

Ion GeneStudio™ S5 Instrument

USER GUIDE

for use with:

Ion GeneStudio™ S5 Instrument

Ion GeneStudio™ S5 Plus Instrument

Ion GeneStudio™ S5 Prime Instrument

Catalog Numbers A38194, A38195, A38196

Publication Number MAN0017528

Revision E



Revision history: MAN0017528 E (English)

Revision	Date	Description
E	5 December 2025	Updated initialization troubleshooting.
D.0	23 October 2025	- Added recommendation for preventive maintenance. - Added symbols statement to copyright page.
C.0	1 February 2023	Removed the following skus from the Ion S5™ Installation Kit contents table. - A28195 - A27897 - A29022
B.0	22 February 2022	Update the quantity of the Ion S5™ Cartridge Tool in the Ion S5™ Installation Kit.
A.0	5 January 2018	New instrument user guide that includes instructions for Ion GeneStudio™ S5 Series instrument operation and maintenance.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF THE DOCUMENT OR THE PRODUCT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

Trademarks: All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. MAXYMum Recovery is a trademark of Axygen, Inc. Tygon is a trademark of Saint-Gobain Performance Plastics Corp.

©2018-2025 Thermo Fisher Scientific Inc. All rights reserved.

Contents

About this guide	6
Purpose of the guide	6
■ CHAPTER 1 Product information	7
Product description	7
System components	7
Required materials and equipment not supplied	8
Instrument installation by only trained personnel	10
Nucleic acid contamination	10
Instrument vibration and clearances	10
■ CHAPTER 2 System components	11
Ion GeneStudio™ S5 System component configuration and layout	11
Internet connectivity	13
Ion GeneStudio™ S5 System component positions	14
Ion GeneStudio™ S5 Sequencer input and output connections	15
■ CHAPTER 3 Instrument operation	16
Precautions	16
Avoid nucleic acid contamination	16
Avoid strong electromagnetic radiation	16
Protection by equipment	16
Power the Ion GeneStudio™ S5 Sequencer on or off	16
Power on	16
Power off	17
Update the Ion GeneStudio™ S5 System software	17
Ion GeneStudio™ S5 instrument run times	18
Maintain the sequencer	19
Materials required	19
Clean or decontaminate the sequencer	19
Perform the sequencer cleaning manually	19

Perform an instrument reset run with an initialized sequencer that is loaded with an unused Reagents cartridge	21
Instrument planned maintenance	21
■ APPENDIX A Touchscreen reference	22
Clean, Initialize, and Run	22
Settings	23
Network Settings	23
System Tools	26
Check for software updates	29
Data management	30
Manually delete run data	31
Perform the instrument Clean operation	32
Instrument settings	33
Touchscreen icons	35
Alarms, Notifications, and Events	37
■ APPENDIX B Troubleshooting	40
Instrument alarms and events	40
Initialization fails	41
Troubleshooting using control libraries	43
Troubleshooting using Control Ion Sphere™ Particles	43
Solid-state buffer leak	44
Pulse clean the sequencer	44
■ APPENDIX C Supplemental procedures	46
Set up and test the Ion Chip™ Minifuge	46
Install the Ion S5™ /Ion Proton™ Rotor and Buckets	46
Test the minifuge	48
■ APPENDIX D Instrument warranty	50
■ APPENDIX E Safety	51
Safety alerts on this instrument	51
Location of safety labels on this instrument	52
Safety information for instruments not manufactured by Thermo Fisher Scientific	52
Instrument safety	53
General	53
Cleaning and decontamination	53
Electrical safety	54

Safety and electromagnetic compatibility (EMC) standards	55
Safety	55
EMC	55
Environmental design	56
Chemical safety	57
Documentation and support	59
Related documentation	59
Customer and technical support	59
Limited product warranty	60

About this guide



CAUTION! ABBREVIATED SAFETY ALERTS. Hazard symbols and hazard types specified in procedures may be abbreviated in this document. For the complete safety information, see the “Safety” appendix in this document.

IMPORTANT! Before using this product, read and understand the information in the “Safety” appendix in this document.

Purpose of the guide

This instrument user guide provides reference information for using and maintaining three Ion GeneStudio™ S5 series sequencers.

- Ion GeneStudio™ S5 System (Cat. No. [A38194](#))
- Ion GeneStudio™ S5 Plus System (Cat. No. [A38195](#))
- Ion GeneStudio™ S5 Prime System (Cat. No. [A38196](#))

Note: In this user guide, Ion GeneStudio™ S5 Sequencer or Ion GeneStudio™ S5 System refers generically to the three Ion GeneStudio™ S5 sequencers or systems, unless otherwise specified.

The discontinued Ion S5™ and Ion S5™ XL instruments continue to be supported by Thermo Fisher Scientific. See the *Ion S5™ and Ion S5™ XL Sequencer User Guide* (Pub. No. [MAN0010811](#)) for reference information for the Ion S5™ and Ion S5™ XL Instruments.



Product information

Product description

The Ion GeneStudio™ S5 series of sequencers are semiconductor-based next-generation sequencing (NGS) systems that enable simple targeted sequencing workflows. With cartridge-based reagents, the systems are simple to use and offers scalability and flexibility. They use Ion 5 Series chips to support a broad range of high-throughput sequencing for clinical research and research applications from microbial genomes and gene panels to exomes and transcriptomes. They share the same foundational technology and fluidic architecture across all three instruments in the series.

- Ion GeneStudio™ S5 System (Cat. No. [A38194](#))—This system is optimized to provide affordability for laboratories focused on smaller panels and lower sequencing throughput requirements.
- Ion GeneStudio™ S5 Plus System (Cat. No. [A38195](#))—From genes to exomes, this system is designed to enable flexibility to help laboratories manage small and large projects across multiple applications.
- Ion GeneStudio™ S5 Prime System (Cat. No. [A38196](#))—For core and service laboratories requiring maximum throughput and rapid turnaround time from a benchtop NGS system, this system enables outstanding breadth and depth, from large panels or exomes to clinical oncology research focused on low allele frequency applications.

System components

The Ion GeneStudio™ S5 series of sequencers includes the following system components.

Ion GeneStudio™ S5 System (Cat. No. [A38194](#))

Components	Part No.
Ion GeneStudio™ S5 Sequencer	A37904
Ion S5™ Installation Kit	A27215

Ion GeneStudio™ S5 Plus System (Cat. No. [A38195](#))

Components	Part No.
Ion GeneStudio™ S5 Plus Sequencer	A37259
Ion S5™ Installation Kit	A27215

Ion GeneStudio™ S5 Prime System (Cat. No. [A38196](#))

Components	Part No.
Ion GeneStudio™ S5 Prime Sequencer	A37260
Ion Torrent™ Server	A36305
Ion S5™ Installation Kit	A27215

Ion S5™ Installation Kit (Part No. A27215; Not available for separate purchase.)

Contents	Part No.	Quantity	Storage
Ion 540™ Chip Kit	A27765	4 pack	15°C to 30°C
Ion S5™ Sequencing Solutions	A27767	1	
Ion S5™ Sequencing Reagents	A27768	1	-30°C to -10°C
Ion S5™ Cartridge Tool	A28308	1	15°C to 30°C

Required materials and equipment not supplied

Unless otherwise indicated, all materials are available through [thermofisher.com](https://www.thermofisher.com). "MLS" indicates that the material is available from [fisherscientific.com](https://www.fisherscientific.com) or another major laboratory supplier.

IMPORTANT! Thermo Fisher Scientific has verified the procedures in this user guide using these specific materials. Substitution can adversely affect system performance.

✓	Item	Source
	Ion Chip™ Minifuge (120 V or 230 V; required for use only with the Ion OneTouch™ 2 System)	4479672 (120V) or 4479673 (230V)
	Minifuge Proton Rotor and Bucket Kit (required for use only with the Ion OneTouch™ 2 System)	4482578
	(Optional) Uninterruptible Power Supply (UPS) ^[1]	MLS
	Thermal cycler with a heated lid	MLS
	Microcentrifuge ^[2]	MLS
	1.5-mL or 1.7-mL microcentrifuge tubes	MLS
	0.2-mL MAXYMum Recovery™ Thin Wall PCR Tubes, Flat Cap (do not use polystyrene tubes)	PCR-02-L-C Axygen™
	Pipettes (P2, P10, P20, P200, P1000) and appropriate low-retention filtered tips	MLS
	Isopropanol (100%)	MLS

(continued)

✓	Item	Source
	Nuclease-free water molecular biology grade	MLS
	Standard laboratory vacuum line or vacuum pump	MLS
	Liquid trap	MLS
	Tygon™ tubing ^[3]	MLS
	Vortex mixer	MLS

^[1] For laboratories that experience frequent power outages or line voltage fluctuations, we recommend that you use an uninterruptible power supply that is compatible with 2500 W output or higher.

^[2] Must fit standard 1.5- and 0.2-mL microcentrifuge tubes and generate $15,500 \times g$.

^[3] As needed to connect laboratory vacuum to liquid trap and liquid trap to P200 pipette tip.

Instrument installation by only trained personnel

IMPORTANT! The Ion GeneStudio™ S5 System is installed only by trained service personnel and must not be relocated without assistance from trained service personnel. See “Customer and technical support” on page 59.

Nucleic acid contamination

IMPORTANT! A primary source of contamination is DNA fragments from previously processed samples. Do not introduce amplified DNA into the library preparation laboratory or work area.

Instrument vibration and clearances

IMPORTANT! Significant vibration during sequencing may add noise and reduce the quality of the measurements. The Ion GeneStudio™ S5 System must be installed on a bench that is free from vibrations or in contact with equipment that can cause vibrations to the bench (freezers, pumps, and other similar equipment).

IMPORTANT! Place the instrument at least 40 in. (1 meter) away from major sources of electronic noise such as refrigerators or microwaves.



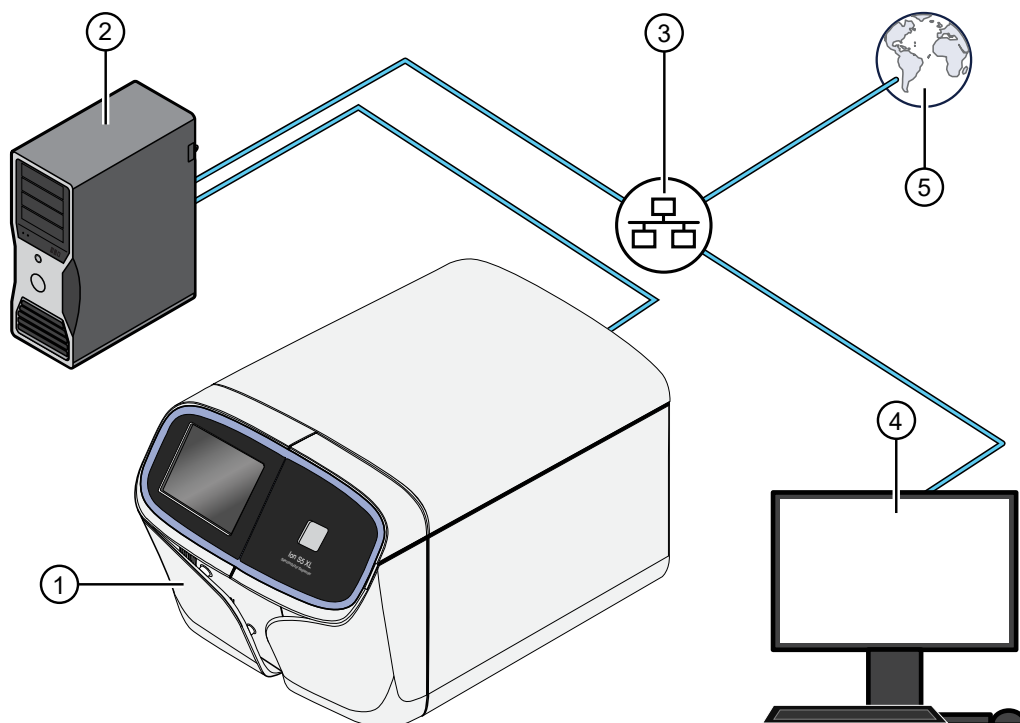
System components

Ion GeneStudio™ S5 System component configuration and layout

The Ion GeneStudio™ S5 Sequencer uses a virtual server, and the Ion GeneStudio™ S5 Plus Sequencer is equipped with an internal Ion Torrent™ Server. These instruments are not connected to an external Ion Torrent™ Server. The Ion GeneStudio™ S5 Prime System includes the Ion GeneStudio™ S5 Prime Sequencer and an external Ion Torrent™ Server.

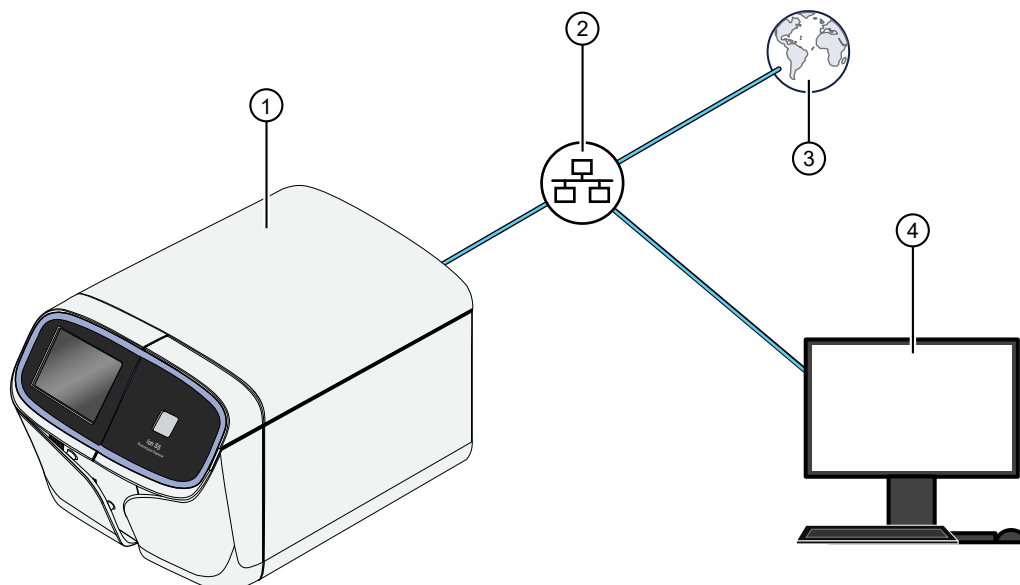
We support the layout in which the external Ion Torrent™ Server is connected directly to the Ion GeneStudio™ S5 Prime Sequencer, instead of through a local area network from a remote location such as a server room. Data are robustly transferred from the Ion GeneStudio™ S5 Prime Sequencer to the Ion Torrent™ Server when they are connected directly by a standard Category 6 Ethernet cable that is provided with the installation materials.

IMPORTANT! The Ion GeneStudio™ S5 Prime Sequencer must be connected to the Ion Torrent™ Server by a standard Category 6 Ethernet cable. We do not troubleshoot data transfer problems that are associated with an indirect connection between the Ion GeneStudio™ S5 Prime Sequencer and the Ion Torrent™ Server.



Ion GeneStudio™ S5 Prime System configuration

- ① Ion GeneStudio™ S5 Prime Sequencer
- ② Ion Torrent™ Server
- ③ Local area network
- ④ Client computer
- ⑤ Internet



Ion GeneStudio™ S5 System or Ion GeneStudio™ S5 Plus System configuration

- ① Ion GeneStudio™ S5 Sequencer or Ion GeneStudio™ S5 Plus Sequencer
- ② Local area network
- ③ Internet
- ④ Client computer

Internet connectivity

The Ion GeneStudio™ S5 Sequencer, the Ion GeneStudio™ S5 Plus Sequencer, or the Ion Torrent™ Server (in an Ion GeneStudio™ S5 Prime System) should be connected to a network with internet access. Connecting to the Internet allows you to update your software and access remote system support easily. Software updates through the network/internet are free. If you do not connect your instrument or server to a network, software updates must be manually installed via USB.

IMPORTANT! Thermo Fisher Scientific does not support the USB method for updating software.

Any problems (for example, file corruption, incomplete updates) resulting from update of the sequencer or server using a USB device that require correction of the faulty update is not covered by your Ion GeneStudio™ S5 Sequencer, Ion GeneStudio™ S5 Plus Sequencer, or Ion GeneStudio™ S5 Prime Sequencer warranty or any service contract you may have bought. You are required to schedule an on-site Time and Materials visit by a field service engineer to correct the problem at your own expense.

In providing outbound access to the Internet from the server, you enable the support team to provide inbound support. The Ion GeneStudio™ S5 Sequencer, the Ion GeneStudio™ S5 Plus Sequencer, and the Ion Torrent™ Server run a remote monitor agent that can provide service personnel with critical system information, such as installed software versions and instrument alarms. With your permission, the agent also allows service personnel to access remotely the Ion GeneStudio™ S5 Sequencer, the Ion GeneStudio™ S5 Plus Sequencer, or the Ion Torrent™ Server, which is required for system support. Without remote access, service personnel cannot access, view, and troubleshoot problems regarding machine performance.

To enable full support, the Ion Torrent™ Server must have outbound internet access (ports 22, 80, and 443) and be behind an appropriately configured firewall. Although not recommended, you can enable access to the Torrent Browser (the web server running on the Ion Torrent™ Server from the Internet). If you provide such access, you must restrict access to the server using HTTP and AUTH firewall rules, or a combination of the two. Implementing and maintaining such restrictions is the responsibility of your server administrator and not of Thermo Fisher Scientific.

Note: For further support and troubleshooting for Ion Torrent™ Server network access requirements, see Software Administration in *Torrent Suite™ Software Help*.

Ion GeneStudio™ S5 System component positions

Note: These positions apply to the Ion S5™ Systems and Ion S5™ XL Systems.

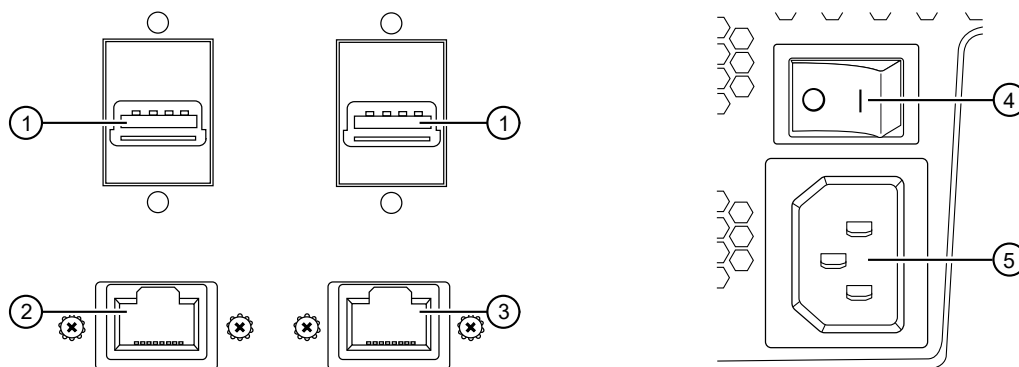


- ① Touchscreen
- ② Power button
- ③ Ion S5™ Sequencing Reagents cartridge
- ④ Chip clamp
- ⑤ Ion S5™ Wash Solution bottle. The waste reservoir is located behind the Ion S5™ Wash Solution bottle (shown on the right).
- ⑥ Ion S5™ Cleaning Solution bottle
- ⑦ Waste reservoir

Note:

- The system uses RFID technology to verify that the proper reagents are loaded in positions 3, 5, and 6. Reagents that exceed their expiration date or usage count generate an error message prompting you to replace the reagent before performing the run.
- RFID regulatory information is on the main screen under **Options ▶ Regulatory info**.

Ion GeneStudio™ S5 Sequencer input and output connections



- ① USB ports—Connects a USB device to the instrument.
- ② Ethernet port—An RJ45 port that provides Ethernet (Gigabit) communication between the sequencer and a local area network.
- ③ Ethernet port—An RJ45 port that provides Ethernet (Gigabit) communication with the Ion Chef™ Instrument.
- ④ On/off switch—Power switch, where the states are on (|) or off (O).
- ⑤ Power port—100–240VAC port that provides power to the instrument.



Instrument operation

Precautions

Avoid nucleic acid contamination

IMPORTANT! A primary source of contamination is spurious nucleic acid fragments from previous sample processing steps. Do not introduce amplified DNA into the work area where the instrument is located.

Avoid strong electromagnetic radiation



WARNING! Do not use the instrument in close proximity to sources of strong electromagnetic radiation (for example, unshielded intentional RF sources), as these sources can interfere with proper operation.

Protection by equipment



WARNING! The protection that is provided by the equipment can be impaired if the instrument is operated outside the environment and use specifications, the user provides inadequate maintenance, or the equipment is used in a manner that is not specified by the manufacturer (Thermo Fisher Scientific).

Power the Ion GeneStudio™ S5 Sequencer on or off

Note: If the Ion GeneStudio™ S5 Sequencer is powered on, and the touchscreen is blank, touch the screen to "wake" the touchscreen.

Power on

If the touchscreen is unresponsive, check the power switch on the back of the instrument to ensure that the switch is in the on (I) position. If the power switch is in the off (O) position, proceed with step 1. If the power switch is already in the on position, proceed to step 2.

1. Turn the power switch on the back of the instrument to the on (I) position.
2. Press the power button on the left side of the instrument.
The button illuminates. When the instrument touchscreen Main Menu appears, the instrument is ready for use.

Power off

It is not necessary to power off the instrument overnight or over the weekend. If the instrument will not be used for more than three days, power off the instrument as follows:

1. In the Main Menu, tap **Settings** ▶ **System Tools** ▶ **Shut Down**.
2. Select either **Shut Down** or **Reboot**.

Note: If you select **Shut Down**, a confirmation message appears. Select **Yes** to power off the instrument.

Note: Do *not* press the power button during a run. Interrupting power to the instrument during a run can result in sequencing run failure and loss of sample.

Update the Ion GeneStudio™ S5 System software

Note: An internet connection is required for the Ion GeneStudio™ S5 System to receive alerts that software updates are available.

If an update to the Ion GeneStudio™ S5 Sequencer software is available, the **Notifications/Alarms** button illuminates red in the touchscreen Main Menu. Press the red **Alarms** button to see the detailed messages. If a message states **New Software Available**, follow these steps to update the software:

1. In the Main Menu, press **Settings** ▶ **Check for Updates**.
2. Press **Update** to download and install the updates automatically.
3. Press **Done**.

The instrument automatically restarts when the software update is complete.

Ion GeneStudio™ S5 instrument run times

Approximate Ion GeneStudio™ S5 instrument run times for a given chip type, read length, and number of reads are shown in the following table. Run times indicate typical time from run start to BAM file delivery (1 run per initialization).

Chip type	Number of reads	Read length (bases) and output (Gb)	Instrument run time		
			Ion GeneStudio™ S5 Sequencer	Ion GeneStudio™ S5 Plus Sequencer	Ion GeneStudio™ S5 Prime Sequencer
Ion 510™ Chip	2–3 M	200 bases 0.3–0.5 Gb	4.5 hours	3 hours	3 hours
		400 bases 0.6–1 Gb	10.5 hours	5 hours	5 hours
Ion 520™ Chip	4–6 M	200 bases 0.6–1 Gb	7.5 hours	3.5 hours	3 hours
		400 bases 1.2–2 Gb	12 hours	5.5 hours	5.5 hours
	3–4 M	600 bases 0.5–1.5 Gb	12 hours	5.5 hours	5.5 hours
Ion 530™ Chip	15–20 M	200 bases 3–4 Gb	10.5 hours	5 hours	4 hours
		400 bases 6–8 Gb	21.5 hours	8 hours	6.5 hours
	9–12 M	600 bases 1.5–4.5 Gb	21 hours	8 hours	7 hours
Ion 540™ Chip	60–80 M	200 bases 10–15 Gb	19 hours	10 hours	6.5 hours
Ion 550™ Chip	100–130 M	200 bases 20–25 Gb	—	11.5 hours	8.5 hours

Maintain the sequencer

Materials required

- Lint-free wipes
- 70% isopropanol
- *(Optional)* 10% bleach solution

Clean or decontaminate the sequencer

In the event of a spill or leak on or inside the instrument, clean and decontaminate the sequencer.

Note: Dispose of all waste in appropriate liquid or solid waste containers.

1. Remove the Ion S5™ Wash Solution bottle, then remove and empty the waste reservoir.
2. Remove the Ion S5™ Sequencing Reagents cartridge.
3. Inspect the waste and nucleotide reagent bays for liquid.
4. Use absorbent paper to soak up as much liquid as possible, then wash the affected area with 10% bleach solution.
5. Wipe the affected surfaces with 70% isopropanol, then allow to air-dry.

Perform the sequencer cleaning manually

The Ion GeneStudio™ S5 Sequencers require cleaning before initialization. Cleaning is normally performed automatically at the completion of the previous sequencing run. To enable two sequencing runs on a single initialization, you must deselect the "Enable post-run clean" checkbox for the first sequencing run. The post-run cleaning is then performed normally after the second sequencing run. However, if the "Enable post-run clean" checkbox is deselected for the second run, the cleaning is not performed after either run. If the second sequencing run is not performed, or the cleaning is not performed automatically after the second run, the instrument will not allow the subsequent initialization to proceed until a manual cleaning has been performed.

If an Ion GeneStudio™ S5 Sequencer is initialized and

- a sequencing run is not started within 24 hours, *or*
- a run is not started or completed due to a power failure or an abort

do not perform a manual cleaning. An instrument reset run is required before reinitialization. See "Perform an instrument reset run with an initialized sequencer that is loaded with an unused Reagents cartridge" on page 21 for more information.

When a cleaning is necessary, use the following procedure:

1. On the home screen, press **Settings ▶ Clean Instrument**.
The instrument door unlocks allowing access to the consumables.
2. Remove the Ion S5™ Wash Solution bottle to access the waste reservoir, then remove and empty the waste reservoir.



3. Reinstall the empty waste reservoir and a *used* Ion S5™ Wash Solution bottle.
4. Ensure the Ion S5™ Sequencing Reagents cartridge and Ion S5™ Wash Solution bottle are properly installed.

IMPORTANT! Perform the cleaning with a used reagent cartridge and wash solution bottle installed. The cleaning procedure pumps cleaning solution into the wash solution bottle and reagent cartridge making them unsuitable for sequencing.

5. Place a used sequencing chip in the chip clamp, then push the chip clamp in all the way to engage.
6. Close the instrument door, then press **Next**.
Cleaning takes ~35 minutes to complete. On completion, the instrument door automatically unlocks and the chip and cartridge clamps disengage.

Perform an instrument reset run with an initialized sequencer that is loaded with an unused Reagents cartridge

Cleaning is normally automatically performed at completion of a sequencing run. If an Ion GeneStudio™ S5 System is initialized, an instrument reset run is required to facilitate proper cleaning before reinitialization in one of the following situations.

- A sequencing run is not started within 24 hours after initialization.
- A sequencing run is not completed due to a power failure or an abort, and <200 flows occurred before the stoppage.

Do NOT perform a manual cleaning with an unused, initialized Ion S5™ Sequencing Reagents cartridge.

Note:

- If a power failure or abort occurs during the second of two runs started after a single initialization, a manual cleaning is sufficient (see [page 19](#)).
 - If the number of flows that occurred before a power failure or abort is unknown, perform an instrument reset run.
-

To perform an instrument reset run, use the following procedure before reinitialization.

1. In the instrument touchscreen main menu, tap **Run**.
The instrument door unlocks and the chip clamp disengages.
2. Ensure that a used sequencing chip is in the chip clamp, then push the chip clamp in all the way to engage.
3. Close the instrument door, then tap **Next**.
4. When prompted, select **Planned Run (none)**. Ensure that the **Enable post-run clean** checkbox is selected, then tap **Review**.
5. In the **Select Run** screen, tap **Edit**, then in the **Detail** screen set the number of flows to **200**. Ensure that the **Post-Run/Clean** checkbox is selected, then tap **Close**.
6. Tap **Start run**, then tap **Accept** to confirm that Post-Run Clean is enabled, and to start the run.

When the instrument reset run completes, the instrument automatically performs the cleaning procedure. After cleaning, the touchscreen returns to the main menu.

Instrument planned maintenance

The Ion GeneStudio™ S5 Sequencer needs annual planned maintenance to keep it operating in peak condition. A notification appears when maintenance is due. To schedule a planned maintenance visit, contact your Thermo Fisher Scientific Field Service Engineer.



Touchscreen reference

Clean, Initialize, and Run

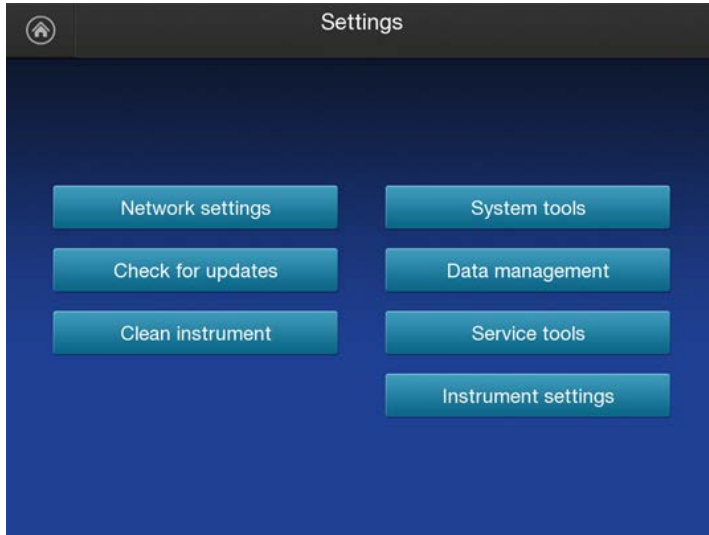


In the **Home** screen, the **Clean**, **Initialize**, and **Run** programs lead you through the necessary steps to prepare the instrument for sequencing and to start a sequencing run. Press the main dial to start a program. Simple, easy to follow instructions are provided on the touchscreen.

- Cleaning must be performed before each initialization to ensure that the reagents from the previous run are cleared from the fluid lines. The **Clean** program is normally performed automatically at the completion of the previous sequencing run. Perform a **Clean** if for any reason the sequencing run was not properly completed. Follow the instructions that are provided on the touchscreen.
- The **Initialize** program must be performed before each run to load and prepare the run reagents. The **Initialize** program walks you through:
 - Emptying the waste reservoir
 - Loading the reagent cartridge, wash solution, and cleaning solution
- The **Run** program walks you through steps leading up to and through sequencing, including:
 - Placing a loaded chip on the instrument
 - Selecting a Planned Run created in the Torrent Suite™ Software
 - Performing sequencing

Settings

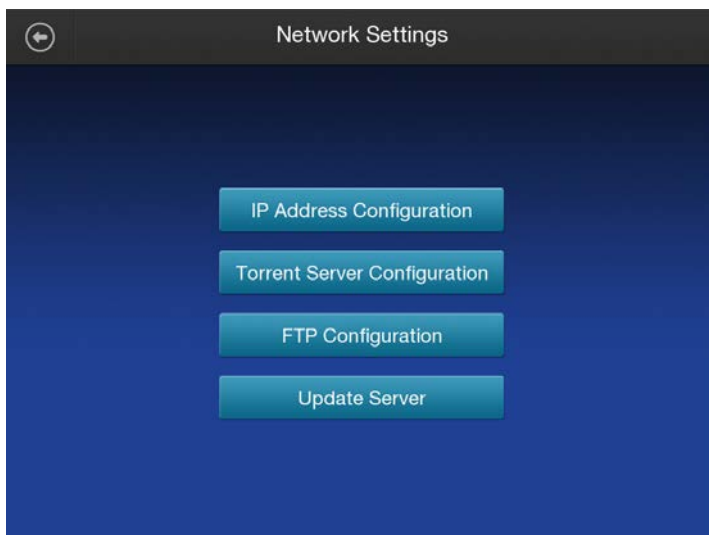
In the **Settings** menu, view or change instrument settings, manage data and network configurations, and update the installed software.

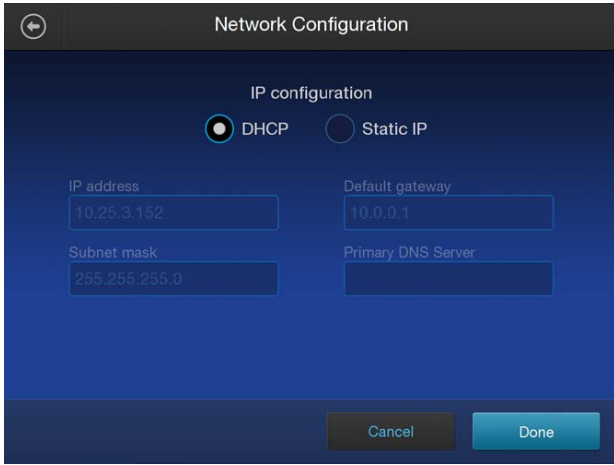
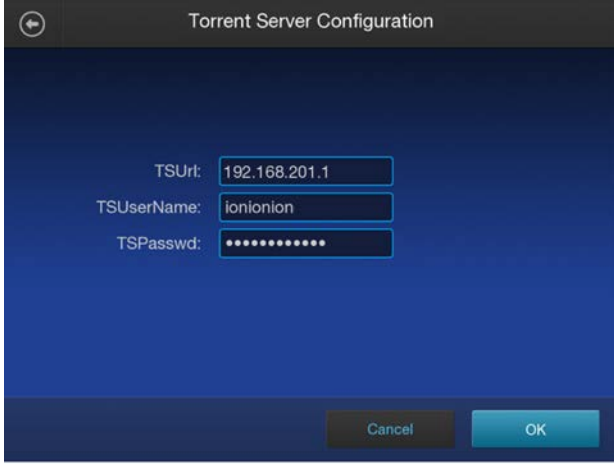


Note: The **Service tools** option is for use only by trained service personnel.

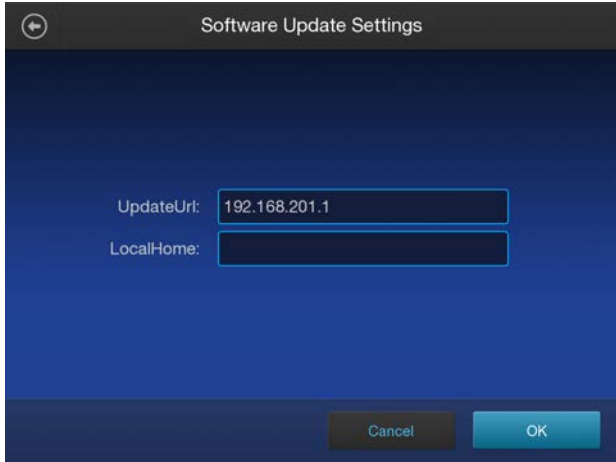
Network Settings

Use the **Network Settings** menu to configure the IP address, configure the Torrent Server, configure FTP settings, and update the server.



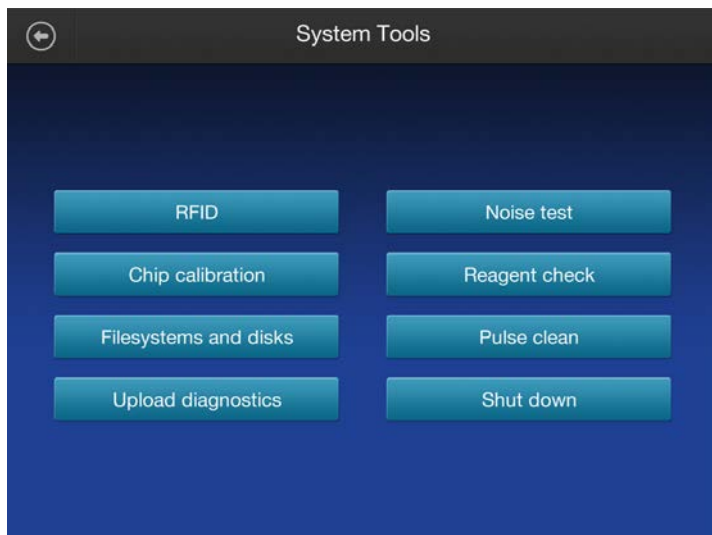
Item	Description	When and how to use
IP Address Configuration	Use this screen to set or change the IP configuration (DHCP or Static IP). 	1. Select either the DHCP or Static IP option. 2. Touch the screen in the field that you want to edit to activate the field. A virtual keyboard appears. 3. Enter the new information, then press Done.
Torrent Server Configuration	Use this screen to change the Ion Torrent™ Server IP address and user information.  Enter one of the following IP addresses in the TSUrl field for your instrument: • Ion GeneStudio™ S5/Ion GeneStudio™ S5 Plus Sequencer: 192.168.122.249 • Ion GeneStudio™ S5 Prime Sequencer: 192.168.201.1	Use when a change to the Ion Torrent™ Server IP address or user information is required.

(continued)

Item	Description	When and how to use
FTP Server Configuration	Use this screen to change the FTP Server IP address and user information.  Enter the following IP address in the FTPserver field for your instrument. - Ion GeneStudio™ S5/Ion GeneStudio™ S5 Plus Sequencer: 192.168.122.249 - Ion GeneStudio™ S5 Prime Sequencer: 192.168.201.1	Use when a change to the FTP Server IP address or user information is required.
Update Server	Use this screen to change the software update server settings.  Enter the following IP address in the UpdateUrl field for your instrument: - Ion GeneStudio™ S5/Ion GeneStudio™ S5 Plus Sequencer: 192.168.122.249 - Ion GeneStudio™ S5 Prime Sequencer: 192.168.201.1	Use when a change to the URL for the software update servers is required. The values on this screen are set at installation. Do not change the settings except at the direction of Thermo Fisher Scientific service staff.

System Tools

Use the **System Tools** menu to upload instrument diagnostics, manage data, and shut down or reboot the instrument.

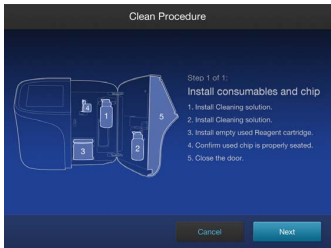


Item	Description	When and how to use
RFID	Lists the product, product expiration date, and remaining uses of the product. 	Use to determine if reagents that were left on an instrument during a period of instrument shutdown have expired. 1. Press the screen within the row of the product to view an RFID Tag Detail page. 2. Press Done to return.
Noise Test	Provides real-time measurement of electrical noise readings on the chip. 	Use for troubleshooting if directed to do so by Technical Support.

(continued)

Item	Description	When and how to use
Chip Calibration	Checks the status of a chip. 	1. Insert a chip into the chip clamp. 2. Press Chip Calibration to start the chip calibration. 3. In the dropdown lists, select the Image to display and the type of Calibration to perform. 4. Press Cancel to quit the chip calibration.
Reagent Check	Performed by the instrument as part of initialization. Confirms that initialization passes by testing reagent flow and signal detection.  Example Reagent Check for an Ion 540™ Chip  Example Reagent Check for an Ion 550™ Chip	Use for troubleshooting if directed to do so by Technical Support. • Press Run to start the reagent check. • Press Last Trace to review the most recent reagent check result trace. Reagent Check confirms that the chip calibration passes, and fails only if the chip is defective, or if the Ion S5™ Wash Solution is not flowing.

(continued)

Item	Description	When and how to use
Filesystems and disks	Provides real-time drive activity status and disk space used. 	Use for troubleshooting if directed to do so by Technical Support. 1. In the System Tools menu, press Filesystems and disks. The Filesystems screen opens. 2. Press the screen within the row of the Filesystem that you want to view, then press Details. The Disk Drives in <Filesystem> screen opens. 3. Press the screen within the row of the Drive to view details of the disk drive. The Details for <Drive> screen opens. 4. Press Done to return to the previous screen.
Pulse clean	Provides step-by-step instructions to perform a pulse clean of the instrument if a solid-state buffer leak occurs. 	Use for troubleshooting if directed to do so by Technical Support. 1. Install consumables and a used sequencing chip. Install cleaning solution in both the wash solution (1) and cleaning solution (2) locations. Use an empty Reagent cartridge. 2. Close the instrument door, then press Next. The pulse cleaning procedure starts automatically. See “Solid-state buffer leak” on page 44 for more information.

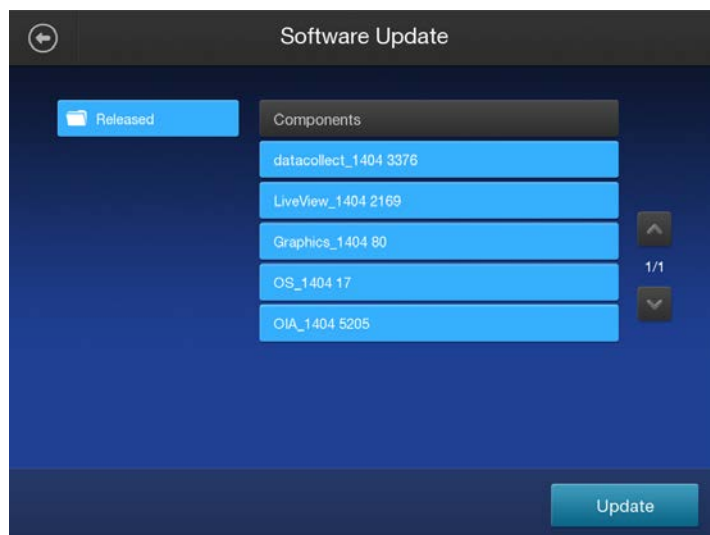
(continued)

Item	Description	When and how to use
Upload Diagnostics	Allows upload of instrument diagnostics files for review by support personnel. 	Use for troubleshooting if directed to do so by Technical Support. The system must be connected to an active network port. If no network is present, the file is pushed to the Torrent Browser. You can then download the file and email it to Technical Support. To retrieve the file using the Torrent Browser, navigate to Help ► Instrument Diagnostics, then follow the on-screen instructions.
Shut Down	Access to Shut Down and Reboot commands. 	Use for troubleshooting If directed to do so by Technical Support. It is not necessary or recommended to shut down the instrument overnight or over the weekend. If needed to shut down the instrument, see “Power off” on page 17.

Check for software updates

When software updates are available, a notification appears on the **Home** screen.

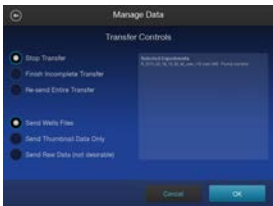
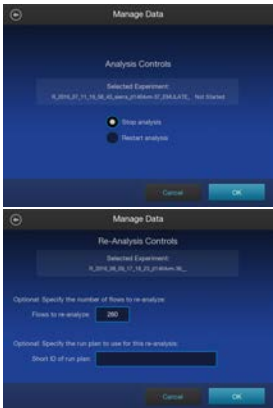
1. In the **Settings** menu, tap **Check for Updates**.
The **Software Update** screen lists the available updates.



2. Tap **Update** to download and install the updates.

Data management

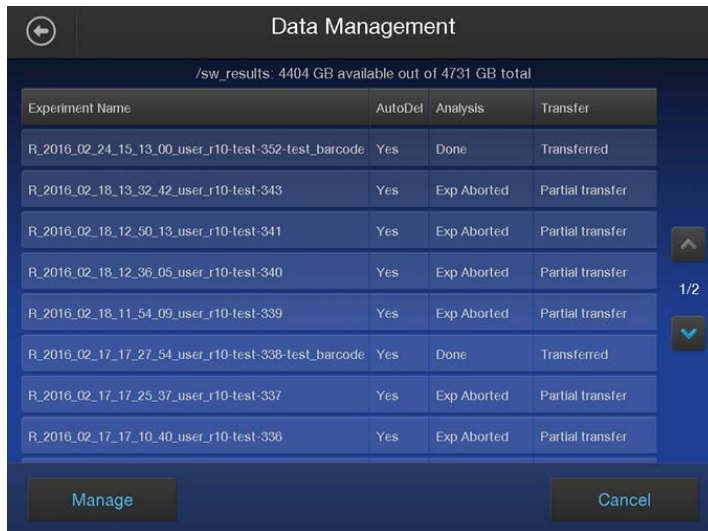
Use the **Data management** function to delete run data manually, analyze the data, or transfer the data to the Ion Torrent™ Server. Under normal conditions, run data are automatically transferred to the server, then deleted from the instrument hard drive.

Item	Description	When and how to use
Delete data	Manually delete run data from the instrument.	If the instrument hard drive becomes full, see “Manually delete run data” on page 31.
Transfer data		Transfer run data from the instrument hard drive to the Ion Torrent™ Server. 1. Tap the screen in the Experiment Name row of the individual experiment to be transferred. 2. Tap Transfer. 3. Select the action to be performed. 4. Select the how you want the files to be transferred. 5. Tap OK.
Analysis Controls	Use this screen to stop or reanalyze the data of individual runs. 	1. Tap the screen in the Experiment Name row of the individual experiment to be analyzed, then press Analysis. 2. Select the Stop Analysis or Restart Analysis option, then tap OK. 3. If performing a reanalysis, follow the on-screen instructions, then tap OK.

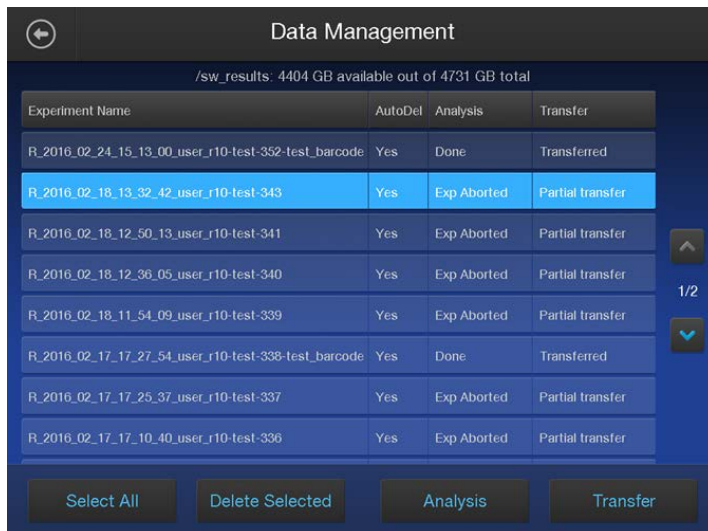
Manually delete run data

To troubleshoot data management problems the **Data Management** function allows users to delete run data manually, or transfer the data to the server. Under normal conditions, run data are automatically transferred to the Ion Torrent™ Server, then deleted from the instrument hard drive.

1. In the **Settings** menu, tap **Data management** to access the **Data Management** screen, then tap **Manage**.



2. Tap **Select All** to select all the available experiments, or tap the screen in the **Experiment Name** row of the individual experiment to be managed.



3. Press **Delete Selected**.

Perform the instrument Clean operation

The **Clean** program is normally performed automatically at the completion of the previous sequencing run. Perform a **Clean** if the sequencing run:

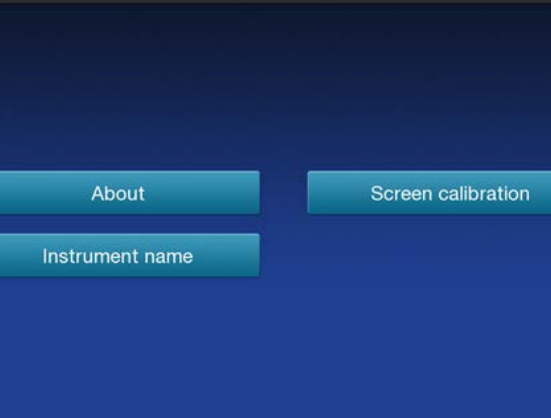
- was aborted or had a power failure during the second of two runs started after a single initialization.
- is not completed and >200 flows occurred before the stoppage.
- the post-sequencing run cleaning was not completed.

IMPORTANT! Do NOT perform a manual cleaning with an unused, initialized Ion S5™ Sequencing Reagents cartridge. Perform an instrument reset run instead. See “Perform an instrument reset run with an initialized sequencer that is loaded with an unused Reagents cartridge” on page 21 for more information.

1. In the **Settings** menu, tap **Clean Instrument**.
The Clean procedure begins automatically.
2. Follow the on-screen instructions (see page 19 for more information), then tap **Next**.



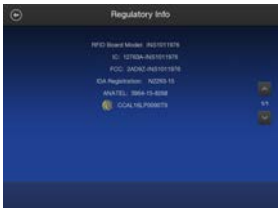
The user interface returns to the Home screen when the cleaning is complete.



The screenshot displays the 'Instrument Settings' interface. At the top, a dark grey header bar contains a circular back arrow icon on the left and the text 'Instrument Settings' in the center. The main content area has a dark blue background. Three light blue rectangular buttons with rounded corners are arranged in the center. The button on the left is labeled 'About'. To its right is a button labeled 'Screen calibration'. Below the 'About' button is a third button labeled 'Instrument name'.

Ion S5™ Sequencer User Guide

(continued)

Item	Description	When and how to use
Regulatory Info	Lists instrument-specific regulatory information. 	Use to view instrument regulatory information.
Instrument Name		Use to change the instrument name.
Screen Calibration		Use for troubleshooting. If directed to do so by Technical Support. Touch the red cross with your finger or a stylus each time it appears. In total, you will touch the screen 4 times, one time in each corner.

Touchscreen icons



Number	Icon	Description
1		Network connectivity – connected
		Network connectivity – not connected
2		Instrument idle
		Sequencing in progress
		Instrument ready
		Error
3		Chip status – Absent
		Chip status – Standby
		Chip status – Connecting
		Chip status – Ready
		Chip status – Imaging
		Chip status – Error

(continued)

Number	Icon	Description
4		Manifold air pressure – Good
		Manifold air pressure – Bad. If the icon reads "0" and there is an alarm, contact Technical Support.
5		Regulator air pressure – Good
		Regulator air pressure – Bad. If the icon reads "0" and there is an alarm, contact Technical Support.
6		Manifold temperature – Good
		Manifold temperature – Bad. Check for related alarms. If the alarm persists contact Technical Support.
7		Chip block temperature – Good
		Chip block temperature – Bad. If the icon reads "0" and there is an alarm, contact Technical Support.
8		Instrument File System Space, the percent of file space used is indicated. ^[1] The instrument checks for sufficient disk space before each run and notifies the user if there is not enough.
9		Torrent Server File System Space, the percent of file space used is indicated. Note: If the indicator turns red, archive data from the server to free up disk space. Refer to the Torrent Suite™ Software help for more information on archiving data.

^[1] Indicator turns yellow when disk space is ≥67% full, indicator turns red when ≥90% full.

Alarms, Notifications, and Events

If the Alarms/Notifications indicator appears, press the indicator to open the **Notifications** screen, then press **Alarms** or **Events** see detailed messages.



① Alarms and Notifications indicator.

Alarm	Description	Recommended Action
Fatal alarms		
Fluidics Calibration Needed.	—	Contact Technical Support.
Chip Cooler can not reach desired temperature.	—	Contact Technical Support.
Drive sdX failed smartctl health check.	Where X is the letter of the drive that failed. Indicates that a drive is failing.	Contact Technical Support.
Failed to mount the results filesystem.	One or more drives are missing from the RAID that makes up sw_results.	Contact Technical Support.
System sensor check failed ipmi -sensors.	A component on the motherboard is failing or has failed.	Contact Technical Support.
System sensor check failed, CC.CC V (LL.LL/HH.HH)	Where CC.CC = current voltage, LL.LL = the low voltage threshold, HH.HH = the high voltage threshold. Indicates a voltage drop on the motherboard.	Contact Technical Support.



(continued)

Alarm	Description	Recommended Action
CPU temperature check failed, Physical Id X.	Where X is the processor number. Indicates that the processor is over heating.	Contact Technical Support.
Front/Rear FanX speed below threshold YYYY.	Where X is the fan number and YYYY is the low threshold speed for the system fan.	Contact Technical Support.
FPGA failed to connect in emulate mode.	Internal error.	Contact Technical Support.
Valve Link Failure.		
OIA (On Instrument Analysis) is not running.		
FPGA temperature is above the threshold of 70°C.	One or more of the system fans has failed.	Contact Technical Support.
Boot drive exceeded 75%.	The boot drive is unexpectedly full.	Contact Technical Support.
RAID degraded one disk has failed.	One of the Ion Torrent™ Server drives has failed. System is still operational but must be serviced soon.	Contact Technical Support.
RAID failure detected, more than one disk failed.	More than one of the Ion Torrent™ Server drives has failed. System is no longer operational.	
/sw_results directory not mounted.	One or more of the drives used during data collection has failed.	
No GPU detected.	GPU has failed.	Contact Technical Support.
No GPU Driver detected.		
Air Compressor failure.	Air compressor is either leaking or has stopped working.	Contact Technical Support.
Non fatal alarms		
Instrument must be cleaned before starting another run.	Maximum number of runs since last clean has been exceeded.	Perform an instrument cleaning. See “Perform the sequencer cleaning manually” on page 19.
Ambient temperature is above the threshold of 35°C.	Verify the temperature in the lab.	Contact Technical Support if the temperature of the lab is within normal operating temperature (20°C to 30°C).
RAID degraded one disk has failed.	One of the Ion Torrent™ Server drives has failed. System is still operational but must be serviced soon.	Contact Technical Support.

(continued)

Alarm	Description	Recommended Action
No Connectivity to Torrent Server.	Check that the Ethernet cable connecting the instrument to the server is properly connected, see page 15 for the Ethernet port location.	Reconfigure the Ion Torrent™ Server as needed, see “Network Settings” on page 23 .
No Connectivity to FTP server.		Reconfigure the FTP server as needed, see “Network Settings” on page 23 .
Network Manager not connected.	Check that the Ethernet cable connecting the instrument to the local area network is properly connected, see page 15 for the Ethernet port location.	If the problem persists, contact Technical Support.
Newer Software Available	Notification that software updates are available.	Update your software. See “Check for software updates” on page 29 .

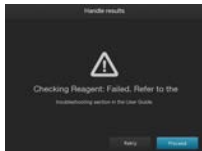


Troubleshooting

Instrument alarms and events

Observation	Possible cause	Recommended action
Red "Alarms" and/or "Events" message in Main Menu 	- Available software updates were detected. - Connectivity issues were detected. - Instrument did not detect required files or hardware.	Tap the red Alarms icon to see detailed messages. - If a message states "Newer Software Available": IMPORTANT! After updates are installed, the instrument must be restarted. - In the main menu, tap Settings ▶ Check for Updates. - Select the Released Updates checkbox, then tap Update. - When installation is complete, follow the onscreen prompts to restart the instrument. Note: In some cases, the instrument restarts automatically after software installation. - If a message states "No Connectivity to Torrent Server", "No Connectivity to ftp server", or "Network Manager not connected", disconnect and re-connect the ethernet cable, confirm that the router is operational, and verify that the network is up and running. - For any other messages: - Power off the instrument: On the home screen, tap Settings ▶ System Tools ▶ Shut Down ▶ Shut Down. - Wait 30 seconds, then press the button on the side of the instrument to power on the instrument. - If the red "Alarms" and/or "Events" message still appears in the main menu, contact Technical Support.

Initialization fails

Observation	Possible cause	Recommended action
Chip Check fails	- The chip clamp was not closed. - The chip was not properly seated. - The chip was damaged.	- Open the chip clamp, remove the chip, and look for signs of water outside the flow cell. - If the chip appears damaged, replace it with a new one. - Close the clamp, then repeat the Chip Check. - If the chip passes, click Next. If the chip fails, replace it with a new chip, then press Chip Check. - If Chip Check continues to fail, there could be a problem with the chip socket. Contact Technical Support.
Reagent Check fails 	A chip failure was detected.	- Replace the used sequencing chip used during initialization with a different used chip. IMPORTANT! If you are using Torrent Suite™ Software version 5.12.1 or earlier and if a used Ion 550™ Chip was used for initialization, replace the used chip with a used Ion 510™ Chip, Ion 520™ Chip, Ion 530™ Chip, or Ion 540™ Chip for the initialization. Using an Ion 550™ Chip causes an error. - Press Retry. - If the initialization completes without failure, touch Home, then continue with your sequencing run. - If the Reagent Check continues to fail, contact Technical Support.
	A wash failure was detected.	- Perform a manual cleaning of the sequencer, see “Perform the sequencer cleaning manually” on page 19. - Replace the Reagents cartridge, Wash Solution bottle, and Cleaning Solution bottle, then repeat initialization of the sequencer. - If the initialization completes without failure, press Home, then continue with your sequencing run. - If the Reagent Check continues to fail, contact Technical Support.
		Ensure that you install the Ion S5™ Wash Solution bottle with the instrument door fully open. If you install the bottle impeded by a partially closed door, the port can be obstructed, and a faulty installation and wash failure can result.
Pressure loss in Wash Solution bottle 	- Improper wash solution bottle installation. - Incomplete thawing of reagent cartridge.	- Tighten wash solution bottle cap. - Reseat wash solution bottle and reagent cartridge. - Repeat the initialization. - If leak check continues to fail, replace wash solution bottle and reagent cartridge, and repeat the initialization. - If leak test continues to fail, contact Technical Support.

Observation	Possible cause	Recommended action
Pressure loss in Reagent Cartridge Chamber 	- Problem with reagent cartridge. - Incomplete thawing of reagent cartridge.	1. Tighten wash solution bottle cap. 2. Reseat wash solution bottle and reagent cartridge. 3. Repeat the initialization. 4. If leak check continues to fail, replace wash solution bottle and reagent cartridge, and repeat the initialization. 5. If leak test continues to fail, contact Technical Support.
Pressure loss in Reagent Cartridge Bags 	- Problem with reagent cartridge. - Incomplete thawing of reagent cartridge.	1. Reseat reagent cartridge. 2. Repeat the initialization. 3. If the initialization continues to fail, replace the reagent cartridge and repeat the initialization. 4. If the initialization continues to fail, contact Technical Support.
Pressure loss in Cleaning Solution bottle 	- Problem with cleaning solution bottle. - Improper cleaning solution bottle installation.	1. Tighten cleaning solution bottle cap. 2. Reseat cleaning solution bottle. 3. Repeat the initialization. 4. If the initialization to fail, replace cleaning solution bottle and repeat the initialization. 5. If the initialization continues to fail, contact Technical Support.
Confirm Blowoff Valve works 	Problem with instrument valve.	1. Restart the instrument and repeat initialization. 2. If the alarm repeats, open a service call.

Troubleshooting using control libraries

Observation	Possible cause	Recommended action
Ion Sphere™ Test Fragments are not present in the Test Fragment Report section of the run report, and library sequencing is poor	- Chip loading was poor. - Control Ion Sphere™ Particles were not added to the sample	1. Confirm that the Control Ion Sphere™ Particles (included in the Ion S5™ Installation Kit) were added. 2. If controls were added, contact Technical Support.
Control Ion Sphere™ Particles are present in the run report, but AQ20 throughput is poor	- The quality of the library was poor. - The quality of the template preparation was poor.	Verify the quality of the library and template preparations using quality assessment procedures recommended in the appropriate library and template preparation user guides. 1. Use the Human CEPH Genomic DNA Control or Human CEPH Control 200 Library, included in the Ion S5™ Controls Kit (Cat. No. A27760), to prepare template-positive ISPs with the Ion OneTouch™ 2 Instrument, or the Ion Chef™ Instrument. 2. Use the ISPs in an Ion GeneStudio™ S5 run. 3. If AQ20 throughput is still below specification, verify the quality of unenriched and enriched ISPs to identify a problem in template preparation. 4. If ISP quality is good, but AQ20 throughput is below specification, contact Technical Support.

Troubleshooting using Control Ion Sphere™ Particles

To prepare Control Ion Sphere™ Particles, provided in the Ion S5™ Installation Kit, for an installation or troubleshooting sequencing run:

1. Create a Planned Run.
2. Clean and initialize the Ion GeneStudio™ S5 Sequencer.
3. Prepare the Control Ion Sphere™ Particles for sequencing:
 - a. Vortex the Control Ion Sphere™ Particles for 5 seconds, then centrifuge for 2 seconds before taking aliquots.
 - b. Add 66 µL of Control Ion Sphere™ Particles to an empty 0.2-mL PCR tube (non-polystyrene).
 - c. Add 150 µL of Ion S5™ Annealing Buffer to the tube.
4. Anneal the sequencing primer to the enriched ISPs, then follow the remaining procedures in your appropriate user guide to load a chip and start the sequencing run.

Solid-state buffer leak

Observation	Possible cause	Recommended action
Two or more of the following observed - Excessive read trimming or the number of short reads is significantly higher than expected. - Percentage of low quality reads is significantly higher than expected. - Poor sequencing performance or no Test Fragments.	A solid-state buffer leak occurred.	Contact Technical Support.
		Perform the pulse cleaning protocol if directed to do so by Technical Support or your Field Service Engineer (FSE).

Pulse clean the sequencer

Pulse clean your Ion GeneStudio™ S5 Sequencer only if directed by Technical Support, or your Field Service Engineer (FSE), to recover instrument performance from a diagnosed solid-state buffer leak.

Contact Technical Support, or your FSE, to diagnose whether a solid-state buffer leak has occurred and obtain the required materials to perform the pulse cleaning.

1. Ensure that the sequencer has completed a normal post-run cleaning. In the main menu:

Display text	Description
Run	Instrument requires cleaning before a pulse cleaning can be performed. See “Perform the sequencer cleaning manually” on page 19.
Clean	Instrument requires cleaning before a pulse cleaning can be performed. See “Perform the sequencer cleaning manually” on page 19.
Initialize	Instrument is ready to be pulse cleaned. Proceed to step 2.

2. Press  (**Settings**) ▶ **System Tools** ▶ **Pulse clean**.

3. Follow the on-screen prompts to prepare the instrument.

- Install new bottles of Ion S5™ Cleaning Solution in both the Wash and Cleaning solution positions.
- Replace the Ion S5™ Sequencing Reagents cartridge with an empty used cartridge.
- Install a used sequencing chip.



4. Close the instrument door, then touch **Next**.

The pulse cleaning procedure begins. The user interface returns to the main menu when the cleaning is complete and the sequencer is ready to be initialized.



Supplemental procedures

Set up and test the Ion Chip™ Minifuge

Note: The Ion Chip™ Minifuge is required only when using the Ion OneTouch™ 2 System for template preparation. Ion Chef™ Instrument users do not require an Ion Chip™ Minifuge.

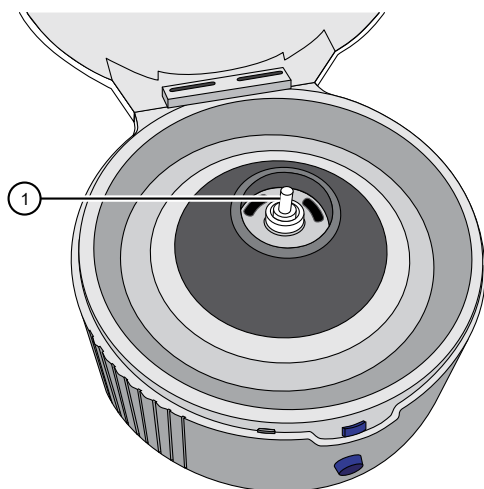
The Ion Chip™ Minifuge (Cat. No. 4479672 or 4479673) is used to load sequencing chips manually for use on Ion PGM™, Ion Proton™, and Ion GeneStudio™ S5 System sequencing platforms. To accommodate the larger chip size, and before using the minifuge to load chips for the first time, Ion Proton™ and Ion GeneStudio™ S5 System sequencer users must:

- Install the Ion S5™ or Ion Proton™ Minifuge Proton Rotor and Bucket (Cat. No. 4482578).
- Test the minifuge to confirm that no liquid is lost during centrifugation.

Note: The protocols in this section may also be used to convert the Ion Chip™ Minifuge back for use with Ion PGM™ sequencing chips.

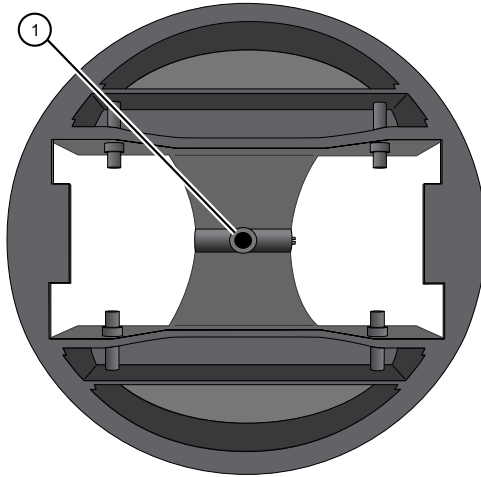
Install the Ion S5™ /Ion Proton™ Rotor and Buckets

1. Grasp the existing rotor and pull straight up to remove the rotor from the motor shaft.



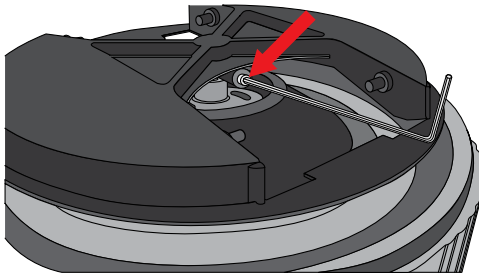
① Motor shaft

2. Press the Ion S5™/Ion Proton™ Rotor down onto the motor shaft to install.



① Insert motor shaft here

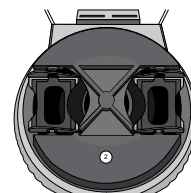
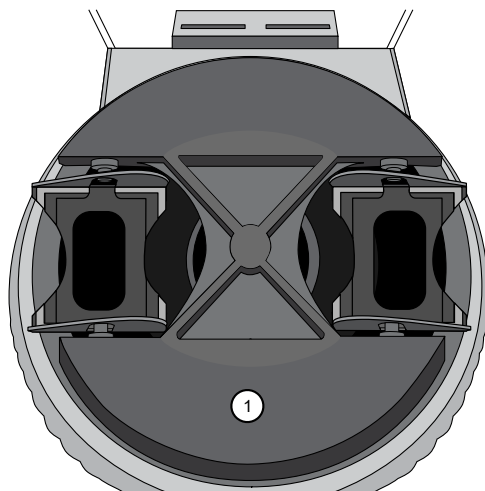
3. Tighten the set screw (arrow) with a 1.5-mm hex wrench.



Note: A newer version of the rotor lacks a set screw. In this case, simply press the rotor firmly onto the motor shaft to install.



4. Install the two buckets. Position the buckets with the larger semi-circular cut-outs facing out, and ensure that the buckets hang freely.



- ① Correct orientation
- ② Incorrect orientation

Test the minifuge

1. Prepare two previously-used chips:
 - a. Inject 100 μL of isopropanol two times into the loading port of each chip. After each injection, remove the expelled liquid from the opposite port.

Note: Use 50 μL volume of isopropanol if testing Ion PGM™ sequencing chips.

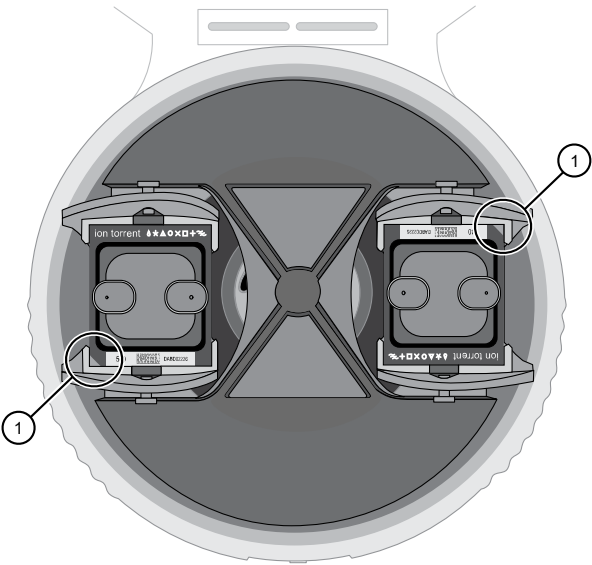
- b. Aspirate the remaining isopropanol from the flow cells for 5–10 seconds. Confirm that the chips are dry.

Note: To aspirate the isopropanol, attach a P200 pipette tip to a vacuum line, then place the pipette tip in the chip loading port.



2. Place the two chips prepared in step 1 in the centrifuge buckets, with the chip notch pointing out. Add 55 µL of nuclease-free water to each chip loading well (do not inject into the chip loading port).

Note: Use 35 µL volume of nuclease-free water if testing Ion PGM™ sequencing chips.



① Chip notch

3. Centrifuge for 5–10 seconds, then examine each chip.
The flow cell in each chip should be completely filled with liquid with no air bubbles. A small volume of liquid will remain in the loading well; this is normal.

Result	Action
The chips are NOT completely filled...	Contact Technical Support.
The chips ARE completely filled...	Centrifuge the chips for an additional 10 minutes, then check the chips again for air bubbles, especially near the inlet and outlet ports.
The chips have air bubbles after the additional 10 minute centrifugation...	Contact Technical Support.
The chips remain completely filled...	The centrifuge is ready to use for chip loading.



Instrument warranty

For new Ion Torrent™ instruments, Life Technologies warrants to and only to buyer for twelve (12) months from the date of shipping, that the Ion Torrent™ software and Ion Torrent™ instruments are free from defects in material and workmanship and conform to Life Technologies' published specifications in all material respects. Where a valid and timely claim in respect of breach of Ion Torrent™ warranty is submitted to Life Technologies, Life Technologies may, at its discretion, replace, repair or modify the Ion Torrent™ instrument. Any agreed replacement shall be at 1:1, like-kind basis, at no cost to the buyer. For Ion Torrent™ chips or reagents reasonably determined by Life Technologies to be defective, independent of user error, shall be replaced by Life Technologies on a 1:1, like-kind basis at no cost to buyer, provided that such defective Ion Torrent™ chips or reagents were used by buyer prior to their expiration date, or if there is no expiration date, the Ion Torrent™ chips or reagents were used within six (6) months of receipt, and the defect was promptly reported with appropriate detail to Life Technologies' technical support.

NO OTHER WARRANTIES SHALL BE APPLICABLE TO ION TORRENT PRODUCTS (WHETHER OR NOT ANY FURTHER WARRANTY DOCUMENTATION MAY BE INCLUDED IN THE SHIPMENT), WITH THE EXCEPTION OF THIRD PARTY WARRANTIES WITH RESPECT TO THIRD PARTY PRODUCT. ANY THIRD PARTY PRODUCTS ARE NOT COVERED BY THIS SECTION AND ANY WARRANTIES FOR THIRD PARTY PRODUCTS ARE PROVIDED BY THE ORIGINAL MANUFACTURER OF THE THIRD PARTY PRODUCT. Warranties are made only to buyer purchasing the Ion Torrent™ Product directly from Life Technologies, are not transferable and do not extend to the benefit of any other person or entity, unless otherwise expressly stated in writing by Life Technologies. ANY PRODUCT NOT COVERED BY AN EXPRESS WRITTEN WARRANTY IS SOLD AND PROVIDED "AS IS," WITHOUT WARRANTY OF ANY KIND, STATUTORY, EXPRESS OR IMPLIED. Any description of Ion Torrent™ Product recited in Life Technologies' quotation is for the sole purpose of identifying Ion Torrent™ Product, and any such description is not part of any contract between Life Technologies and buyer and does not constitute a warranty that Ion Torrent™ Product shall conform to that description. Any sample or model used in connection with Life Technologies' quotation is for illustrative purposes only, and is not part of any contract between Life Technologies and buyer and does not constitute a warranty that Ion Torrent™ Product will conform to the sample or model. No affirmation of fact or promise made by Life Technologies, whether or not in Life Technologies' quotation, shall constitute a warranty that Ion Torrent™ Product will conform to the affirmation or promise. Unless otherwise specified in writing in documentation shipped with Ion Torrent™ Product or otherwise agreed by Life Technologies in writing. Life Technologies does not provide service or support for custom products or other products made to buyer's specifications. THE WARRANTIES IDENTIFIED IN THIS CLAUSE ARE LIFE TECHNOLOGIES' SOLE AND EXCLUSIVE WARRANTIES WITH RESPECT TO Ion Torrent™ PRODUCT AND ARE IN LIEU OF ALL OTHER WARRANTIES, STATUTORY, EXPRESS OR IMPLIED, ALL OF WHICH OTHER WARRANTIES ARE EXPRESSLY DISCLAIMED, INCLUDING WITHOUT LIMITATION ANY IMPLIED WARRANTY OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, NON-INFRINGEMENT, OR REGARDING RESULTS OBTAINED THROUGH THE USE OF ANY PRODUCT (INCLUDING, WITHOUT LIMITATION, ANY CLAIM OF INACCURATE, INVALID OR INCOMPLETE RESULTS), WHETHER ARISING FROM A STATUTE OR OTHERWISE IN LAW OR FROM A COURSE OF PERFORMANCE, DEALING OR USAGE OF TRADE.



Safety



WARNING! GENERAL SAFETY. Using this product in a manner not specified in the user documentation may result in personal injury or damage to the instrument or device. Ensure that anyone using this product has received instructions in general safety practices for laboratories and the safety information provided in this document.

- Before using an instrument or device, read and understand the safety information provided in the user documentation provided by the manufacturer of the instrument or device.
- Before handling chemicals, read and understand all applicable Safety Data Sheets (SDSs) and use appropriate personal protective equipment (gloves, gowns, eye protection, and so on). To obtain SDSs, visit thermofisher.com/support.

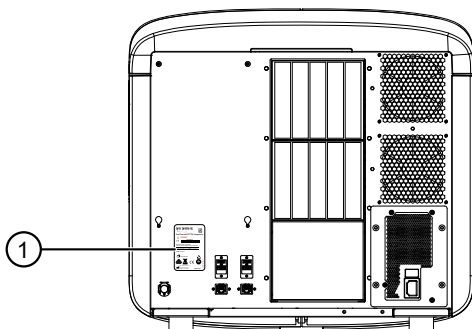
Safety alerts on this instrument

Additional text may be used with one of the symbols described above when more specific information is needed to avoid exposure to a hazard. See the following table for safety alerts found on the instrument.

English	Français
 CAUTION! Hazardous chemicals. Read the Safety Data Sheets (SDSs) before handling.	 MISE EN GARDE ! Produits chimiques dangereux. Lire les fiches signalétiques (FS) avant de manipuler les produits.
 CAUTION! Hazardous waste. Refer to SDS(s) and local regulations for handling and disposal.	 MISE EN GARDE ! Déchets dangereux. Lire les fiches signalétiques (FS) et la réglementation locale associées à la manipulation et à l'élimination des déchets.

Location of safety labels on this instrument

Ion GeneStudio™ S5 Series sequencers have warnings at the location shown:



Ion GeneStudio™ S5 Sequencer labels

Safety information for instruments not manufactured by Thermo Fisher Scientific

Some of the accessories provided as part of the instrument system are not designed or built by Thermo Fisher Scientific. Consult the manufacturer's documentation for the information needed for the safe use of these products.

Instrument safety

General



CAUTION! Do not remove instrument protective covers. If you remove the protective instrument panels or disable interlock devices, you may be exposed to serious hazards including, but not limited to, severe electrical shock, laser exposure, crushing, or chemical exposure.

Cleaning and decontamination



CAUTION! Cleaning and Decontamination. Use only the cleaning and decontamination methods specified in the manufacturer's user documentation. It is the responsibility of the operator (or other responsible person) to ensure the following requirements are met:

- No decontamination or cleaning agents are used that could cause a HAZARD as a result of a reaction with parts of the equipment or with material contained in the equipment.
- The instrument is properly decontaminated a) if hazardous material is spilled onto or into the equipment, and/or b) prior to having the instrument serviced at your facility or sending the instrument for repair, maintenance, trade-in, disposal, or termination of a loan (decontamination forms may be requested from customer service).
- Before using any cleaning or decontamination methods (except those recommended by the manufacturer), users should confirm with the manufacturer that the proposed method will not damage the equipment.



MISE EN GARDE ! Nettoyage et décontamination. Utiliser uniquement les méthodes de nettoyage et de décontamination indiquées dans la documentation du fabricant destinée aux utilisateurs. L'opérateur (ou toute autre personne responsable) est tenu d'assurer le respect des exigences suivantes:

- Ne pas utiliser d'agents de nettoyage ou de décontamination susceptibles de réagir avec certaines parties de l'appareil ou avec les matières qu'il contient et de constituer, de ce fait, un DANGER.
- L'instrument doit être correctement décontaminé a) si des substances dangereuses sont renversées sur ou à l'intérieur de l'équipement, et/ou b) avant de le faire réviser sur site ou de l'envoyer à des fins de réparation, de maintenance, de revente, d'élimination ou à l'expiration d'une période de prêt (des informations sur les formes de décontamination peuvent être demandées auprès du Service clientèle).
- Avant d'utiliser une méthode de nettoyage ou de décontamination (autre que celles recommandées par le fabricant), les utilisateurs doivent vérifier auprès de celui-ci qu'elle ne risque pas d'endommager l'appareil.

Electrical safety



WARNING! Ensure appropriate electrical supply. For safe operation of the instrument:

- Plug the system into a properly grounded receptacle with adequate current capacity.
- Ensure the electrical supply is of suitable voltage.
- Never operate the instrument with the ground disconnected. Grounding continuity is required for safe operation of the instrument.



AVERTISSEMENT ! Veiller à utiliser une alimentation électrique appropriée. Pour garantir le fonctionnement de l'instrument en toute sécurité :

- Brancher le système sur une prise électrique correctement mise à la terre et de puissance adéquate.
- S'assurer que la tension électrique est convenable.
- Ne jamais utiliser l'instrument alors que le dispositif de mise à la terre est déconnecté. La continuité de la mise à la terre est impérative pour le fonctionnement de l'instrument en toute sécurité.



WARNING! Power Supply Line Cords. Use properly configured and approved line cords for the power supply in your facility.



AVERTISSEMENT ! Cordons d'alimentation électrique. Utiliser des cordons d'alimentation adaptés et approuvés pour raccorder l'instrument au circuit électrique du site.



WARNING! Disconnecting Power. To fully disconnect power either detach or unplug the power cord, positioning the instrument such that the power cord is accessible.



AVERTISSEMENT ! Déconnecter l'alimentation. Pour déconnecter entièrement l'alimentation, détacher ou débrancher le cordon d'alimentation. Placer l'instrument de manière à ce que le cordon d'alimentation soit accessible.

Safety and electromagnetic compatibility (EMC) standards

The instrument design and manufacture complies with the following standards and requirements for safety and electromagnetic compatibility.

Safety

Reference	Description
EU Directive 2014/35/EU	European Union “Low Voltage Directive”
IEC 61010-1 UL 61010-1 CSA C22.2 No. 61010-1	<i>Safety requirements for electrical equipment for measurement, control, and laboratory use – Part 1: General requirements</i>
IEC 61010-2-010	<i>Safety requirements for electrical equipment for measurement, control and laboratory use – Part 2-010: Particular requirements for laboratory equipment for the heating of materials</i>

EMC

Reference	Description
Directive 2014/30/EU	European Union “EMC Directive”
EN 61326-1	<i>Electrical Equipment for Measurement, Control and Laboratory Use – EMC Requirements – Part 1: General Requirements</i>
AS/NZS CISPR 22 2009+A1 2010	<i>Limits and Methods of Measurement of Electromagnetic Disturbance Characteristics of Industrial, Scientific, and Medical (ISM) Radiofrequency Equipment</i>
ICES-003, Issue 5	<i>Industrial, Scientific and Medical (ISM) Radio Frequency Generators</i>
FCC Part 15 Subpart B (47 CFR)	<i>U.S. Standard Radio Frequency Devices</i>

Environmental design

Waste and hazardous substances

Reference	Description
Directive 2012/19/EU	European Union “WEEE Directive” – Waste electrical and electronic equipment
Directive 2011/65/EU	European Union “RoHS Directive” – Restriction of hazardous substances in electrical and electronic equipment

Radio compliance

Reference	Description
Directive 2014/53/EU	European Radio Equipment Directive (RED)
RFID	FCC Notice (for U.S. Customers): This device complies with Part 15 of the FCC Rules. Operation is subject to the following conditions: 1. This device may not cause harmful interference, and 2. This device must accept any interference received, including interference that may cause undesired operation Changes and Modifications not expressly approved by Thermo Fisher Scientific can void your authority to operate this equipment under Federal Communications Commissions rules. Canada: <i>This device complies with Industry Canada licence-exempt RSS standard(s).</i> Operation is subject to the following two conditions: (1) this device may not cause interference, and (2) this device must accept any interference, including interference that may cause undesired operation of the device. Le présent appareil est conforme aux CNR d'Industrie Canada applicables aux appareils radio exempts de licence. L'exploitation est autorisée aux deux conditions suivantes : (1) l'appareil ne doit pas produire de brouillage, et (2) l'utilisateur de l'appareil doit accepter tout brouillage radioélectrique subi, même si le brouillage est susceptible d'en compromettre le fonctionnement.

Chemical safety



WARNING! GENERAL CHEMICAL HANDLING. To minimize hazards, ensure laboratory personnel read and practice the general safety guidelines for chemical usage, storage, and waste provided below. Consult the relevant SDS for specific precautions and instructions:

- Read and understand the Safety Data Sheets (SDSs) provided by the chemical manufacturer before you store, handle, or work with any chemicals or hazardous materials. To obtain SDSs, see the “Documentation and Support” section in this document.
- Minimize contact with chemicals. Wear appropriate personal protective equipment when handling chemicals (for example, safety glasses, gloves, or protective clothing).
- Minimize the inhalation of chemicals. Do not leave chemical containers open. Use only with adequate ventilation (for example, fume hood).
- Check regularly for chemical leaks or spills. If a leak or spill occurs, follow the manufacturer's cleanup procedures as recommended in the SDS.
- Handle chemical wastes in a fume hood.
- Ensure use of primary and secondary waste containers. (A primary waste container holds the immediate waste. A secondary container contains spills or leaks from the primary container. Both containers must be compatible with the waste material and meet federal, state, and local requirements for container storage.)
- After emptying a waste container, seal it with the cap provided.
- Characterize (by analysis if necessary) the waste generated by the particular applications, reagents, and substrates used in your laboratory.
- Ensure that the waste is stored, transferred, transported, and disposed of according to all local, state/provincial, and/or national regulations.
- **IMPORTANT!** Radioactive or biohazardous materials may require special handling, and disposal limitations may apply.



AVERTISSEMENT ! PRÉCAUTIONS GÉNÉRALES EN CAS DE MANIPULATION DE PRODUITS CHIMIQUES. Pour minimiser les risques, veiller à ce que le personnel du laboratoire lise attentivement et mette en œuvre les consignes de sécurité générales relatives à l'utilisation et au stockage des produits chimiques et à la gestion des déchets qui en découlent, décrites ci-dessous. Consulter également la FDS appropriée pour connaître les précautions et instructions particulières à respecter :

- Lire et comprendre les fiches de données de sécurité (FDS) fournies par le fabricant avant de stocker, de manipuler ou d'utiliser les matériaux dangereux ou les produits chimiques. Pour obtenir les FDS, se reporter à la section « Documentation et support » du présent document.
- Limiter les contacts avec les produits chimiques. Porter des équipements de protection appropriés lors de la manipulation des produits chimiques (par exemple : lunettes de sûreté, gants ou vêtements de protection).
- Limiter l'inhalation des produits chimiques. Ne pas laisser les récipients de produits chimiques ouverts. Ils ne doivent être utilisés qu'avec une ventilation adéquate (par exemple, sorbonne).
- Vérifier régulièrement l'absence de fuite ou d'écoulement des produits chimiques. En cas de fuite ou d'écoulement d'un produit, respecter les directives de nettoyage du fabricant recommandées dans la FDS.
- Manipuler les déchets chimiques dans une sorbonne.

- Veiller à utiliser des récipients à déchets primaire et secondaire. (Le récipient primaire contient les déchets immédiats, le récipient secondaire contient les fuites et les écoulements du récipient primaire. Les deux récipients doivent être compatibles avec les matériaux mis au rebut et conformes aux exigences locales, nationales et communautaires en matière de confinement des récipients.)
- Une fois le récipient à déchets vidé, il doit être refermé hermétiquement avec le couvercle fourni.
- Caractériser (par une analyse si nécessaire) les déchets générés par les applications, les réactifs et les substrats particuliers utilisés dans le laboratoire.
- Vérifier que les déchets sont convenablement stockés, transférés, transportés et éliminés en respectant toutes les réglementations locales, nationales et/ou communautaires en vigueur.
- **IMPORTANT !** Les matériaux représentant un danger biologique ou radioactif exigent parfois une manipulation spéciale, et des limitations peuvent s'appliquer à leur élimination.

Documentation and support

Related documentation

Title	Pub. No.
<i>Ion 510™ & Ion 520™ & Ion 530™ Kit – Chef User Guide</i>	MAN0016854
<i>Ion 540™ Kit – Chef User Guide</i>	MAN0010851
<i>Ion 550™ Kit – Chef User Guide</i>	MAN0017275
<i>Ion 520™ & Ion 530™ ExT Kit – Chef User Guide</i>	MAN0015805
<i>Ion 520™ & Ion 530™ Kit – OT2 User Guide</i>	MAN0010844
<i>Ion 540™ Kit – OT2 User Guide</i>	MAN0010850
<i>Ion GeneStudio™ S5 System Site Preparation Guide</i>	MAN0017529

Customer and technical support

Visit thermofisher.com/support for the latest service and support information.

- Worldwide contact telephone numbers
- Product support information
 - Product FAQs
 - Software, patches, and updates
 - Training for many applications and instruments
- Order and web support
- Product documentation
 - User guides, manuals, and protocols
 - Certificates of Analysis
 - Safety Data Sheets (SDSs; also known as MSDSs)

Note: For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

Limited product warranty

Life Technologies Corporation and its affiliates warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have questions, contact Life Technologies at www.thermofisher.com/support.

