

Agilent 8697 Headspace Samplers

Safety Manual



Notices

© Agilent Technologies, Inc. 2023

No part of this manual may be reproduced in any form or by any means (including electronic storage and retrieval or translation into a foreign language) without prior agreement and written consent from Agilent Technologies, Inc. as governed by United States and international copyright laws.

Manual Part Number

G4511-90000

Edition

Second edition, April 2023

First edition, February 2021

Printed in USA

Agilent Technologies, Inc.
2850 Centerville Road
Wilmington, DE 19808-1610 USA

安捷伦科技（上海）有限公司
上海市浦东新区外高桥保税区
英伦路 412 号
联系电话：（800） 820 3278

Warranty

The material contained in this document is provided “as is,” and is subject to being changed, without notice, in future editions. Further, to the maximum extent permitted by applicable law, Agilent disclaims all warranties, either express or implied, with regard to this manual and any information contained herein, including but not limited to the implied warranties of merchantability and fitness for a particular purpose. Agilent shall not be liable for errors or for incidental or consequential damages in connection with the furnishing, use, or performance of this document or of any information contained herein. Should Agilent and the user have a separate written agreement with warranty terms covering the material in this document that conflict with these terms, the warranty terms in the separate agreement shall control.

Technology Licenses

The hardware and/or software described in this document are furnished under a license and may be used or copied only in accordance with the terms of such license.

Restricted Rights Legend

U.S. Government Restricted Rights. Software and technical data rights granted to the federal government include only those rights customarily provided to end user customers. Agilent provides this customary commercial license in Software and technical data pursuant to FAR 12.211 (Technical Data) and 12.212 (Computer Software) and, for the Department of Defense, DFARS 252.227-7015 (Technical Data -Commercial Items) and DFARS 227.7202-3 (Rights in Commercial Computer Software or Computer Software Documentation).

Safety Notices

CAUTION

A CAUTION notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in damage to the product or loss of important data. Do not proceed beyond a CAUTION notice until the indicated conditions are fully understood and met.

WARNING

A WARNING notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in personal injury or death. Do not proceed beyond a WARNING notice until the indicated conditions are fully understood and met.

Introduction

Important Safety Warnings	4
Hydrogen Safety	8
Fume Hood	9
Disconnect Device	9
Magnets	10
Safety and Regulatory Certifications	11
Intended Use	16
Cleaning and Disinfecting	16
Recycling the Product	18

Important Safety Warnings

Before moving on, there are several important safety notices that you should always keep in mind when using the Agilent 8697.

WARNING

When handling/using chemicals for preparation or use within the headspace sampler, all applicable local and national laboratory safety practices must be followed. This would include, but is not limited to, correct use of Personal Protective Equipment (PPE), correct use of storage vials, using vent hoods, and correct handling of chemicals, as defined in the laboratory's internal safety analysis and standard operating procedures. Failure to adhere to laboratory safety practices could lead to injury or death.

Many internal parts of the headspace sampler carry dangerous voltages

If the headspace sampler is connected to a power source, even if the power switch is set to standby, potentially dangerous voltages exist on:

- The wiring between the Headspace Sampler power cord and the AC power supply, the AC power supply itself, and the wiring from the AC power supply to the power switch.

With the power switch on, potentially dangerous voltages also exist on:

- All electronics boards in the instrument.
- The internal wires and cables connected to these boards.
- The wires for the oven heater.

WARNING

All these parts are shielded by covers. With the covers in place, it is difficult to accidentally make contact with dangerous voltages. Unless specifically instructed to, never remove a cover unless the instrument is unplugged.

WARNING

If the power cord insulation is frayed or worn, the cord must be replaced. Contact your Agilent service representative.

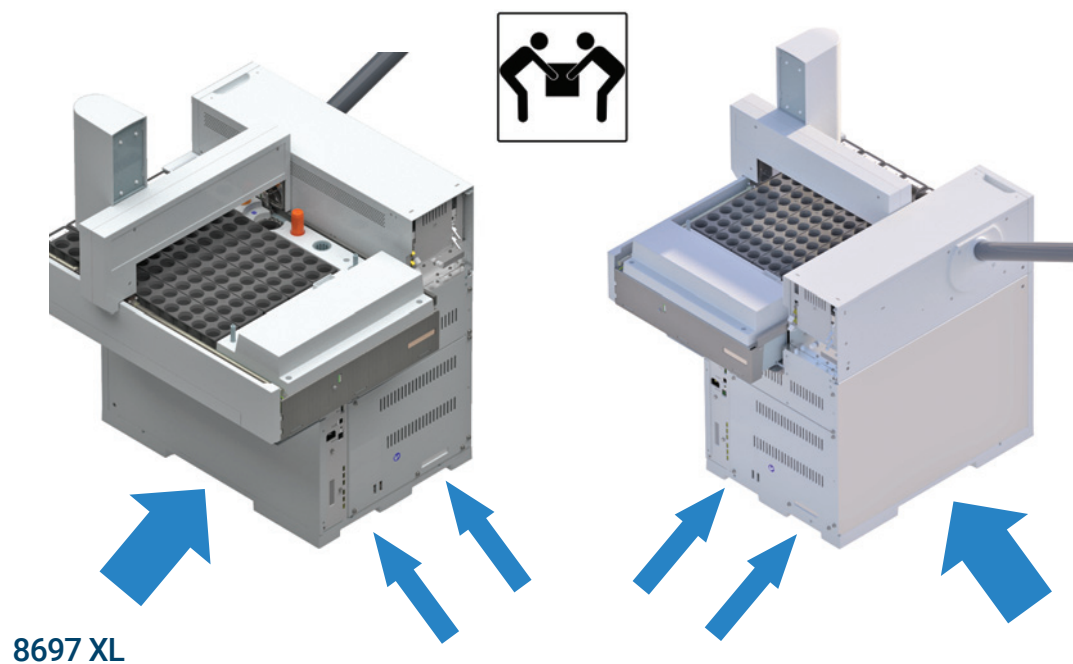
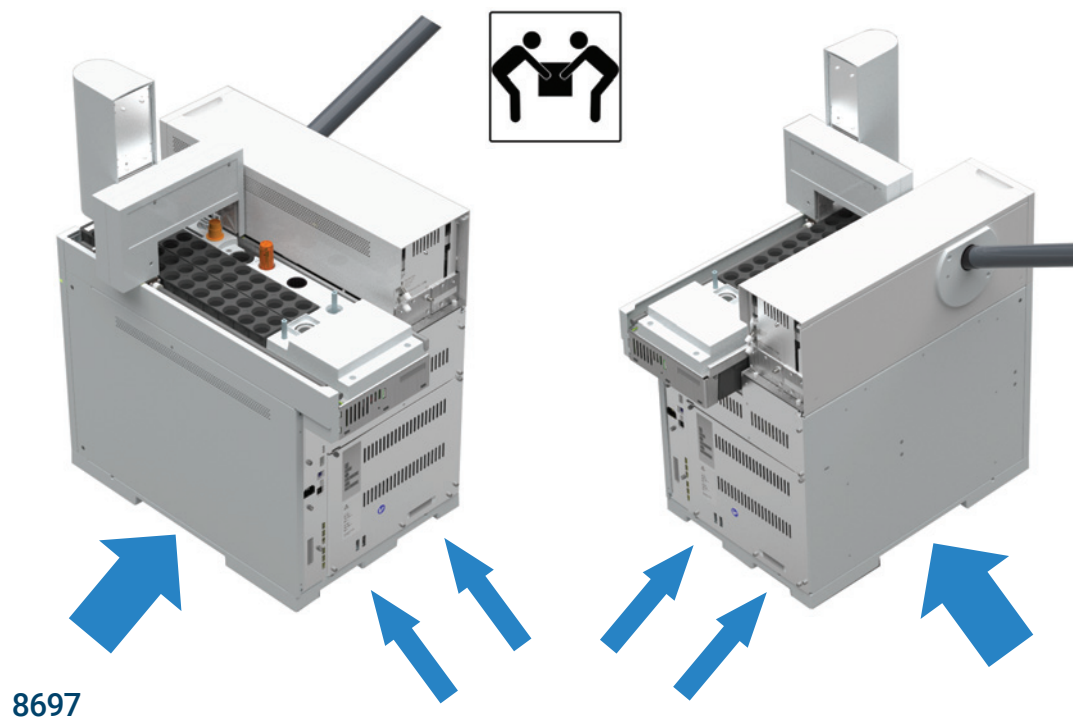
WARNING

This is a Safety Class 1 Product (provided with a protective earthing ground, incorporated in the power cord). The mains plug shall only be inserted in a socket outlet provided with a protective earth contact. Any interruption of the protective conductor inside or outside of the instrument is likely to make the instrument dangerous. Intentional interruption is prohibited.

Lifting

WARNING

The headspace sampler is heavy. To avoid injury, use a two-person lift.



CAUTION

Make sure that all samples and trays have been removed prior to moving or relocating the headspace sampler. Prior to moving, all covers and enclosures must also be seated properly. Make sure all covers and enclosures are secure.

CAUTION

Make sure your headspace sampler has adequate distance from the wall and other instruments.

Electrostatic discharge is a threat to headspace sampler electronics

The printed circuit (PC) boards in the headspace sampler can be damaged by electrostatic discharge. Do not touch any of the boards unless it is absolutely necessary. If you must handle them, wear a grounded wrist strap and take other antistatic precautions. Wear a grounded wrist strap any time you must remove the side cover.

Many parts are dangerously hot

Many parts of the headspace sampler operate at temperatures high enough to cause serious burns. These parts include but are not limited to:

- The oven carousel/tray and its contents
- The probe and heated blocks
- The six port valve and heated blocks
- Vials recently removed from the heated oven

You should always cool these areas of the headspace sampler to room temperature before working on them. They will cool faster if you first set the temperature of the heated zone to room temperature. Turn the zone off after it has reached the setpoint. If you must perform maintenance on hot parts, use a wrench and wear thermally protective gloves.

Whenever possible, cool the instrument part that you will be maintaining before you begin working on it.

Oven thermal leaks

WARNING

Objects passing through the oven lid seal can cause thermal leaks which create hazardous hot spots which cause burns and melt equipment.

Do not allow wiring or temperature probes to pass through the oven lid seal.

Never use a flammable gas for vial pressurization

Flammable gases, such as hydrogen and argon/methane, create a potential explosion hazard during vial pressurization and in venting. The Agilent 8897 Headspace Sampler cannot be configured to use a flammable gas for vial pressurization.

General gas safety

- Wear eye protection when using compressed gas to avoid eye injury.
- Fasten all compressed gas cylinders securely to an immovable structure or permanent wall.
- Store and handle compressed gases in accordance with relevant safety codes.
- Do not put gas cylinders in the path of a hot air vent (including a GC oven exhaust).

General warnings

- Perform periodic leak checks on supply lines, fittings, and pneumatic plumbing to prevent a potentially hazardous condition.
- To avoid a potential shock hazard when using liquid solution to locate leaks, turn the main power switch to standby and disconnect the main power cord. Be careful not to spill leak solution on electrical leads.
- Turn off all HS and GC gas supplies at their sources before performing maintenance.

Use with chemicals

When handling/using chemicals for preparation or use within the instrument, follow all applicable local and national laboratory safety practices. This includes, but is not limited to, correct use of PPE, correct use of storage vials, and correct handling of chemicals, as defined in the laboratory's internal safety analysis and standard operating procedures. Failure to adhere to laboratory safety practices could lead to injury or death.

Hydrogen Safety

Hydrogen gas may be used as carrier gas. When mixed with air, hydrogen can form explosive mixtures.

WARNING

When using hydrogen (H₂) as the carrier gas, be aware that hydrogen gas can flow into the headspace instrument or GC oven and create an explosion hazard. Therefore, be sure that the hydrogen supply is turned off until all connections are made. When using H₂, make sure the sample loop is properly installed and that the transfer line is connected to the GC before supplying hydrogen gas to the instrument. Hydrogen is flammable. Leaks, when confined in an enclosed space, may create a fire or explosion hazard. In any application using hydrogen, leak test all connections, lines, and valves before operating the instrument. Always turn off the hydrogen supply at its source before working on the instrument.

WARNING

Never use flammable gas for vial pressurization. Flammable gases, such as hydrogen and argon/methane, can create a potential fire hazard or explosion hazard when used for vial pressurization. The 8897 Headspace Sampler does not support use of flammable gases for vial pressurization.

Hydrogen is a commonly used GC carrier gas. Hydrogen is potentially explosive and has other dangerous characteristics.

- Hydrogen is combustible over a wide range of concentrations. At atmospheric pressure, hydrogen is combustible at concentrations from 4% to 74.2% by volume.
- Hydrogen has the highest burning velocity of any gas.
- Hydrogen has a very low ignition energy.
- Hydrogen that is allowed to expand rapidly from high pressure into the atmosphere can self-ignite due to an electrostatic spark.
- Hydrogen burns with a nonluminous flame which can be invisible under bright light.

Fume Hood

During normal operation of the headspace sampler, some sample can exhaust from the sample vial out the headspace sampler vent. If any sample components are toxic or noxious, or if hydrogen is used as the carrier gas, these exhausts must be vented to a fume hood. If the Headspace is in an enclosed, small space then a fume hood should be used regardless of the Headspace gases connected. Place the headspace sampler in the hood or attach a large diameter venting tube to the outlet for proper ventilation.

Disconnect Device

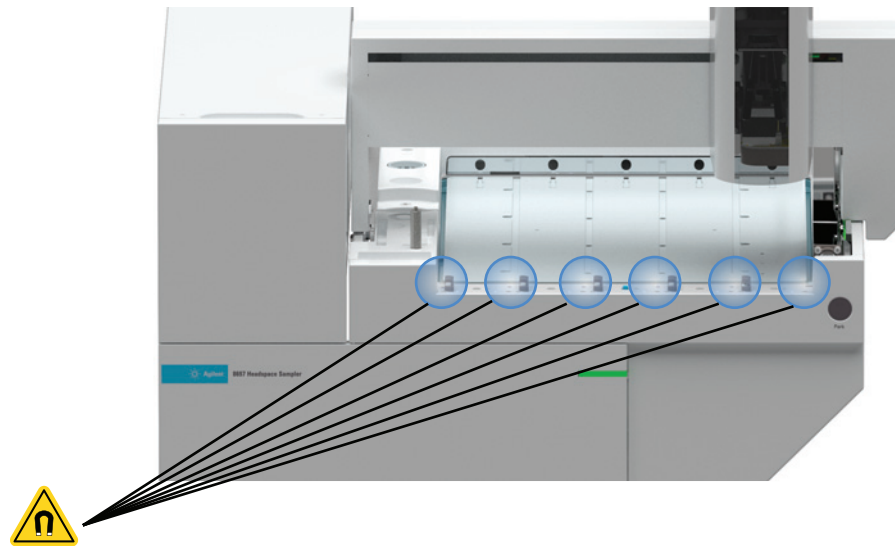
WARNING

The appliance inlet coupler (main input power cord) is the power disconnect device. Do not position the instrument such that access to the coupler or plug is impaired.

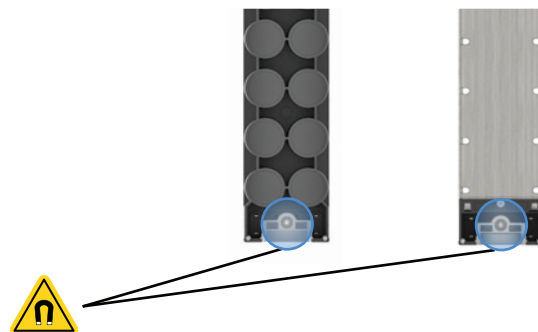
WARNING

The front power switch does not disconnect power for the instrument. Pressing the disconnect switch places the instrument in a standby mode. To avoid an electric shock hazard, disconnect the device by unplugging the power cord before performing instrument service.

Magnets



Magnets in vial tray base (8697 XL shown, 8697 is similar)



Magnets in bottom front of vial racks



Magnetic Field. This product contains magnets, which may affect the functioning of pacemakers, implanted heart defibrillators, or other active implanted medical devices (AIMD). If you wear an AIMD keep at least a 30 mm (1.2 inch) distance from the magnets and follow the AIMD manufacturer guidance when using this product.

Safety and Regulatory Certifications

The Agilent 8697 Headspace Sampler conforms to the following safety standards:

- Canadian Standards Association (CSA): C22.2 No. 61010-1
- CSA/Nationally Recognized Test Laboratory (NRTL): ANSI/UL 61010-1
- International Electrotechnical Commission (IEC): 61010-1, 61010-2-010, 61010-2-081
- EuroNorm (EN): 61010-1

The Agilent 8697 Headspace Sampler conforms to the following regulations on Electromagnetic Compatibility (EMC) and Radio Frequency Interference (RFI):

- CISPR 11/EN 55011: Group 1, Class A

Group 1 ISM equipment: group 1 contains all Industrial, Scientific and Medical (ISM) equipment in which there is intentionally generated and/or used conductively coupled radio-frequency energy which is necessary for the internal functioning of the equipment itself.

Class A equipment is equipment suitable for use in all establishments other than domestic and those directly connected to a low voltage power supply network which supplies buildings used for domestic purposes.

- IEC/EN 61326-1
-  AUS/NZ

This ISM device complies with Canadian ICES-001. Cet appareil ISM est conforme a la norme NMB-001 du Canada.

- CAN ICES-1/NMB-1, ISM 1-A

The Agilent 8697 Headspace Sampler is designed and manufactured under a quality system registered to ISO 9001. Declaration of Conformity available.

This product complies with the EU RoHS Directive 2011/65/EU and conforms to EN 50581.



CE Compliance

Your Agilent instrument has been designed to comply with the requirements of the applicable directives of the European Union, such as Electromagnetic Compatibility (EMC) Directive, Low Voltage Directive (LVD), Machinery Directive (MD), RoHS Directive, etc. Agilent has confirmed that each product complies with the relevant Directives by testing samples against the harmonized EN (European Norm) standards published on the Official Journal of the European Union (OJEU).

Proof that a product complies with these directives is indicated by:

- The CE Marking appearing on the rear of the product, and

- The documentation package that accompanies the product containing a copy of the Declaration of Conformity. The Declaration of Conformity is the legal declaration by Agilent that the product complies with the relevant directives listed above, and it shows the EN standards to which the product was tested to demonstrate compliance.

UK Compliance



Your Agilent instrument has been designed to comply with the requirements of the applicable regulations of the United Kingdom, such as The Electromagnetic Compatibility Regulations 2016, The Electrical Equipment (Safety) Regulations 2016, The Supply of Machinery (Safety) Regulations 2008, The Restriction of the Use of Certain Hazardous Substances in Electrical and Electronic Equipment Regulations 2012, etc.

Agilent has confirmed that each product complies with the relevant Regulations by testing samples against the designated standards published on GOV.UK.

Proof that a product complies with these regulations is indicated by:

- the UK CA Marking appearing on the rear of the product, and
- the documentation package that accompanies the product containing a copy of the Declaration of Conformity. The Declaration of Conformity is the legal declaration by Agilent that the product complies with the relevant regulations listed above, and shows the designated standards to which the product was tested to demonstrate compliance.

Electromagnetic compatibility

This device complies with the requirements of CISPR 11, Group 1, Class A. Operation is subject to the following two conditions:

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

If this equipment does cause harmful interference to radio or television reception, which can be determined by turning the equipment off and on, the user is encouraged to try one or more of the following measures:

- 1 Relocate the radio or antenna.
- 2 Move the device away from the radio or television.
- 3 Plug the device into a different electrical outlet, so that the device and the radio or television are on separate electrical circuits.
- 4 Make sure that all peripheral devices are also certified.
- 5 Make sure that appropriate cables are used to connect the device to peripheral equipment.
- 6 Consult your equipment dealer, Agilent Technologies, or an experienced technician for assistance.

Changes or modifications not expressly approved by Agilent Technologies could void the user's authority to operate the equipment.

EMC Declaration for South Korea

사용자안내문

이 기기는 업무용 환경에서 사용할 목적으로 적합성평가를 받은 기기로서 가정용 환경에서 사용하는 경우 전파간섭의 우려가 있습니다 .

※ 사용자 안내문은 " 업무용 방송통신기자재 " 에만 적용한다 .

This equipment has been evaluated for its suitability for use in a commercial environment. When used in a domestic environment, there is a risk of radio interference.

Detachable Power Cord Declaration for Japan

(弊社提供の電源コードセットが汎用性がない旨を示す)

電源コードセットの取扱いについて (日本国内向け)

製品には、同梱された電源コードセットをお使いください。同梱された電源コードセットは、他の製品では使用できません。

Notice - Power Cords for Japanese Market

Your product must only use the power cord that was shipped with this product. Do not use this power cord with any other product.

Information

The Agilent 8697 Headspace Sampler meets the following IEC (International Electro-technical Commission) classifications: Safety Class 1, Transient Overvoltage Category II, Pollution Degree 2.

This unit has been designed and tested in accordance with recognized safety standards and is designed for use indoors in non-classified locations. If the instrument is used in a manner not specified by the manufacturer, the protection provided by the instrument may be impaired. Whenever the safety protection of the Agilent 8697 Headspace Sampler has been compromised, disconnect the unit from all power sources and secure the unit against unintended operation.

Refer servicing to qualified service personnel. Substituting parts or performing any unauthorized modification to the instrument may result in a safety hazard.

Symbols

The apparatus is marked with this symbol when the user should refer to the instruction manual in order to protect risk of harm to the operator and to protect the apparatus against damage.



Indicates a hot surface. Keep hands, electrical cables, gas lines, chart paper, combustible material and other items safely clear of the hot surface.



Indicates hazardous voltages.



Indicates earth (ground) terminal.



Indicates potential explosion hazard.



Indicates radioactivity hazard.



Indicates electrostatic discharge hazard.



Indicates a hazard. See the Agilent Headspace Sampler user documentation for the item labeled.



or



Indicates that you must not discard this electrical/electronic product in domestic household waste



To prevent personal injury, two person lift recommended



Manufacturing date.



Power symbol indicates **On/Standby**. The apparatus is not completely disconnected from the main supply when the power switch is in the **Standby** position.



Warnings in the manual or on the instrument must be observed during all phases of operation, service, and repair of this instrument. Failure to comply with these precautions violates safety standards of design and the intended use of the instrument. Agilent Technologies assumes no liability for the customer's failure to comply with these requirements.

Sound Emission Certification for Federal Republic of Germany

Sound pressure

Sound pressure $L_p < 70 \text{ dB(A)}$ according to DIN-EN 7779.

Schalldruckpegel

Schalldruckpegel $LP < 70 \text{ dB(A)}$ nach DIN-EN 7779.

Intended Use

Agilent products must only be used in the manner described in the Agilent product user guides. Any other use may result in damage to the products or personal injury. Agilent is not responsible for any damages caused, in whole or in part, by improper use of the products, unauthorized alterations, adjustments or modifications to the products, failure to comply with procedures in Agilent product user guides, or use of the products in violation of applicable laws, rules or regulations.

Cleaning and Disinfecting

The United States Centers for Disease Control and Prevention (CDC) lists a variety of chemicals and commercial products for disinfection, however, some of these are not suitable for cleaning Agilent instruments. A compatible, CDC-recommended cleaning solution for external instrument surfaces is a solution of 70% isopropanol (isopropyl alcohol) and 30% water. This solution is commonly sold in stores as rubbing alcohol.

Concentrations of isopropanol above 70% may damage the instrument.

WARNING

When working with solvents observe appropriate safety procedures (for example, goggles, safety gloves and protective clothing) as described in the safety data sheet supplied by the solvent vendor, especially when toxic or hazardous solvents and flammable liquids are used. For this cleaning procedure, the appropriate personal protective equipment (PPE) consists of nitrile or similar gloves, safety glasses, and a laboratory coat.

General procedure

Use the following procedure to clean often-touched external surfaces of Agilent instruments.

NOTE

Any thermal zone on an instrument is likely to already be decontaminated, provided it was operated above 121 °C for at least one hour. For example, on a gas chromatograph, the inlets, oven, columns, transfer lines, and detectors can be decontaminated by setting the temperature zones to 200 °C for at least one hour. To avoid burns, cool these temperature zones to room temperature before accessing.

- 1 Before beginning, establish the current baseline performance as appropriate. For example, if using an MS perform an autotune and generate a report.
- 2 Put on the appropriate PPE.
- 3 Cool all thermal zones and allow them to reach room temperature.
- 4 Wait until the instrument is cool, then turn off the instrument and disconnect the power cord. Never clean an instrument while it is powered on or plugged in. Never clean an instrument while it is hot.
- 5 Remove any external devices, if necessary, to allow cleaning access to all potentially contaminated surfaces. For example, remove an ALS and clean separately. (If desired you can disconnect a GC or LC from a mass spectrometer or other connected instrument as

described in its user documentation to allow cleaning between the instruments. Be sure to use all of the recommended procedures and two persons if it is necessary to move the instruments.)

- 6 If needed, create an amount of 70% isopropanol/30% water solution that is sufficient to clean the instruments.

WARNING

Never spray any liquids directly onto the product.

- 7 Moisten a clean, lint-free cloth with the mixture of 70% isopropanol/30% water.
 - Do not use a paper towel or other fibrous material.
 - The cloth should be moist, not dripping wet.
- 8 Gently wipe external surfaces to be cleaned using the moistened cloth. Do not allow any liquid to drip into the instrument. Many areas inside Agilent instruments contain electrical wiring, printed circuit boards, and other parts that can be damaged by liquids or potentially create an electrical safety hazard when wet.
 - Start with the top and side surfaces, and end with external cabling, external gas supply lines, and transfer lines.
 - Wipe the outer surface of cables and their connectors but do not touch the electrical connections (for example, pins and plugs inside the cable or connection port).
 - When cleaning a display or touchscreen, carefully wipe in one direction, moving from the top to the bottom.
- 9 Allow all surfaces to completely air-dry. No moisture should be present on the instrument surfaces.
- 10 Plug in the instruments. If separated, return instruments to their original locations on the lab bench, and restore to normal operating conditions as described in the instrument user documentation.
- 11 Verify performance after reassembly, as appropriate.
- 12 Discard the gloves and other personal protective equipment or clean them in an approved process, then wash your hands.

Refer to the instrument documentation for stabilization times after turning everything back on.

You may notice some visible changes to cosmetic finishes over time as a result of the cleaning techniques mentioned in this document.

Other cleaners

While there are many other common household cleaners and disinfectants recommended by the CDC for disinfecting surfaces, many of these can damage painted surfaces and electronic products. Do not use: bleach (including sodium hypochlorite), peroxides (including hydrogen peroxide), acetone, ammonia, quaternary ammonium-based cleaners, ethyl alcohol, methylene chloride, sodium dichloroisocyanurate (sodium dichloro-S-triazinetriene), or any petroleum-based solvents (for example gasoline, paint thinner, benzene, or toluene). This includes wipes, liquids, and sprays. Read the product labels.

Recycling the Product



Instructions for Disposal of Waste Equipment by Users in the European Union. This symbol on the product or its packaging indicates that this product must not be disposed of with other waste. Instead, it is your responsibility to dispose of your waste equipment by handing it over to a designated collection point for the recycling of waste electrical and electronic equipment. The separate collection and recycling of your waste equipment at the time of disposal will help conserve natural resources and ensure that it is recycled in a manner that protects human health and the environment. For more information about where you can drop off your waste equipment for recycling, please contact your local city recycling office or the dealer from whom you originally purchased the product.

For recycling, contact your local Agilent sales office.