

1. IDENTIFICATION OF THE SUBSTANCE/PREPARATION AND THE COMPANY/UNDERTAKING

Product Code 401446
Product Name Premix Buffer Concentrate
Trade Name(s) Not Applicable

Recommended Use For research use only. Not for use in diagnostic procedures.

Contact Manufacturer
APPLIED BIOSYSTEMS
7 KINGSLAND GRANGE, WOOLSTON
WARRINGTON, CHESHIRE WA1 4SR
Telephone: +44 1925 82 56 50 (UK)
Fax: +44 1925 28 27 02

Emergency Telephone Number CHEMTREC: +1-703-527-3887 (INTERNATIONAL)
1-800-424-9300 (NORTH AMERICA)

2. COMPOSITION/INFORMATION ON INGREDIENTS

Chemical nature of the preparation Mixture of organic compounds.

Haz/Non-haz

Chemical Name	EINECS-No.	CAS-No	Weight %	Classification
Acetic acid	200-580-7	64-19-7	10-30	R10 C;R35

For the full text of the R phrases mentioned in this Section, see Section 16.

3. HAZARDS IDENTIFICATION

The preparation is classified as dangerous in accordance with Directive 1999/45/EC.

Classification

Xi;R36/38
Xi - Irritant.

Most Important Hazards

Irritating to eyes and skin..

4. FIRST AID MEASURES

General advice If symptoms persist, call a physician. Show this safety data sheet to the doctor in attendance.

Eye contact Rinse immediately with plenty of water, also under the eyelids, for at least 15 minutes. If symptoms persist, call a physician.

Skin contact Wash off immediately with soap and plenty of water for at least 15 minutes while removing all contaminated clothing and shoes. If symptoms persist, call a physician.

Ingestion	Rinse mouth. Drink plenty of water. Do NOT induce vomiting. Never give anything by mouth to an unconscious person. Consult a physician.
Inhalation	Move to fresh air. If symptoms persist, call a physician. If breathing is irregular or stopped, administer artificial respiration.
Notes to physician	Treat symptomatically.
Protection of first-aiders	Use personal protective equipment.

5. FIRE-FIGHTING MEASURES

Suitable extinguishing media

Water spray. Carbon dioxide (CO₂). Foam. Dry powder.

Extinguishing media which must not be used for safety reasons

No information available.

Specific hazards

Keep product and empty container away from heat and sources of ignition.

Special protective equipment for firefighters

Wear self-contained breathing apparatus and protective suit.

6. ACCIDENTAL RELEASE MEASURES

Environmental precautions	Do not flush into surface water or sanitary sewer system. Prevent further leakage or spillage if safe to do so.
Personal precautions	Use personal protective equipment. Avoid contact with the skin and the eyes. Evacuate personnel to safe areas. Keep people away from and upwind of spill/leak.
Methods for cleaning up	Dam up. Soak up with inert absorbent material (e.g. sand, silica gel, acid binder, universal binder, sawdust). Sweep up and shovel into suitable containers for disposal. After cleaning, flush away traces with water.
Methods for Containment	Prevent further leakage or spillage if safe to do so.
Other information	Refer to protective measures listed in Sections 7 and 8.

7. HANDLING AND STORAGE

Handling	Ensure adequate ventilation. Wear personal protective equipment. Do not breathe vapours or spray mist. Avoid contact with skin, eyes and clothing. Remove and wash contaminated clothing before re-use. Handle in accordance with good industrial hygiene and safety practice.
Storage	Keep container tightly closed. Keep out of the reach of children. Keep in properly labelled containers. Keep away from heat and sources of ignition. Do not store near combustible materials.

8. EXPOSURE CONTROLS / PERSONAL PROTECTION

Exposure Limits

Chemical Name	EU	The United Kingdom	France	Spain	Germany
Acetic acid 64-19-7	TWA: 10 ppm TWA: 25 mg/m ³	TWA: 10 ppm STEL: 15 ppm TWA: 25 mg/m ³ STEL: 37 mg/m ³	VLCT: 10 ppm VLCT: 25 mg/m ³	VLA-EC: 15 ppm VLA-EC; 37 mg/m ³ VLA-EC VLA-ED: 10 ppm VLA-ED; 25 mg/m ³ VLA-ED	

Chemical Name	Italy	Portugal	The Netherlands	Finland	Denmark
Acetic acid 64-19-7		STEL: 15 ppm TWA: 10 ppm		TWA: 13 mg/m ³ TWA: 5 ppm STEL: 10 ppm STEL: 25 mg/m ³	TWA: 10 ppm TWA: 25 mg/m ³

Chemical Name	Austria	Switzerland	Poland	Norway	Ireland
Acetic acid 64-19-7	STEL: 20 ppm STEL; 50 mg/m ³ STEL MAK: 10 ppm MAK; 25 mg/m ³ MAK	STEL: 20 ppm STEL; 50 mg/m ³ STEL MAK: 10 ppm MAK; 25 mg/m ³ MAK	NDSch: 30 mg/m ³ NDS: 15 mg/m ³	TWA: 10 ppm	TWA: 10 ppm TWA: 25 mg/m ³ STEL: 15 ppm STEL: 37 mg/m ³

Occupational Exposure Controls

Engineering measures

Showers, eyewash stations, and ventilation systems.

Personal protective equipment

Respiratory protection

In case of inadequate ventilation wear respiratory protection.

Eye protection

Face-shield.

Skin and body protection

Wear protective gloves/clothing.

Hand protection

Protective gloves.

Other protective equipment

Eye wash bottle with pure water

Hygiene measures

Remove and wash contaminated clothing before re-use. Regular cleaning of equipment, work area and clothing. Wash hands before breaks and immediately after handling the product.

Environmental exposure controls

Do not allow material to contaminate ground water system. Prevent product from entering drains.

9. PHYSICAL AND CHEMICAL PROPERTIES

Appearance	Colourless.	Odour	Irritating, Pungent.
Physical state	Liquid.	pH	No information available
Flash point	Not flammable.	Autoignition temperature	No information available.
Boiling point/range	No information available.	Melting point/range	No information available.
Flammability Limits in Air	No information available.	Water solubility	Soluble in water.
Specific Gravity	No information available.	Evaporation rate	No information available.
Solubility	No information available	Vapour density	No information available.
Vapour pressure	No information available.	Viscosity	No information available.
VOC content	40		

10. STABILITY AND REACTIVITY

10. STABILITY AND REACTIVITY

Stability	Stable under normal conditions.
Conditions to avoid	Heat, flames and sparks.
Materials to avoid	Incompatible with oxidizing agents. Acids. Combustible materials.
Hazardous decomposition products	Sulfur oxides.

11. TOXICOLOGICAL INFORMATION

Acute Toxicity

Product Information Product does not present an acute toxicity hazard based on known or supplied information..

Chronic Toxicity

Irritation	Irritating to eyes and skin.
Reproductive Toxicity	No information available.
Developmental Toxicity	No information available.
Target Organ Effects	Eyes. Respiratory system. Skin. Teeth.

12. ECOLOGICAL INFORMATION

Ecotoxicity

Ecotoxicity effects of component substances.

Chemical Name	Toxicity to algae	Toxicity to fish	Microtox Data	Daphnia magna (Water flea)
Acetic acid		LC50= 75 mg/L Lepomis macrochirus 96 h LC50= 88 mg/L Pimephales promelas 96 h	EC50 = 8.8 mg/L 15 min EC50 = 8.8 mg/L 25 min EC50 = 8.8 mg/L 5 min	EC50 = 95 mg/L 24 h

Persistence and degradability	No information available.
Bioaccumulative potential	Product has a low potential to bioconcentrate.
Mobility	The product contains volatile organic compounds (VOC) which will evaporate easily from all surfaces. The product is water soluble, and may spread in water systems.

Chemical Name	log Pow
Acetic acid	-0.31

13. DISPOSAL CONSIDERATIONS

Waste from residues / unused products	Dispose of as hazardous waste in compliance with local and national regulations.
Contaminated packaging	Empty containers should be taken for local recycling, recovery or waste disposal.
EWC waste disposal No	07 07 04 16 05 06

Other information According to the European Waste Catalogue, Waste Codes are not product specific, but application specific. Waste codes should be assigned by the user based on the application for which the product was used.

14. TRANSPORT INFORMATION

IMDG/IMO

Proper shipping name Acetic acid, solution (Mixture)
Hazard Class 8
UN-No UN2790
Packing group III
EmS F-A, S-B
Description UN2790, Acetic acid, solution,8,PG III,Mixture

RID

Proper shipping name Acetic acid, solution (Mixture)
Hazard Class 8
UN-No UN2790
Packing group III
Classification Code C3
Description UN2790 Acetic acid, solution,8,III,RID,Mixture
ADR/RID-Labels 8

ADR

Proper shipping name Acetic acid, solution (Mixture)
Hazard Class 8
UN-No UN2790
Packing group III
Classification Code C3
Description UN2790 Acetic acid, solution,8,III,ADR,Mixture
ADR/RID-Labels 8

ICAO

UN-No UN2790
Proper shipping name Acetic acid solution (Mixture)
Hazard Class 8
Packing group III
Description Acetic acid solution,8,UN2790,PG III,Mixture

ICAO/IATA

UN-No UN2790
Proper shipping name Acetic acid solution (Mixture)
Hazard Class 8
Packing group III
ERG Code 8L
Description UN2790,Acetic acid solution,8,PG III,Mixture

15. REGULATORY INFORMATION

The preparation is classified as dangerous in accordance with Directive 1999/45/EC.

Labelling

Symbol(s) Xi - Irritant.

Xi

**R-phrase(s)**

R36/38 - Irritating to eyes and skin.

S-phrase(s)

S26 - In case of contact with eyes, rinse immediately with plenty of water and seek medical advice.

S37 - Wear suitable gloves.

S23 - Do not breathe gas/fumes/vapour/spray.

S60 - This material and its container must be disposed of as hazardous waste.

International Inventories

Chemical Name	EINECS	ELINCS	TSCA	DSL	NDSL	PICCS	ENCS	China	AICS	KECL
Acetic acid	200-580-7	-	Present	X	-	X	X	X	X	KE-00013

16. OTHER INFORMATION

Text of R phrases mentioned in Section 2

R35 - Causes severe burns.

R10 - Flammable.

Issuing date	23-Apr-2007
---------------------	-------------

Revision Date	29-Jun-2007
----------------------	-------------

Revision Note	No information available
----------------------	--------------------------

Disclaimer

The information provided on this SDS is correct to the best of our knowledge, information and belief at the date of its publication. The information given is designed only as a guide for safe handling, use, processing, storage, transportation, disposal and release and is not to be considered as a warranty or quality specification. The information relates only to the specific material designated and may not be valid for such material used in combination with any other material or in any process, unless specified in the text.

End of Safety Data Sheet