer to be zeroed.

Authorised distributor

In Australia:

For customer service, call 1300-735-292

To email an order, ordersau@thermofisher.comTo order online: thermofisher.com**In New Zealand:**

For customer service, call 0800-933-966

To email an order, ordersnz@thermofisher.comTo order online: thermofisher.com**ThermoFisher**
S C I E N T I F I C