

## **Agilent GC, MSD, and ALS**

### **Site Preparation Checklist**

Dimensions and Weight	2
Power Consumption	4
Heat Dissipation	6
Environmental Conditions	7
Gas Selection	8
Gas Supply	10
Other Considerations	12
Exhaust venting requirements for the GC/MSD	12
Basic tools	13

This checklist outlines the space and resource requirements for a GC, MSD, and automatic liquid sampler (ALS) installation. For a successful and timely installation of the instrument, the site must meet these requirements before beginning installation. Necessary supplies (gases, tubing, operating supplies, consumables, and other usage-dependent items such as columns, vials, syringes, and solvents) must also be available. Note that performance verification requires the use of helium carrier gas and, for models using chemical ionization, methane reagent gas. Refer to the Agilent Web site at [www.agilent.com/chem](http://www.agilent.com/chem) for the most up-to-date listing of GC, MSD, and ALS supplies and consumables.



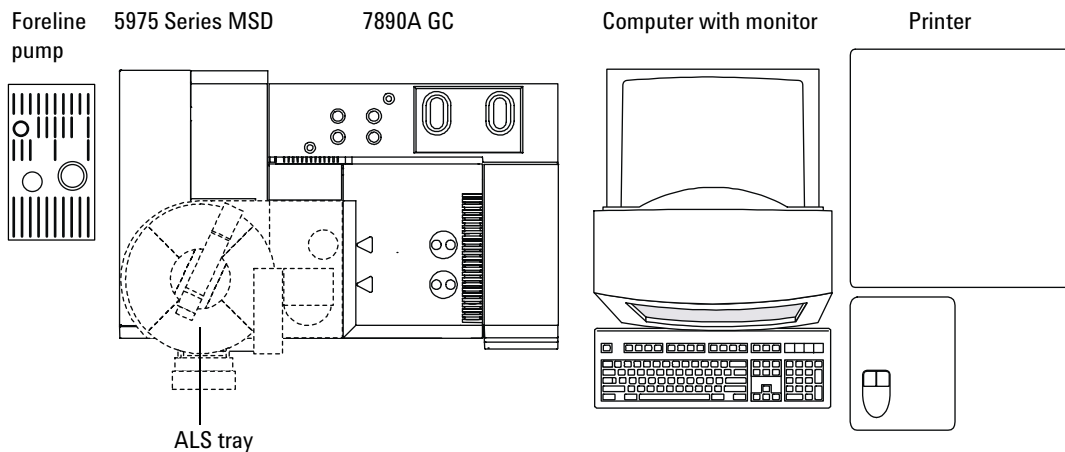
## Dimensions and Weight

Select the laboratory bench space before the system arrives. Pay special attention to the total height requirements. Avoid bench space with overhanging shelves. See Table 1.

Allow at least 20-cm clearance between back of GC and wall to dissipate air.

**Table 1** Required height, width, depth, and weight

	Height	Width	Depth	Weight
5975 Series MSDs				
• Diffusion pump	41 cm (16 in)	30 cm (12 in)	54 cm (22 in)	39 kg (85 lb)
• Standard turbo pump	41 cm (16 in)	30 cm (12 in)	54 cm (22 in)	39 kg (85 lb)
• Performance turbo pump	41 cm (16 in)	30 cm (12 in)	54 cm (22 in)	41 kg (90 lb)
• Performance CI/EI turbo pump	41 cm (16 in)	30 cm (12 in)	54 cm (22 in)	46 kg (100 lb)
Foreline pump				
Standard	21 cm (8 in)	13 cm (5 in)	31 cm (12 in)	11 kg (23.1 lb)
Oil-free	19 cm (7.5 in)	32 cm (13 in)	28 cm (11 in)	16 kg (35.2 lb)
7890A Series GCs	54 cm (22 in)	59 cm (23 in)	54 cm (22 in)	50 kg (112 lb)
6890 Series GCs	54 cm (22 in)	55 cm (22 in)	54 cm (22 in)	50 kg (112 lb)
6850 Series GCs	51 cm (20 in)	29 cm (12 in) 34 cm (14 in) CO <sub>2</sub> 37 cm (15 in) 6850 ALS	57 cm (23 in)	< 23 kg (51 lb)
Additional space requirements				
• MSD operational and maintenance access	Requires 30 cm (1 ft) to its left			
• Typical laser printer	Requires 41 cm (16 in)			
• GC operational oven access	Requires ≥ 30 cm (12 in) open space above			
• GC with ALS	Requires 44 cm (17.3 in) above the GC			
• GC with CTC PAL Autosampler	Requires 66 cm (26 in) above the GC and 4 to 20 cm (1.5 to 8 in) to the left or right, depending on configuration			



**Figure 1** Top view of typical installation (GC/MSD system with ALS)

A simple system that includes a GC, an ALS, and a computer would require about 153 cm (5 ft) of bench space. Allowing for operational access and a printer, a total of 244 cm (8 ft) of bench space should be available for a full GC/MSD system. Some repairs to the MSD or to the GC will also require access to the back of the instrument(s).

Note that the length of the vacuum hose is 130 cm (4 ft 3 in) from the high vacuum pump to the foreline pump, and the length of the foreline pump power cord is 2 m (6 ft 6 in).

## Power Consumption

Table 2 lists site power requirements.

- The number and type of electrical outlets depend on the size and complexity of the system.
- Power consumption and requirements depend on the country the unit ships to.
- The electrical outlet for the unit should have a dedicated ground.

**Table 2** Power requirements

Product	Line voltage (VAC)	Frequency (Hz)	Current rating (amps)	Maximum continuous power consumption (VA)	Outlets required
5975 Series MSD	120 (–10% / +5%)	50/60 ± 5%	8	1100 (400 for foreline pump only)	1
	220–240 (–10% / +5%)	50/60 ± 5%	8	1100 (400 for foreline pump only)	1
	200 (–10% / +5%)	50/60 ± 5%	8	1100 (400 for foreline pump only)	1
ChemStation PC system (monitor, CPU, printer)	120 (–10% / +5%)	50/60 ± 5%	15	1000	3–5
	200–240 (–10% / +5%)	50/60 ± 5%	15	1000	3–5
Agilent 7890A and 6890	Standard Americas: 120* single phase (–10% / +5%)	48–66	18.8	2250	1
	220/230/240 single/split phase (–10% / +5%)	48–66	10.2/9.8/9.4	2250	1
	Fast Japan: 200 split phase (–10% / +5%)	48–66	14.8	2950	1
	220/230/240 <sup>†‡</sup> single/split phase (–10% / +5%)	48–66	13.4/12.8/12.3	2950	1
6850	Standard Japan: 100 single phase (–10% / +10%) <sup>**</sup>	48–66	15	1440	1

**Table 2** Power requirements (continued)

Product	Line voltage (VAC)	Frequency (Hz)	Current rating (amps)	Maximum continuous power consumption (VA)	Outlets required
	Americas: 120 single phase (-10% / +10%)**	48-66	12	1440	1
	230 single/split phase (-10% / +10%)**	48-66	9	2000	1
Fast	120 single phase (-10% / +10%)**	48-66	20	2400	1
	220/230/240 single/split phase (-10% / +10%)**	48-66	11	2400	1
	200/208 single/split phase (-10% / +10%)**	48-66	12	2400	1

\* Americas 120 VAC requires 20 amp dedicated line. Americas 240 VAC requires 15 amp dedicated line.

† Option 003, 208 VAC fast oven, uses a 220 VAC unit with operating range of 193 to 231 VAC. Most labs have 4-wire service resulting in 208 VAC at the wall receptacle. It is important to measure the line voltage at the receptacle for the GC.

‡ Power line conditioners should not be used with 6890 or 7890A GCs.

\*\* Requires an isolated ground and dedicated outlet.

## Heat Dissipation

Use Table 3 to estimate the additional BTUs of heat dissipated from this equipment. Maximums represent the heat given off when heated zones are set for maximum temperatures.

**Table 3** Heat dissipation

Oven	Agilent 7890A and 6890 Series	6850 Series	5975 Series
Standard oven ramp	7681 BTU/hour maximum	< 4800 BTU/hour maximum	3000 BTU/hour including GC/MSD interface
Fast oven ramp (options 002 and 003)	10,071 BTU/hour maximum	< 4800 BTU/hour maximum	3000 BTU/hour including GC/MSD interface

## Environmental Conditions

Operating the GC/MSD within the recommended ranges optimizes instrument performance and lifetime. The instrument needs space for proper convection of heat and ventilation. Performance can be affected by sources of heat and cold from heating, air conditioning systems, or drafts. See Table 4.

**Table 4** Environmental conditions for operation and storage

Product	Conditions	Operating temp range	Operating humidity range	Maximum altitude
Agilent 7890A and 6890 Series	Standard oven ramp	20 to 27 °C	50 to 60%	4,615 m
	Fast oven ramp (options 002 and 003)	20 to 27 °C	50 to 60%	4,615 m
	Storage	5 to 40 °C	5 to 95%	
6850 Series	Standard oven ramp	15 to 35 °C	5 to 95%	4,615 m
	Fast oven ramp (options 002 and 003)	15 to 35 °C	5 to 95%	4,615 m
	Storage	5 to 40 °C	5 to 95%	
5975 Series	Operation	15 to 35 °C (59 to 95 °F)	40 to 80%	4,615 m*
	Storage	-20 to 70 °C (-4 to 158 °F)	0 to 95%	

\* 5975B VL MSD: 2,300 m

## Gas Selection

Table 5 lists gases usable with Agilent GCs and capillary columns.

**Table 5** Gases usable with Agilent GCs

Detector type	Carrier	Preferred makeup	Alternate choice	Detector, anode purge, or reference
Electron capture (ECD)	Hydrogen	Argon/Methane	Nitrogen	Anode purge must be same as makeup
	Helium	Argon/Methane	Nitrogen	
	Nitrogen*	Nitrogen	Argon/Methane	
	Argon/Methane*	Argon/Methane	Nitrogen	
Flame ionization (FID)	Hydrogen	Nitrogen	Helium	Hydrogen and air for detector
	Helium	Nitrogen	Helium	
	Nitrogen*	Nitrogen	Helium	
Flame photometric (FPD)	Hydrogen	Nitrogen		Hydrogen and air for detector
	Helium	Nitrogen		
	Nitrogen*	Nitrogen		
	Argon*	Nitrogen		
Nitrogen-Phosphorus (NPD)	Helium	Nitrogen	Helium	Hydrogen and air for detector
	Nitrogen*	Nitrogen	Helium	
Thermal conductivity (TCD)	Hydrogen	Must be same as carrier and reference	Must be same as carrier and reference	Reference must be same as carrier and makeup
	Helium			
	Nitrogen*			

\* Not generally suitable for MSD carrier gas.



Agilent recommends that carrier and detector gases be 99.9995% pure. See Table 6. Air needs to be zero grade or better. Agilent also recommends using high quality traps to remove hydrocarbons, water, and oxygen.

**Table 6** Carrier and reagent gases purity

<b>Carrier and reagent gas requirements</b>	<b>Purity</b>	<b>Notes</b>
Helium (carrier)	99.9995%	Hydrocarbon free
Hydrogen (carrier)	99.9995%	SFC grade
Methane reagent gas <sup>*</sup>	99.999%	Research or SFC grade
Isobutane reagent gas <sup>†</sup>	99.99%	Instrument grade
Ammonia reagent gas <sup>†</sup>	99.9995%	Research or SFC grade
Carbon dioxide reagent gas <sup>†</sup>	99.995%	SFC grade

\* Required reagent gas for installation and performance verification, CI MSDs only.

† Optional reagent gases, CI MSDs only

## Gas Supply

Supply instrument gases using tanks, an internal distribution system, or gas generators. If used, tanks require two-stage pressure regulators with packless, stainless steel diaphragms. The GC/MSD requires 1/8-inch Swagelok connections to its gas supplies. Plumb the gas supply tubing/regulators so that one 1/8-inch Swagelok female connector is available for each gas.

Table 7 lists minimum and maximum delivery pressures for inlets and detectors, measured at the bulkhead fittings on the back of the instrument.

**Table 7** Delivery pressures required at the GC/MSD, in kPa (psig)

	Detector type					Inlet type				
	FID	NPD	TCD	ECD	FPD	Split/Splitless 150 psi	Split/Splitless 100 psi	On-column	Purged packed	PTV
Hydrogen	240–690 (35–100)	240–690 (35–100)			310–690 (45–100)					
Air	380–690 (55–100)	380–690 (55–100)			690–827 (100–120)					
Makeup	380–690 (55–100)	380–690 (55–100)	380–690 (55–100)	380–690 (55–100)	380–690 (55–100)					
Reference			380–690 (55–100)							
Carrier (max)						1,172 (170)	827 (120)	827 (120)	827 (120)	827 (120)
Carrier (min)						(20 psi) above pressure used in method				

Conversions: 1 psi = 6.8947 kPa = 0.068947 Bar = 0.068 ATM

Notes:

- If you have not requested option 305, you must supply pre-cleaned, 1/8-inch copper tubing and a variety of 1/8-inch Swagelok fittings to connect the GC to inlet and detector gas supplies.
- Cryogenic cooling with Liquid N<sub>2</sub> requires 1/4-inch insulated copper tubing.

- Cryogenic cooling with Liquid CO<sub>2</sub> requires 1/8-inch heavy-walled, stainless steel tubing.
- Valve actuation requires a separate pressurized, dry air supply at 380 kPa (55 psig).
- Never use liquid thread sealer to connect fittings. Never use chlorinated solvents to clean tubing or fittings.

Table 8 lists the limits on total gas flow into the 5975 Series MSD.

**Table 8** 5975 Series total gas flow limitations

Feature	G3170A	G3171A	G3172A	G3174A
High vacuum pump	Diffusion	Standard turbo	Performance turbo	Performance turbo, EI/PCI/NCI
Optimal gas flow mL/min <sup>*</sup>	1.0	1.0	1.0 to 2.0	1.0 to 2.0
Maximum recommended gas flow, mL/min	1.5	2.0	4.0	4.0
Maximum gas flow, mL/min <sup>†</sup>	2.0	2.4	6.5	4.0
Max column id	0.25 mm (30 m)	0.32 mm (30 m)	0.53 mm (30 m)	0.53 mm (30 m)

\* Total gas flow into the MSD: column flow plus reagent gas flow (if applicable).

† Expect degradation of spectral performance and sensitivity.

Table 9 lists typical flows resulting from selected carrier and reagent gas source pressures.

**Table 9** 5975 Series carrier and reagent gases

Carrier and reagent gas requirements	Typical pressure range	Typical flow (mL/min)
Helium (required) (column and split flow)	345 to 552 kPa (50 to 80 psi)	20 to 50
Hydrogen (optional) <sup>*</sup> (column and split flow)	345 to 552 kPa (50 to 80 psi)	20 to 50

**Table 9** 5975 Series carrier and reagent gases (continued)

Carrier and reagent gas requirements	Typical pressure range	Typical flow (mL/min)
Methane reagent gas (required for CI operation)	103 to 172 kPa (15 to 25 psi)	1 to 2
Isobutane reagent gas (optional)	103 to 172 kPa (15 to 25 psi)	1 to 2
Ammonia reagent gas (optional)	34 to 55 kPa (5 to 8 psi)	1 to 2
Carbon dioxide reagent gas (optional)	103 to 138 kPa (15 to 20 psi)	1 to 2

\* Hydrogen gas can be used for the carrier gas but specifications are based on helium as the carrier gas. Please observe all hydrogen gas safety cautions.

## Other Considerations

### Exhaust venting requirements for the GC/MSD

Vent the MSD externally to the building via an ambient-pressure vent system, within 460 cm (15 ft) of both the GC split vent and MSD foreline pump, or vent to a fume hood. Note that an exhaust vent system is not part of the building environmental control system, which recirculates air. Exhaust venting must comply with all local environmental and safety codes. Contact your Environmental Health & Safety (EHS) specialist.

For GCs with the exhaust deflector option installed, the exhaust is about 65 ft<sup>3</sup>/min (1.840 m<sup>3</sup>/min). Without the deflector, the exhaust rate is about 99 ft<sup>3</sup>/min (2.8 m<sup>3</sup>/min). The deflector outlet diameter is 10 cm (4 in).

## Basic tools

The GC/MSD comes with a few basic tools and consumables depending on the specific inlet and detector that you ordered. Below is a general list of what comes with the instrument.

<b>Tool or consumable</b>	<b>Used for</b>
T10 and T20 Torx wrenches	Removing tray. Removing covers to access gas control modules, traps, and pneumatic connections.
1/4-inch nut driver	FID jet replacement.
FID flow measuring insert	FID troubleshooting.
Column cutter, ceramic or diamond	Column installation.
1/8-inch Tee, Swagelok, brass	Connect gas supplies.
1/8-inch nuts & ferrules, Swagelok, brass	Connect gas supplies.
Inlet septa appropriate for type	Inlet seal.
Inlet insert or liner	Contains sample during vaporization in inlet.
1.5-mm and 2.0-mm hex driver	Source maintenance (disassembly).
Tool bag	Holding GC and MS tools.
Q-Tips	Cleaning source parts.
Cloths	Keeping surfaces and parts clean.
Gloves	Reducing contamination on GC and MSD parts.
Funnel	Changing oil.
Hex key, 5 mm	Removing oil plug and screws in safety shield handle.

## 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.



**Agilent Technologies**

First Edition March 2007

This information is subject to change without notice.

© Agilent Technologies, Inc. 2007

Printed in U.S.A., March 2007

G3170-90025



G3170-90025