

# **5973N and 5973 inert Mass Selective Detectors**

## **Site Preparation Guide**



**Agilent Technologies**

## Notices

© Agilent Technologies, Inc. 2003

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

G2589-90070

### Edition

First edition, August 2003

Replaces part number G2589-90010

Printed in USA

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

### Acknowledgements

Microsoft® and Windows® are U.S. registered trademarks of Microsoft Corporation

## 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.**

## 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.

---

# Contents

## 1 General Information

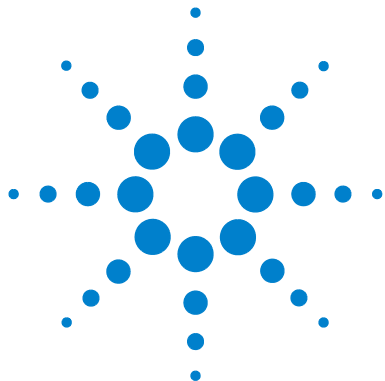
Instrument Identification	6
5973N MSD version	6
5973 inert MSD version	6
Important Safety Warnings	7
Many internal parts of the MSD carry dangerous voltages	7
Electrostatic discharge is a threat to MSD electronics	8
Many parts are dangerously hot	8
Hydrogen	9
Chemical safety	10
Safety and Regulatory Certifications	11
Information	11
Symbols	12
Electromagnetic compatibility	13
Sound emission certification for Federal Republic of Germany	13
Cleaning	14
Recycling the Product	14

## 2 Site Preparation

Overview	16
Customer responsibility	16
Agilent responsibility	17
Other documentation	19
Space and Weight Requirements	20

Communications Requirements	22
Telephone	22
Building network	22
Electrical Requirements	23
Voltage ranges of major components	23
Power configurations	24
Power requirements	25
Power plugs and cords	26
Other electrical considerations	27
Air Conditioning Requirements	29
Temperature and humidity	29
Airborne dust	30
Exhaust venting	30
Fume (exhaust) hood	31
Carrier and Reagent Gas Requirements	32
Regulators, tubing, and fittings	33
Laboratory Supply Requirements	34
Cleaning solvents	34
Data system supplies	34
Spare parts and consumables	35
Receiving the System	38
Delivery and unloading	38
Inspecting for damage	38
Storage	39
Unpacking	39
Installation and Verification	40
Installation	40
Verification	40
Sensitivity specifications	41

## A Power Cords



# 1 General Information

Instrument Identification	6
Important Safety Warnings	7
Safety and Regulatory Certifications	11
Cleaning	14
Recycling the Product	14

This section provides information on how to identify specifically what type of MSD you are working with as well as how to properly clean and dispose of the instrument. Also included in this section are the “[Important Safety Warnings](#)” which list critical safety precautions for all users.



## Instrument Identification

Each 5973N and 5973 inert Mass Selective Detector (MSD) is identified by a unique 10-character serial number. This serial number is located on a label on the lower left side near the front of the instrument. On CI upgrades, a serial number is located on the flow module.

When corresponding with Agilent Technologies about your instrument, be sure to include the model number and both full 10-character serial numbers.

Write the serial number of your 5973 MSD here for reference:

### 5973N MSD version

The 5973N MSD is equipped with a diffusion pump. The serial number label displays a product number that tells what kind of MSD you have:

Product number	MSD type
G2577A	Diffusion Pump, EI

### 5973 inert MSD version

The 5973 inert MSD is equipped with either a standard turbomolecular pump or a performance turbomolecular pump. Chemical Ionization is available for the turbo MSDs only. The serial number label displays a product number that tells what kind of MSD you have:

Product number	MSD type
G2578A	Standard Turbo Pump, EI
G2579A	Performance Turbo Pump, EI
G2589A	Performance Turbo Pump, EI, NCI, PCI

This site preparation guide will refer to both 5973N and 5973 inert MSDs as 5973 Series MSDs unless noted otherwise.

## Important Safety Warnings

Before moving on, there are several important safety notices that you should always keep in mind when using the 5973 Series Mass Selective Detector.

### Many internal parts of the MSD carry dangerous voltages

If the MSD is connected to a power source, even if the power switch is off, potentially dangerous voltages exist on:

- The wiring between the MSD 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 any heater.

#### **WARNING**

**All these parts are shielded by covers. With the covers in place, it should be difficult to accidentally make contact with dangerous voltages. Unless specifically instructed to, never remove a cover unless the detector, inlet, or oven are turned off.**

---

#### **WARNING**

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

---

Connecting the MSD to power sources that are not equipped with protective earth contacts creates a shock hazard for the operator and can damage the instrument.

Interrupting the protective conductor inside or outside the MSD or disconnecting the protective earth terminal creates a shock hazard for the operator and can damage the instrument.

Make sure the power cords supplied with the MSD are appropriate for your country and site before using them. Maintain easy access to the power cords so they can be disconnected during maintenance.

The use of incorrect or makeshift fuses or the short-circuiting of fuse holders creates a shock hazard for the operator and can damage the instrument. Replace fuses only with fuses of identical current rating and type.

Excessive fluctuations in the line voltage can create a shock hazard and can damage the instrument. Make sure the supply voltage does not fluctuate more than +5% or -10% from the rated voltage. This equipment must be installed in a Category II environment as defined in IEC 664.

### **Electrostatic discharge is a threat to MSD electronics**

The printed circuit (PC) boards in the MSD 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 MSD covers.

### **Many parts are dangerously hot**

Many parts of the MSD operate at temperatures high enough to cause serious burns.

You should always cool heated areas of the MSD 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 gloves. Whenever possible, cool the part of the instrument that you will be maintaining before you begin working on it.



**WARNING**

Be careful when working behind the gas chromatograph (GC). During cool-down cycles, the GC emits hot exhaust which can cause burns.

---

**WARNING**

The insulation around the GC inlets, detectors, valve box, and the insulation cups is made of refractory ceramic fibers. To avoid inhaling fiber particles, we recommend the following safety procedures: ventilate your work area; wear long sleeves, gloves, safety glasses, and a disposable dust/mist respirator; dispose of insulation in a sealed plastic bag; wash your hands with mild soap and cold water after handling the insulation.

---

## Hydrogen

Hydrogen gas may be used as carrier gas, and/or as fuel for an FID. When mixed with air, hydrogen can form explosive mixtures.

**WARNING**

When using hydrogen (H<sub>2</sub>) as the carrier gas or fuel gas, be aware that hydrogen gas can flow into the oven and create an explosion hazard. Therefore, be sure that the supply is off until all connections are made, and ensure that the GC inlet and detector column fittings are either connected to a column or capped at all times when hydrogen gas is supplied 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**

The MSD cannot detect leaks in gas streams. For this reason, it is vital that column fittings should always be either connected to a column, or have a cap or plug installed.

---

## General Information

When using hydrogen gas, check the system for leaks as described by your local Environmental Health and Safety (EHS) requirements to prevent possible fire and explosion hazards. Always check for leaks after changing a tank or servicing the gas lines. Always make sure the vent line is vented into a fume hood.

### Chemical safety

#### **WARNING**

**The foreline pump exhaust and split vent exhaust will contain traces of the chemicals you are analyzing. These could potentially be toxic. Vent the foreline pump exhaust and split vent exhaust outside your laboratory or into a fume hood. Be sure to comply with all local environmental regulations.**


---

## Safety and Regulatory Certifications

The 5973 Series Mass Selective Detector conforms to the following safety standards:

- Canadian Standards Association (CSA): C22.2 No. 1010.1
- CSA/Nationally Recognized Test Laboratory (NRTL): UL 61010A-1
- International Electrotechnical Commission (IEC): 61010-1
- EuroNorm (EN): 61010-1

The 5973 Series Mass Selective Detector conforms to the following regulations on Electromagnetic Compatibility (EMC) and Radio Frequency Interference (RFI):

- CISPR 11/EN 55011: Group 1, Class A
- IEC/EN 61326
- AUS/NZ 

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



The 5973 Series Mass Selective Detector is designed and manufactured under a quality system registered to ISO 9001.

### Information

The Agilent Technologies 5973 Series Mass Selective Detector meets the following IEC (International Electro-technical Commission) classifications: Safety Class I, 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. 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 MSD 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

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.

See accompanying instructions for more information.



Indicates a hot surface.



Indicates hazardous voltages.



Indicates earth (ground) terminal.



Indicates explosion hazard.



Indicates radioactivity hazard.



Indicates electrostatic discharge hazard.



## Electromagnetic compatibility

This device complies with the requirements of CISPR 11.

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

## Sound emission certification for Federal Republic of Germany

### Sound pressure

Sound pressure  $L_p < 70$  dB according to EN 27779:1991.

### Schalldruckpegel

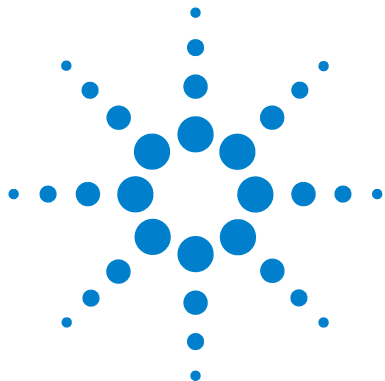
Schalldruckpegel  $L_P < 70$  dB am nach EN 27779:1991.

## **Cleaning**

To clean the unit, disconnect the power and wipe down with a damp, lint-free cloth.

## **Recycling the Product**

For recycling, contact your local Agilent sales office.



## 2 Site Preparation

Overview	16
Space and Weight Requirements	20
Communications Requirements	22
Electrical Requirements	23
Air Conditioning Requirements	29
Carrier and Reagent Gas Requirements	32
Laboratory Supply Requirements	34
Receiving the System	38
Installation and Verification	40

This section describes how to properly prepare your site for a new MSD. Follow these instructions carefully as delays due to improper site preparation may result in loss of instrument use during the warranty period.



### Overview

Before the 5973 Series Mass Selective Detector system can be installed, the site must be properly prepared. Site preparation includes, but is not limited to, ensuring that adequate facilities are available. Among the site requirements are:

- Adequate space is available for the MSD system.
- A suitable supporting bench is available.
- Adequate electrical power is available at the correct voltages and frequencies.
- Environmental control systems are adequate to maintain a correct, stable operating environment.
- Preparations for safe exhaust venting are adequate.
- Supplies necessary for instrument operation are available, including solvents, carrier and reagent gasses, and printer paper.

#### NOTE

Installation and verification will **only** be performed using helium carrier gas.

Delays due to inadequate site preparation could cause loss of instrument use during the warranty period. In extreme cases, Agilent Technologies may ask to be reimbursed for the additional time required to complete the installation. Agilent Technologies provides service during the warranty period and under maintenance agreements only if the specified site requirements are met.

### Customer responsibility

Unless previous arrangements have been made with Agilent Technologies, site preparation is the customer's responsibility. Customer responsibilities include, but are not limited to:

- Planning, scheduling, and preparing the site according to the specifications in this manual



- Verifying that the electrical environment is safe and adequate for the MSD system installation and operation
- Complying with all local laws (codes, ordinances, and regulations) for mechanical, building, and electrical distribution systems, hazardous waste disposal, and chemical storage.  
*Compliance must exist prior to installation.*
- Providing lifting equipment adequate to unload the system from the delivery vehicle and transport it to the site where it will be installed
- Providing someone to help lift the MSD onto the laboratory bench
- Providing adequate secure storage space for the system until it can be installed by an Agilent Technologies representative

### Agilent responsibility

An Agilent Technologies service representative will install the MSD and verify its performance. The service representative's responsibilities are limited to:

- Unpacking the MSD system and verifying that all components are present and undamaged
- Connecting the carrier gas line to the instrument from the tank, regulators, and lines previously installed by the customer
- Installing, connecting, and turning on MSD system components
- Verifying that the system meets Agilent Technologies performance standards
- Providing **basic** user familiarization for system hardware and software
- Signing the customer up through the Response Centers for instrument and software support

Agilent Technologies is not responsible for:

- Any task not listed in the *5973 Series MSD Hardware Installation Manual* (G2589-90073) or the installation guides for the GC, data system, and other accessories
- Installing helium carrier gas or reagent gas tanks and regulators
- Connecting or verifying the performance of hardware and software not provided by Agilent Technologies

---

**NOTE**

Items that do not include installation by Agilent Technologies must be installed by the customer.

- 
- Connecting the MSD/ChemStation system to the building network
  - Testing the MSD against customer standards or samples
  - Providing detailed instruction in the operation of the computer operating system and MSD software

---

**NOTE**

Contact Agilent Technologies for information concerning training classes.

- 
- Setting up laboratory procedures

---

**NOTE**

Contact Agilent Technologies for information concerning Application Assist or Chemical Analysis Consulting services.

- 
- Operating the MSD system following installation

## Other documentation

Additional information is contained in the following documentation:

- *5973N or 5973 inert Mass Selective Detector Hardware Manual*, which is located on the *5973N and 5973 inert Mass Selective Detector User Information* CD-ROM
- *5973 Network Mass Selective Detector Site Preparation manual* or *5973 inert Mass Selective Detector Site Preparation Guide*
- 6890 Series GC manuals
- GC accessories (autosampler, etc.) manuals
- G1701DA MSD Productivity ChemStations software manuals and online help
- The appropriate sensitivity specifications for your instrument:  
5973N: Agilent publication 5968-7358E  
5973 inert: Agilent publication 5988-9991EN
- *Hydrogen Carrier Gas Safety Guide* (5955-5398)

For updated information, see the Agilent Technologies website at <http://www.agilent.com/chem>.

## Space and Weight Requirements

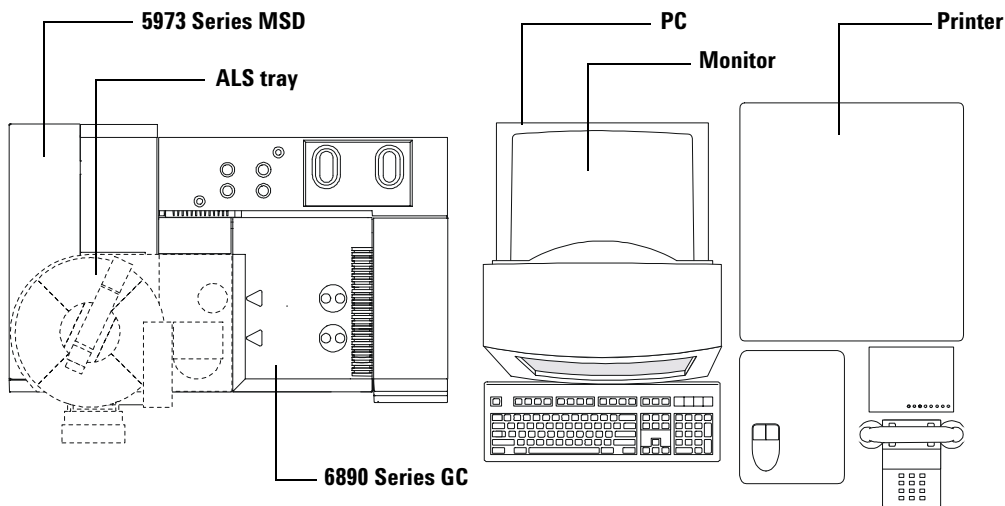
Table 1 lists dimension and weight information for the MSD and related components. Your site must have enough bench space for the MSD, GC, data system, and accessories. In addition, there must be sufficient space around the system for ventilation and maintenance access. At least 30 cm (12 inches) to the left of the MSD and at least 20 cm (8 1/4 inches) behind the MSD must be kept clear.

Benches must be sturdy enough to support the weight of the entire system.

The foreline pump can be located on the laboratory bench or on the floor. It must be close to the MSD because it is connected by a 200-cm (79-inch) hose. The hose is stiff and cannot be bent sharply.

### CAUTION

Do not put the foreline pump on your laboratory bench if vibration-sensitive equipment is located on the bench.



**Table 1** Product dimensions\*

Product	Dimensions, cm (in)			
	Height	Width	Depth	Weight, kg (lb)
5973 Series MSD <sup>†</sup>				
G2577A Diffusion pump EI MSD	40.6 (16)	29.8 (11.75)	54 (21.25)	36 (80)
G2578A Standard turbo EI MSD	40.6 (16)	29.8 (11.75)	54 (21.25)	39 (85)
G2579A Performance turbo EI MSD	40.6 (16)	29.8 (11.75)	54 (21.25)	44 (95)
G2589A EI/PCI/NCI MSD	40.6 (16)	30.3 (12.01)	54 (21.25)	46 (100)
6890 Series GC	50.0 (21)	58.5 (23)	50.0 (21)	50.0 (110)
Data System	Data system size and weight depend on the components included in the data system. Reserve at least 100 cm (39 in) of bench space for the data system. A typical data system weight is 34 kg (75 lb).			

\* All dimensions are approximate

† Be sure to plan for the space for your carrier and reagent gas source(s).

**NOTE**

The specifications in [Table 1](#) are for a single MSD ChemStation configuration. Two-MSD configurations require additional table space. The second GC/MSD system can be located on a different table, connected to the ChemStation with a network cable.

## Communications Requirements

### Telephone

A telephone with a cord long enough to be used at the computer will allow the MSD operator to communicate with Agilent Technologies support personnel. In addition, a second analog phone line is strongly recommended to allow remote control and diagnosis of the MSD through a modem. This is **required** for some service contracts.

### Building network

If you intend to connect your MSD system to your building's network, you must have an additional shielded twisted pair (STP) network cable.

#### NOTE

Agilent Technologies is not responsible for connecting to or establishing communication with your building network. The representative will test the 5973 Series MSD's ability to communicate on a mini-hub only.

---

#### NOTE

The IP addresses assigned to the instrument(s) and the ChemStation PC must be **fixed** (permanently assigned) addresses. If you intend to connect your MSD system to your building's network, each piece of equipment (the MSD, the GC, and the PC) must have a unique IP address assigned to it.

---

## Electrical Requirements

You are responsible for providing appropriate electrical power and power outlets for all of the components in your 5973 Series MSD system. Power considerations include:

- Voltage ranges of major components
- Power configurations
- Power requirements
- Power plugs and cords

### Voltage ranges of major components

The 5973 Series MSD includes a full-range power supply that can operate without reconfiguration on either of two wide ranges of single-phase alternating current (AC) electrical power:

- 120 VAC +5/-10%, 50/60 Hz  $\pm$  5% (typical for North America)
- 200-240 VAC +5/-10%, 50/60 Hz  $\pm$  5% (typical for Europe and Japan)

The foreline pump also draws its power from the MSD. However, a different foreline pump is supplied depending on the voltage range on which the MSD will be operating. The pump is supplied according to the standard voltage in the country from which the order originates. For example, if an order originates from an Agilent Technologies sales office in Germany, the foreline pump supplied will be configured to operated on the standard voltage and frequency of electrical power in Germany.

#### NOTE

The MSD does not support 100 VAC input power. In countries that have only 100 VAC power, a step-up transformer must be used to provide the appropriate voltage to the MSD.

### CAUTION

If an instrument is being ordered from one location, but is to be installed in another location with different electrical power characteristics, this must be noted on the order. A special note must also be made if the electrical power at the site is different from the standard electrical power in that country.

---

### Power configurations

Electrical power for the MSD may be delivered in either single-phase or 208-Wye configuration. Correct grounding for the 208-Wye configuration must be verified by an electrician. The neutral wire **cannot** be used for safety grounding. The ground wire should carry zero current except for ground-fault current or static electric discharge. The entire system should share an isolated, noise-free electrical ground. This system ground should be electrically separate from the ground for the rest of the building, back to the main ground for the facility. Power configurations are provided in [Table 2](#).

### WARNING

**Connecting an MSD to a power source which is not equipped with a protective earth contact (ground) creates a shock hazard for the operator and can damage the instrument.**

---

### WARNING

**Interrupting the protective conductor inside or outside the MSD or disconnecting the protective earth terminal (ground) creates a shock hazard for the operator and can damage the instrument.**

---



**Table 2** Power configurations

Configuration	Measurement	Nominal voltage
Single phase, 120 VAC	Line to neutral	120 VAC*
	Line to ground	120 VAC*
	Ground to neutral	< 0.5 V rms
Single phase, 200-240 VAC	Line to neutral	200, 220, or 240 VAC*
	Line to ground	200, 220, or 240 VAC*
	Ground to neutral	< 0.5 V rms
208-Wye	Line to line (phase A to phase B)	208 VAC
	Line to ground (phase A to ground)	120 VAC
	Line to ground (phase B to ground)	120 VAC

\* Varies with country and/or region

## Power requirements

Table 3 lists the power requirements for the 5973 Series MSD and related equipment. Extra power capacity for future additions is a very good idea.

Each product listed requires a dedicated circuit. The MSD, GC, and data system must each have a separate circuit breaker. All of the equipment **must** share a common ground.

Power must meet the stability specifications listed in Table 3. Use a line monitor to check power stability. If your line power is unstable, you may need to install a line conditioner.

**Table 3** Power requirements

Product	Line voltage	Maximum continuous power consumption	Supply circuit rating	Outlets required
5973 Series MSD <sup>*</sup>	120 VAC +5/-10%, 50/60 Hz ± 5%	900 VA (400 VA for foreline pump only)	15 A	1
	200-240 VAC +5/-10%, 50/60 Hz ± 5%	900 VA	15 A	1
6890 Series GC and ALS	120 VAC +5/-10%, 50/60 Hz ± 5%	1000 VA <sup>2</sup>	20 A	1 or 2
	200-240 VAC +5/-10%, 50/60 Hz ± 5%	1000 VA <sup>†</sup>	15 A	1 or 2
Gauge controller	100-240 VAC			1
ChemStation PC	100-120 or 220-240 VAC +5/-10%, 50/60 Hz ± 5%	1000 VA <sup>2</sup>	15 A	3 to 5

\* The MSD operates on either voltage range. The foreline pump and the diffusion pump are supplied for one voltage range or the other based on the standard voltage in the country where the order originates.

† Depends on product configuration.

**WARNING**

**Excessive fluctuations in the voltage of the power supply can create a shock hazard and can damage the instrument. Make sure the supply voltage does not fluctuate more than +5% or –10% from the rated voltage. This equipment must be installed in a Category II environment as defined in IEC 664.**

### Power plugs and cords

The MSD is supplied with a power cord and plug appropriate for the country from which the order originates. For example, if an order originates from a Agilent Technologies sales office in Germany, the power cord and plug supplied will be compatible with the standard voltage and outlet configuration in Germany. See [“Power Cords”](#) on page 43 for illustrations of the power cords available.

Data system components also include power cords with plugs appropriate for the country where the order was placed.

Power cord lengths for the MSD and the data system components and accessories are approximately 2.3 m (7.5 ft).

**CAUTION**

If an instrument is being ordered from one location, but is to be installed in another location with different electrical power characteristics, this must be noted on the order. A special note must also be made if the electrical power at the site is different from the standard electrical power in that country.

---

**WARNING**

**Make sure the power cords supplied with the MSD are appropriate for your country and site before installing the instrument. See “Power Cords” on page 43.**

---

**WARNING**

**Do not use extension cords with the MSD or GC. Extension cords normally do not provide enough power and can be a safety hazard.**

---

**NOTE**

Maintain easy access to the power cords so they can be disconnected for maintenance.

---

## Other electrical considerations

Additional electrical considerations include:

- Electromagnetic interference (EMI), such as is generated by NMRs, radio transmitters, and microwave links, may interfere with system performance.
- Protect the system from static electricity by observing humidity and temperature requirements. Minimize the presence of non-conductive products such as carpets and vinyl floor tiles.
- Install emergency-off pushbuttons that can disconnect power to the ventilation system and all electric equipment in the room except overhead lighting.

## Site Preparation

- Provide separate convenience outlets for building maintenance and other appliances. Convenience outlets must be on circuits separate from the MSD system. Convenience outlets must share the normal building distribution ground, **not** the MSD system ground.
- In some geographical areas it may be advisable to install lightning protection for personnel and equipment.

## Air Conditioning Requirements

Air conditioning considerations include temperature, humidity, airborne dust, and exhaust venting. Each of these is considered in more detail in the following material.

### Temperature and humidity

The 5973 Series MSD is specified for operation under the following conditions:

- Operation requires constant temperature (variations < 2 °C/hr)
- Operation and storage require a non-condensing, non-corrosive atmosphere
- The temperature and humidity limits in [Table 4](#)

**Table 4** Temperature and humidity limits

	Temperature	Humidity (relative)
Operation	15 °C to 35 °C (59 °F to 95 °F)	40% to 80%
Storage	-20 °C to 70 °C (-4 °F to 158 °F)	0% to 95%

Environmental control systems must maintain these temperature and humidity ranges.

The MSD is rated for 900 Watts (3000 BTU/hr). The GC and the data system also contribute significantly to the cooling load although the exact amounts depend on their configurations. Additional allowances should be made for other heat sources such as heat from other equipment, heat from adjacent rooms, and heat from laboratory personnel.

### Airborne dust

Agilent Technologies recommends a maximum airborne particle density of  $55 \mu\text{g}/\text{m}^3$ . If you suspect your site exceeds this limit, contact your local Agilent Technologies Customer Service Organization. Customer Engineers with special training and equipment can test for airborne particle density. They can offer suggestions for reducing airborne dust.

### Exhaust venting

There are two sources of exhaust on the GC/MSD system: the foreline pump and the GC split vent. The foreline pump outputs gas removed from the vacuum manifold by the high vacuum pumps. The foreline pump exhaust will also contain traces of solvent and sample.

#### **WARNING**

**User safety requires that the exhaust gases from the MSD be vented externally to the building and not recirculated by the environmental control system. Health hazards include chemical toxicity of solvents, samples, derivitizing agents, pump fluid vapor, and aerosolized biological samples.**

---

#### **WARNING**

**The pump exhaust contains carrier gas and traces of solvents, analytes, and foreline pump oil. The supplied oil trap stops only pump oil. It does *not* trap or filter out toxic chemicals. If you are using toxic solvents or toxic or flammable carrier gas, or analyzing toxic chemicals, do not install the oil trap. Install a hose to take the pump exhaust to a fume hood. Do not use the oil trap on any CI MSD.**

---

The foreline pump exhaust **cannot** be vented into the laboratory if any hazardous materials will be introduced into the GC/MS system, including samples, solvents, carrier gases, and reagent gases. It must be vented external to the building or vacuum exhausted to a fume hood.

If a negative pressure vent is not available, the length of tubing from the foreline pump to an ambient pressure vent should not exceed 460 cm (15 ft). The exhaust can **not** be connected to a positive pressure vent.

Exhaust gas venting must comply with all local environmental and safety codes.

### Fume (exhaust) hood

An auxiliary work space and fume hood are needed for some maintenance procedures.

#### **WARNING**

**Hydrogen is potentially explosive. Take extreme care when using hydrogen as the GC carrier gas in a GC/MS system. Read the *Hydrogen Carrier Gas Safety Guide* (5955-5398) before operating the MSD with hydrogen carrier gas.**

---

## Carrier and Reagent Gas Requirements

Table 5 lists the specifications for the carrier and reagent gases.

**Table 5** Carrier and reagent gas requirements

	Purity	Typical pressure range (psi)	Typical flow (mL/min)
Helium (required)	99.999% or better and hydrocarbon free	50 to 80	20 to 50 (column and split flow)
Methane reagent gas (required for CI operation)	Research or SFC grade, 99.999%	15 to 25	1 to 2
Isobutane reagent gas (optional)	Instrument grade, 99.999%	15 to 25	1 to 2
Ammonia reagent gas (optional)	Research or SFC grade, 99.9995%	5 to 8	1 to 2
Carbon dioxide reagent gas (optional)	SFC Grade, 99.995%	15 to 25	1 to 2

**WARNING**

**Methane is flammable and ammonia is corrosive and toxic. Read and follow the safety information that comes with the gases.**



## Regulators, tubing, and fittings

You must supply an appropriate regulator for your carrier and reagent gas. The regulator must be able to supply gas in the specified pressure range. It must have one outlet with 1/4-inch Swagelok fittings. See the Agilent Technologies Chemical Analysis Columns and Supplies Catalog or visit the Agilent Technologies Chemical Analysis website at <http://www.agilent.com/chem> for dual-stage regulators available from Agilent Technologies.

You must supply fittings, ferrules, and connectors of a Swagelok design for the 1/8-inch tubing.

## Laboratory Supply Requirements

Laboratory supply requirements vary with applications.

### Cleaning solvents

Cleaning tasks for the MSD require the following HPLC-grade (or better) solvents:

- Methylene chloride
- Isopropyl alcohol
- Methanol
- Acetone
- Water

Proper storage, handling, and disposal of these chemicals is required for personal and environmental safety.

#### CAUTION

Carbon disulfide used as a sample solvent is corrosive and will damage the MSD analyzer. Using carbon disulfide will void the warranty on the analyzer.

---

#### WARNING

**Chemical solvents should be considered hazardous and must be handled with care. Contact your chemical supplier for solvent handling and safety information, preferably a material safety data sheet (MSDS).**

---

### Data system supplies

You will need paper for printing the results of the testing done during installation and later for printing reports of your analyses. You will also need appropriate hardware and media (flexible disks, tape cartridges, CD writers, etc.) for making backup copies of your data files.

**NOTE**

Agilent Technologies does not include CD writers with MSD systems. These are available from many suppliers.

## Spare parts and consumables

The supplies and parts listed in [Table 6](#) are used in the operation and maintenance of a 5973 Series MSD system. Keeping these parts on hand can reduce system downtime related to instrument maintenance and repair.

**Table 6** Maintenance supplies and consumables

Description	Part number
<b>Miscellaneous consumables</b>	
One-year maintenance kit	5183-0296
He/H <sub>2</sub> gas filter for GC	5182-3468
Abrasive sheets (5/pk)	5061-5896
Cloths, lint-free (15/pk)	05980-60051
Cotton swabs (100/pk)	5080-5400
Diffusion pump fluid, 18.5 mL, approximately 30 mL needed (diffusion pump MSDs only)	6040-0809
Foreline pump oil, Inland 45, 1 liter	6040-0834
El filament (2/pk)	05972-60053
El high temp filament	G2590-60053
Octafluoronaphthalene (OFN)	8500-5441
<b>Gloves, clean</b>	
large	8650-0030
small	8650-0029
Nitrile rubber gloves, medium	9300-1751
<b>Filaments</b>	
CI (2/pk)	G1099-80053

**Table 6** Maintenance supplies and consumables (continued)

El filament (2/pk)	05972-60053
El high temp filament	G2590-60053
<b>Chemical Consumables</b>	
PFTBA (El calibration fluid)	05971-60571
PFDTD (Cl calibration fluid)	8500-8130
EVAL A (evaluation sample)	05990-60045
OFN (Octafluoronaphthalene)	8500-5441
Benzophenone	8500-5440
Diffusion pump fluid (2 required)	6040-0809
Foreline pump oil, Inland 45, 1 liter	6040-0834
<b>Gas Filters</b>	
He/H <sub>2</sub> gas purifier for GC	RMSH-2
Methane/Isobutane gas purifier for Cl MSD	G1999-80410
<b>Commonly used tools</b>	
Safety glasses	9300-1159
Wrenches	
1/2-inch x 9/16-inch, open end	8710-0877
<b>Ferrules</b>	
Blank, graphite-vespel	5181-3308
GC/MSD interface	
0.3-mm id, 85% Vespel 15% graphite, for 0.10-mm id columns	5062-3507

**Table 6** Maintenance supplies and consumables (continued)

0.4-mm id, 85% Vespel 15% graphite, for 0.20-mm id and 0.25-mm id columns	5062-3508
0.5-mm id, 85% Vespel 15% graphite, for 0.32-mm id columns	5062-3506
0.8-mm id, 85% Vespel 15% graphite, for 0.53-mm id columns	5062-3538
<b>Injection port</b>	
0.27-mm id, 90% Vespel 10% graphite, for 0.10-mm id columns	5062-3518
0.37-mm id, 90% Vespel 10% graphite, for 0.20-mm id columns	5062-3516
0.40-mm id, 90% Vespel 10% graphite, for 0.25-mm id columns	5181-3323
0.47-mm id, 90% Vespel 10% graphite, for 0.32-mm id columns	5062-3514
0.74-mm id, 90% Vespel 10% graphite, for 0.53-mm id columns	5062-3512

# Receiving the System

When your MSD system is delivered, it is your responsibility to provide for removal of the shipping containers from the truck and their storage until installation. Contact your Agilent Technologies service representative as soon as your shipment arrives to arrange an installation date.

## Delivery and unloading

The shipping containers are large and heavy. The largest container in the shipment is 121.9 cm (48 inches) x 121.9 cm x 121.9 cm. It contains several smaller cartons that can be removed from the large container after delivery to facilitate moving them to the location where the instrument is to be installed. After removing the smaller cartons and discarding the outer container, the largest remaining carton, which contains the GC, is approximately:

- 88.9 cm (35 inches) high
- 81.3 cm (32 inches) wide
- 73.7 cm (29 inches) deep
- 60 kg (132 lbs)

All doorways, hallways, floors, and elevators must be able to accommodate the largest, heaviest container.

## Inspecting for damage

Once the shipping containers are unloaded, examine them for any obvious **external** damage. If any of the containers appear damaged, note on the carrier's bill of lading that there is: *Apparent damage - subject to inspection and test*. Arrange for both the carrier's claims representative and your Agilent Technologies service representative to be present when the containers are unpacked.

Do not open any of the shipping containers unless a representative of Agilent Technologies is present. **Opening any of the containers without an Agilent Technologies representative being present will void the receiving warranty on the instrument.**

## Storage

It is your responsibility to store the containers until installation. If your site does not have adequate storage space, the containers may be stored at your expense in a bonded warehouse. Allow space for data system and accessory containers too.

The environment in the storage area should be between 5 °C and 50 °C (41 °F and 122 °F), 20% to 80% relative humidity, non-condensing and non-corrosive.

## Unpacking

Do not open any shipping containers until an Agilent Technologies representative is present. Warranty claims for missing items will not be honored unless an Agilent Technologies representative is present to verify the contents of each container as it is unpacked.

The actual shipping containers become your property and should not be returned to Agilent Technologies.

# Installation and Verification

## Installation

Once the installation has begun, it should progress in a timely manner to completion. Delays due to inadequate site preparation could cause loss of instrument use during the warranty period. In extreme cases, Agilent Technologies may ask to be reimbursed for the additional time required to complete the installation.

The primary user and, if possible, all other users of the MSD should be present during installation to receive familiarization instruction from the Agilent Technologies service representative.

## Verification

The final step in the installation process is system verification. Your Agilent Technologies service representative will test the system against Agilent Technologies specifications as documented for the product(s) or system(s) you have purchased. (Not all tests are performed for every system.)

**The Agilent Technologies service representative will not test your system against your standards or samples.**

Further, the Agilent Technologies service representative will not set up your laboratory procedures. Assistance with laboratory procedures can be obtained from your local Agilent Technologies Applications Engineer (AE) on a consulting basis at additional cost.



## Sensitivity specifications

See the appropriate sensitivity specifications for your instrument type:

5973N: Request Agilent publication 5968-7358E

5973 inert: Request Agilent publication 5988-9991EN

### NOTE

SIM performance verification is not performed at installation and must be purchased if desired.

---

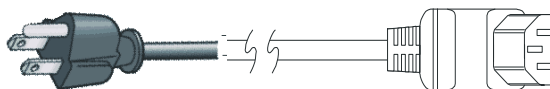
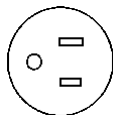




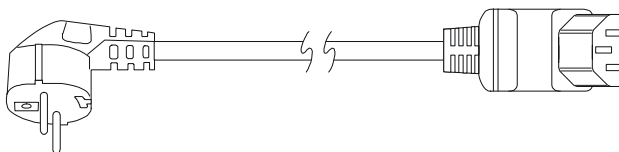
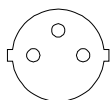
## A Power Cords

This appendix shows the power cords available for the 5973 Series MSD. See [page 26](#) for information about ensuring that the correct power cord is supplied with your system.

US and Canada, NEMA 6-15P (Part number 8120-6825)

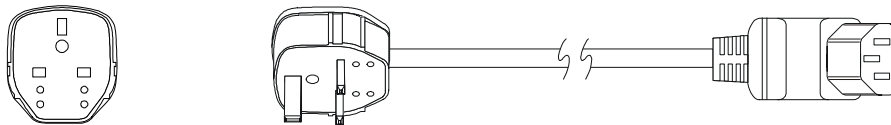


European Power, CEE 7/7 (Part number 8120-1689)

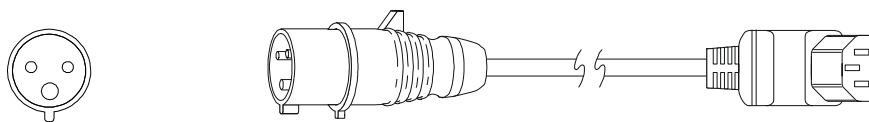


## Power Cords

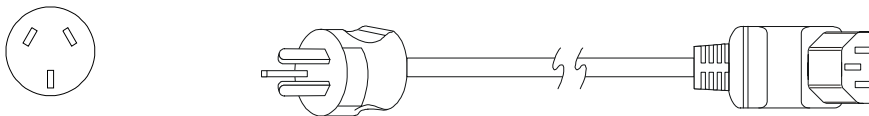
UK / Hong Kong, BS 1363 (Part number 8120-1351)



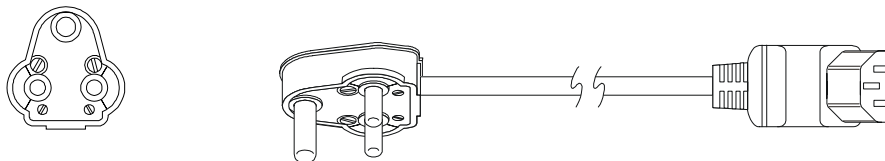
Denmark/Greenland, IEC 309 (Part number 8120-3997)



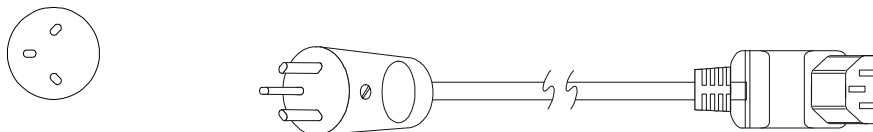
Australia / New Zealand, AS 3112-1981 (Part number 8120-1369)



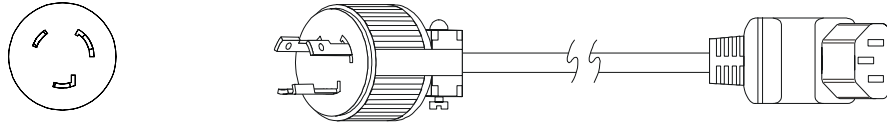
India / South Africa, BS 546 (Part number 8120-4211)



Israel, SI 32 (Part number 8120-5182)



Japan, NEMA L6-20P (Part number G2025-60189)



## Power Cords





© Agilent Technologies, Inc.  
Printed in USA, August 2003



G2589-90070