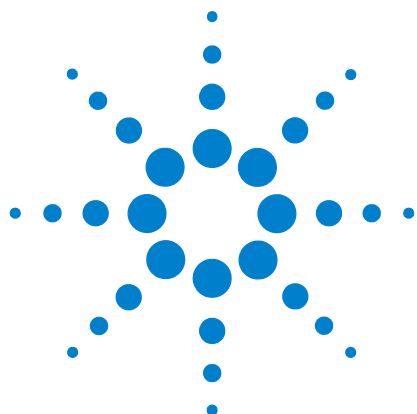




Agilent 7696A Sample Prep WorkBench

Site Preparation



Agilent Technologies

Notices

© Agilent Technologies, Inc. 2010

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

G4529-90016

Edition

First edition, December 2010

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.

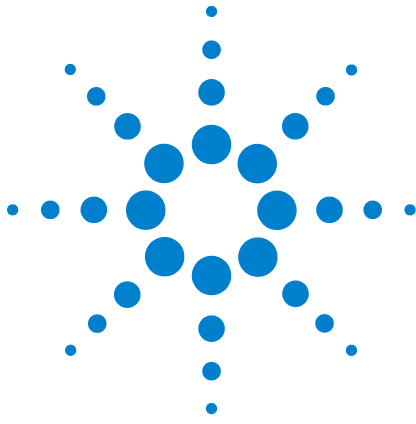
Safety Notices

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.



Agilent 7696A Sample Prep WorkBench Site Preparation

Customer Responsibilities 4

Dimensions and Weight 5

Power Consumption 6

Environmental Conditions 8

Maximum Length of Cables 9

Site LAN Network 10

Basic Tools 12

This section outlines the space and resource requirements for Agilent 7696A Sample Prep WorkBench (WorkBench) installation. For a successful and timely installation of the instrument, the site must meet these requirements before beginning installation. Necessary supplies (operating supplies, consumables, and other usage-dependent items such as vials and solvents) must also be available. Refer to the Agilent website at www.agilent.com/chem for the most up-to-date listing of supplies and consumables.



Customer Responsibilities

The specifications in this manual outline the necessary space, electrical outlets, operating supplies, consumables, and other usage-dependent items such as vials and solvents required for the successful installation of instruments and systems.

If Agilent is delivering installation and familiarization services, users of the instrument should be present throughout these services; otherwise, they will miss important operational, maintenance, and safety information.

If Agilent is delivering installation and familiarization services, 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.

Dimensions and Weight

Prepare the laboratory bench space before the system arrives. Pay special attention to the total height requirements. It is recommended that you leave at least 10 cm (4 in) of clearance for tower removal/installation. Avoid bench space with overhanging shelves.

See [Figure 1](#) for an example of total dimensions when used.

Table 1 Required height, width, depth, and weight

Product	Height	Width	Depth	Weight
Agilent 7696A Sample Prep WorkBench	70.4 cm (27.72 in)	67.6 cm (26.62 in)	49.3 cm (19.41 in)	40.8 kg (90 lbs)

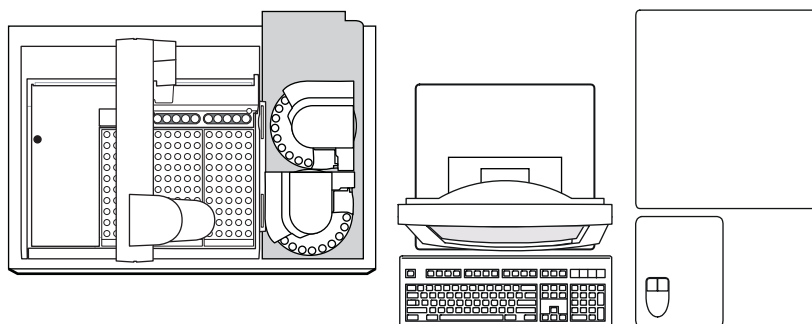


Figure 1 Example bench space needed for a 7696A WorkBench system with computer and printer.

A system that includes a 7696A WorkBench and computer would require about 130 cm (52 in) of bench space. Allowing for operational access and a printer, a total of 174 cm (68 in) of bench space should be available. Some repairs to the system will also require access to the back of the instrument.

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 to which the unit ships.
- The voltage requirements for your instrument are printed near the power cord attachment.
- The electrical outlet for the unit should have a dedicated ground.
- All instruments should be on a dedicated circuit.
- Power line conditioners should not be used with Agilent instruments.

Table 2 Power consumption

Product	Line voltage (VAC)	Frequency (Hz)	Maximum continuous power consumption (VA)	Current rating (amps)
7696A Sample Prep WorkBench	Americas: 120 single phase (–10% / +10%)	50–60	800	7.4
	220/230/240 single/split phase (–10% / +10%)	50–60	800	4.0/3.9/3.7

WARNING

Do not use extension cords with Agilent instruments. Extension cords normally are not rated to carry enough power and can be a safety hazard.

Although your instrument should arrive with the parts needed for operation in your country, compare its voltage requirements with those listed in Table 2. If the voltage option you ordered is not suitable for your installation, contact Agilent Technologies.

CAUTION

A proper earth ground is required for instrument operations. Any interruption of the grounding conductor or disconnection of the power cord could cause a shock that could result in personal injury.

To protect users, the metal instrument panels and cabinet are grounded through the three-conductor power line cord in accordance with International Electrotechnical Commission (IEC) Requirements.

The three-conductor power line cord, when plugged into a properly grounded receptacle, grounds the instrument and minimizes shock hazard. A properly grounded receptacle is one that is connected to a suitable earth ground. Be sure to verify proper receptacle grounding. The instrument requires an isolated ground.

Environmental Conditions

Operating the instrument within the recommended ranges optimizes instrument performance and lifetime. Performance can be affected by sources of heat and cold from heating, air conditioning systems, or drafts. See [Table 3](#). The conditions assume a noncondensing, noncorrosive atmosphere.

Table 3 Environmental conditions for operation and storage

Product	Conditions	Operating temp range	Operating humidity range	Maximum altitude
Agilent 7696A WorkBench	Operation	15 to 35 °C (59 to 95 °F)	5 to 95%	4,300 m
	Storage	–20 to 70 °C (–4 to 158 °F)	0 to 95%	

Maximum Length of Cables

The distance between system modules may be limited by some of the cabling and the vent or vacuum hoses. Plan to set up the work space as shown in [“Dimensions and Weight”](#).

- The length of the Agilent-supplied remote cable is 2 meters (6.6 feet).
- The length of the Agilent-supplied LAN cable is 10 meters (32.8 feet).
- The lengths of the power cords are 2 meters (6.6 feet).

Site LAN Network

If you intend to connect your system to your site's LAN network, you must have an additional shielded twisted pair network cable.

NOTE

Agilent Technologies is not responsible for connecting to or establishing communication with your site LAN network. The representative will test the system's ability to communicate on a mini-hub or LAN switch only.

NOTE

The IP addresses assigned to the instrument(s) must be fixed (permanently assigned) addresses. If you intend to connect your system to your site's network, each piece of equipment must have a unique, fixed (static) IP address assigned to it.

Laboratory Safety Practices

When handling/using chemicals for preparation or use within the Workbench unit, 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, 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.

WARNING

Use extreme caution when handling heavy parts. A two-person lift is recommended. Failure to perform a two-person lift may result in personal injury.

Basic Tools

The WorkBench comes with a few basic tools and consumables. Below is a general list of what comes with the instrument.

Table 4 Basic tools

Tool or consumable	Used for
Crimper, 20 mm seals	Crimp-top sample vials
T-10 Torx key driver	Maintenance
Short Arm L Key T-20 Torx	Maintenance
Cross Head Screw Driver	Maintenance
Calibration Vial	Calibration procedure
Screw Cap Vials Clear (100)	
Blue Screw Caps	