

Agilent 8697 Headspace Sampler

## **Firmware Bulletin – Revision 1.6.0.5**



# Notices

## Copyright

© Agilent Technologies, Inc. 2025

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., governed by United States and international copyright laws.

### Agilent Technologies, Inc.

5301 Stevens Creek Blvd.  
Santa Clara, CA 95051  
USA

## Firmware Revision

This guide is valid for the 1.6.0.5 of  
[Agilent 8697 HS firmware](#)

## 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 conflicts 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.

## 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 damage the product or lose 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 could result in personal injury or death if not correctly performed or adhered to. Do not proceed beyond a **WARNING** notice until the indicated conditions are fully understood and met.

|   |   |    |
|---|---|----|
| 1 | Introduction  | 4  |
| 2 | Agilent 8697 HS Firmware – What's new?              | 5  |
| 3 | Compatibility                                       | 6  |
| 4 | Support Information                                 | 7  |
| 5 | Resolved Issues                                     | 8  |
| 6 | Agilent 8697 HS Firmware - Previous Version Changes | 9  |
| 7 | Changelog   | 11 |

## Introduction

The purpose of this document is to provide users of the Agilent 8697 Headspace Sampler guidance to relevant support documentation that contains further levels of information and detail.

**Table 1** Terms and abbreviations used in this document

| Terms | Description                               |
|-------|---|
| SSB   | Software Status Bulletin                  |
| CDS   | Chromatographic Data System               |
| U     | Update                                    |
| SP    | Service Pack                              |
| HF    | Hotfix                                    |
| SR    | Service Release                           |
| [xx]  | Final Build Numbers of a released product |
| DSA   | Data System Adapter                       |

### For our Regulated Customers

When any change is made to Agilent software/firmware, the validation status of the software needs to be re-established by the user/customer.

Whenever software/firmware changes, a validation analysis should be conducted to validate an individual change and determine the extent and impact of that change on the entire system.

### Trademarks

Microsoft, Windows, Windows Server, and Microsoft .NET are registered trademarks owned by Microsoft Corporation.

## 2

## Agilent 8697 HS Firmware – What's new?

### What's new in 8697 HS Firmware?

- *Defects fixes and Enhancements.*

### Supported Agilent Data Systems

- OpenLab CDS (2.8, 2.7)
- OpenLab CDS ChemStation Edition (LTS 01.11, C.01.11)
- GC/MS MassHunter Acquisition (13.1, 10.2 SR3, 10.2)
- OpenLab CDS EZChrom Edition (A.04.10 (Update 3))

### Supported Agilent GC Driver

- Agilent GC Driver (4.4.92 or later)

### Supported Agilent HS Driver

- Agilent HS Driver (4.4.28 or later)

### Supported Agilent GC Firmware

- Agilent GC Firmware (3.2.0.130 or later) for the 8890, 8860, and Intuvo 9000 GCs

#### Important Information:

1. Please use Firmware Update Utility 3.2.8.0 or newer.

### Non-Agilent Chromatography Data System (CDS)

- The 3rd-party CDS software vendor is responsible for compatibility testing with the respective CDS revision.
- The 3rd-party CDS software vendor defines the minimum firmware revision required for CDS compatibility.
- The 3rd-party CDS release notes issued by the respective CDS vendor may use different terminology for the firmware requirements, such as "tested firmware," "supported firmware," "firmware requirements," "minimum tested firmware," etc.

Agilent recommends using the latest firmware revision to avoid interoperability issues.

### **OQ/PV - Validation Information**

- If a firmware upgrade has been performed, no revalidation of the module/system is normally required. This is stated in the released Service Note applicable to the release.
- In addition, this information is stated further in this document.

### **Agilent on the web**

Visit [Agilent's website](#) for up-to-date information on Agilent Systems.

## 5 Resolved Issues

The current firmware for the 8697 HS is 1.6.0.5.

*Released in Nov 2025*

[https://download.chem.agilent.com/firmware/8697\\_FW\\_1.6.0.5.tgz.gpg](https://download.chem.agilent.com/firmware/8697_FW_1.6.0.5.tgz.gpg)

To download this version, use GC Firmware Update Utility 3.2.8.0 or newer.

*Updating to this firmware does not require instrument revalidation.*

Service Note Number: D0139638

**Table 2** Resolved Issues (8697 HS Firmware)

| Issue ID | Current Behavior   | Resolution |
|----------|--|------------|
| 046503   | 8697A does not have Argon as vial pressure and carrier gas | Resolved   |
| 046725   | Headspace EPC Module FW Update Support                     | Resolved   |



## 6

## Agilent 8697 Firmware - Previous Version Changes

### Previous Firmware Versions.

This section details previous versions of 8697 firmware giving release notes.

#### **1.5.0.4 8697 HS Firmware.**

***Released in Jan 2025***

*Updating to this firmware does not require instrument revalidation.*

Changes/Fixes made since the last version. (1.4.0.10):

| Issue ID | Current Behavior  | Resolution |
|----------|---|------------|
| 044302   | Aborted headspace injection leads to unnecessary "Headspace pressure pre-shutdown" warning                                    | Resolved   |
| 045911   | HS FW needs to change to support later versions of EPCs   | Resolved   |
| 046379   | System report "Headspace: Pressure pre-shutdown" and GC beep for few minutes during purge time with no impact to sequence run | Resolved   |

#### **1.4.0.10 8697 HS Firmware.**

***Released in Jul 2023***

*Updating to this firmware does not require instrument revalidation.*

Changes/Fixes made since the last version. (1.3.0.59):

| Issue ID | Current Behavior   | Resolution |
|----------|--|------------|
|          | Initial support for 8697-XL  | Resolved   |
| 041998   | Use new board ID bits to implement a faster glitch filter on overcurrent detection | Resolved   |
| 044042   | Tray calibration with broken axis corrupts RAM                                     | Resolved   |
| 044071   | To add 8697 oven fan check   | Resolved   |
| 037180   | Fan speed dependent on settings and conditions.                                    | Resolved   |
| 043747   | EPC Validation   | Resolved   |

## Agilent 8697 Firmware - Previous Version Changes

|        |  |          |
|--------|--|----------|
| 045128 | Enable dynamic oven fan error checking | Resolved |
|--------|--|----------|

### 1.3.0.59 8697 HS Firmware.

*Released in Feb 2023*

*Updating to this firmware does not require instrument revalidation.*

Changes/Fixes made since the last version. (1.2.0.36):

| Issue ID | Current Behavior  | Resolution |
|----------|---|------------|
| 041844   | Unable to move the shutter from half open to full open without having an error                  | Resolved   |
| 042198   | If y-axis cannot reach the right rear calibration post it now return a calibration failed error | Resolved   |
| 042248   | The quick calibration check at power on should return an error if one of the 3 axes is broken   | Resolved   |
| 042551   | The calibration status command 00AS does not work   | Resolved   |
| 042587   | Bryce pneumatic restriction pressure has issue  | Resolved   |
| 042630   | The Home check for z-axis always fails  | Resolved   |
| 042631   | The z-axis motor grinds at the top of the travel at power-on                                    | Resolved   |
| 042709   | Time off in Bryce tray  | Resolved   |
| 042722   | Tray: Calibration commands not respecting stop  | Resolved   |
| 042756   | Headspace tray extended calibration   | Resolved   |
| 042776   | Oven calibration resets the shutter offset value  | Resolved   |
| 042939   | The PRT being used for the chiller plate has a wide 700 deg C temperature span                  | Resolved   |
| 043015   | The carousel power during lifter initialization was inadvertently reduced                       | Resolved   |
| 043053   | Gripper pad maintenance flag warning came up  | Resolved   |
| 043246   | There is a sudden drop in the vial pressure actual while filling a vial in constant volume mode | Resolved   |
| 043311   | Calibration post fail   | Resolved   |

## Agilent 8697 Firmware - Previous Version Changes

|        |   |          |
|--------|---|----------|
| 043315 | Incorrect "tray type mismatch" error after stalling both x and z axes                                     | Resolved |
| 043511 | A DEFOA board from ATS had incorrect EEPROM contents  | Resolved |
| 043671 | In final test, we rely on the oven fan spinning by default to measure whether the fan is working properly | Resolved |
| 043675 | The Diag and ELVDS status LEDs are much brighter than the other status LEDs                               | Resolved |
| 041624 | Speed up tray calibration   | Resolved |

### 1.2.0.36 8697 HS Firmware.

*Released November 2021*

*Updating to this firmware does not require instrument revalidation.*

Changes/Fixes made since the last version. (Initial version):

| Issue ID | Current Behavior  | Resolution |
|----------|---|------------|
| 042333   | Left rear post calibration fails when recalibration if gripper offsets change in X or Y by large amount | Resolved   |
| 039896   | Insufficient vial pressurization gas pressure fails to make Bryce "Not ready"                           | Resolved   |
| 040114   | Park button does not work after disconnecting / reconnecting tray cable                                 | Resolved   |
| 041035   | System calibration fails with wrong vial type error during shutter offset measurement                   | Resolved   |
| 041045   | Bryce EPC unable to reach 200ml/min of N2 below 60psi input pressure                                    | Resolved   |
| 041070   | Pressure pre-shutdown warning occurs during loop fill for constant volume of 0 ml is active             | Resolved   |
| 041631   | 8697 Transfer Line has "Heating too slowly" error reported  | Resolved   |

## Revision History

Table 19 Revision History

| Revision | Release Date | What's New/What Changed |
|----------|--------------|-------------------------|
| Rev 1.00 | 05 Nov 2025  | Initial release         |

[www.agilent.com](http://www.agilent.com)

© Agilent Technologies, Inc. 2025

Edition Nov/2025

D0139638