

**Site Preparation Specification (Preliminary rev. 20060310)**

**Preliminary—the following information is believed to be correct but may change prior to shipment.**

**Purpose of Procedure**

To ensure that the installation site is properly evaluated and prepared with the appropriate utilities, consumables and supplies for the successful installation of Agilent instruments and systems.

**Customer Responsibilities**

Customers should ensure that all necessary operating supplies, consumables and usage dependent items such as columns, vials, syringes, solvents and buffers required for the successful installation of instruments and systems are available. Installation sites should be prepared in accordance with the following specifications.

**Important Information**

This checklist is designed to be used in conjunction with the Agilent 6410A Triple Quadrupole LC/MS Site Preparation Manual. If you have problems providing any of the following, please contact your local Agilent sales office for assistance. Assistance with user specific applications may be provided but should be contracted separately. Users of the instrument should be present throughout the installation and familiarization otherwise important operational, maintenance and safety information may be missed.

**Procedure Checklist**



**Agilent G6410A Triple Quad Mainframe:**  
*footprint:\**

Depth: 66 cm      Width: 111 cm  
          26 in            43.5 in

**Tick Boxes**

*Maximum cabinet dimensions:\*\**

Weight: 110 kg      Height: 47 cm  
          242 lb            18.5 in  
Depth: 79 cm      Width: 128 cm  
          31 in              50.5 in

**E2M28 Mechanical Roughing Pump:**

Weight: 40 kg      Height: 28 cm  
          88 lb              11 in  
Depth: 58 cm      Width: 18 cm  
          23 in              7.1 in

**Agilent G1947B APCI Source:**

Weight: 1.7 kg      Height: 23 cm  
          3.75 lb            9.2 in  
Depth: 13.0 cm      Width: 18 cm  
          5.1 in            7.1 in

**Agilent G1948A/B API-ES Source:**

Weight: 1.7 kg      Height: 17 cm  
          3.75 lb            6.8 in  
Depth: 9.5 cm      Width: 18.0 cm  
          3.7 in            7.1 in

**Agilent G1978B Multimode Source:**

Weight: 2.29 kg      Height: 17 cm  
          5.05 lb            6.8 in  
Depth: 13.0 cm      Width: 18.0 cm  
          5.1 in            7.1 in

\* The footprint dimensions represent the minimum dimensions of the supporting surface. This surface must also be relatively vibration-free and capable of supporting at least 114 kg (250 lbs).

\*\* Maximum cabinet dimensions are for an Agilent G6410A with an Agilent G1947B APCI, or G1978B Multimode source installed. At least 30 cm (1 ft) to the left of the cabinet and at least 55 cm (1.8 ft) above the cabinet must be added to these dimensions to provide adequate instrument access. In addition, the Agilent G6410A has a 183 cm (6 ft) metal hose connected from the back of the instrument to its mechanical roughing pump.

**Site Preparation Specification (Preliminary rev. 20060310)**

**Power**



**Americas & Japan:** 200 to 220 VAC; 2500 VA max, 50/60 Hz, 15 A circuit<sup>1</sup>  
**Europe & Asia Pacific:** 220 to 240 VAC; 2500 VA max, 50 Hz, 15 A circuit<sup>1</sup>

<sup>1</sup>Single outlet for G6410A LC/MS.

**Tick Boxes**



**Heat Dissipation**



**Output:** 1100 Watts (3700 BTU / hour)<sup>1</sup>

<sup>1</sup>Approximately 600 Watts are removed with the source exhaust.



**Nitrogen Gas Supply**



**The G6410A LC/MS will require two sources of nitrogen—one for the Nebulizer and Drying Gas, and the other for the Collision Cell.**

**Nebulizer and Drying Gas:**

**Purity:** 99.5% or better - Gas cylinder  
 98.0% or better - N<sub>2</sub> gas generator or liquid N<sub>2</sub> Dewar.

Balance of impurity should consist of oxygen and/or argon. Gas must be hydrocarbon free (< 0.1 ppm).

**Outlet Pressure:** 80-100 psi. A 1/4" Swagelok outlet (male) fitting is required to connect the LC/MS.

**Volume:** Up to 15 liters/min.



**Collision Cell:**

**Purity:** 99.999% or better  
**Outlet Pressure:** 10-30 psi. A 1/8" Swagelok outlet (male) fitting is required to connect to the LC/MS.  
**Volume:** Up to 2 sccm (2 cc per min at atmospheric pressure)



**Laboratory Supply Requirements**



**Mobile Phases:** Water, Methanol, Isopropanol, Acetonitrile<sup>1</sup>

**Purity:** HPLC-grade or better

**Buffers:** Ammonium Formate

**Acids:** Acetic or Formic Acid

**Purity:** Ammonium formate, 97% or better  
 Acetic acid, 99.7% or better  
 Formic acid, 96% or better

<sup>1</sup>Methanol/water required for G6410A installation.





Site Preparation Specification (Preliminary rev. 20060310)

Tick Boxes

*Exhaust Venting Requirements*



**Capacity:** Up to 15 liters/min. total.

**Connections:** Separate 1/2" hose barbs required for rough pump and ion source (ES, APCI, or MM).

<sup>1</sup>A 7.6m (20 ft.) length of 1/2 inch i.d. Tygon™ tubing is included for venting source exhaust (drain bottle) and rough pump. (Sufficient for two 3.8 m (10 ft) lengths.)

*Environmental Requirements*



**Temperature:** 15 to 35 °C (59 to 95 °F)  
at constant temperature (variations < 3 °C/hr).

**Humidity:** 20 to 85 %