



Agilent 8850 GC Firmware

Release Notes – Revision 3.2.0.130

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. 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 3.2.0.130
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 3.2.0.130	9
7	Previous Versions	11
8	Firmware Update Program	15

1

Document History

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

Date	Version	Description
11 th September 2024	Rev A.01	Initial Release
11 th February 2025	Rev A.02	Added 3.0.3.21 Release Note
20 th June 2025	Rev A.03	Added 3.1.0.380 Release Note
20 th September 2025	Rev A.04	Added 3.2.0.130 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 (2.8, 2.7, 2.6, 2.5 or higher)

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

GC/MS MassHunter Acquisition (13.0)

OpenLab CDS EZChrom Edition (A.04.10 Update 3)

Supported Agilent GC Driver

Agilent GC Driver (4.4.92, 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 3.2.0.130?

- Enhancements
 - Support new 8850 features (MSD, et al)
- Bug fixes as detailed below.

6

Resolved Issues in revision 3.2.0.130

The current shipping firmware for the 8850 is 3.2.0.130

Released on 20th September 2025

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 2 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

7

Previous Versions

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

3.1.0.380 Firmware

Released on 20th June 2025

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 3 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 6th February 2025

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 4 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 10th September 2024

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 5 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.

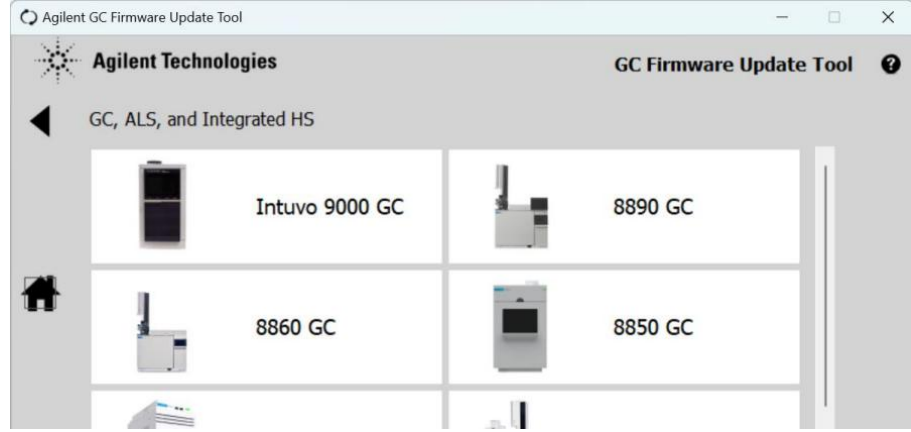
8

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.



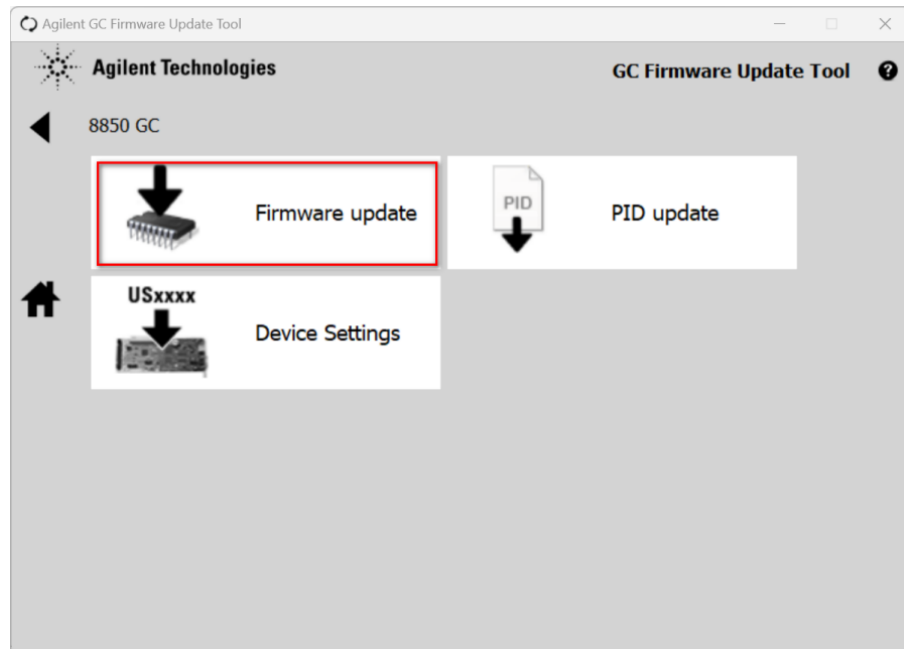
This enables the user to select which instrument type needs to have its firmware downloaded. In this case we select the GC and ALS.



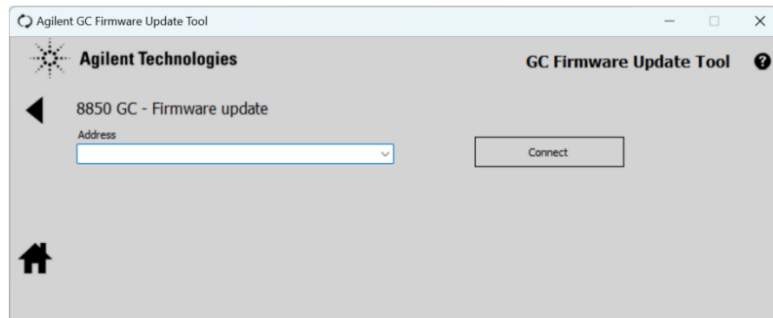
The **8850 GC** is chosen from the Product menu.

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

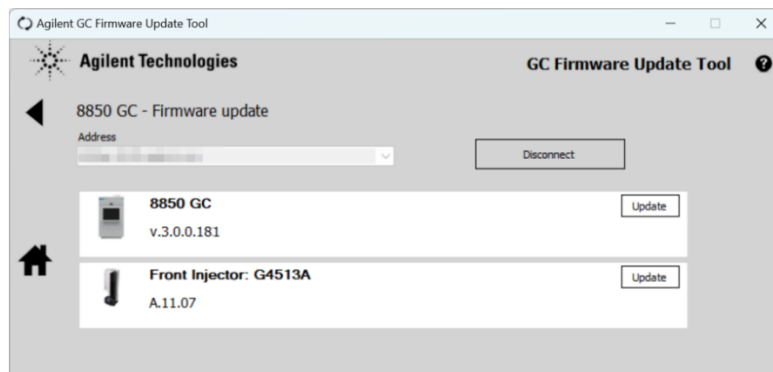
Firmware update is selected.



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

Name	Status	Date modified	Type	Size
8850_FW_3.0.0.181.tgz.gpg		8/29/2024 4:08 PM	GPG File	1,318,269 KB



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

© Agilent Technologies, Inc. 2025

Edition 09/2025

D0120223

