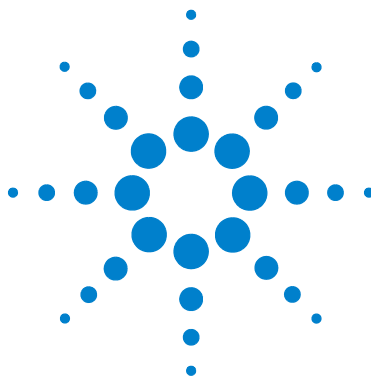




Agilent 7697A Headspace Sampler

Getting Started



Agilent Technologies

Notices

© Agilent Technologies, Inc. 2011

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

G4556-90010

Edition

First edition, January 2011

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

If software is for use in the performance of a U.S. Government prime contract or subcontract, Software is delivered and licensed as “Commercial computer software” as defined in DFAR 252.227-7014 (June 1995), or as a “commercial item” as defined in FAR 2.101(a) or as “Restricted computer software” as defined in FAR 52.227-19 (June 1987) or any equivalent agency regulation or contract clause. Use, duplication or disclosure of Software is subject to Agilent Technologies’ standard commercial license terms, and non-DOD Departments and Agencies of the U.S. Government will receive no greater than Restricted Rights as defined in FAR 52.227-19(c)(1-2) (June 1987). U.S. Government users will receive no greater than Limited Rights as defined in FAR 52.227-14 (June 1987) or DFAR 252.227-7015 (b)(2) (November 1995), as applicable in any technical data.

1 Introduction

This document provides an overview of the Agilent Technologies 7697A Headspace Sampler system, available user information materials, and installation steps.



Figure 1 Agilent 7697A Headspace Sampler with 7890A GC and 7693A ALS



Where to Find Information



In addition to this document, Agilent provides several learning products that document how to install, operate, maintain, and troubleshoot the 7697A HS. This information can be found on the Agilent Technologies GC and GC/MS Hardware User Information & Instrument Utilities Software DVDs that ship with your instrument.



The Agilent Technologies GC and GC/MS Hardware User Information & Instrument Utilities Software DVDs that ship with your instrument provides an extensive collection of online help, videos, and books for current Agilent gas chromatographs, mass selective detectors, and GC samplers. Included are localized versions of the information you need most, such as:

- Familiarization information
- Safety and regulatory guides
- Site preparation information
- Installation information
- Operating guides
- Maintenance information
- Troubleshooting details

Site Preparation

Start here. The *Site Preparation* manual lists all of the parts and requirements needed to install and set up the headspace sampler.

Installation and First Startup

The *Installation and First Startup* manual provides all of the steps required to successfully install and check out the headspace sampler in your system.

Safety

Read the *Safety* manual before using the headspace sampler, and whenever needed. The *Safety* manual describes the important safety precautions to follow when installing and using the headspace sampler, as well as important regulatory information.

Familiarization

The *Familiarization* guide explains how to use the headspace sampler using the Agilent GC ChemStation software. It shows you how to set up a method and run sequences of samples. Agilent shares its instrument control technology between its data systems, so users of other Agilent data systems will still find this manual useful.

Operation

The *Operation* guide describes the typical hardware and software tasks needed to operate the headspace sampler on a daily basis. Use this manual to familiarize yourself with the instrument and software, and to find part numbers for consumables such as vials, vial labels, caps, and so forth.

Advanced Operation

Method developers should use the *Advanced Operation* guide to learn about the full capabilities of the headspace sampler. Standalone users (users without Agilent data system control) will find all keypad functionality here. Also included are procedures and concepts for configuration, energy conservation, advanced method development, converting methods made for other headspace sampler models, and using the 7697A early maintenance feedback (EMF) features.

Maintenance

When needed, refer to the *Maintenance* manual for routine maintenance schedule and all related user maintenance procedures. The *Maintenance* manual also includes a complete list of user-replaceable parts.

Troubleshooting

When needed, refer to the *Troubleshooting* manual for hints and procedures to improve performance and to correct sampler-related chromatographic problems.

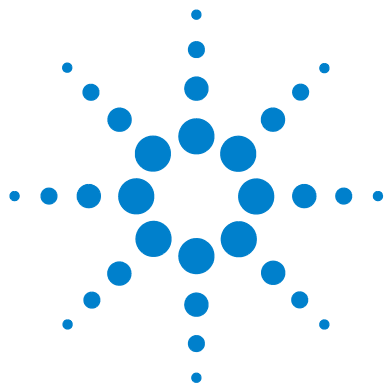
The Agilent 7697A Headspace Sampler

Front view, G4557A (111 vial model)



Back view, G4557A (111 vial model)





2 Installation Summary

Following is a brief overview of the Agilent 7697A Headspace Sampler installation steps. For detailed instructions, see the *Agilent 7697A Installation and First Startup* guide on the Agilent GC and GC/MS Hardware User Information & Instrument Utilities Software DVD.

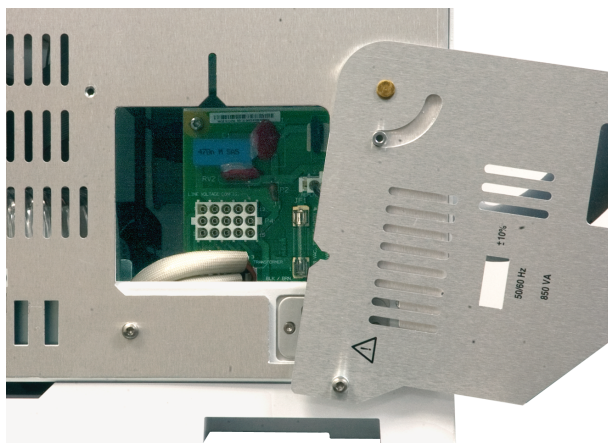


Installation Summary

- 1 Unpack.
- 2 Place on bench.



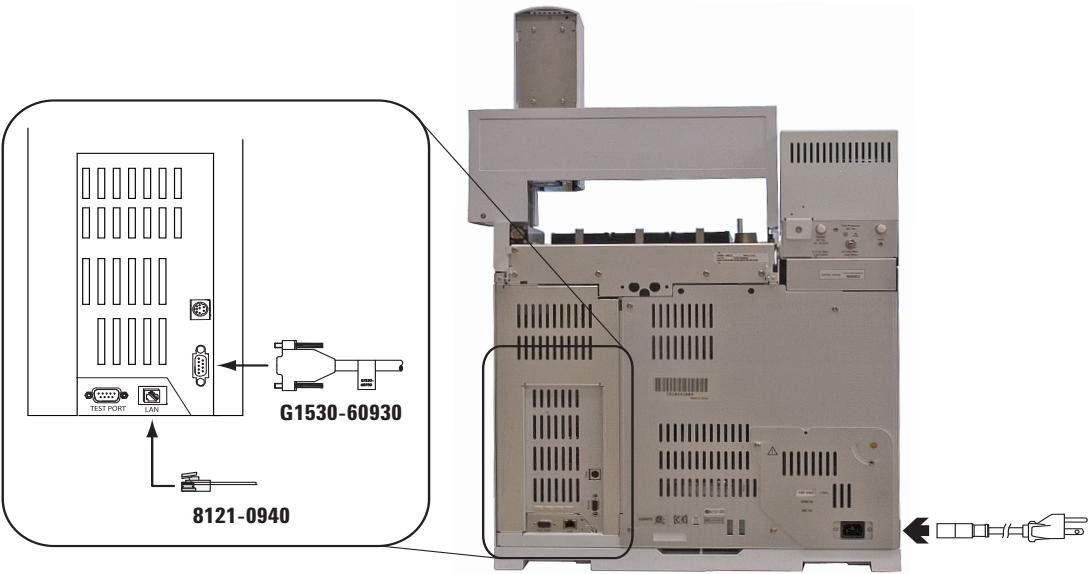
3 Install power configuration plug.



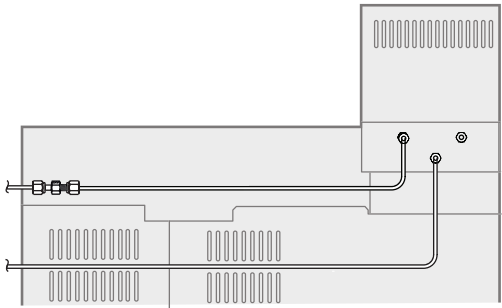
4 Install vial tray (111 vial model) and vial racks.



5 Connect cabling.

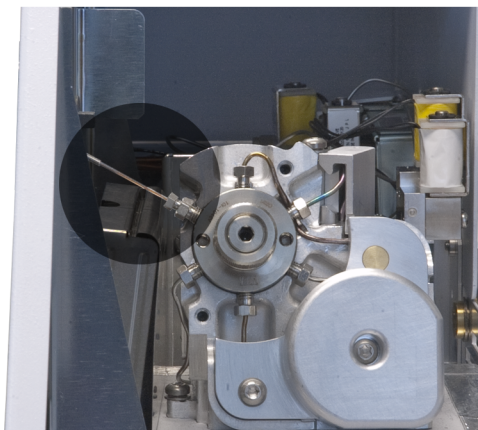


6 Connect gases.



Gas stream	Recommended starting supply pressure
Vial pressurization	410 kPa (80 psi)
Carrier (if using optional G4562A 7697A Carrier Gas EPC Module Accessory)	410 kPa (80 psi)

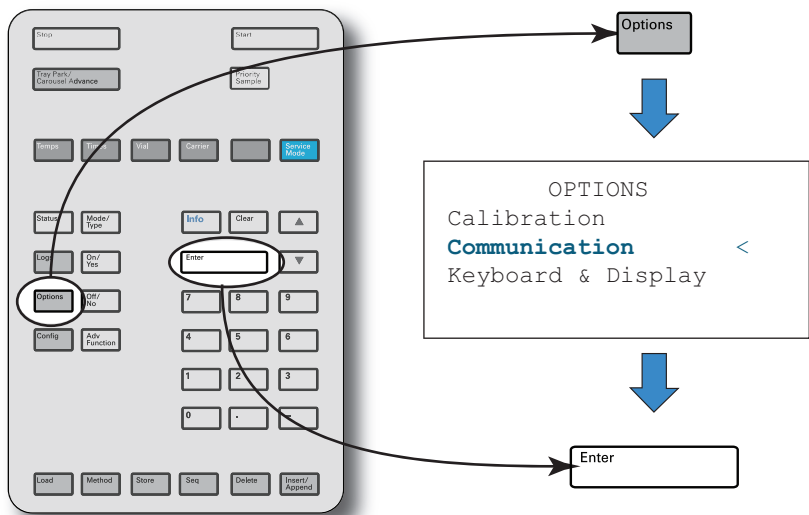
7 Install and connect the transfer line.



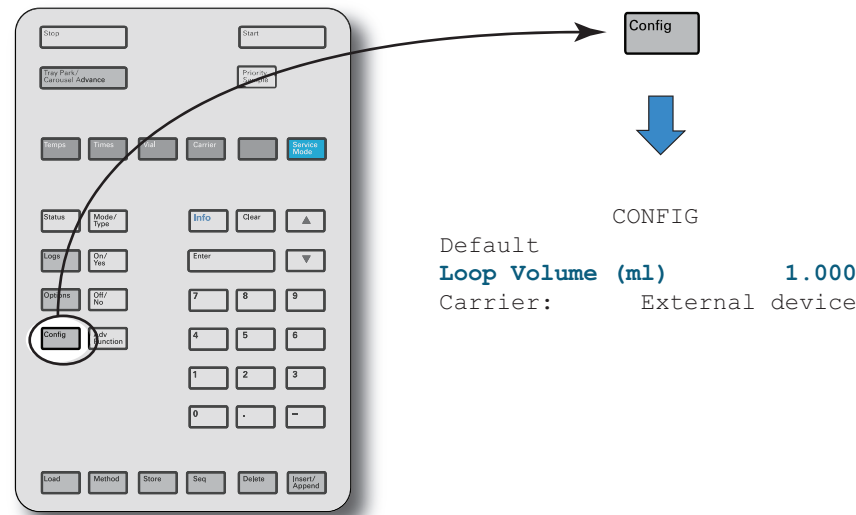
8 Turn on.



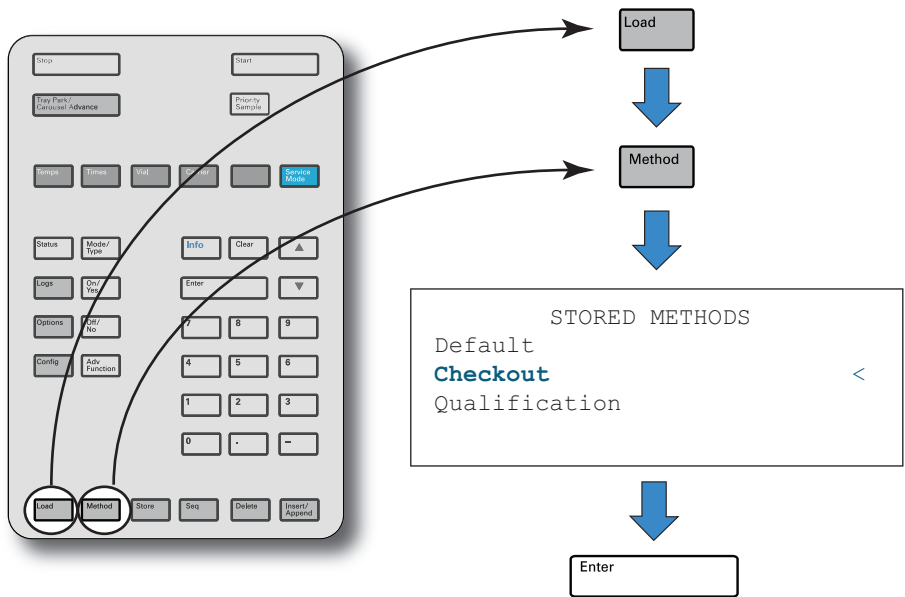
9 Set IP address.



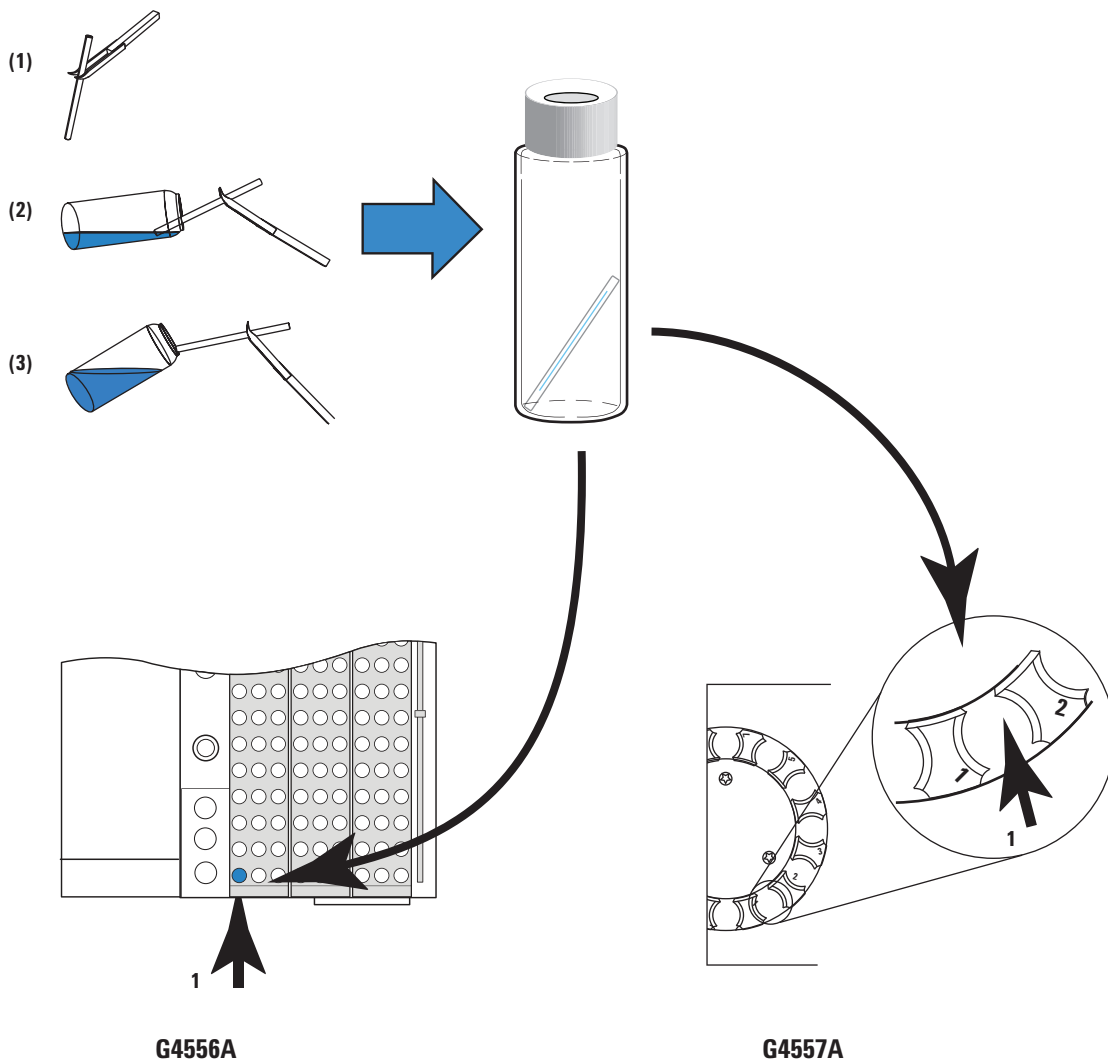
10 Configure.



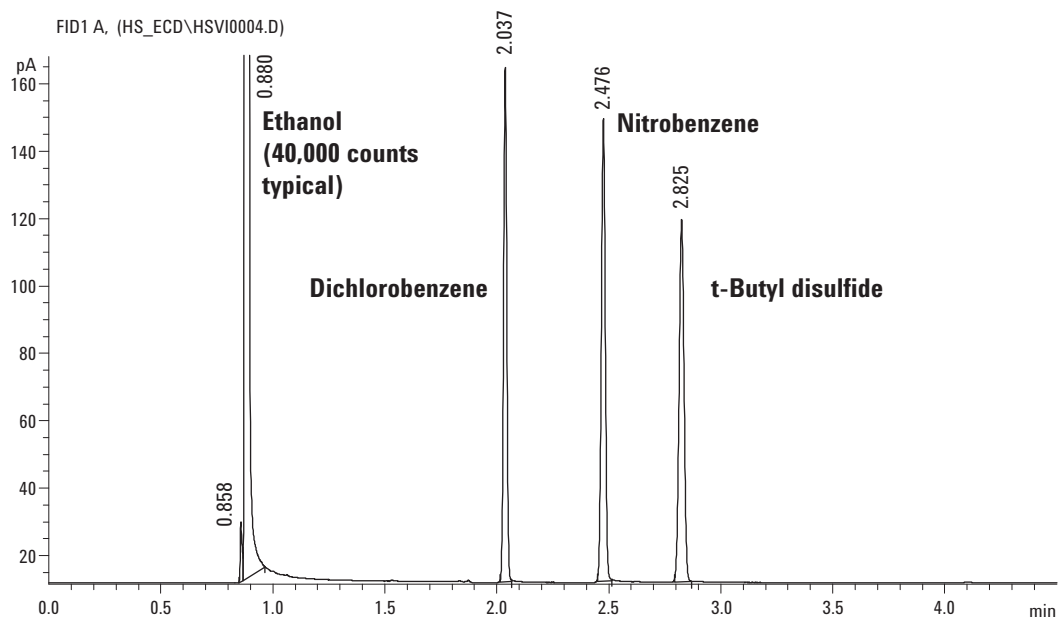
11 Load checkout method.



12 Prepare sample and load sample vial into tray.



13 Run. Compare results.





Agilent Technologies

© Agilent Technologies, Inc.

Printed in USA or China

First edition, January, 2011



G4556-90010