



Agilent 8850 GC Firmware

## **Release Notes – Revision 3.2.5.4**

# Notices

## Copyright

© Agilent Technologies, Inc. 2026

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.

### Agilent Technologies, Inc.

5301 Stevens Creek Blvd.  
Santa Clara, CA 95051  
USA

## Firmware/Software Revision

This guide is valid for the 3.2.5.4  
revision of *Agilent 8850 GC 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 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.

## 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.

# Table of Contents

|   |                                      |    |
|---|--------------------------------------|----|
| 1 | Document History                     | 4  |
| 2 | Introduction                         | 5  |
| 3 | Compatibility                        | 6  |
| 4 | Support Information                  | 7  |
| 5 | Agilent 8850 GC System – What's new? | 8  |
| 6 | Resolved Issues in revision 3.2.5.4  | 9  |
| 7 | Previous Versions                    | 10 |
| 8 | Firmware Update Program              | 16 |

# 1

## Document History

The table below lists all changes that have been made to this document.

| Date                            | Version  | Description                  |
|---------------------------------|----------|------------------------------|
| 11 <sup>th</sup> September 2024 | Rev A.01 | Initial Release              |
| 11 <sup>th</sup> February 2025  | Rev A.02 | Added 3.0.3.21 Release Note  |
| 20 <sup>th</sup> June 2025      | Rev A.03 | Added 3.1.0.380 Release Note |
| 20 <sup>th</sup> September 2025 | Rev A.04 | Added 3.2.0.130 Release Note |
| 7 <sup>th</sup> January 2026    | Rev A.05 | Added 3.2.2.3 Release Note   |
| 25 <sup>th</sup> March 2026     | Rev A.06 | Added 3.2.3.3 Release Note   |
| 31 <sup>st</sup> July 2026      | Rev A.07 | Added 3.2.5.4 Release Note   |

## 2

## Introduction

The purpose of this document is to provide users of the Agilent [8850 GC](#) system 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                       |

# 3

## Compatibility

### Supported Agilent Data Systems

OpenLab CDS (3.0, 2.8, 2.7, 2.6, 2.5, 2.4 or higher)

OpenLab CDS ChemStation Edition (C.01.10, C.01.09, LTS.01.11)

GC/MS MassHunter Acquisition (13.1, 13.0 or higher)

OpenLab CDS EZChrom Edition (A.04.10 Update 3)

### Supported Agilent GC Driver

Agilent GC Driver (4.4.117, 4.5.228 or higher)

## 4

## Support Information

- Agilent recommends using always the latest firmware revision of a firmware in order to avoid interoperability issues.
- Generally, Agilent recommends keeping the GC instrument firmware always current.

### Notes for Agilent 8850 controlled by non-Agilent Chromatography Data Systems (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.

### OO/PV - Validation Information

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



## 5

# Agilent 8850 GC System – What's new?

### What's new in 8850 GC firmware 3.2.5.4?

- Bug fixes as detailed below.

## 6

## Resolved Issues in revision 3.2.5.4

The current shipping firmware for the 8850 is 3.2.5.4

Released on 31<sup>st</sup> March 2026

[https://download.chem.agilent.com/firmware/8850\\_FW\\_3.2.5.4.tgz.gpg](https://download.chem.agilent.com/firmware/8850_FW_3.2.5.4.tgz.gpg)

To download this version, use GC Firmware Update Tool 3.2.8.0 or greater.

Updating to this firmware does not require instrument requalification.

Service Note Number: D0151332

### Resolved Issues:

**Table 2 Resolved Issues (Agilent 8850 GC System 3.2.5.4)**

| Issue ID | Current Behavior                                                                                                       | Resolution |
|----------|------------------------------------------------------------------------------------------------------------------------|------------|
| 047112   | Headspace MHC Didn't Move on from Second Extraction                                                                    | Resolved   |
| 047873   | Align the 8697 sample overlap behavior with 7697                                                                       | Resolved   |
| 048093   | Gather Logs Should Include MF Command Error Info When Headspace Configured                                             | Resolved   |
| 048182   | Uploading New PID File Using GC FW Update Utility Fails                                                                | Resolved   |
| 048561   | Developer Exceptions shown in user logs                                                                                | Resolved   |
| 048668   | Method parameters are not actually downloaded and applied to GC (with 8697 HS connection but no 8697 HS configuration) | Resolved   |
| 048779   | Headspace MHC Didn't Move on from Second Extraction                                                                    | Resolved   |
| 048873   | New style EPC is missing the SF09 Reference rawSTP value in EEPROM, causing Gas Identification error                   | Resolved   |
| 048911   | New style EPC N2 SSL w/HS High Flow Noise                                                                              | Resolved   |
| 048912   | New style EPC N2 MMI w/HS High Flow Noise                                                                              | Resolved   |
| 048920   | PID Updating Incorrectly                                                                                               | Resolved   |

# 7

## Previous Versions

This section details previous versions of 8850 firmware giving release notes.

### 3.2.3.3 Firmware

Released on 25<sup>th</sup> March 2026

[https://download.chem.agilent.com/firmware/8850\\_FW\\_3.2.3.3.tgz.gpg](https://download.chem.agilent.com/firmware/8850_FW_3.2.3.3.tgz.gpg)

To download this version, use GC Firmware Update Tool 3.2.8.0 or greater.

[Updating to this firmware does not require instrument requalification.](#)

Service Note Number: D0145780

#### Resolved Issues:

**Table 3 Resolved Issues (Agilent 8850 GC System 3.2.3.3)**

| Issue ID | Current Behavior                                                                                                                | Resolution |
|----------|---------------------------------------------------------------------------------------------------------------------------------|------------|
| 048135   | New design EPC module sensor error "Pressure Sensor Fault" with vacuum detectors connected when the pressure goes below -6.5psi | Resolved   |
| 048182   | Uploading New PID File Using GC FW Update Utility Fails                                                                         | Resolved   |

### 3.2.2.3 Firmware

Released on 7th January 2026

[https://download.chem.agilent.com/firmware/8850\\_FW\\_3.2.2.3.tgz.gpg](https://download.chem.agilent.com/firmware/8850_FW_3.2.2.3.tgz.gpg)

To download this version, use GC Firmware Update Tool 3.2.8.0 or greater.

[Updating to this firmware does not require instrument requalification.](#)

Service Note Number: D0142569

#### Resolved Issues:

**Table 4 Resolved Issues (Agilent 8850 GC System 3.2.2.3)**

| Issue ID | Current Behavior                                                           | Resolution |
|----------|----------------------------------------------------------------------------|------------|
| 047366   | Gather Logs Does not Work from BrowserUI                                   | Resolved   |
| 047371   | System Checkout is listed under Diagnostic Tests for the 3.2.0.130 release | Resolved   |

|        |                                                                                                        |          |
|--------|--------------------------------------------------------------------------------------------------------|----------|
| 047495 | Mainframe Test Broken, doesn't turn on relay or flaps                                                  | Resolved |
| 047502 | The Max Event number is inconsistent between Datasheet and DSA                                         | Resolved |
| 047505 | Need to confirm and align on power requirements for 8850 2XX for "UK type" labs, may want to update FW | Resolved |
| 047518 | System Fault in 3.2.0.133 when running sequence                                                        | Resolved |
| 047524 | "New Style" EPC pressure sensor fault under vacuum                                                     | Resolved |
| 047525 | Delete invalid HSS messages in Empower Run Log                                                         | Resolved |
| 047526 | GC System Fault - Code: 0x00000200                                                                     | Resolved |
| 047530 | H2 Sensor Calibration check not functioning correctly                                                  | Added    |
| 047619 | The Max Event number is inconsistent between Datasheet, LUI & BUI                                      | Added    |

### 3.2.0.130 Firmware

Released on 20<sup>th</sup> September 2025

[https://download.chem.agilent.com/firmware/8850\\_FW\\_3.2.0.130.tgz.gpg](https://download.chem.agilent.com/firmware/8850_FW_3.2.0.130.tgz.gpg)

To download this version, use GC Firmware Update Tool 3.2.8.0 or greater.

[Updating to this firmware does not require instrument requalification.](#)

Service Note Number: D0137759

#### Resolved Issues:

**Table 5 Resolved Issues (Agilent 8850 GC System 3.2.0.130)**

| Issue ID | Current Behavior                                                               | Resolution |
|----------|--------------------------------------------------------------------------------|------------|
| 046828   | Allow users to disable SMB Share                                               | Added      |
| 046910   | Touch Screen - Disable start button when instrument is in sleep mode           | Resolved   |
| 046918   | Browser Interface - One event type of "Detector Negative Polarity" not display | Resolved   |
| 046919   | Browser Interface - Untranslation happens on the list of signal source         | Resolved   |

|        |                                                                                                                                       |          |
|--------|---------------------------------------------------------------------------------------------------------------------------------------|----------|
| 046920 | Browser Interface - Untranslation happens on Browser Interface method report                                                          | Resolved |
| 046943 | Browser Interface - In Method\Events, the comp and column comp detector items are removed                                             | Resolved |
| 046957 | Untranslation: "Cancelled while waiting for ready" in logs                                                                            | Resolved |
| 046959 | The terminology of the message should be changed from setpoint/actual to expected/target in PreRun Flow Test Log Message              | Resolved |
| 047037 | Detector Evaluation Report not present - undefined error FW 3.0.3.21                                                                  | Resolved |
| 047089 | Request to show inlet pressure during column bakeout maintenance procedure                                                            | Added    |
| 047107 | Enhancement: Include instrument IP in Browser title bar                                                                               | Added    |
| 047251 | GC Firmware update tool revision 3.1.30.0 report error "Nothing to update ...." when upgrade 7697 firmware                            | Resolved |
| 047253 | 8697 pressure unit is always Psi even after changing GC unit                                                                          | Resolved |
| 047262 | False displayed message during Inlet Pressure zeroing                                                                                 | Resolved |
| 047283 | Improvement to the auto-zero process within the diagnostic test to properly account for the septum purge valve being not bubble tight | Resolved |

### 3.1.0.380 Firmware

Released on 20<sup>th</sup> June 2025

[https://download.chem.agilent.com/firmware/8850\\_FW\\_3.1.0.380.tgz.gpg](https://download.chem.agilent.com/firmware/8850_FW_3.1.0.380.tgz.gpg)

To download this version, use GC Firmware Update Tool 3.1.30.0 or greater.

[Updating to this firmware does not require instrument requalification.](#)

Service Note Number: D0133025

#### Resolved Issues:

**Table 6 Resolved Issues (Agilent 8850 GC System 3.1.0.380)**

| Issue ID | Current Behavior                                             | Resolution |
|----------|--------------------------------------------------------------|------------|
| 045867   | Browser Interface Help 8850 Miscellaneous topic forgets Oven | Resolved   |

|        |                                                                                                           |          |
|--------|-----------------------------------------------------------------------------------------------------------|----------|
| 045974 | Browser Interface: Un-translation happens on signal selection window                                      | Resolved |
| 046154 | 150 psi Split/Splitless inlet type is not identified in the Service Mode listing for the Pneumatic Module | Resolved |
| 045271 | Oven lid sensor is not responsive                                                                         | Resolved |
| 046345 | 8850_Untranslation happens on instrument maintenance and gas identification                               | Resolved |
| 046353 | Calibrating TCD when it's in a fault state                                                                | Resolved |
| 046505 | Text in the Touch Screen: History Logs is cut off                                                         | Resolved |
| 046593 | Invalid Syringe Volume should abort Run                                                                   | Resolved |
| 046639 | 8850 G3940A#001 Oven setting for 100V(JP)                                                                 | Resolved |

### 3.0.3.21 Firmware

Released on 6<sup>th</sup> February 2025

[https://download.chem.agilent.com/firmware/8860\\_FW\\_3.0.3.21.tgz.gpg](https://download.chem.agilent.com/firmware/8860_FW_3.0.3.21.tgz.gpg)

To download this version, use GC Firmware Update Tool 3.0.1.0 or greater.

Updating to this firmware does not require instrument requalification.

Service Note Number: D0127638

#### Resolved Issues:

**Table 7 Resolved Issues (Agilent 8850 GC System 3.0.3.21)**

| Issue ID | Current Behavior                                                                 | Resolution |
|----------|----------------------------------------------------------------------------------|------------|
| 046366   | GC system fault-Fault code, 0x00000200, Error code, 1000600a, sequence abort     | Resolved   |
| 046425   | PP Inlet - post run defect in flow control mode                                  | Resolved   |
| 046522   | GC Fault 1000600a//0x00000200 with FW 2.8                                        | Resolved   |
| 046544   | H2 Sensor appears to be zeroing after each run. Adding time to the GC Cycle time | Resolved   |

|        |                                                                                                                         |          |
|--------|-------------------------------------------------------------------------------------------------------------------------|----------|
| 046579 | Waters: They report a need for 500 mL flow rate                                                                         | Resolved |
| 046586 | Latest VICI actuator requires GC FW changes to pass EFT                                                                 | Resolved |
| 046595 | 8850 Channel Partner and AIB detector configuration support                                                             | Resolved |
| 046597 | Allow AIB compatibility with the 8850                                                                                   | Resolved |
| 046606 | Phone Browser UI Stuck on Loading Wheel Upon Startup                                                                    | Resolved |
| 046607 | H2 Sensor giving error messages reporting "Hydrogen Sensor: Not Functioning" and "Hydrogen Sensor: Invalid Environment" | Resolved |
| 046633 | Can't Download Diagnostic Data Collection CSV file                                                                      | Resolved |
| 046716 | Phone UI: Instrument firmware revision is empty                                                                         |          |
| 046717 | Phone UI: Attempting to change language does not bring up language selection menu                                       |          |

### 3.0.0.181 Firmware

Released on 10<sup>th</sup> September 2024

[http://download.chem.agilent.com/firmware/8850\\_FW\\_3.0.0.181.tgz.gpg](http://download.chem.agilent.com/firmware/8850_FW_3.0.0.181.tgz.gpg)

To download this version, use GC Firmware Update Tool 3.0.1.0 or greater.

Updating to this firmware does not require instrument requalification.

Service Note Number: D0120220

#### Resolved Issues:

**Table 8 Resolved Issues (Agilent 8850 GC System 3.0.0.181)**

| Issue ID | Current Behavior                                                                                   | Resolution |
|----------|----------------------------------------------------------------------------------------------------|------------|
| 037012   | Browser Interface should disable 'Enable Blank Evaluation' when detector signals are not selected. | Resolved   |
| 037227   | Browser UI: Oven Ramp setpoints not invalidated after changing Max Temp                            | Resolved   |
| 041981   | Detector Evaluation initial and final baseline should not fail because user selects Zero Signal    | Resolved   |

|        |                                                                                                                                          |          |
|--------|------------------------------------------------------------------------------------------------------------------------------------------|----------|
| 044226 | Add Gas Supply Pressure check back - in addition to method supply pressure check                                                         | Resolved |
| 044756 | Enhancement request to print the serial number of the main parts of the GC.                                                              | Added    |
| 045387 | 8697: System reported "flow not ready" and "Pressure not ready" after every injection when purge time starts                             | Resolved |
| 045496 | 8850 negative temperature ramps not working correctly                                                                                    | Resolved |
| 045541 | Maintenance Mode Options: remain cool or return to previous operating conditions                                                         | Resolved |
| 045561 | Detector Gas type labels do not agree between interfaces                                                                                 | Resolved |
| 045577 | Touchscreen and browser do not show any diagnostic progress for Browser sequence option 'Perform Pre-sequence Gas Supply Pressure check' | Resolved |
| 045633 | Browser UI column configuration doesn't correctly handle decreasing temperature limits                                                   | Resolved |
| 045719 | 8850: Touchscreen does not show Carrier Gas Flow Correction after download from browser.                                                 | Resolved |
| 045800 | 8850 Reset and Factory Reset offers 2 sets of Valves in Browser UI.                                                                      | Resolved |
| 045807 | 8850 Browser does not display or print Total Flow (for COC); inconsistent with touchscreen and GC Driver                                 | Resolved |
| 045845 | Incorrect 8697 logs after gripper failure happens                                                                                        | Resolved |
| 045918 | Real Time Plot for Duty Cycle is incorrectly scaled                                                                                      | Added    |
| 045925 | 8697: Appended line on a sequence doesn't start the run                                                                                  | Resolved |
| 045939 | 8850 COC Browser method print forgets Inlet Control Mode and Septum Purge Flow                                                           | Resolved |
| 045943 | Browser UI: GC Model issues when selecting certain events in the Method\Events\Run Table Events Setpoints dialog                         | Resolved |
| 045963 | Each advance from System Setup Date/Time screen changes the clock regardless of going back or abort & exit                               | Resolved |
| 045975 | 8850 Browser UI separated when loading reference chromatogram.                                                                           | Resolved |
| 046107 | 8850: Oven ramp upper limit in method parameters edit allows setting of 400C/min which exceeds the data sheet limit                      | Resolved |

|        |                                                                                                        |          |
|--------|--------------------------------------------------------------------------------------------------------|----------|
|        |                                                                                                        |          |
| 046215 | H2 Sensor purged into the oven during a run.                                                           | Resolved |
| 046250 | GC aborted sequence mid-run and recovered without notifying user                                       | Resolved |
| 046307 | GC Driver Early Maintenance Feedback counters: Linked field for User Defined counters is not available | Resolved |

**3.0.0.181 8850 GC Firmware**

*First manufacturing release September 2024.*

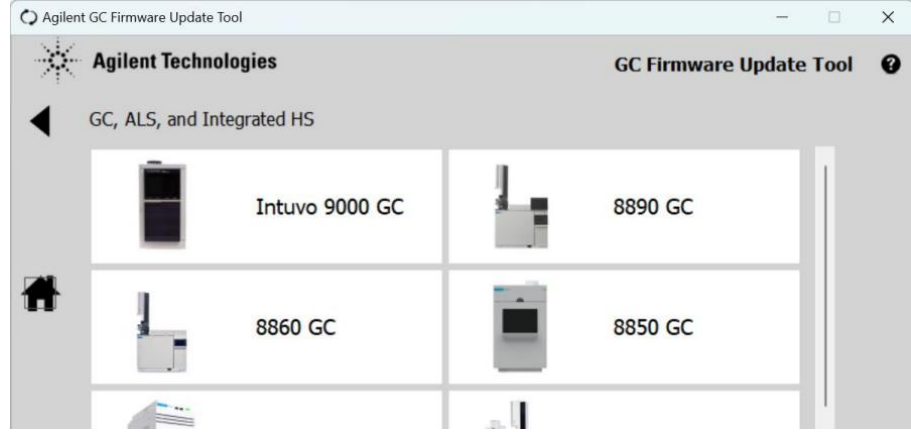
## 8

## Firmware Update Program

The GC firmware update program is supplied as part of the Firmware Update Package. Always use the latest version of the update tool. After the program has been loaded and executed the following screen appears.



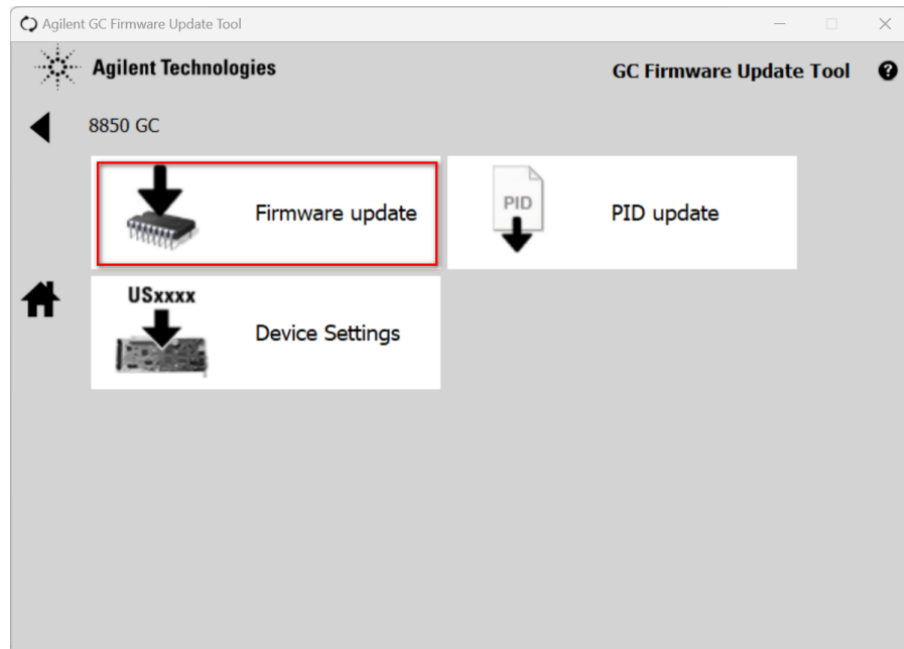
This enables the user to select which instrument type needs to have its firmware downloaded. In this case we select the GC and ALS.



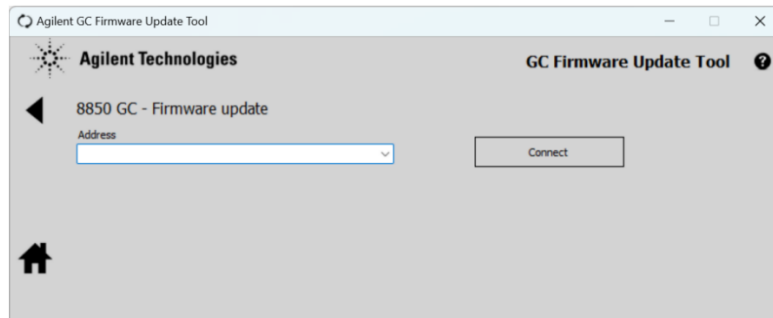
The **8850 GC** is chosen from the Product menu.

The next screen allows us to either modify firmware or PID values for the EPC modules.

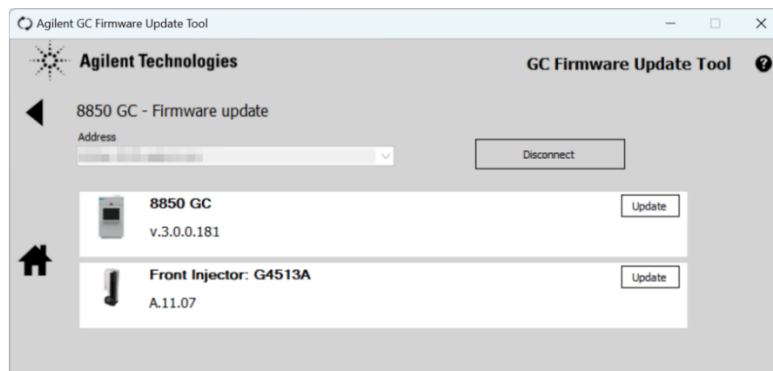
**Firmware update** is selected.



After choosing to update firmware, you must then enter the IP address of the 8850 GC that you wish to connect to and then press **Connect**.



Information such as current installed firmware will then be displayed.

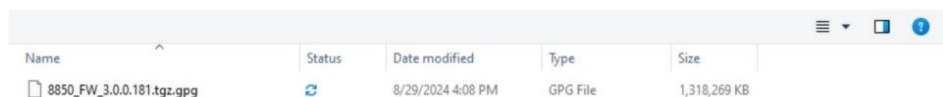


Press **Update**.

A Windows Explorer directory will then appear, and you must select the firmware revision that needs to be downloaded.

If the required firmware version is not displayed then you can "browse" your hard drive to find it.

The firmware file has the structure 8850-FW-X.X.X.X.tgz.gpg





After you have selected the correct file the firmware will automatically be downloaded.

After the download is finished the GC will reboot and the firmware update is complete.



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

© Agilent Technologies, Inc. 2026

Edition 07/2026

D0120223

