

Agilent 8860B Gas Chromatograph

Operation Manual



Notices

Manual Part Number

G2790-90022

Edition

First edition May 2026

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.
2850 Centerville Road
Wilmington, DE 19808-1610 USA
www.agilent.com

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 Legend

U.S. Government Restricted Rights. Software and technical data rights granted to the federal government include only those rights customarily provided to end user customers. Agilent provides this customary commercial license in Software and technical data pursuant to FAR 12.211 (Technical Data) and 12.212 (Computer Software) and, for the Department of Defense, DFARS 252.227-7015 (Technical Data - Commercial Items) and DFARS 227.7202-3 (Rights in Commercial Computer Software or Computer Software Documentation).

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.

Contents

1	Getting Started	8
	The 8860B Gas Chromatograph	9
	Before Operating Your GC	10
	Chromatography Using a GC	11
	Inlets	12
	Automatic injectors	12
	Automatic sampling valves	12
	GC Columns and Oven	13
	Detectors	14
	Touchscreen and Control Buttons	15
	System Operation	16
	Touchscreen and Control Buttons	16
	Browser Interface	16
	Data System	17
	Browser Interface	18
	Status Indicator	22
	GC Status	23
	Status Colors and details	23
	Alert tones	24
	Error conditions	24
	Clearing a shutdown condition	24
	Overview of Running a Sample	25
	Correcting Problems	26
2	Help and Information	27
	Where to Find Information	28
	Help from a Browser	29
	Context-Sensitive Help	32
3	Startup and Shutdown	34
	To Start Up the GC	35
	To Put the GC into Standby Mode	36
	To Shut Down the GC	37
4	Touchscreen and Controls Operation	38
	Navigation	39
	Run controls	40
	Status/control tray	40
	Entering Data	43
	Home View	44
	Status page	44
	Plot page	45
	Shortcuts	46
5	Methods	48
	What is a Method?	49
	What Is Saved in a Method?	50
	What Happens When You Load a Method?	51

Create a Method	52
Load a Method	53
Running Methods	54
Pre Run and Prep Run	54
Manually prepare for the run	54
To manually inject a sample with a syringe and start a run	54
To run a method to process a single ALS sample	55
To abort a method	55
Events	56
Using run time events	56
Programming run time events	57
The run table	57
Editing events in the run table	57
Deleting run time events	57
Inlets	58
Carrier gas flow rates	58
About gas saver	58
About Oven Temperature Programming	60
Oven ramp rates	60
Columns	62
Flow modes	62
Pressure modes	62
Nickel Catalyst Tube	63
About the nickel catalyst tube	63
Detectors	64
FID	64
FPD+	65
NPD	66
TCD	67
ECD	69
Valves	70
The 8860B valve box	70
Valve control	70
Valve types	71
Controlling a valve	72
GC Output Signals	74
Analog signals	74
Digital signals	76
6 Sequences	79
What is a Sequence?	80
Recoverable Errors	81
7 Diagnostics	82
About Diagnostics	83
Automated testing	83
The system health report	84
Self-guided diagnostics	84
Gather Log Files	85
Warnings and Errors	85
Using the Diagnostics View	86

	Performing Diagnostic Tests	87
8	Maintenance	88
	Automated Maintenance	89
	Early Maintenance Feedback (EMF)	90
	Counter types	90
	Default Thresholds	90
	Available Counters	91
	Viewing Maintenance Counters	93
	To Enable, Reset, or Change a Limit for an EMF Counter	94
	EMF Counters for Columns	95
	EMF Counters for Autosamplers	96
	Counters for 7693A and 7650 ALS with EMF-enabled firmware	96
	Counters for ALS with earlier firmware	96
	EMF Counters for MS Instruments	97
9	Logs	98
	Logs View	99
	Maintenance logs	99
	Run log	99
	System log	99
	Sequence log	99
	Run history log	99
10	Settings	100
	About Settings	101
	Service Mode	102
	About	103
	Calibration	104
	Maintaining EPC calibration—inlets, detectors, PCM, and AUX	104
	To zero a specific flow or pressure sensor	105
	System Settings	106
	Configuring the IP address for the GC	106
	To change the system locale	106
	To access stored run data	107
	To control Browser Interface access	108
	Tools (Browser Interface)	109
	Performing a column compensation run	109
11	Configuration	110
	About Configuration	111
	Configuration Changes	112
	Valve Configuration	113
	To configure valves	113
	Inlet Configuration	114
	Columns	115
	To configure a single column	115
	To configure a packed column	116
	Additional notes on column configuration	116
	Oven	118
	To configure the oven	118

	Detector Configuration	119
	Analog Output Settings	120
	MSD Configuration	121
	MSD configuration	121
	To enable or disable MS communications	121
	To use the GC when the MS is shut down	121
	Headspace Sampler (8697)	123
	Readiness	124
	Valve Box	125
	PCMs	126
	Aux EPCs	127
12	Resource Conservation	128
	Resource Conservation	129
	Sleep methods	129
	Wake and condition methods	130
	To Set the GC to Conserve Resources	132
13	Programming	133
	Clock Time Programming	134
	Using clock time events	134
	Adding events to the clock table	134
	Deleting clock time events	134
14	Flow and Pressure Modules	135
	About Flow and Pressure Control	136
	Maximum Operating Pressure	137
	Auxiliary Pressure Controllers	138
	Restrictors	139
	Selecting a frit	139
	Example: Using the PCM Channels	141
	Channel 1: Forward-pressure	141
	Channel 2: Two-way channel	141
	PIDs	142
15	Inlets	143
	Inlet Overview	144
	About the Split/Splitless Inlet	145
	Selecting the correct S/SL inlet liner	145
	About the Purged Packed Column Inlet	147
	About the Packed Column Inlet	148
	About the Cool On-Column Inlet	149
	Setup modes of the COC inlet	149
	Retention gaps	149
16	Columns and Oven	150
	Hydrogen Sensor	151
	Usage	151
	Audible Alarm	151
	Instrument logs	151

	Calibration	151
	Status information	152
	Operation with an Agilent data system	152
17	Chromatographic Checkout	154
	About Chromatographic Checkout	155
	To Prepare for Chromatographic Checkout	156
	To check FID performance with a purged packed, split splitless or cool-on column inlet	158
	To check FID performance with a packed column inlet (PCI)	161
	To check TCD performance with a purged packed, split splitless or cool-on column inlet	164
	To check TCD performance with a packed column inlet (PCI)	167
	To Check NPD Performance	170
	To Check ECD Performance	173
	To Check FPD+ Performance (Sample 5188-5953)	176
	Preparation	176
	Phosphorus performance	176
	Sulfur performance	179
	To Check FPD+ Performance (Sample 5188-5245, Japan)	180
	Preparation	180
	Phosphorus performance	180
	Sulfur performance	183
18	Chinese Metrology Testing	185
	FPD+ and ECD Unit Conversion Factors	186
	Conversion factors for the FPD+	186
	Conversion factor for the ECD	186
	Using the Conversion Factors	187
	References	188
19	Glossary	189

Getting Started

The 8860B Gas Chromatograph

Before Operating Your GC

Chromatography Using a GC

Inlets

Automatic injectors

Automatic sampling valves

GC Columns and Oven

Detectors

Touchscreen and Control Buttons

System Operation

Browser Interface

Status Indicator

GC Status

Overview of Running a Sample

Correcting Problems

This document provides an overview of the Agilent 8860B Gas Chromatograph (GC) along with detailed operating instructions.

Some graphics may not represent your GC. Not all features listed may be available at introduction of product.

The 8860B Gas Chromatograph



Figure 1. The 8860B GC

Before Operating Your GC

Before operating your GC, be sure to read the safety and regulatory information available on the Agilent website or the Browser Interface with a connected web browser. The most common safety hazards when working on the GC are:

- Burns caused by touching heated areas on or in the GC and MSD.
- Release of pressurized gas containing hazardous chemical compounds caused by opening inlets.
- Glass cuts or puncture wounds caused by sharp capillary column ends.
- Use of hydrogen as a GC carrier gas.

Magnetic Fields:

Magnets may affect the functioning of pacemakers, implanted heart defibrillators, or other active implanted medical devices (AIMD).



Figure 2. 8860B GC

The 8860 GC contains magnets on the inside of the oven door. Users that wear an AIMD are advised to avoid these magnets when the oven door is open.

Chromatography Using a GC

Chromatography is the separation of a mixture of compounds into individual components.

There are three major steps involved with separating and identifying components of a mixture using a GC. They are:

1. **Injecting** a sample into the GC. (This takes place at the inlet.)
2. **Separating** the sample into individual components. (This takes place inside the column in the oven.)
3. **Detecting** what compounds were in the sample. (This is done in the detector.)

During this process, status messages from the GC are displayed, and user changes to parameter settings can be made through the Browser Interface or the data system.

Inlets

Inlets are where samples are injected into the GC. The Agilent 8860B GC can have a maximum of two inlets, identified as Front Inlet and Back Inlet.

The following inlet types are available:

- Split/splitless inlet (SSL)
- Purged packed inlet (PP)
- Packed column inlet (PCI)
- Cool on column inlet (COC)

The type of inlet chosen is based on the type of analysis being done, the type of sample being analyzed, and the column being used.

Samples can be injected into the inlets by hand using a syringe, or an automatic sampling device (such as an Agilent Automatic Liquid Sampler or Agilent Headspace Sampler).

Automatic injectors

The Agilent 8860B GC can accommodate up to two autoinjectors, identified as Front Injector and Back Injector.

Automatic sampling valves

The optional sampling valves are simple mechanical devices that introduce a sample of fixed size into the carrier gas stream. Valves are most frequently used to sample gases in constantly flowing streams.

The Agilent 8860B GC can accommodate up to four total valves. Up to two non-heated liquid sampling valves, or up to two gas sampling valves may be used in any combination.

The valves are located inside the sampling valve box.

GC Columns and Oven

GC columns are located inside a temperature-controlled oven. Generally, one end of the column is attached to the inlet, while the other end is attached to the detector.

Columns vary in length, diameter, and internal coating. Each column is designed for use with different compounds.

The purpose of the column and the oven is to separate the injected sample into individual compounds as it travels through the column. To aid this process, the GC oven can be programmed to speed the sample flow through the column.

The Agilent 8860B GC can accommodate up to three columns, identified as Column #1, Column #2, and Column #3

To open the GC 8860 column oven, press the oven door button.

Detectors

Detectors identify the presence of compounds as they exit the column.

As each compound enters the detector, an electrical signal proportional to the amount of compound detected is generated. This signal is generally sent to a data analysis system—such as Agilent OpenLab CDS ChemStation edition—where it shows up as a peak on a chromatogram.

The 8860 GC can accommodate up to three detectors (two mounted on top of the GC, and one mounted in a side carrier, or in the aux position on the top), identified as Front Detector, Back Detector, and Aux 2 Detector.

A complete selection of detectors (FID, TCD, NPD, FPD+, ECD, and MSD) are available. The type of detector chosen is based on the type of analysis required.



Figure 3. Detector locations

NOTE

A side mounted detector may only be a TCD, ECD, or FID. Other detector types are not supported in this location.

An Aux Detector mounted on top may only be a FPD+ or a TCD.

Touchscreen and Control Buttons

The touchscreen shows GC status, activity information, shortcuts and settings, while the control buttons allow you to start, stop, and prepare the GC to run a sample.

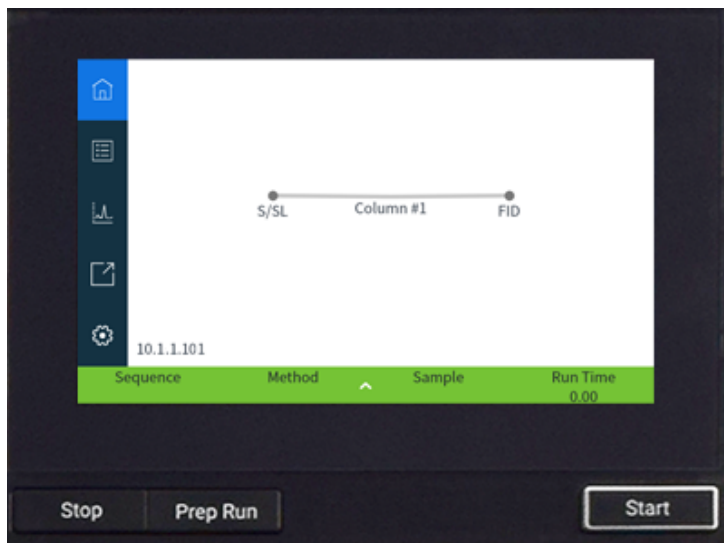


Figure 4. View status, status listing, live plots, shortcuts, settings, and control runs.

See [Touchscreen and Controls Operation](#) for a detailed description of the touchscreen and control button functions and capabilities.

System Operation

The GC can be controlled using the touchscreen and control buttons, the Browser Interface, and an Agilent data system.

Touchscreen and Control Buttons

The touchscreen provides controls for:

- Setting the GC IP address
- Selecting the touchscreen language
- Viewing status data and real-time plots
- Monitor GC and run status
- Monitor system health
- Make system settings, for example system locale, IP address setup, and so forth

Control button	Function
Prep Run	Typically required before manual injections to exit any gas saver mode and to prepare the inlet flows for injection.
Start	Starts a run.
Stop	Stops the current run.

Some tasks, such as setting the IP address, can only be performed using the touchscreen.

Browser Interface

The Browser Interface provides many of the same functions as the touchscreen and control buttons. The Browser Interface provides for instrument setup and control, as well as the ability to run the instrument stand-alone (without a connected data system). The Browser Interface can be viewed using any typical web browsing device, such as a computer or tablet, provided the device is connected to the same gateway as the GC. Use the Browser Interface to:

- Create methods
- Run samples and sequences of samples
- Run diagnostic tests
- Check status and performance of the GC
- Perform maintenance procedures
- Check maintenance details and reset counters
- Change language settings
- Access on-line help

To connect to the Browser Interface:

1. On a tablet or computer connected to the same gateway as the GC, open a browser window.
2. Navigate to **http://GC Name or IP Address**. For example, if the GC IP address is 10.1.1.100, navigate to **http://10.1.1.100**.

3. If prompted, enter the access code. (The access code is available at the GC touchscreen.)

For more information regarding the Browser Interface, see **Browser Interface**.

If multiple users connect to a GC using the Browser Interface, only the first connection has normal control.

When a data system is controlling the GC, setpoint changes and other features may be disabled in the Browser Interface.

Data System

The data system adapter provided with Agilent data systems, such as OpenLab CDS, provides full control of the GC for creating methods, running samples, and so forth. Use the data system to:

- Create methods
- Run samples and sequences of samples
- Monitor GC and run status
- Monitor system health
- Track consumables usage and usable life remaining
- Make some system settings
- View system logs
- Set GC clock table and resource conservation events

In addition, the data system adapter provides access to the complete help and user information. From anywhere within the method editor for the GC, select **Help and Information Browser** from the navigation tree.

Also note:

- The data system adapter does not provide direct access to all of the GC's diagnostic and maintenance features.
- The data system cannot directly use any methods or sequences created in the Browser Interface.
- When connected to the GC, the data system adapter can be set to restrict anyone from making certain changes at the touchscreen or with a Browser Interface.

Browser Interface

You can control and monitor the GC using a web browser that is on the same gateway as the GC. An internet connection is not required. This Browser Interface can be accessed by using computer browser clients and mobile device client browsers, such as tablets. The Browser Interface provides complete control of the GC. Use it to perform tasks such as:

- Configuring GC gas types and flows
- Performing automated maintenance procedures
- Running diagnostic tests
- Creating, editing, and loading methods
- Running samples
- Creating, editing, and loading sequences of samples
- Monitoring GC performance (viewing logs, current statuses, and plots)

While connected to your GC with a data system, you cannot edit methods, edit sequences, and start or stop runs using the Browser Interface. See [What is a Method?](#) for more details.

The screenshot displays the Browser Interface with the **Method** tab selected. The interface includes a top navigation bar with tabs: **Method**, **Sequences**, **Diagnostics**, **Maintenance**, **Logs**, **Settings**, and **Help**. Below the navigation bar is a toolbar with icons for navigation and actions. The main content area shows a table of components with columns: **Name**, **Type**, **On**, **Load Time (min)**, and **Inject Time (min)**. The table lists 10 items, all marked as **Not Installed**. On the left, a sidebar shows a tree view of components: **ALS/Tray**, **Front Injector**, **Other**, **Valves**, **Inlets** (with sub-items **Front Inlet - SS** and **Back Inlet - MMI**), **Columns**, **Oven**, **Detectors** (with sub-items **Front Detector - FID**, **Back Detector - TCD**, and **Aux Detector 2 - TCD**), **Analog Out**, **Analog Out 1**, and **Analog Out 2**. At the bottom, a status bar shows **STATUS: READY** and a table with columns **Sequence**, **Method**, **Sample Name**, and **Est. Remaining**. The **Method** column shows **AddedCol3** and the **Est. Remaining** column shows **1**. On the right side of the interface, there are three callout boxes: **Edit GC Methods, Sequences, Settings, and more** (pointing to the top navigation bar), **Access your Help & Information Suite** (pointing to the Help tab), and **View context sensitive help about your current screen** (pointing to the sidebar).

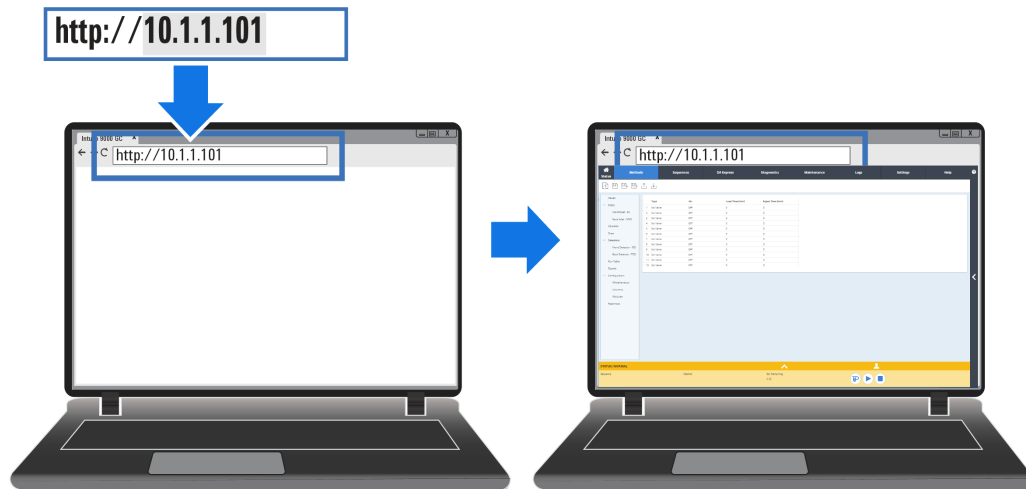
When using the Browser Interface, the GC stores run data internally. Access can be granted to allow a connected computer on the same gateway to access this data and copy it for analysis. Note that run data can only be deleted by using the Browser Interface directly, or by the GC according to either date or disk space settings.

To connect to the GC using a browser:

1. If you do not know the GC's IP address or host name, use the touchscreen to find it.

NETWORK	ACCESS	STORAGE	POWER	MISC	ABOUT
DHCP	☐ Enabled				
IP Address	10.1.1.101				
Netmask	255.255.255.0				
Gateway	10.1.1.100				
Mac Address	00:00:00:00:00:00				
					Apply
Sequence		Method	Sample	Run Time 0.00	

- Open a web browser. Supported browsers include Chrome, Safari (on a tablet), and Edge. Make sure the browser version is up to date.
- Enter **http://xxx.xx.xx.xxx**, where **xxx.xx.xx.xxx** is the GC's IP address. (If using a host name, enter it instead.) In this example, the GC IP address is 10.1.1.101. Accessing the Browser Interface only requires that the tablet or computer be connected to the same gateway as the GC; no internet connection is required.



For more information on how to utilize the Browser Interface, click the Help tab to access the Help & Information suite, or click the < on the right side of the screen to access the context sensitive help. See [Help from a Browser](#) and [Context-Sensitive Help](#) for more information.

Getting Started

Browser Interface

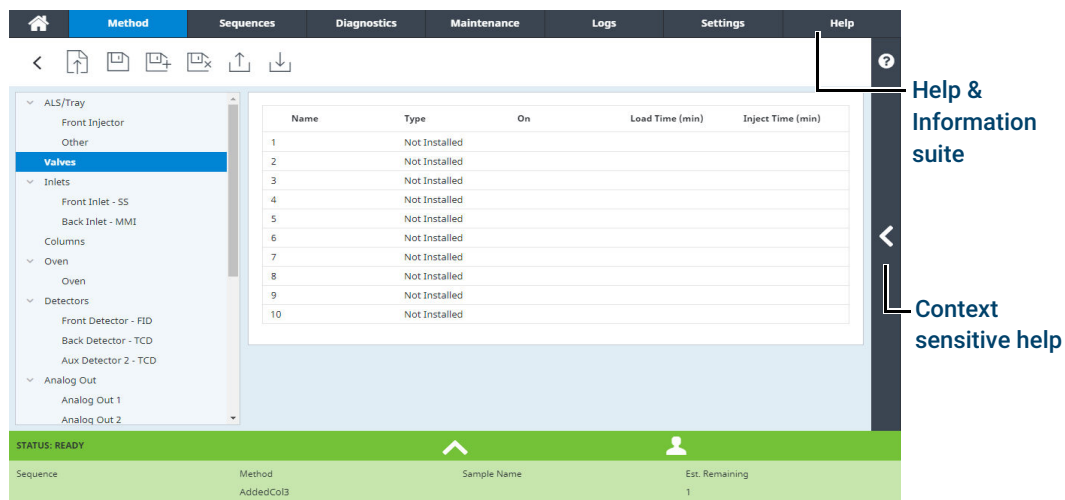


Figure 5. Accessing help from the Browser Interface

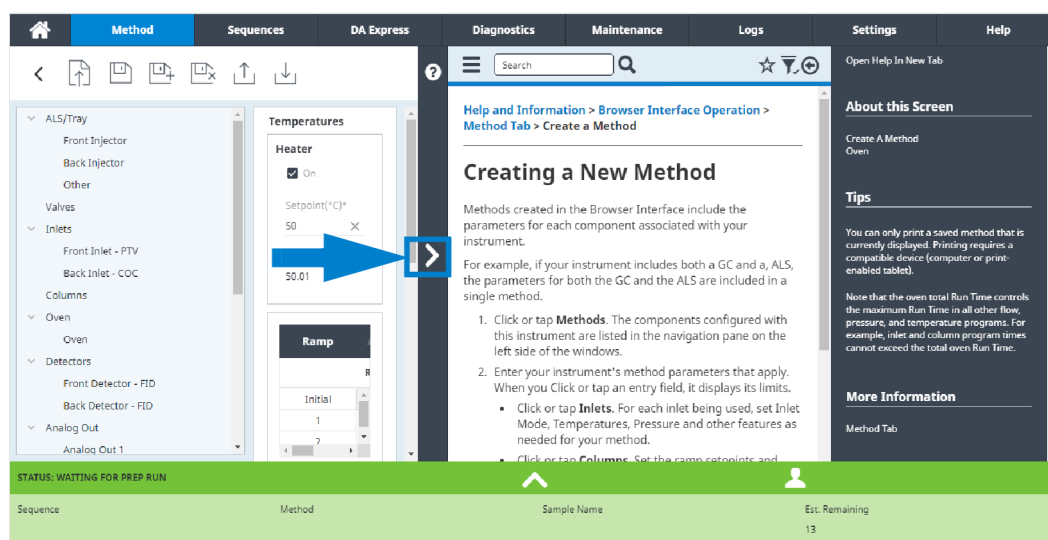


Figure 6. Context sensitive help

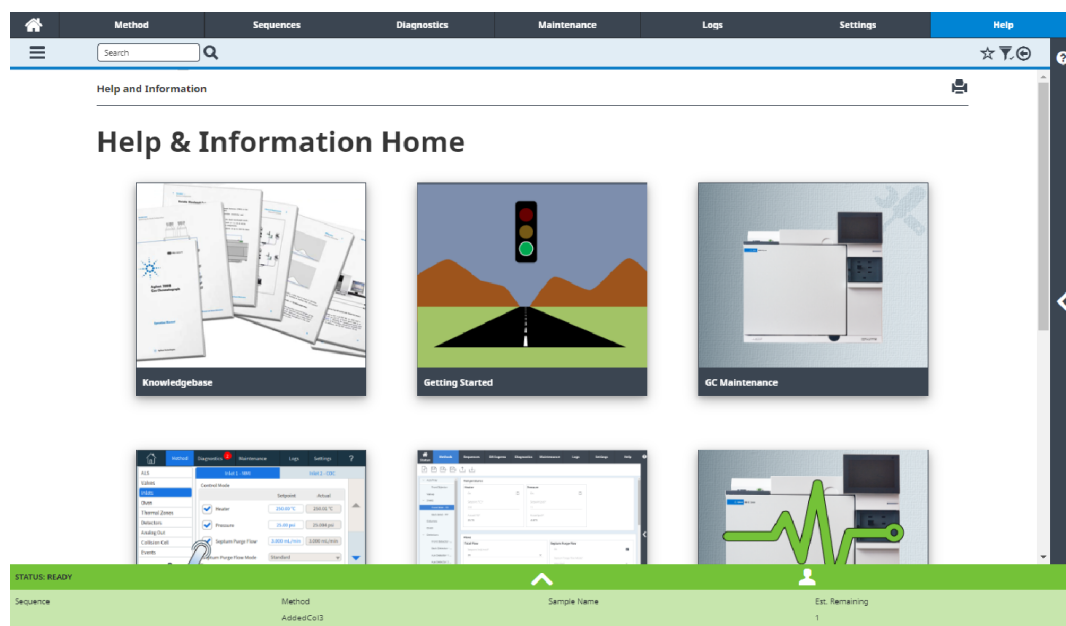


Figure 7. Complete Help and Information content

Status Indicator

The GC includes a status indicator on the front panel to allow you to quickly determine the status and readiness of the GC. The status indicator changes color depending on the current state of the GC.

-  (Green): Indicates that GC is ready for operation.
-  (Yellow): Indicates that the GC is not ready for operation. Power is on and available, but not all parameters have reached operating setpoints, or a maintenance or diagnostic task is in progress. A warning or other message may exist. Check the GC touchscreen for additional information.
-  (Red): Indicates a fault or other serious condition. A fault or other message may exist. Check the GC touchscreen for additional information. The GC cannot be used until the fault condition is resolved.



Figure 8. Status Indicator

GC Status

When the GC is ready to begin a run, the Browser Interface shows **STATUS: READY FOR INJECTION**. Alternately, when a component of the GC is not ready to begin a run, the Browser Interface shows **STATUS: NOT READY** and the status indicator on the GC front panel is yellow. Clicking the **Diagnostics** tab displays indications of why the GC is not ready.

Status Colors and details

Colors provide quick GC status:

Ready (all settings at setpoint)

STATUS: NORMAL

Run in progress

STATUS: RUN

Not ready

STATUS: NOT READY

Maintenance mode (an automated maintenance task is running)

STATUS: MAINTENANCE MODE

Preparing sample

STATUS: PREPARING SAMPLE

GC is in sleep mode

STATUS: SLEEP

GC is in an error state

STATUS: ERROR

Status text displayed in the status tray can include:

Status text	Comment
CONDITIONING RUN	GC is performing a conditioning run after waking from Sleep.
COOLDOWN	The GC is performing a cooldown between runs.
DIAGNOSTIC MODE	The GC is currently performing a diagnostic test.
ERROR	An error is preventing normal operation. Check the Diagnostics and Logs tabs for more information.
H2 SHUTDOWN	The GC is shut down: heaters are turned off, the oven flaps are open, and gas flows are off.
MAINTENANCE MODE	The GC is running a user maintenance procedure.
MSD VENTING	The GC is set to provide flows appropriate for MSD venting.
MSD VENTED	The MSD is vented.
NORMAL	No conditions prevent operation. The GC may or may not be ready to begin a run.
POST RUN	The GC is executing a method post run program.
PREPARING SAMPLE	During a run, a connected ALS is performing sample preparation.
PREP RUN	The GC is in the Prep Run state, preparing for an injection.
RUN	The GC is performing a run.
SLEEP	The GC is in sleep mode. To wake the GC, apply a setpoint change, download a method, or from an Agilent data system select "Wake Up."

Status text	Comment
SLEEP DEFERRED	Due to ongoing activity, the GC did not enter sleep mode when scheduled. The GC will enter sleep mode after it is has been idle for at least 15 minutes.
SHUTDOWN	A GC component is preventing the GC from operating normally. Check for a diagnostic condition or message at the Diagnostics tab.

Alert tones

The GC provides information via beeps.

A series of warning beeps sounds before a shutdown occurs. The GC starts with one beep. The longer the problem persists, the more the GC beeps. After a short time the component with the problem shuts down, the GC emits one beep, and a brief message is displayed. For example, a series of beeps sounds if the inlet gas flow cannot reach setpoint. The message **Inlet flow shutdown** is briefly displayed. The flow shuts down after 2 minutes.

WARNING

Before resuming GC operations, investigate and resolve the cause of the hydrogen shutdown.

One beep sounds when a problem exists, but the problem will not prevent the GC from executing the run. The GC will emit one beep and display a message. The GC can start the run and the warning will disappear when a run starts.

Fault messages indicate hardware problems that require user intervention. Depending on the type of error, the GC emits no beep or a single beep.

Error conditions

If a problem occurs, the GC status bar changes to Not Ready, the GC status indicator turns yellow, a number appears next to the Diagnostics tab, and the status tray and Diagnostics tab list the condition that is causing the Not Ready condition. Select **Diagnostics** to view the issue and resolve it.

Clearing a shutdown condition

When a component is shut down, the GC becomes Not Ready, the status indicator and status bar turn yellow, and the Diagnostics tab displays a number in a red circle to indicate that there are a number of warnings or other issues that require attention. **Diagnostics > Warnings and Errors** will display any current conditions that require attention.

To clear a shutdown condition, select **Diagnostics > Warnings and Errors**, then select a condition to view its possible causes and resolutions. Choose a **Setpoint Action**. Select **Keep Shutdown Zone On** to clear the alert and turn on all zones. Select **Turn Shutdown Zone Off** to clear the alert and turn on all zones except the shutdown zone.

Note that Hydrogen shutdowns initiated from the hydrogen sensor can only be cleared by power cycling (rebooting) the GC. After the GC restarts, the affected hydrogen flows and heaters will remain turned off. (Note that the GC can only shut down hydrogen flows for devices under its control.) Determine the cause of the shutdown and resolve it. See the Troubleshooting manual.

Overview of Running a Sample

Operating the GC involves the following tasks:

- Setting up the GC hardware for an analytical method.
- Starting up the GC. See **To Start Up the GC**.
- Preparing any attached sampler. Install the method-defined syringe; configure solvent and waste bottle usage and syringe size; and prepare and load solvent, waste, and sample vials, as applicable. Refer to the documentation provided with your ALS or Headspace Sampler (HS) for details on its installation, operation, and maintenance.
- Loading the analytical method or sequence into the GC control system.
 - See the Agilent data system documentation.
 - For standalone GC operation see **Load a Method**.
- Running the method or sequence.
 - See the Agilent data system documentation.
 - For standalone GC operation, see **To manually inject a sample with a syringe and start a run** and **To run a method to process a single ALS sample**.
- Monitoring sample runs from the GC touchscreen, Browser Interface, or Agilent data system. See **Home View**, or the Agilent data system documentation.
- Shutting down the GC. See **To Put the GC into Standby Mode** or **To Shut Down the GC**.

Correcting Problems

If the GC stops operation because of a fault, check the touchscreen or Browser Interface for any messages. The GC includes diagnostics functions to help you determine the cause of a fault.

1. Use the touchscreen, Browser Interface, or data system to view the alert. (See **Home View** and **Diagnostics** for details.)
2. Press the "Stop" button located beneath the touchscreen, click the stop button on the Browser Interface, or turn off the offending component in the data system.
3. Diagnose the problem using the built in diagnostics tools on the GC. See **Diagnostics**.
4. Resolve the problem by, for example, changing gas cylinders or fixing the leak.

Once the problem is fixed, you may need to power cycle the instrument. Most shutdown errors can be cleared from the LUI, but some may require a power cycle.



2 Help and Information

Where to Find Information

Help from a Browser

Context-Sensitive Help

Where to Find Information

Agilent provides an extensive amount of documentation for installing, operating, and maintaining the instrument directly from the 8860B GC. In addition, there are multiple ways to access the comprehensive Help and Information suite either through a browser, or from your Agilent data system.

- **“Help from a Browser”.**

The complete set of user information is available directly from the GC by using a connected web browser.

- **“Context-Sensitive Help”.**

In addition to the complete set of user documentation, context-sensitive information is also available from the Browser Interface.

- **The Agilent web site** (www.agilent.com).

The Agilent web site provides information about the complete Agilent line of products, including samplers and mass selective detectors. The available information includes manuals, applications, technical briefs, and more.

When unpacking the instrument, make sure you take a look at the provided Agilent 8860B GC *Quick Start* Poster to help you quickly get familiar with your GC, as well as setup and configure your instrument. This poster can also be found in your browser help, in the familiarization section.

Agilent
8860 Gas Chromatograph
Quick Start

1 8860 GC System Ship Kit (83490-6088)

Agilent GC and GC/MS User Manuals and Tools (44000-64000)

2 Remove Protective Packaging

Remove a Embalagem de Proteção

Удалите защитную упаковку

移除保护包装

3

LAN

4

5 Select Settings tab to enter Network Information

Selecione a guia Parâmetros para inserir informações da rede

Выберите вкладки параметров и введите информацию о сети

选择“设置”选项卡输入网络信息

6 Access the Browser Interface to configure the GC and get detailed installation instructions by connecting to the GC via your web browser (no internet required).

The GC and PC must be configured on the same gateway (for example, isolated LAN).

Acesso a interface do navegador para configurar o GC e obter instruções detalhadas de instalação conectando-se ao GC através do seu navegador da Web (não é necessário usar a Internet).

O GC e o PC devem ser configurados no mesmo gateway (por exemplo, uma LAN isolada).

Доступ к интерфейсу веб-браузера для настройки прибора и получения подробных инструкций по установке, подключению к ПК через веб-браузер (подключение к интернету не требуется).

Пк и ПК должны быть настроены на одном и том же адресе веб-браузера (например, изолированная LAN).

ウェブブラウザを通じてGCのブラウザインターフェイスに接続し、詳細なインストール手順を閲覧してください（インターネットに接続する必要はありません）。

GCとPCは同じゲートウェイ（例えばLAN）に接続してください。

访问浏览器界面配置 GC 并浏览详细安装说明（不需要互联网连接）。连接至 GC 获取详细安装说明。GC 和 PC 必须配置在相同网关（例如，独立网络环境中）。

Browser Interface

Provides instrument control

<http://10.1.1.101/>

Installation instructions

GC and GC/MS User Manuals & Tools

<http://10.1.1.101/instel/>

Where to find additional information

Onde encontrar informação adicional

Где найти дополнительную информацию

添加情報を検索する方法

如何获得相关信息

Help and Information

<http://10.1.1.101/info>

Accessing Help and Information from Agilent Data Systems

Acessando Ajuda & Informação através do Sistema de Dados da Agilent

Доступ к справке и информации по системе данных Agilent

通过 Agilent 数据系统访问 帮助和信息

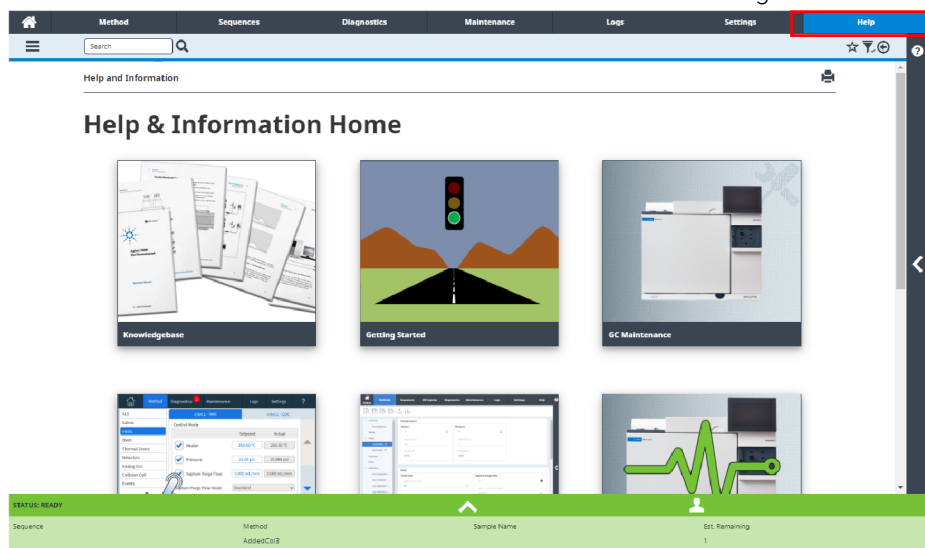
© Agilent Technologies, Inc., 2015
Printed in USA or China

Help from a Browser

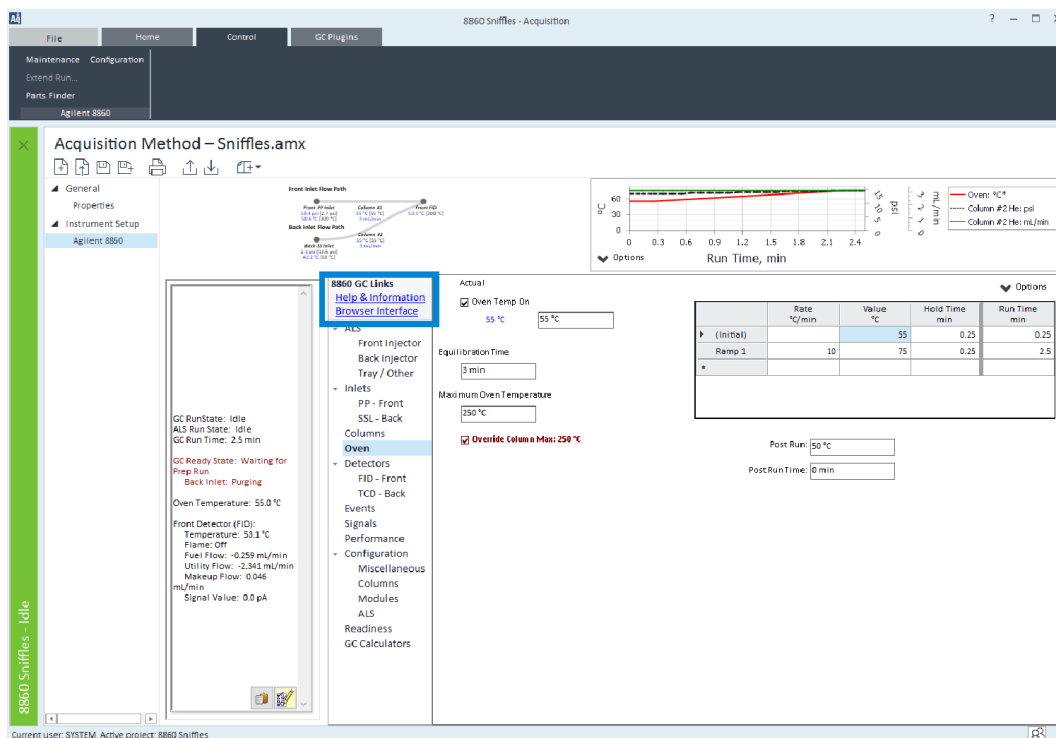
You can access an extensive amount of on-board documentation that is designed to assist with topics such as getting started, familiarization, installation, operation, maintenance, troubleshooting, and other useful information. **Access to the internet is not required to use this enhanced help package.** Accessing this help only requires that your PC or tablet be connected to the same gateway as your GC.

This enhanced version of **Help & Information** can be easily accessed through:

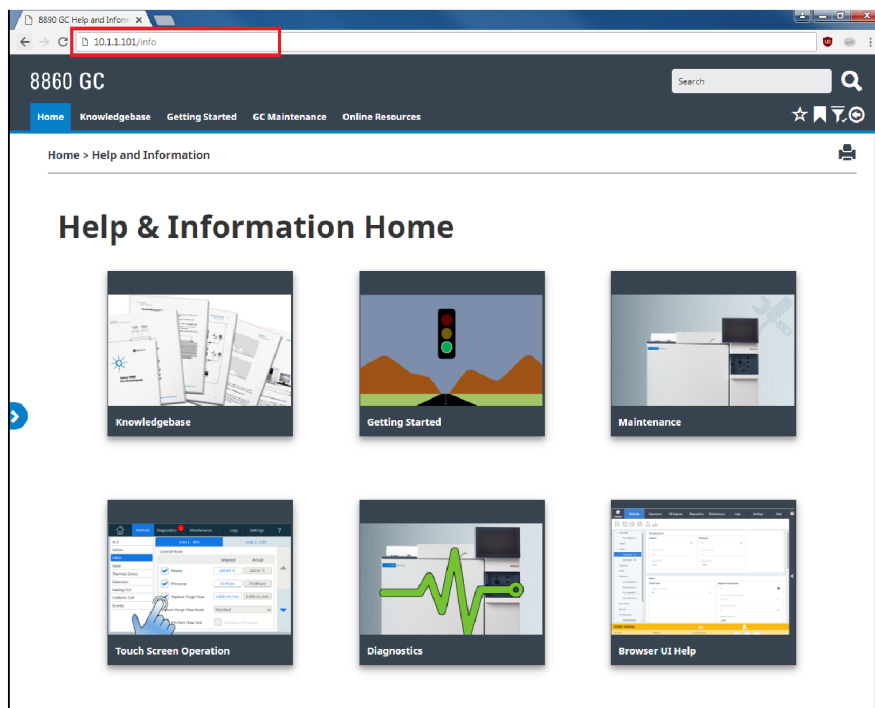
- The Browser Interface. Access the Help & Information suite by clicking the **Help** tab in the Browser Interface. See [Browser Interface](#) for instructions on connecting to the Browser Interface.



- Your Agilent data system. Access the Help & Information suite by clicking the **Help & Information Browser Interface** link at the top of the tree view.



- A web browser on any device that is on the same gateway as the GC. Access the Help & Information suite by typing **http://xxx.xx.xx.xxx/info** into your web browser of choice, where **xxx.xx.xx.xxx** is the IP address or host name of your GC.



Accessing the Help & Information suite provides you with the following information:

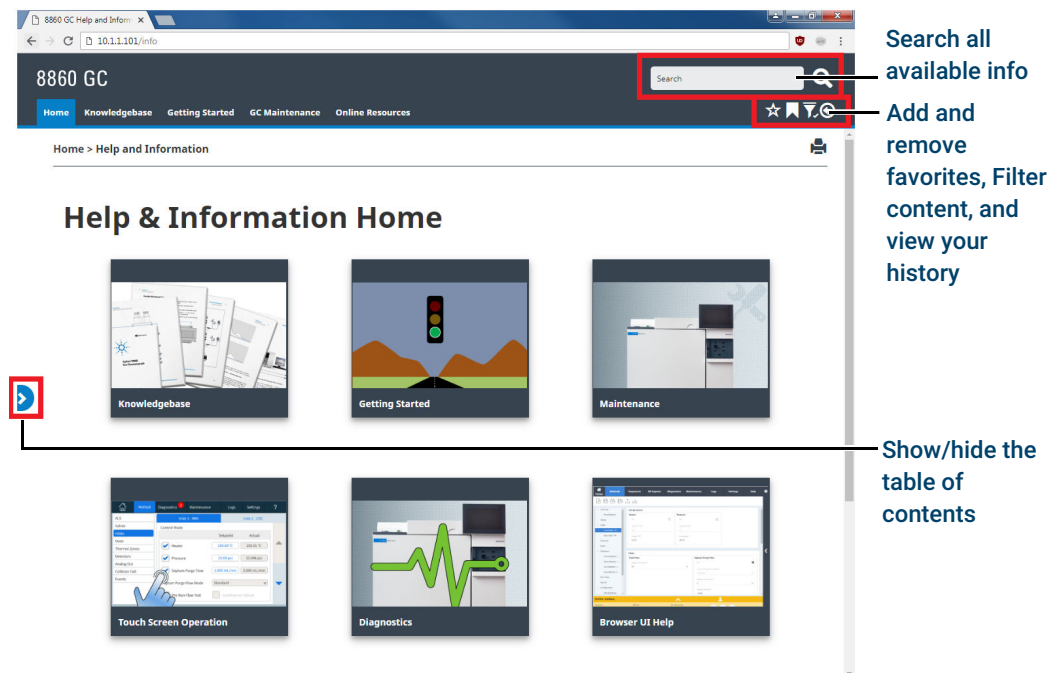
- Knowledgebase:** manuals, installation info, safety info, and site prep.

Help and Information

Help from a Browser

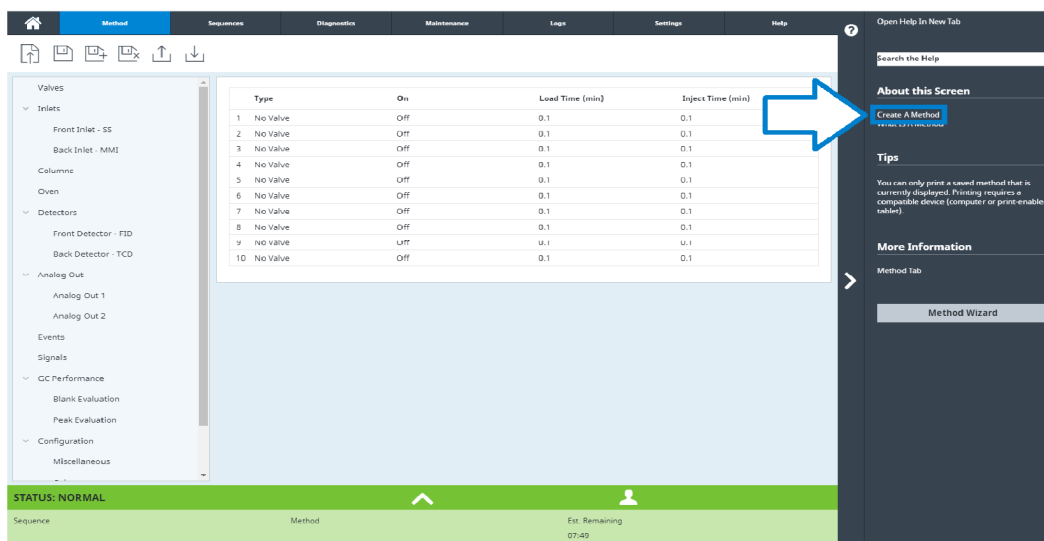
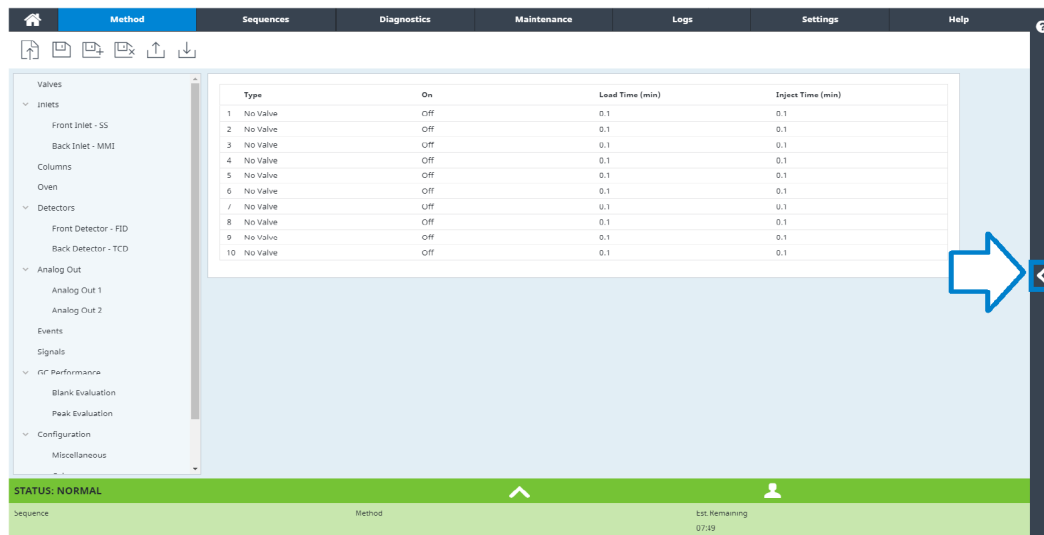
- **Getting Started:** BrowserUI tutorial, Quick Start Poster, eIntroduction, a GC Feature Tour, the System Setup Wizard, and instructional videos.
- **Maintenance:** information for performing maintenance on your configuration (inlets, detectors, etc.).
- **System Operation:** information on operating your GC, such as its settings, or EMF counters.
- **Diagnostics:** information on the diagnostic tests, self-guided diagnostics, and tasks available on your GC.
- **Browser Interface Help:** help and instructions for using the Browser Interface.
- **Online Resources:** links to Agilent University, Agilent YouTube, Agilent Community, services, and more.

By default, the help is filtered to only show information pertaining to your GC's configuration. If information does not appear to be available, be sure to enable all help content. To check your active filters, click the  icon in the top right corner of the Help & Information screen. These filters allow you to show or hide information pertaining to the various components of your GC, such as information about specific detectors or inlets.

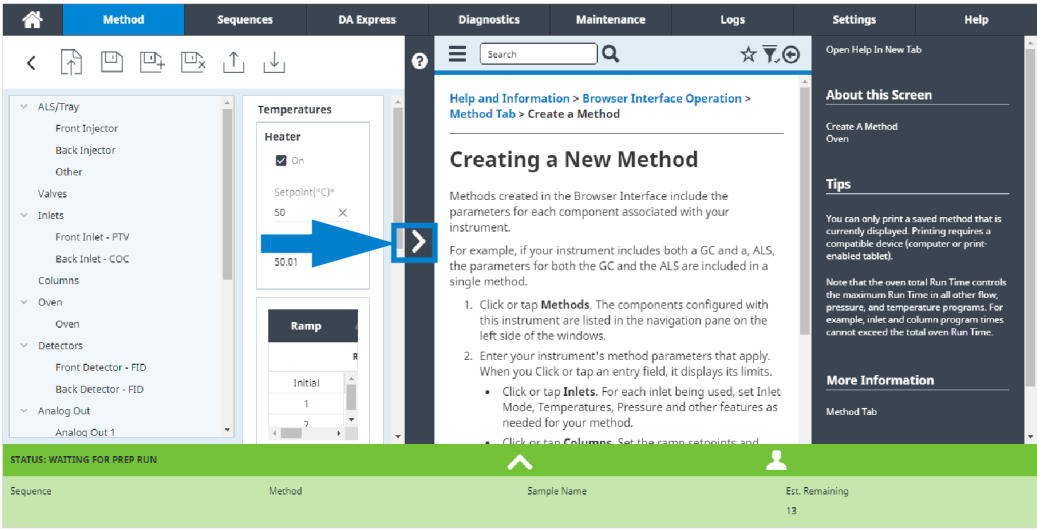


Context-Sensitive Help

Available from each of the Browser Interface pages is the ability to access context-sensitive information or help. Select the **<** from the right side of the screen to access information and tips relevant to the page you are viewing. **About this Screen** provides several links to relevant help topics. When clicked, these links open their corresponding topics within the help tray. Under **Tips**, you will see snippets of useful information about the current page. Additionally, some pages will provide links to additional documentation, such as the link to the Method Wizard in the example below.



To minimize the topic you are viewing, click the **>** between the context sensitive help and the page you are viewing. Click the **>** again to minimize the context sensitive help tray.





3 Startup and Shutdown

To Start Up the GC

To Put the GC into Standby Mode

To Shut Down the GC

To Start Up the GC

Successful operation begins with a properly installed and maintained GC. The utility requirements for gases, power supply, venting of hazardous chemicals, and required operational clearances around the GC are detailed in the Agilent *8860B Gas Chromatograph Site Preparation Guide*.

1. Check gas source pressures. For required pressures, see the Agilent 8860B Gas Chromatograph Site Preparation Guide.
2. Turn on the carrier and detector gases at their sources and open the local shutoff valves.
3. Turn on the GC power. Wait for **Power on successful** to display on the touchscreen.
4. Install the column.
5. Check that the column fittings are leak free.
6. Set a method for the GC to use.
If using a data system, download the method to the GC.
If using the Browser Interface, load a method.
7. Wait for the detector(s) to stabilize before acquiring data. The time required for the detector to reach a stable condition depends on whether the detector was turned off or its temperature was reduced while the detector remained powered.

Table 1. Detector stabilization times

Detector type	Stabilization time starting from a reduced temperature (hours)	Stabilization time starting from detector off (hours)
FID	2	4
TCD	2	4
ECD	4	18 to 24
FPD+	2	12
NPD	4	18 to 24

To Put the GC into Standby Mode

1. Wait for the current run to finish.
2. If the active method has been modified, save the changes.

WARNING

Never leave flammable gas flows on if the GC will be unmonitored. If a leak develops, the gas could create a fire or explosion hazard.

3. Reduce oven temperature to 50 °C or lower. Reduce detector and inlet temperatures to between 150 and 200 °C. If desired, the detector can be turned off. See [To Start Up the GC](#) to determine if it is advantageous to shut down the detector for a short time period. The time required to return the detector to a stable condition is a factor.

To Shut Down the GC

1. Put the GC into general maintenance mode by selecting **Maintenance > Instrument > Perform Maintenance > Start Maintenance**, and wait for the GC to become ready.
2. Turn off the main power switch.

WARNING

Be careful! The oven, inlet, and/or detector may be hot enough to cause burns. If they are hot, wear heat-resistant gloves to protect your hands.



4 Touchscreen and Controls Operation

Navigation

Run controls

Status/control tray

Home View

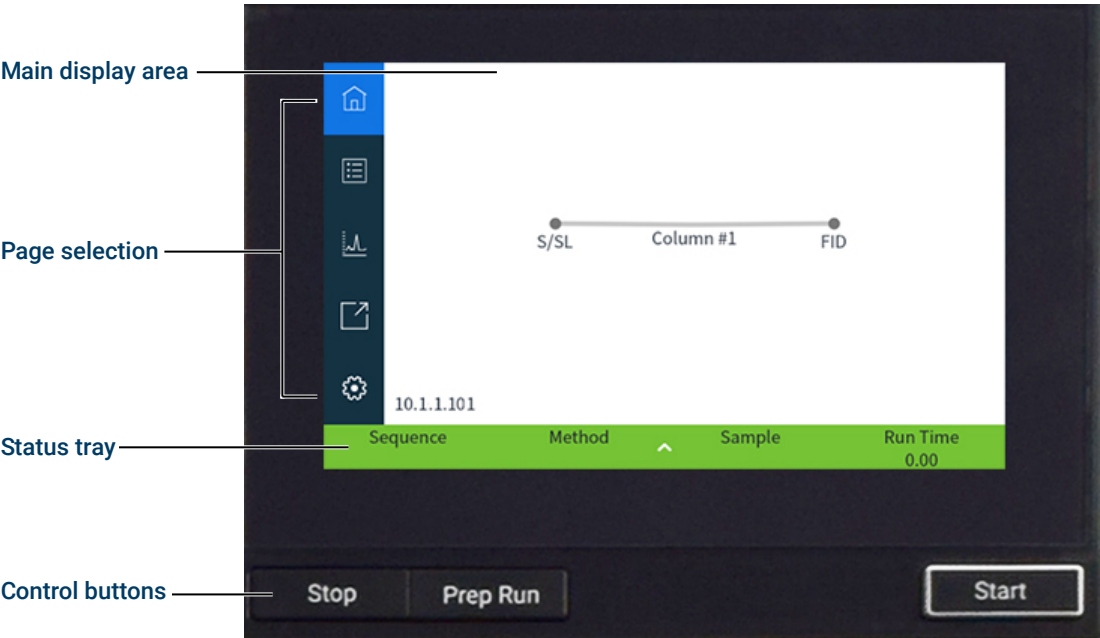
Status page

Plot page

This section describes the basic operation of the Agilent GC touchscreen and control button interface.

Navigation

The touchscreen shows GC status and activity information (current temperatures, flows, pressures, and information about GC readiness). The control buttons allow you to start, stop, and prepare the GC to run a sample.



On the home page, selecting one of the page selection buttons loads the corresponding page.

The main display area provides information related to the selected functional area/page. This area contains status displays, controls, settable parameters, and so on.

Depending on which page is selected, additional controls may appear. This can include page selection buttons, selectable tabs, back and next buttons, scroll buttons, and so on. See [Figure 9](#).

	On/Off	Actual	Setpoint	
Oven Temperature	☐	29.1	50.0	
Front Inlet Temperature	☐	26.1	250.0	
Front Inlet Pressure(psi)	☐	0.06	11.00	^
Front Inlet Total Flow		-4.5	10.0	
Back Inlet Temperature	☐	26.4	250.0	
Back Inlet Pressure(psi)	☐	0.02		
Back Inlet Total Flow		-3.9	25.0	v
Front Detector Temperature	☐	27.5	275.0	
Sequence	Method	Sample	Run Time	
			0.00	

Figure 9. Additional controls

Scroll buttons are enabled if additional information or settings are available via scrolling.

Run controls

Runs are controlled by the buttons (Stop, Prep Run, and Start) located beneath the touchscreen. The run controls are used to start, stop, and prepare the GC to run a sample.

Prep Run	The Prep Run control activates processes required to bring the GC to the starting condition for a run (such as turning off the inlet purge flow for a splitless injection). This is typically required before manual injections to exit any gas saver mode and to prepare the inlet flows for injection.
Start	The Start control starts a run after manually injecting a sample. (When you are using an automatic liquid sampler or gas sampling valve, the run is automatically activated at the appropriate time.)
Stop	The Stop control immediately terminates the run. If the GC is in the middle of a run, the data from that run may be lost.

For details on running methods, see [Running Methods](#).

Status/control tray

The status/control tray provides details on the current status of the GC, the current sequence and method (if connected to an Agilent data system), the time remaining for the current operation being performed by the GC, run controls, and so on.

The status/control tray is color coded to reflect the run or ready status of the GC, see [Status Colors and details](#) for more information.

Any Early Maintenance Feedback (EMF) flags are also displayed. See [Early Maintenance Feedback \(EMF\)](#).

The tray can be expanded by selecting the up arrow on the tray. See [Figure 10](#).

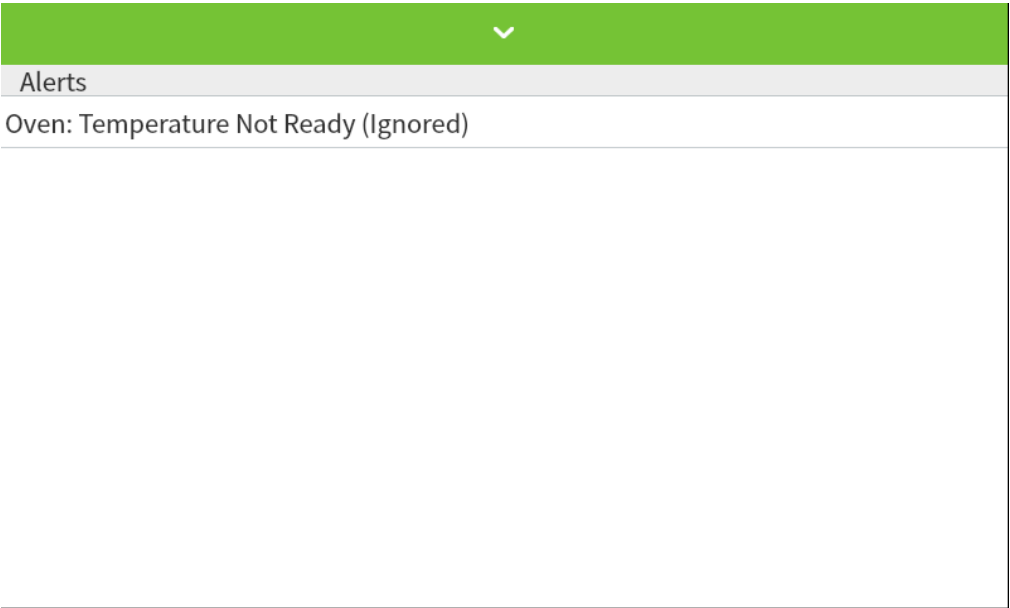


Figure 10. Status/control tray - expanded

The tray can be minimized by selecting the arrow on the tray.

Status Colors and details

Colors provide quick GC status:

Ready (all settings at setpoint)	STATUS: NORMAL
Run in progress	STATUS: RUN
Not ready	STATUS: NOT READY
Maintenance mode (an automated maintenance task is running)	STATUS: MAINTENANCE MODE
Preparing sample	STATUS: PREPARING SAMPLE
GC is in sleep mode	STATUS: SLEEP
GC is in an error state	STATUS: ERROR

Status text displayed in the status tray can include:

Status text	Comment
CONDITIONING RUN	GC is performing a conditioning run after waking from Sleep.
COOLDOWN	The GC is performing a cooldown between runs.
DIAGNOSTIC MODE	The GC is currently performing a diagnostic test.
ERROR	An error is preventing normal operation. Check the Diagnostics and Logs tabs for more information.

Status text	Comment
H2 SHUTDOWN	The GC is shut down: heaters are turned off, the oven flaps are open, and gas flows are off.
MAINTENANCE MODE	The GC is running a user maintenance procedure.
MSD VENTING	The GC is set to provide flows appropriate for MSD venting.
MSD VENTED	The MSD is vented.
NORMAL	No conditions prevent operation. The GC may or may not be ready to begin a run.
POST RUN	The GC is executing a method post run program.
PREPARING SAMPLE	During a run, a connected ALS is performing sample preparation.
PREP RUN	The GC is in the Prep Run state, preparing for an injection.
RUN	The GC is performing a run.
SLEEP	The GC is in sleep mode. To wake the GC, apply a setpoint change, download a method, or from an Agilent data system select "Wake Up."
SLEEP DEFERRED	Due to ongoing activity, the GC did not enter sleep mode when scheduled. The GC will enter sleep mode after it is has been idle for at least 15 minutes.
SHUTDOWN	A GC component is preventing the GC from operating normally. Check for a diagnostic condition or message at the Diagnostics tab.

Entering Data

When you touch a data entry field, a touch keyboard or keypad appears, as applicable. See **Figure 11**.

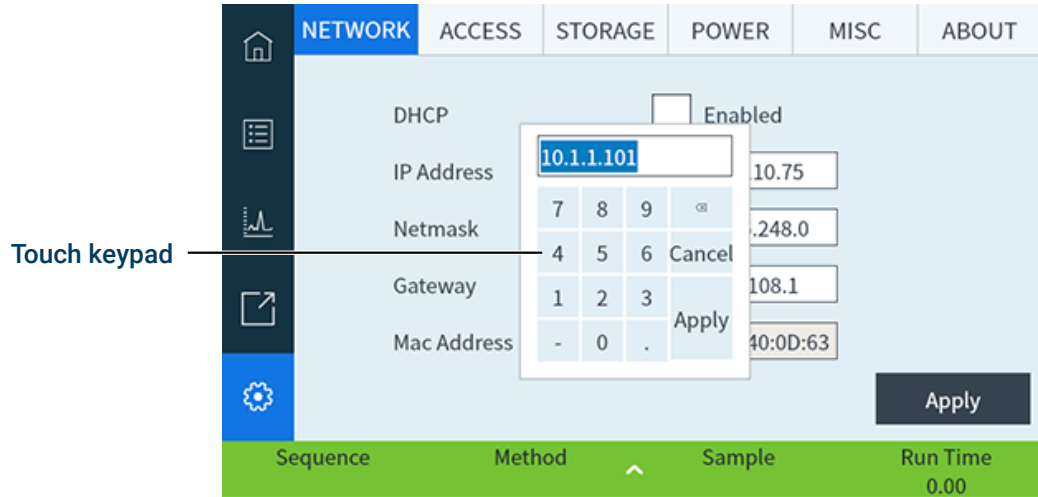


Figure 11. Touch keypad for data entry

If you enter an out of range entry, it is highlighted in a different color.

If the field is a drop-down list box (indicated by a down arrow to the right of the displayed contents of the field), select it to open the list, and then select the desired entry.

Home View

The home view's Flow Path page shows flow path information (including current temperatures and flow rates), run status (including user selectable status items), access to shortcuts to allow quick access to common tasks, a real-time plot of the current chromatogram, and related information. See **Figure 12**.

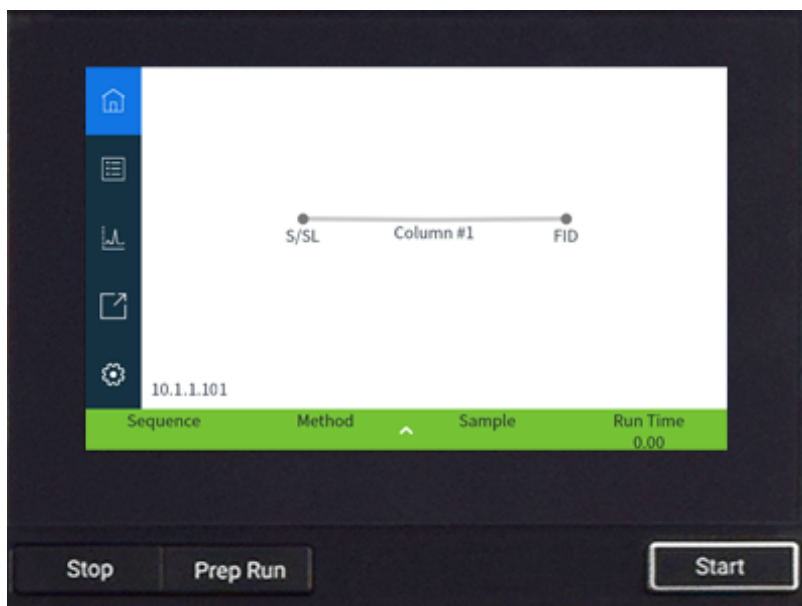


Figure 12. Home view

The other pages available from the home view are:

- Flow path
- Status
- Plot
- Shortcuts
- Settings

These pages are displayed by selecting the corresponding page selection button on the left side of the home view.

The flow path page provides details on sample flow through the GC. This includes visual indications of whether an ALS is installed on the GC, the inlet type(s), column setup, and detector type(s), along with corresponding setpoints.

Each page is described below.

Status page

The status page displays a user-selectable list of parameters, along with their setpoints and actual values. See **Figure 13**.

		On/Off	Actual	Setpoint	
	Oven Temperature	☐	29.1	50.0	^
	Front Inlet Temperature	☐	26.1	250.0	
	Front Inlet Pressure(psi)	☐	0.06	11.00	
	Front Inlet Total Flow		-4.5	10.0	v
	Back Inlet Temperature	☐	26.4	250.0	
	Back Inlet Pressure(psi)	☐	0.02		
	Back Inlet Total Flow		-3.9	25.0	
	Front Detector Temperature	☐	27.5	275.0	
Sequence		Method	Sample	Run Time	
				0.00	

Figure 13. Home view - status page

Plot page

The plot page displays a plot of the currently selected signal. See [Figure 14](#).

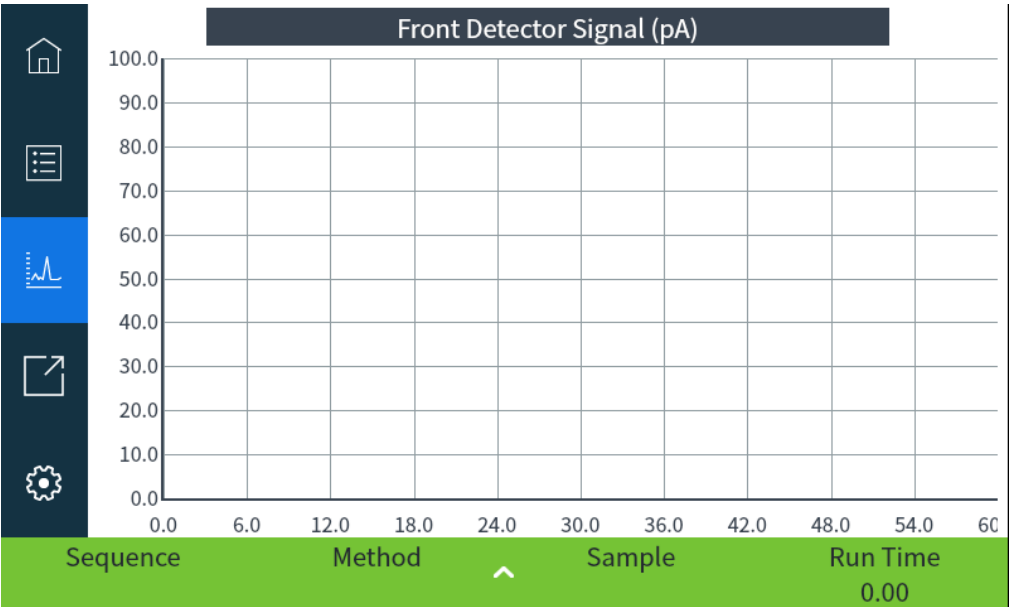


Figure 14. Home view - plot page

Selecting the displayed signal name opens a Plot Options dialog box. This allows you to select which signal to display. See [Figure 15](#).

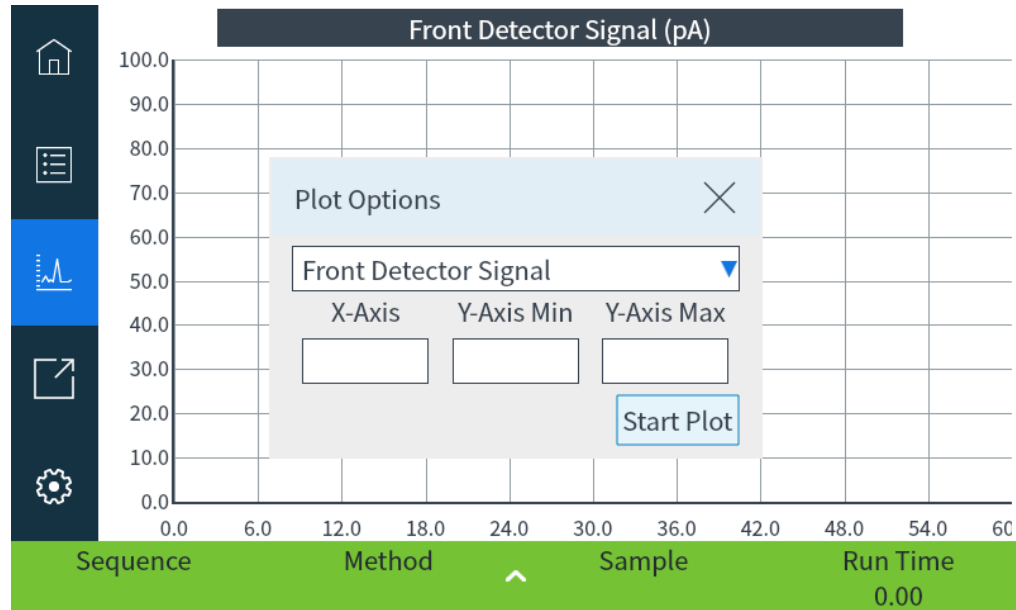


Figure 15. Plot Options dialog box

Use the drop down list box to select which parameter to display on the plot.

The displayed **X-Axis** interval is 1 to 60 minutes. The **Y-Axis Range** is negative infinity to infinity. Selecting either field brings up a keypad which allows you to set the corresponding value.

If the plot is not currently running, selecting **Start Plot** starts it. If the plot is currently running, selecting **Stop Plot** halts data collection and display. (When changing the Signal Name, it may be necessary to select **Stop Plot** and then **Start Plot** to display the signal.)

To zoom the plot, tap it. Tapping the plot cycles through different levels of zoom.

Shortcuts

Selecting the shortcuts button opens the shortcuts menu and shows the user their selected list of shortcuts.

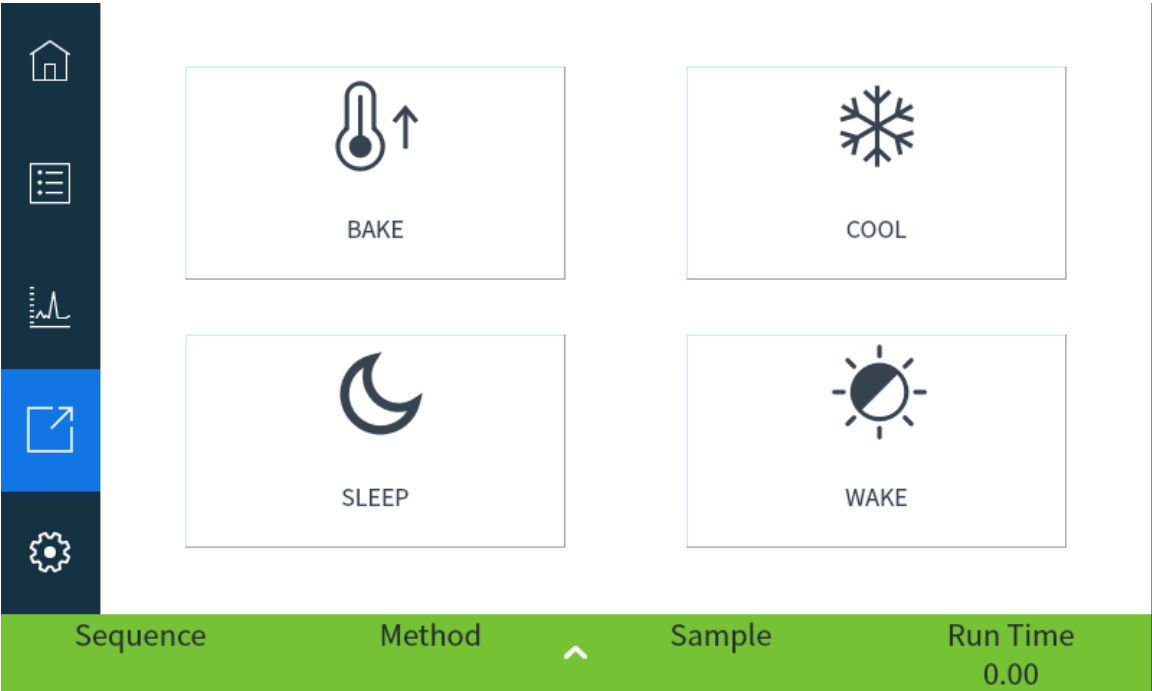


Figure 16. Shortcuts

The shortcuts menu will be pre-populated with standard shortcuts:

Shortcut	Function
Sleep	Invoke the sleep method
Wake	Invoke the wake method
Cool	Invoke the Cool down maintenance procedure
Bake	Start the column Bake maintenance procedure

5

Methods

What is a Method?

What Is Saved in a Method?

What Happens When You Load a Method?

Create a Method

Load a Method

Running Methods

Events

Inlets

About gas saver

About Oven Temperature Programming

Columns

Detectors

FID

FPD+

NPD

TCD

ECD

Valves

The 8860B valve box

Valve control

Valve types

Controlling a valve

GC Output Signals

Analog signals

Digital signals

What is a Method?

A method is the group of settings required to analyze a specific sample.

Since every type of sample reacts differently in the GC—some samples require a higher oven temperature, others require a lower gas pressure or a different detector—a unique method must be created for each specific type of analysis.

The GC also can store several specialized methods. The GC stores three methods used for resource conservation, called **SLEEP**, **CONDITION**, and **WAKE**. When configured for use with an attached MS, the GC also provides a method called **MS VENT**, used to change GC setpoints to values appropriate for a safe MS venting process. Create these methods using an Agilent data system. See [Resource Conservation](#) for more information about these specialized methods.

Methods can be created and edited using the Browser Interface. Methods can also be created, edited, and stored on a connected data system. While a data system is connected to your GC, any attempt to connect to your GC via the Browser Interface will result in the Browser Interface having limited functionality. You will not be able to edit methods, edit sequences, and start or stop runs until the data system is disconnected. Similarly, once a Browser Interface is connected to your GC, any subsequent Browser Interfaces attempting to connect will not be able to edit methods, edit sequences, and start or stop runs until the first session is disconnected.

Methods and sequences created using the Browser Interface are not directly accessible from a connected data system. Methods and sequences created using a connected data system are not directly accessible using the Browser Interface.

What Is Saved in a Method?

Some of the settings saved in a method define how the sample will be processed when the method is used. Examples of method settings include:

- The oven temperature program
- The type of carrier gas and flows
- The type of detector and flows
- The type of inlet and flows
- The type of column
- The length of time to process a sample

NOTE

For GCs equipped with EPR (electronic pneumatics regulation), detector, column, and inlet flows are not saved with the method.

Data analysis and reporting parameters are also stored in a method when it is created on an Agilent data system, for example OpenLab CDS or MassHunter software. These parameters describe how to interpret the chromatogram generated by the sample and what type of report to print.

The GC method also includes sampler setpoints. Refer to the sampler documentation for details on setpoints for the supported device:

- For the 7650A ALS, see its Installation, Operation, and Maintenance manual.
- For the 7693A ALS, see its Installation, Operation, and Maintenance manual.
- For the 7697A HS, see its Installation and First Startup manual and Operation manual.
- For the CTC PAL3 AS, see its Installation and First Startup manual and Operation manual.

Current setpoint parameters are saved when the GC is turned off, and loaded when you turn the instrument back on.

What Happens When You Load a Method?

There are two kinds of methods:

- The active method—This is sometimes referred to as the current method. The settings defined in this method are the settings the GC is currently maintaining.
- Stored methods—User-created methods can be stored in the GC, along with one SLEEP method, one WAKE method, one CONDITION method, an MS VENT method, and a default method.

For EPC (electronic pneumatics control) equipped GCs, when a method is loaded from the GC or Agilent data system, the setpoints of the active method are immediately replaced with the setpoints of the method loaded.

For EPR (electronic pneumatics regulation) equipped GCs, when a method is loaded from the GC or Agilent data system, any manually set parameters are not replaced with the setpoints of the method loaded.

- The method loaded becomes the active (current) method.
- The GC front panel status indicator will be yellow (not ready) and will stay yellow until the GC reaches all of the settings specified by the method that was just loaded.

Create a Method

Methods may be created using your Agilent data system, or via the Browser Interface. For more information on creating a method using your data system, refer to the documentation provided with your data system.

From the Browser Interface:

1. Click **Method** on the control ribbon.
2. Click the **Create New** button . You are prompted to name your method and save.
3. From the navigation tree, select each instrument device, and set its method parameters to the desired values.
4. Click **Configuration > Modules**. Check the gas configuration for each inlet and detector. (Gas configuration was initially set when the Setup Wizard was run at installation.) Make any changes if needed.
Click **Configuration > Columns**. If the installed columns include Smart ID Keys, the column configurations are set by the Smart ID Keys. If not, edit the column configuration as needed. Double-click a column to edit it.
5. Click **Columns**. For each column, set the control mode, select the **On** checkbox, and set the flow on the column.
6. Click **Inlets**, and then select either the front or back inlet. Set the inlet mode, temperatures, and other parameters needed for your method. Repeat for the other inlet, if present.
7. Click **Detectors**. For each detector being used, set the detector temperature and gas flows. Select all check boxes to turn the detector on.

NOTE

Carrier Gas Flow Correction setting recommendations

- **Column + Fuel = Constant** (H₂ carrier in Constant Pressure mode)
 - **Column + Makeup = Constant** (He/N₂ carrier Constant Pressure mode)
 - **Constant Makeup and Fuel Flow** (any carrier in Constant Flow mode)
8. Click **Oven**. Set the initial oven temperature, ramp temperatures, and times for your method. Select the **Heater On** checkbox. For an isothermal run, do not create any ramps.
 9. Click **Signals**. Choose the signals to be included in the data file when a run is initiated. Normally this is **Front Signal** or **Back Signal** for detector output. Select the **Save** check box and select a data rate that meets the needs of your chromatography.
 10. Click **ALS/Tray**, then select either the front or back injector. Set the **Injection Volume**, washes, and pumps.

NOTE

It is important to do both sample and solvent washes to eliminate carryover and keep the syringe clean. Pumping the plunger multiple times eliminates air bubbles in the syringe for better reproducibility.

11. If using a tray, click **ALS/Tray > Other**. Set **Sample Overlap** if desired.
12. Click the **Save** button  to save your method.

Load a Method

1. Connect to the GC using the Browser Interface. See [Browser Interface](#).
2. Open or create the desired method. See [Create a Method](#).
3. If needed, click the **Save** button to save the method .
4. Click the **Download** button  to load the method to your GC.

Running Methods

Pre Run and Prep Run

With some inlets and operating modes, certain instrument setpoints are different between runs than during an analysis. To restore the setpoints for injection, you must place the GC into the Pre Run state.

You must use the Pre Run state when:

- Using gas saver with any inlet.
- Using splitless mode with any inlet.
- Using a pressure pulse mode with any inlet.

NOTE

Gas saver and pressure pulse mode are not supported on EPR (electronic pneumatics regulation) equipped GCs.

There are three ways to begin Pre Run—manually (press the "Prep Run" button beneath the touchscreen before each run), automatically (for Agilent samplers), or using **Auto Prep Run** (for non-Agilent samplers). The three methods are discussed below.

During the Pre Run state:

- The status on the touchscreen changes to indicate the GC is preparing for the injection.
- Setpoints change to the correct values for injection.
- Inlet, detector, and oven equilibration times begin.

When all criteria for a run are met, the Status/Control tray indicates that the GC is ready for sample injection.

Manually prepare for the run

Press the "Prep Run" button before you inject a sample manually. The GC enters the Pre Run state. When the touchscreen indicates that the GC is ready, begin the analysis.

If you are using an Agilent automatic sampling system, the **Prep Run** function is automatic.

To manually inject a sample with a syringe and start a run

1. Prepare the sample syringe for injection.
2. Load the desired method. See [Load a Method](#).
3. Press the "Prep Run" button. See [Run controls](#) for more information.
4. Wait for status **Ready** to display.
5. Insert the syringe needle through the septum and all the way into the inlet.
6. Simultaneously depress the syringe plunger to inject the sample and press the "Start" button.

To run a method to process a single ALS sample

1. Prepare the sample for injection.
2. Load the sample vial into the assigned location in the ALS tray or turret.
3. Load the desired method. (See [Load a Method](#).)
4. Press the "Start" button to initiate the ALS syringe cleaning, sample loading, and sample injection method. After the sample is loaded into the syringe, the sample is automatically injected when the GC reaches the ready state. See [Run controls](#) for more information.

To abort a method

1. Press the "Stop" button.
2. When you are ready to resume running analyses, load the appropriate sequence or method. (See [Load a Method](#).)

Events

Run time programming during a method allows certain setpoints to be changed during a run as a function of the chromatographic run time. Thus an event that is programmed to occur at 2 minutes will occur 2 minutes after every injection.

- Controlling column switching or other valves
- Changing analog signal definition, zero, or range
- Controlling an auxiliary pressure channel
- Changing polarity of a thermal conductivity detector (TCD)
- Make sample flow bypass TCD filament
- Turning the hydrogen flow to a nitrogen-phosphorus detector (NPD) on or off
- Switching digital signal output (requires an Agilent data system)
- Pausing ("freezing") and resuming digital signal output (requires an Agilent data system)
- Performing signal math on the front and back detectors.

The changes are entered into a run table that specifies the setpoint to be changed, the time for the change, and the new value. At the end of the chromatographic run, most setpoints changed by a run time table are returned to their original values.

Valves can be run time programmed but are not restored to their starting position at the end of the run. You must program the reset operation in the run table if this action is desired.

Using run time events

The **Events** page in the **Method** tab is used to program the following timed events.

- Valves (1-10)
- Multiposition valve
- Signal type
- Analog signal definition, zero, and range
- Auxiliary pressures (1 through 9)
- TCD negative polarity (on/off)
- Detector gas flow (on/off), including NPD H₂ fuel gas
- Inlet septum purge flow
- AO signal range
- AO signal source
- AO signal zero
- Bypass TCD Filament

Programming run time events

1. Select **Method**.
2. Select **Events**.
3. Input the time at which you want the event to occur, select the event you want to program, select the position of the hardware that is being controlled, and set the desired setpoint.

The run table

The programmed events are arranged in order of execution time in the Run Table.

Editing events in the run table

1. Select the event you wish to edit. If the desired event is not immediately visible, use the up and down arrows to the right to scroll up and down the table until it is.
2. Select the parameter you want to change.
3. Input the new value

Deleting run time events

1. Select the event you wish to delete.
2. Select **Delete**.

Inlets

The Inlets page in the Method tab is used to modify the method parameters and each inlet connected to your GC. Common parameters include the heater temperature and inlet pressure. To modify your inlets' method parameters:

1. Select **Method > Inlets**.
2. Select the inlet you want to modify.
3. Scroll to the desired setting and edit as necessary.

Carrier gas flow rates

The flow rates in [Table 2](#) are recommended for all column temperatures.

Table 2. Column size and carrier flow rate

Column type	Column size	Carrier flow rate, mL/min		
		Hydrogen	Helium	Nitrogen
Packed	1/8-inch		30	20
	1/4-inch		60	40
Capillary	0.05 mm id	0.5	0.4	n/a
	0.10 mm id	1.0	0.8	n/a
	0.20 mm id	2.0	1.6	0.25
	0.25 mm id	2.5	2.0	0.5
	0.32 mm id	3.2	2.6	0.75
	0.53 mm id	5.3	4.2	1.5

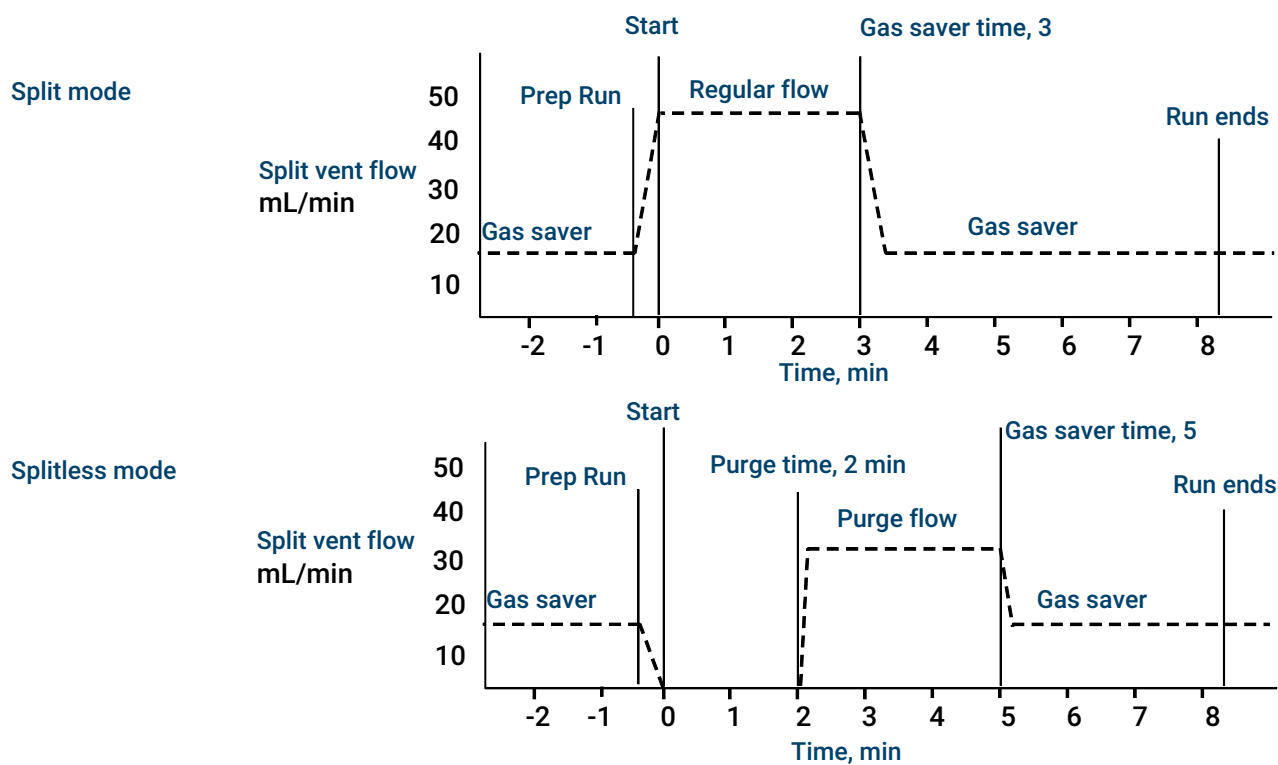
About gas saver

NOTE

Gas saver is not supported on EPR (electronic pneumatics regulation) equipped GCs.

Gas saver reduces carrier flow from the split vent after the sample is on the column. It applies to the Split/Splitless inlet (all modes). It is most useful in split applications.

Column head pressure and flow rate are maintained, while split vent flows decrease. Flows—except column flow—remain at the reduced level until you press the "Prep Run" button.



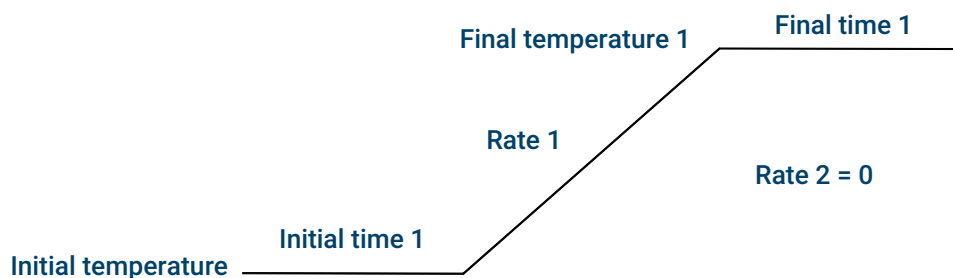
To use gas saver

1. Select **Method > Inlets**.
2. Scroll down until the Gas Saver option is visible.
3. Select the checkbox next to **Gas Saver** to enable it.
4. Input the **Setpoint**. The setpoint must be at least 15 mL/min greater than the column flow.
5. Input the **Time**.

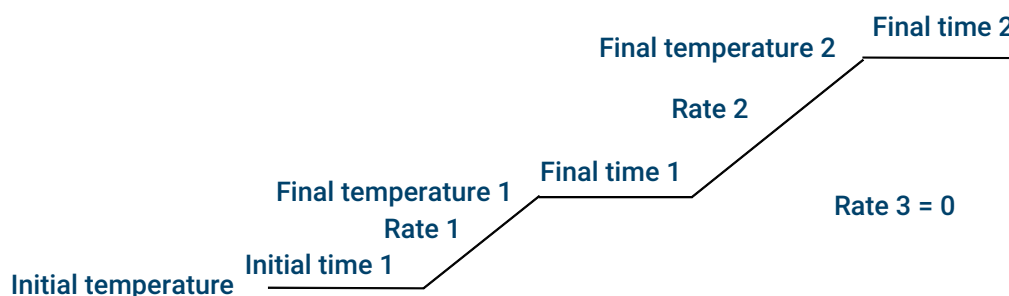
About Oven Temperature Programming

You can program the oven temperature from an initial temperature to a final temperature using up to 5 ramps during a run.

A single ramp temperature program raises the initial oven temperature to a specified final temperature at a specified rate and holds at the final temperature for a specified period of time.



The multiple-ramp temperature program is similar. You can program the oven from an initial temperature to a final temperature, but with various rates, times, and temperatures in between. Multiple ramps can also be programmed for temperature decreases as well as increases.



Oven ramp rates

The highest rate that you can achieve depends on many factors, including the room temperature, temperatures of the inlets and detectors, the amount of material inside the oven (columns, valves, etc.), and whether or not this is the first run of the day.

Table 3 lists typical oven ramp rates.

Table 3. Oven ramp rates

Temperature range (°C)	100 V oven ramp rate (°C/minute)	200/220/230/240 V oven ramp rate (°C/minute)
50 to 70	30	75
70 to 115	30	45
115 to 175	30	40
175 to 300	30	30
300 to 425	20	20

Isothermal runs

An isothermal run is one in which the oven is maintained at a constant temperature. For an isothermal run, set Rate 1 to zero.

1. Select **Oven** to open the oven parameter list.
2. Enter the oven temperature for the isothermal run.
3. Enter the number of minutes (Initial time) that you want the oven to stay at this temperature. This time is the duration of the run.
4. If Rate 1 is not already 0, enter zero for an isothermal run.

Columns

The column flow mode determines whether pressure or flows are available as setpoints for the GC inlets. When all columns in the flow path are defined, you can enter pressures or flows. If any column in the flow path is not defined, then the settings for the inlet may be restricted based on the inlet type and whether the column is in flow or pressure mode.

There are two different flow modes that control the mass flow rate through the column: **Constant flow** and **Ramped flow**. The flow rates are corrected to NTP (normal temperature and pressure, 25 °C and 1 atmosphere).

There are two different pressure modes that control the pressure at the head of the column: **Constant pressure** and **Ramped pressure**. These pressures are gauge pressures—the difference between the absolute pressure and the local atmospheric pressure.

Flow modes

- **Ramped flow** — Increases the mass flow rate in the column during the run according to a program you enter. A column flow profile can have up to three ramps, each consisting of a programmed increase followed by a hold period.
- **Constant flow** — Maintains a constant mass flow rate of carrier gas in the column throughout the run. If the column resistance changes due to a temperature program, the column head pressure is adjusted to keep the flow rate constant. This can shorten runs significantly.

Pressure modes

The pressure modes are not available if the column is not defined or the inlet's mode is set to Flow control. Pressures are gauge pressures—the difference between the absolute pressure and the local atmospheric pressure.

Because most detectors present little resistance to the column flow, the gauge pressure at the column head is usually the same as the pressure difference between column inlet and exit. The mass selective detector and the atomic emission detector are the exceptions.

- **Constant pressure** — Maintains a constant gauge pressure at the head of the column throughout the run. If the column resistance and gas density changes during a run, the gauge pressure does not change but the mass flow rate does.
- **Ramped pressure** — Increases the column head gauge pressure during the run according to a program you enter. A column pressure profile can have up to three ramps, each consisting of a programmed increase followed by a hold period.

NOTE

Flow modes and ramped pressure mode are not supported on SSL EPR (electronic pneumatics regulation) equipped GCs.

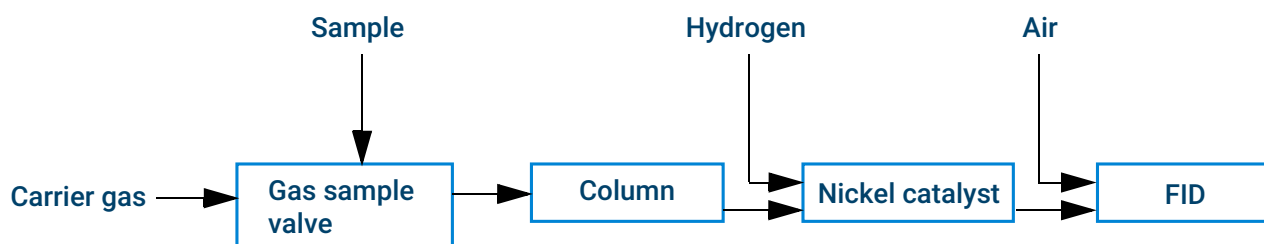
Pressure modes are not supported on PPIP EPC equipped GCs.

Pressure modes and ramped flow mode are not supported on PCI EPR equipped GCs.

Nickel Catalyst Tube

About the nickel catalyst tube

The Nickel Catalyst Tube accessory, G3534A, is used for trace analysis of CO and CO₂ with a flame ionization detector. The gas sample is separated on the column and passed over a heated catalyst in the presence of hydrogen, which converts the CO and CO₂ peaks to CH₄.



Detectors

For help creating a new method, or troubleshooting detector problems, see the recommended starting conditions for each detector.

The makeup gas line of your detector parameter list changes depending on your instrument configuration. If you have an inlet with the column not defined, the makeup flow is constant.

FID

FID automatic reignition (Lit offset)

Lit offset is the expected minimum difference between the FID output with the flame lit and the output with the flame off. The GC checks this value during runs and when loading a method.

During a run, if the output falls below the **Lit offset** value, the FID will attempt to reignite three times. If after the third attempt the output does not increase by at least this value, the detector shuts down all functions except temperature and makeup gas flow.

When loading a method that includes a **Flame On** setting, the GC performs a similar check. If the detector output is less than the **Lit offset**, it will attempt reignition after reaching method setpoints.

The default setting for **Lit offset** is 2.0 picoamps. This is a good working value for all but very clean gases and systems. You may want to lower this setpoint if the detector attempts to reignite when the flame is still on, thus producing a shutdown.

To change **Lit offset**:

1. In the browser interface, go to **Method > Configuration > Modules**.
2. Enter a new **Lit Offset** value for the desired detector.
3. Save the method.

NOTE

The newly entered **Lit offset** is not applied/referenced against the detector output until the next ignition cycle.

Recommended starting conditions

See [Table 4](#) for guidelines and rules to select initial detector settings for new methods.

Table 4. Recommended starting conditions

Gas type	Suggested flow rate
Carrier gas (hydrogen, helium, nitrogen)	
Packed columns	10 to 60 mL/min
Capillary columns	1 to 5 mL/min
Detector gases	
Hydrogen	40 mL/min ¹
Air	450 mL/min ¹
Column plus capillary makeup (N ₂ is recommended, or He as an alternative)	50 mL.min ¹
Detector Temperature	
If < 150 °C, the flame will not light. Agilent recommends a temperature ≥ 300 °C to prevent condensation damage. The detector temperature should be approximately 20 °C greater than highest oven ramp temperature.	

¹The hydrogen / air ratio should be 8% to 12% to keep the flame lit.

FPD+

The sample burns in a hydrogen- rich flame, where some species are reduced and excited. The gas flow moves the excited species to a cooler emission zone above the flame where they decay and emit light. A narrow bandpass filter selects light unique to one species, while a shield prevents intense carbon emission from reaching the photomultiplier tube (PMT).

The light strikes a photosensitive surface in the PMT where a light photon knocks loose an electron. The electron is amplified inside the PMT for an overall gain of up to a million.

Storage

The FPD+ should not be stored at temperatures above 50 °C.

FPD+ linearity

Several mechanisms produce sulfur emission. The excited species is diatomic, so that emission intensity is approximately proportional to the square of the sulfur atom concentration.

The excited species in the phosphorus mode is monatomic, leading to a linear relationship between emission intensity and atom concentration.

Inlet liners for use with the FPD+

Compounds containing sulfur may adsorb on an inlet liner and degrade the GC's performance. Use deactivated, clean liners or a cool on-column inlet, which injects directly onto the column.

For best results use Agilent Ultra Inert liners:

- Splitless 5190-2293
- Split 5190-2295

FPD+ temperature considerations

The FPD+ provides two temperatures zones, one for the transfer line (the main detector temperature) and one for the emission block. For the transfer line temperature, we recommend a temperature that is 25 °C higher than the highest column temperature.

The emission block temperature ranges from 125–175 °C. Typically, the default temperature of 150 °C is sufficient for most applications. However, when setting the emission block temperature, consider the following:

- If using the GC oven at high temperature (> 325 °C) with the transfer line set to 400 °C, set the emission block temperature to 165 °C to avoid a system Not Ready if the emission block temperature cannot be maintained.
- If using the transfer line at 400 °C, set the emission block temperature to at least 150 °C to avoid a system Not Ready.
- For sulfur analyses, the highest area response will be realized with the lowest possible emission block temperature.
- For phosphorus analyses, the area response is independent of the emission block temperature.

Recommended starting conditions

The FPD+ provides two temperatures zones, one for the transfer line (the main detector temperature) and one for the emission block. For the transfer line temperature, we recommend a temperature that is 25 °C higher than the highest column temperature. The emission block temperature ranges from 125–175 °C. Typically, the default temperature of 150 °C is sufficient for most applications.

Flows for the maximum sensitivity FPD+ flame, which is hydrogen-rich and oxygen-poor. Helium, either as carrier or makeup gas, may cool the detector gases below the ignition temperature.

Table 5. Recommended starting conditions

Gas type	Suggested flow rate
Carrier gas (hydrogen, helium, nitrogen)	
Packed columns	10 to 60 mL/min
Capillary columns	1 to 5 mL/min
Detector gases	
Hydrogen	60 mL/min
Air	60 mL/min
Column plus capillary makeup	60 mL.min

As with the FID, the FPD has a Lit Offset associated with it. The default Lit Offset for the FPD is 2.0 pA.

NPD

Setting NPD adjust offset on the clock table

You can use the **Clock table** feature to begin **Adjust offset** at a specified time.

Extending the NPD bead life

These actions, together with the automated heatup and adjust procedures, can extend bead life considerably.

- Use the lowest practical **Adjust Offset** value. This will result in a lower applied bead current during operation.
- Run clean samples.
- Turn the bead off when not in use.
- Keep the detector temperature high (320 to 335 °C).
- Turn the hydrogen flow off during solvent peaks and between runs.

Turning hydrogen off during a solvent peak

When using the NPD, the baseline shifts after a solvent peak and can take some time to stabilize, especially with chlorinated solvents. To minimize this effect, turn off the hydrogen flow during the solvent peak and turn it back on after the solvent elutes. With this technique, the baseline recovers to its original value in less than 30 seconds. This also extends the life of the bead. The hydrogen can be turned on and off automatically as part of a Run Table. See [Events](#).

Turning hydrogen off between runs

To extend bead life, turn off the hydrogen flow between runs. Leave all other flows and the detector temperature on. Turn on the hydrogen flow for the next run; the bead will ignite almost immediately. The process can be automated with Run Table entries.

Recommended Starting conditions

Table 6. Recommended starting conditions

Gas type	Suggested flow rates
Carrier gas (helium, hydrogen, nitrogen ¹)	Capillary: choose an optimal flow rate based on column dimensions.
Detector gases	
Hydrogen	1 to 3 mL/min
Air	60 mL/min
Makeup Flow (He, N ₂ ²)	1 to 20 mL/min, less is recommended
Bead voltage	
• Use Adjust Bead and let the GC set the applied bead current for you. • Bead voltage is not settable as it is calculated based on applied current. The bead voltage can be viewed from the Browser Interface: Settings > Service mode > Detectors > Signals.	

¹Flow rate must be less than 3 mL/min when using hydrogen as carrier gas.

²Helium is recommended for best peak shapes.

TCD

Chemically active compounds reduce TCD filament life

The tungsten-rhenium TCD filament has been chemically passivated to protect against oxygen damage. However, chemically active compounds such as acids and halogenated compounds may attack the filament. The immediate symptom is a permanent change in detector sensitivity due to a change in filament resistance.

If possible, such compounds should be avoided. If this is not possible, the TCD cell may have to be replaced frequently.

Changing the TCD polarity during a run

Negative polarity On inverts the peak so the integrator or Agilent data system can measure it.

Negative Polarity can be a run table entry; see [Events](#).

Detecting hydrogen with the TCD using helium carrier gas

Hydrogen is the only element with thermal conductivity greater than helium, and mixtures of small amounts of hydrogen (<20%) in helium at moderate temperatures exhibit thermal conductivities less than either component alone. If you are analyzing for hydrogen with helium carrier gas, a hydrogen peak may appear as positive, negative, or as a split peak.

There are two solutions to this problem:

- Use nitrogen or argon-methane as carrier gas. This eliminates problems inherent with using helium as carrier, but causes reduced sensitivity to components other than hydrogen.
- Operate the detector at higher temperatures—from 200 °C to 300 °C.

You can find the correct detector operating temperature by analyzing a known range of hydrogen concentrations, increasing the operating temperature until the hydrogen peak exhibits normal shape and is always in the same direction (negative relative to normal response to air or propane) regardless of concentration. This temperature also ensures high sensitivity and linear dynamic range.

Because hydrogen peaks are negative, you must turn negative polarity on at appropriate times so the peak appears positive.

Recommended Starting conditions

Table 7. Recommended starting conditions

Gas type	Suggested flowrates
Carrier gas	Packed: 10 to 60 mL/min
(hydrogen, helium, nitrogen)	Capillary: 1 to 5 mL/min
Reference	15 to 60 mL/min
(same gas type as carrier)	
Capillary makeup	Packed: 2 to 3 mL/min
(same gas type as carrier)	Capillary: 5 to 15 mL/min
Detector temperature	
< 135 °C, cannot turn on filament	
If detector temperature goes below 125 °C, the filament turns off.	
Detector temperature should be 30 °C to 50 °C greater than the highest oven ramp temperature.	

ECD

ECD linearity

The ECD response factor versus concentration curve is linear for four orders of magnitude or more (linear dynamic range = 10^4 or higher) for a broad range of compounds. You should still run a calibration curve on your samples to find the limits of the linear range for your materials.

ECD makeup gas notes

If the carrier gas type is different from the makeup gas type, the makeup gas flow rate must be at least three times the carrier gas flow rate.

ECD sensitivity can be increased by reducing the makeup gas flow rate.

ECD chromatographic speed (for fast peaks) can be increased by increasing the makeup gas flow rate.

ECD temperature programming

The ECD is flow sensitive. If you are using temperature programming, in which the column flow resistance changes with temperature, set up the instrument as follows:

- Set the carrier gas in the **Constant flow** mode. Set detector makeup gas to **Constant makeup**.
- If you choose to work in the constant pressure mode, the makeup gas should be set in the **Column + makeup = constant** mode.

Recommended starting conditions

Use the following information when selecting temperatures and flows. Maximum source pressure must not exceed 100 psi. Use the maximum source pressure to achieve maximum makeup flow rate.

Table 8. Recommended starting conditions

Gas type	Suggested flow rates
Carrier gas	30 to 60 mL/min
Packed columns (nitrogen or argon-methane)	0.1 to 20 mL/min, depending on diameter
Capillary columns (hydrogen, nitrogen, or argon-methane)	
Capillary makeup (nitrogen or argon-methane)	10 to 150 mL/min (30 to 60 mL/min typical)
Temperature	
250 °C to 400 °C	
Detector temperature is typically set 25 °C greater than the highest oven ramp temperature.	

Valves

The 8860B valve box

The GC holds up to four valves in a heated valve box on top of the oven.

The valve box is the preferred location for valves because it is a stable temperature zone, isolated from the column oven.

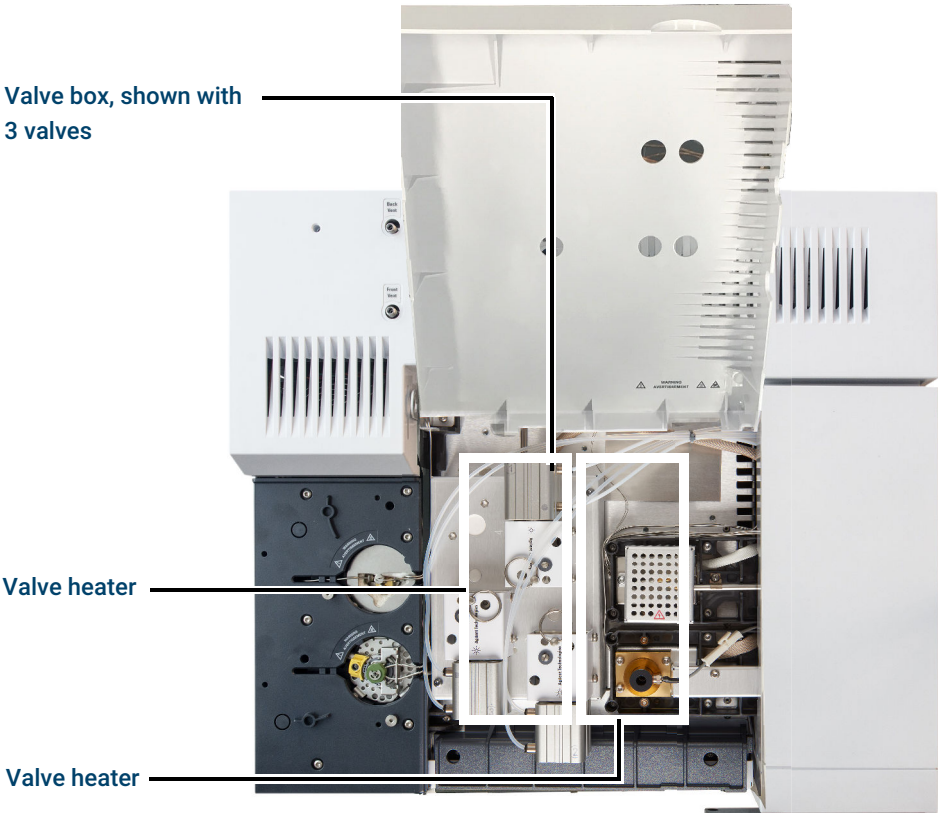


Figure 17. Diagram of valve locations on GC

Valve control

Valves can be controlled manually from the Browser Interface or as part of a clock or run time program. Note that sampling valves automatically reset at the end of a run.

The valve drivers

A valve driver is the software and circuitry in the GC that controls a valve or related function. There are eight drivers, known as Valves 1, 2, 3, 4, 7, 8, 9, and 10. Each valve is independently controlled.

Table 9. Valve drivers

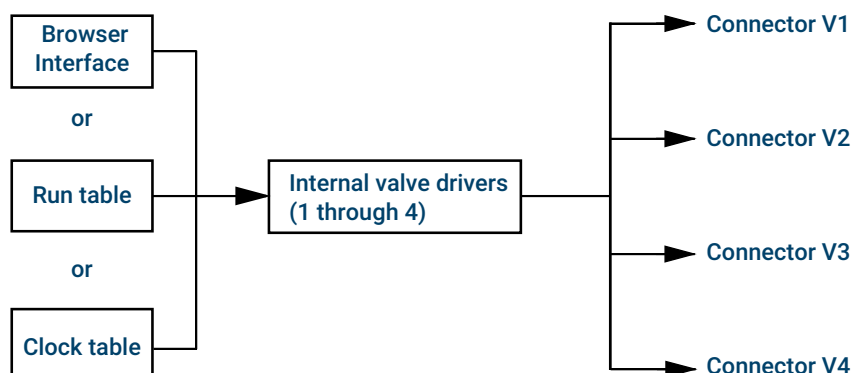
Valve number	Type	Volts	Power or current	Use
1, 2, 3 and 4	Current source	24 VDC	13 watts	Pneumatic valve control

Valve number	Type	Volts	Power or current	Use
9 and 10	Current source	24 VDC	100 mA	Relays and low-power devices
7 and 8	Contact closure	48 VDC or 48 VAC RMS		Control an external current source

The internal valve drivers

Valve drivers 1 through 6 are usually used to control pneumatically operated valves mounted in the valve box. The wiring for these appears at a set of connectors inside the right cover of the GC.

Pneumatically driven valves are controlled by solenoids mounted near the connectors that control the flow of air to the valve actuators.



There is no direct relationship between the location of a valve in the valve box and the driver that controls it. This depends on how the solenoids are wired and the actuators are plumbed.

Manual valves must be switched by hand, and are heated or unheated.

Valve types

The possible valve types are:

Sampling

A two-position (load and inject) valve. In load position, an external sample stream flows through an attached (gas sampling) or internal (liquid sampling) loop and out to waste. In inject position, the filled sampling loop is inserted into the carrier gas stream. When the valve switches from **Load** to **Inject**, it starts a run if one is not already in progress. See the example on [Controlling a valve](#).

Switching

A two-position valve with four, six, or more ports. These are general-purpose valves used for such tasks as column selection, column isolation, and many others.

Other

Could be anything.

Not installed

There is no valve installed in this position.

Controlling a valve

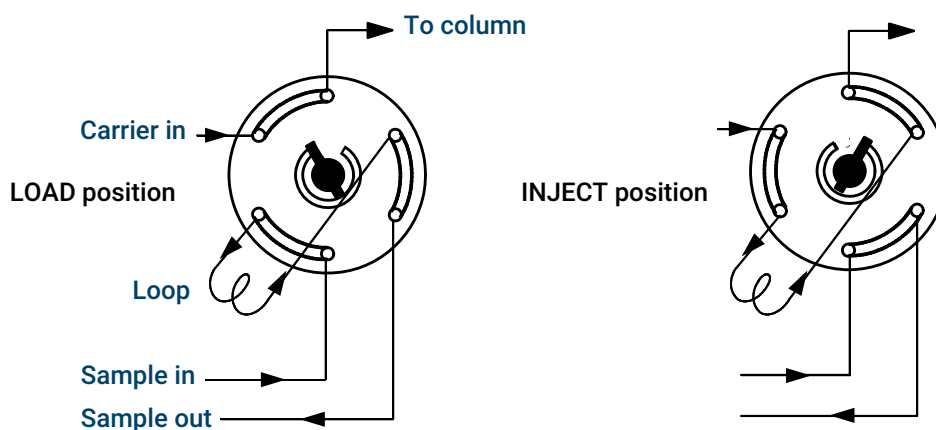
From the run or clock time tables

The **Valve On** and **Valve Off** commands can be run time or clock time programmed. See **“Events”** and **“Clock Time Programming”**.

If a valve is rotated by a run time program, it is not automatically returned to its initial position at the end of the run. You must program this reset operation yourself.

Sampling valve

If a valve is configured as a sampling valve, it starts a run automatically when it is switched to the Inject position. This can be done by a subsequence or clock table entry. You may have two gas sampling valves installed.



Sampling valves have two positions:

Load position

The loop (external for gas sampling, internal for liquid sampling) is flushed with a stream of the sample gas. The column is flushed with carrier gas.

Inject position

The filled loop is inserted into the carrier gas stream. The sample is flushed onto the column. The run starts automatically.

Carrier gas may be provided by an (optional) PCM channel. To do this, configure the column and specify the PCM channel as the inlet. The channel then becomes programmable with four operating modes.

The gas sampling valve control parameters are:

Load time

Time in minutes that the valve remains in the Load position before becoming ready.

Inject time

Time in minutes that the valve remains in the Inject position before returning to the Load position.

The sampling valve cycle is:

1. The sampling valve rotates to the Load position. **Load time** begins. Valve is not ready.
2. **Load time** ends. The valve becomes ready.
3. If everything else is ready, the GC becomes ready. If anything is not ready:
 - If you are using Clock Table or sequence control, the GC waits until everything is ready, then executes the valve inject command.
 - If you are not using Clock Table or sequence control, the valve injection can be made at any time from the keyboard.
4. The sampling valve rotates (keyboard command or sequence control) to the Inject position. **Inject time** begins. The run begins.
5. **Inject time** ends. Return to step 1.

GC Output Signals

Signal is the GC output to a data handling device, analog or digital. It can be a detector output or the output from flow, temperature, or pressure sensors. One signal output channel is provided.

Signal output can be either analog or digital, depending on your data handling device. Analog output is available at either of two speeds, suitable to peaks with minimum widths of 0.004 minutes (fast data rate) or 0.01 minutes (normal rate). Analog output ranges are 0 to 1 V, 0 to 10 V.

Digital output rates are set by your Agilent data system, such as OpenLab CDS or MassHunter Workstation.

See **Table 10** for the conversions from units shown on the GC display to units as shown in Agilent data systems and integrators.

Table 10. Signal conversions

Signal type	1 display unit is equivalent to:
Detector:	
FID, NPD	1.0 pA (1.0×10^{-12} A)
FPD+	150 pA (150×10^{-12} A)
TCD	25 μ V (2.5×10^{-5} V)
ECD	1 Hz
Analog input board (use to connect the GC to non-Agilent detector)	15 μ V
Nondetector:	
Thermal	1 °C
Pneumatic:	
Flow	1 mL/min
Pressure	1 pressure unit (psi, bar, or kPa)
Diagnostic	Mixed, some unscaled

When outputting a column pressure signal, the GC reports the pressure in absolute units. For example, an inlet pressure of 68.9 kPa would be reported as 170.2 kPa.

Analog signals

If you use an analog recorder, you may need to adjust the signal to make it more usable. Zero and Range in the Signal parameter list do this.

Analog zero

Zero

Subtracts value entered from baseline. Either select **On** to set Zero at the current signal value, or enter a number between -500,000 and +500,000 as the setpoint to subtract from the baseline.

This is used to correct baseline elevation or offsets. A common application is to correct a baseline shift that occurs as the result of a valve operation. After zeroing, the analog output signal is equal to the **Value** line of the parameter list minus the **Zero** setpoint.

Zero can be programmed as a run time event. For details, see [Events](#).

Analog range

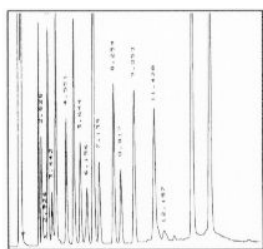
Range

Scales data coming from the detector

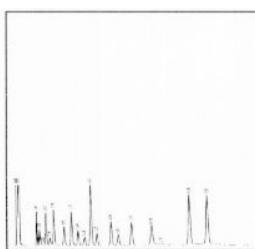
Range is also referred to as gain, scaling, or sizing. It sizes the data coming from the detector to the analog signal circuits to avoid overloading the circuits (clamping). **Range** scales all analog signals.

If a chromatogram looks like A or B in the next figure, the data needs to be scaled (as in C) so that all peaks are visible on the paper.

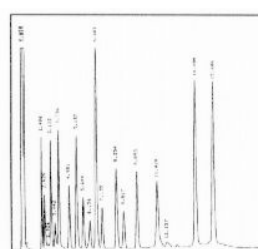
Valid setpoints are from 0 to 13 and represent 2^0 (=1) to 2^{13} (=8192). Changing a setpoint by 1 changes the height of the chromatogram by a factor of 2. The following chromatograms illustrate this. Use the smallest possible value to minimize integration error.



A: Range = 0



B: Range = 3



C: Range = 1

There are limits to usable range settings for some detectors. The table lists the valid range setpoints by detector.

Table 11. Range limits

Detector	Usable range settings (2x)
FID	0 to 13
NPD	0 to 13
FPD+	0 to 13
TCD	0 to 6
ECD	0 to 6
Analog input	0 to 7

Range may be run time programmed. See [Events](#) for details.

Analog data rates

Your integrator or recorder must be fast enough to process data coming from the GC. If it cannot keep up with the GC, the data may be damaged. This usually shows up as broadened peaks and loss of resolution.

Speed is measured in terms of bandwidth. Your recorder or integrator should have a bandwidth twice that of the signal you are measuring.

The GC allows you to operate at two speeds. The faster speed allows minimum peak widths of 0.004 minutes (8 Hz bandwidth), while the standard speed allows minimum peak widths of 0.01 minutes (1.6 Hz bandwidth).

If you use the fast peaks feature, your integrator should operate at around 15 Hz.

Selecting fast peaks (analog output)

1. Select **Settings > Configuration**.
2. Select **Analog Out**.
3. Select the checkbox next to **Fast Peaks**.

Agilent does not recommend using **Fast Peaks** with a thermal conductivity detector. Since the gas streams switch at 5 Hz, the gain in peak width is offset by increased noise.

Digital signals

The GC outputs digital signals only to an Agilent data system. The following discussions describe features that impact the data sent to data systems, not the analog data available to integrators. Access these features from the data system. These features are not accessible from the GC touchscreen or Browser Interface.

Zero signal

Available only from an Agilent data system.

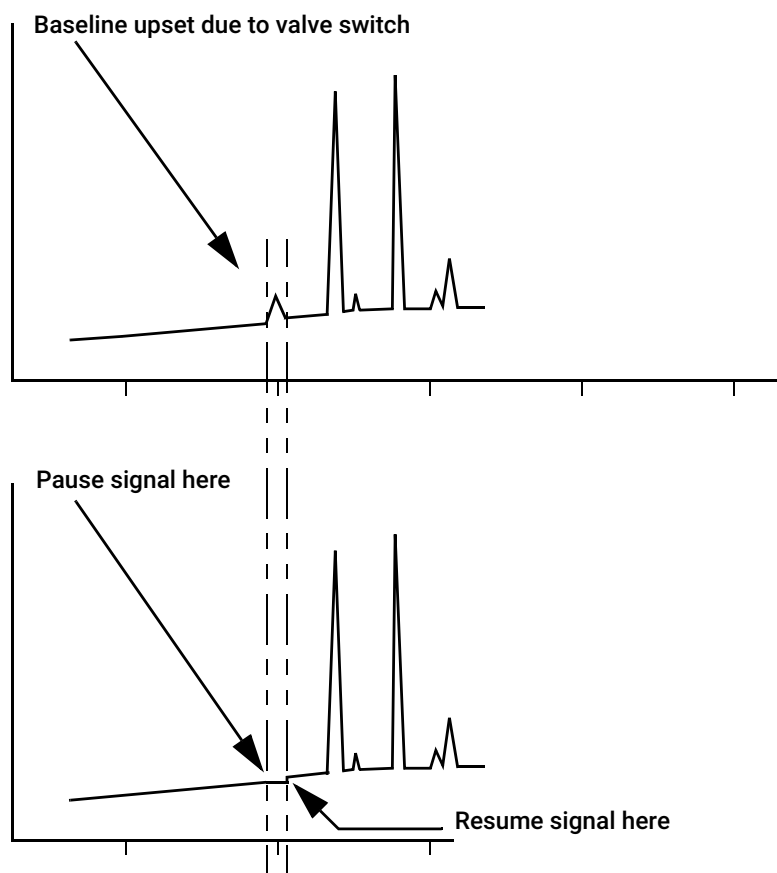
Digital signal outputs respond to a zero command by subtracting the signal level at the time of the command from all future values.

Signal Freeze and Resume

Available only from an Agilent data system.

Some run time operations, such as changing signal assignments or switching a valve, can cause baseline upsets. Other factors can cause baseline upsets also. The GC can compensate for this by pausing (freezing) the signal at a particular value, using that signal value for a specified duration, and then resuming normal signal output.

Consider a system that uses a switching valve. When the valve switches, an anomaly occurs in the baseline. By freezing and resuming the signal, the anomaly can be removed so that the peak identification and integration software operates more smoothly.



Data rates with Agilent data systems

The GC can process data at various data rates, each corresponding to a minimum peak width. The table shows the effect of data rate selection.

Table 12. Agilent data system data processing

Data rate, Hz	Minimum peak width, minutes	Relative noise	Detector	Column type
1000	0.0002	6.96	NPD	Narrow-bore, 0.05 mm
500	0.0004	5	FID/NPD	Narrow-bore, 0.05 mm
200	0.001	3.1	FID/FPD+/NPD	Narrow-bore, 0.05 mm
100	0.002	2.2	FID/FPD+/NPD	Capillary
50	0.004	1.6	ECD/FID/FPD+/NPD	Capillary
20	0.01	1	ECD/FID/FPD+/NPD	Capillary
10	0.02	0.7	ECD/FID/FPD+/NPD	Capillary
5	0.04	0.5	ECD/FID/FPD+/NPD/TCD	Capillary
2	0.1	0.3	ECD	Capillary
1	0.2	0.22	ECD	Capillary
0.5	0.4	0.16	ECD	Capillary
0.2	1.0	0.10	ECD	Capillary
0.1	2.0	0.07	ECD	Slow packed

Methods

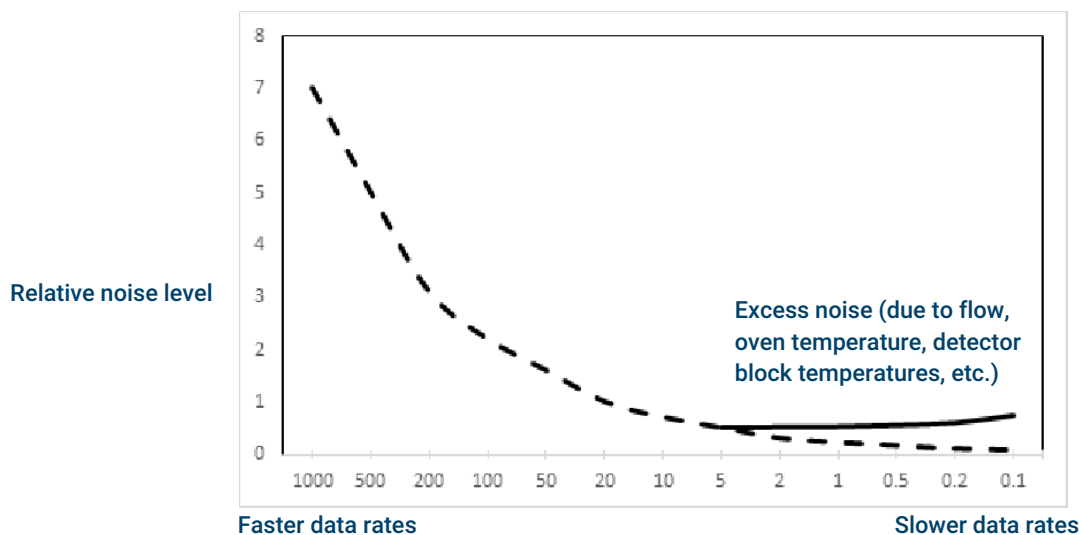
GC Output Signals

You cannot change the data rate during a run.

You will see higher relative noise at the faster sampling rates. Doubling the data rate can double peak height while the relative noise increases by 40%. Although noise increases, the signal-to-noise ratio is better at the faster rates.

This benefit only occurs if the original rate was too low, leading to peak broadening and reduced resolution. We suggest that rates be chosen so that the product of data rate and peak width in seconds is about 10 to 20.

The figure shows the relationship between relative noise and data rates. Noise decreases as the data rate decreases until you get to data rates of around 5 Hz. As the sampling rate slows, other factors such as thermal noise increase noise levels.





6

Sequences

What is a Sequence?

Recoverable Errors

What is a Sequence?

A sequence is a list of samples to be analyzed along with the method to be used for each analysis. Sequences can be set up from the Browser Interface, or from the connected Agilent data system. Refer to the help provided in the Browser Interface or data system for details.

Recoverable Errors

For details on how this feature works in your data system, refer to its help and documentation.

Some types of errors, such as an ALS missing vial error or a headspace sampler vial size mismatch, may not always justify stopping an entire sequence. These errors are called recoverable errors, since you may be able to recover from them and continue running a sequence, if desired. Agilent data systems now provide features to allow you to control how the system will react to these types of errors. When using an Agilent data system, the data system will now control whether or not the sequence pauses, aborts completely, continues with the next sample, and so on, for each type of recoverable error.

Note that the data system only controls what happens to the next run in the sequence, not the *current* run, except when set to immediately abort. (In that case, the data system typically aborts the current run and the sequence.)

For example, pressing the "Stop" button on the GC always halts the current run. However, the data systems can allow you to choose whether to continue with the next run or to pause or to abort the whole sequence.

7

Diagnostics

About Diagnostics

Automated testing

The system health report

Self-guided diagnostics

Gather Log Files

Warnings and Errors

Using the Diagnostics View

Performing Diagnostic Tests

About Diagnostics

The GC provides diagnostics features for inlets, detectors, and other installed components. This includes tests that are performed by the operator and automated testing that is performed by the GC without operator intervention.

The Diagnostics view provides access to current warnings and errors, diagnostic tests, the system health report, blank evaluation reports, system checkout reports, and gas and power usage statistics.

Select a link to access the desired reports or tests.

To view active warning and error conditions, select **Warnings and Errors**. This panel displays the list of conditions that require attention. If the GC has not identified any conditions that require attention, the panel will be blank. Touch a condition to see a list of possible causes and resolutions.

Automated testing

The GC performs continuous, automated testing of the following items. If a failure occurs, an alert appears on the diagnostic tab. In addition, an entry is made in the appropriate log.

See below for a list of the continuously monitored signals and states:

Detector:

- Supply voltage
- ADC (analog to digital converter) references
- FID flameout
- NPD bead open/short
- Igniter open/short
- Collector short
- Electrometer not plugged in
- TCD filament open or short
- TCD switching valve open or short
- FPD flameout

EPC (Electronic Pneumatic Control) - Inlets, detectors, and other modules.

ADC (analog to digital converter) references

Actuator movements

Thermal:

- Sensor open/short
- Missing heater
- Wrong heater

Diagnostics

About Diagnostics

- Heater current:
 - Quiescent
 - Leakage

Configuration Mismatch

The system health report

To access the system health report, select **System Health Report**. The system health report appears.

The system health report includes the following types of information:

- System information
- System Configuration details
- Active instrument conditions
- Column details
- Early Maintenance Feedback details
- Diagnostic test results
- Network information
- Status snapshot information

Self-guided diagnostics

The GC offers several useful diagnostic checks to use when troubleshooting issues with the GC or method. See below for a list of self-guided diagnostic tests available:

Inlet tests:

- Leak & Restriction test
- Gas Supply Pressure test
- Split Vent Restriction test
- Pressure Decay test
- Septum Purge test
- Gas Identification test

Detector tests:

- FID Jet Restriction test
- FID Leakage Current test
- FPD Dark Current check

Other tests:

- Hydrogen sensor calibration check. If the optional hydrogen sensor is installed, this test verifies sensor operation.

Gather Log Files

The GC provides many diagnostic tools to automatically diagnose itself. If you need to contact Agilent for support, please use the convenient Gather Logs feature (found in the browser interface on the Diagnostics tab, and also available in the data system) and provide the collected information when you contact Agilent support. Gather Logs will bundle all relevant logs files and data in a convenient package file. This information will be used to diagnose and resolve any issues. Information gathered includes GC and PC time information, the system logs, and firmware revisions. You can also include setpoints, instrument configuration, the last five runs of sample information, and the last five data runs.

From an Agilent data system, find Gather Logs on the Maintenance Actions tab when accessing the GC method parameters. The data system command will also include the data system version and related basic details.

Log options

In addition to the normal logs (run log, system log, maintenance log, and so forth), optionally include the active method, run information log, run data logs, and the trend analysis database.

When including the trend analysis database, the log file will include EMF, diagnostic, system check-out, and blank evaluation data. Collecting this information will increase the time needed to generate the log file.

Available Logs

Use the selection list in Available Logs to search for, load, and delete existing log files as desired.

Warnings and Errors

When the GC determines a condition that requires attention, it lists the condition in **Diagnostics > Warnings and Errors**. Select a condition to see a list of possible causes and resolutions. A blank screen means that the GC has not identified any issues that require attention.

< Diagnostics		Warnings and Errors	?
	MPV: Missing	2023-02-02 23:23:08	
	Front Inlet: Invalid PIDs	2023-02-02 23:23:08	
	Instrument: Unsupported Electronics Detected	2023-02-02 23:23:08	
	Instrument: Power Supply Communication Fault	2023-02-02 23:23:08	

Figure 18. Example listing of warning and error conditions

Using the Diagnostics View

To use the Diagnostics view:

1. Select **Diagnostics**. The Diagnostics view appears.
2. Select **Diagnostic Tests**. The Diagnostic Tests page appears.
3. Select the desired test from the list.

Performing Diagnostic Tests

To perform a diagnostic test:

1. Access the desired test from the Diagnostics view. See [Using the Diagnostics View](#).
2. Select the desired test. The corresponding test page appears. The test page includes a test description and an indication of the parameter being tested.
3. Select **Start Test**. Testing is initiated. Test details are displayed along with test results.

The currently running test can be aborted by selecting **Cancel**. This brings up a dialog box which allows you to confirm that you want to cancel the test.

8 Maintenance

Automated Maintenance

Early Maintenance Feedback (EMF)

Counter types

Default Thresholds

Available Counters

Viewing Maintenance Counters

To Enable, Reset, or Change a Limit for an EMF Counter

EMF Counters for Columns

EMF Counters for Autosamplers

Counters for 7693A and 7650 ALS with EMF-enabled firmware

Counters for ALS with earlier firmware

EMF Counters for MS Instruments

Selecting the maintenance navigation tab shows buttons for counters by component, lets you view the maintenance log, and lets you start automated, step-by-step maintenance procedures for various components. These procedures also include automated tests as needed to help- be sure the GC is ready when maintenance is complete. See the GC's *Maintaining Your GC* manual for more information.

Automated Maintenance

For many common maintenance procedures, your GC contains a detailed walkthrough of each procedure. For all GC models, these procedures may be accessed by selecting **Maintenance > Perform Maintenance**, or by selecting **Maintenance > your desired component > Perform Maintenance**. Once here, select your desired maintenance procedure, and select **Start Maintenance** to begin the walkthrough.

Automated maintenance procedures not only step you through performing the maintenance task, but also prepare the GC, perform checks at appropriate times, and provide videos to assist you in performing hardware tasks.

Many non-automated maintenance procedures require you to put the GC in maintenance mode prior to performing any maintenance. To do so, select **Maintenance > Instrument > Perform Maintenance**, select the checkbox next to **Maintenance Mode**, and then select **Start Maintenance**.

Putting the GC into a maintenance mode can include:

- Setting low temperatures to avoid burns and other injuries
- Setting reduced flows to avoid safety hazards and to prevent damage to the instrument
- Venting a mass selective detector (MSD)
- Making other settings to prevent damage to the instrument (electronics, columns, and so forth) or to connected instruments (MSD)

For example, to change the gas clean filter, select **Maintenance > Instrument > Perform Maintenance**, select the checkbox next to Gas Clean Filter Maintenance, and then select **Start Maintenance**. This cools down the required components of the GC, and then walks you through the procedure for replacing the gas clean filter.

When performing an automated maintenance procedure:

- The GC status indicator turns yellow
- The touchscreen status bar color changes
- The GC is Not Ready

STATUS: MAINTENANCE MODE

To exit maintenance mode, either complete the running maintenance procedure or exit/abort it. Maintenance procedures also end when the GC is turned off. If aborting a running maintenance procedure, make sure that all GC parts and consumables have been installed, configured, and so forth.

Early Maintenance Feedback (EMF)

The GC provides injection- and time-based counters for various consumable and maintenance parts. Use these counters to track usage and replace or recondition these items before potential degradation impacts the chromatographic results.

If using an Agilent data system, these counters can be set and reset from within the data system.

Counter types

Counters are provided for injections, runs, and time. Each type is described below.

Injection counters increment whenever an injection occurs on the GC via an ALS injector, head-space sampler, or sampling valve. Manual injections do not increment counters. The GC differentiates between front and back injections, and only increments counters associated with the configured injection flow path.

For example, consider the following GC:

Configured front flow path	Configured back flow path
Front injector	Back injector
Front inlet	Back inlet
Column 1 (GC oven)	Column 2 (GC oven)
Purged Union / Aux EPC 1	Back detector
Column 3 (GC oven)	
Front detector	

In this example, for a front ALS injection, the GC will increment counters for the front injector, front inlet, and the front detector, but it would not increment the back injector, back inlet, or back detector counters. For the columns, the GC would increment the injection counters for columns 1 and 3, and the oven cycles counter for all 3 columns.

Run counters increment against the number of runs performed on the GC.

Time counters increment against the GC clock. Changing the GC clock changes the age of tracked consumables.

Default Thresholds

Selected counters have default thresholds to use as a starting point.

If you want to change a default limit, enter a conservative limit based on your experience. Use the warning feature to alert you when service is approaching, then track performance to determine if the **Service Due** threshold is too high or too low.

For any EMF counter, you may need to adjust the threshold values based on the demands of your applications.

Available Counters

Table 13 lists the most common counters available. The available counters will vary based on the installed GC options, consumables, and future updates.

Table 13. Common EMF counters

GC Component	Parts with a counter	Type	Default value
Detectors			
FID	Collector	Number of injections	
	Jet	Number of injections	
	Ignitor	Number of ignition attempts	
TCD	Switching solenoid	On time	
	Filament on time	On time	
ECD	Insert liner	Number of injections	
	Time since wipe test	On time	6 months
NPD	Bead	Number of injections	
	Collector	Number of injections	
	Bead baseline offset	pA Value	
	Bead baseline voltage	Voltage value	Blos bead: 1.045
	Bead current integral	pA-sec Value	
FPD+	Bead on time	On time	Blos bead: 2400 h
	Ignitor	Number of ignition attempts	
	PMT	Number of injections	
	PMT	On time	6 months
Inlets			
SSL	Gold seal	Number of injections	5000
	Gold seal	Time	90 days
	Liner	Number of injections	200
	Liner	Time	30 days
	Liner O-ring	Number of injections	1000
	Liner O-ring	Time	60 days
	Septum	Number of injections	200
	Split vent trap	Number of injections	10,000
	Split vent trap	Time	6 months
PP	Liner	Number of injections	200
	Liner	Time	30 days
	Septum	Number of injections	200
	Top weldment O-ring	Number of injections	10,000
	Top weldment O-ring	Time	1 year
COC	Septum	Number of injections	200
Columns			
Column	Injections onto column	Number of injections	
	Oven cycles	Number of injections	
	Length	Value	
	Run count	Number of runs	

Maintenance

Available Counters

GC Component	Parts with a counter	Type	Default value
Valves			
Valve	Rotor	Activations (number of injections)	
	Maximum temperature	Value	
Instrument			
Instrument	On time	Time	
	Run count	Number of runs	
	Filters	Time	
Hydrogen sensor (optional)			
Hydrogen sensor	On time since calibration	Time	6 months
ALS Injectors			
ALS	Syringe	Number of injections	800
	Syringe	Time	2 months
	Needle	Number of injections	800
	Plunger moves	Value	6000
Mass spectrometers			
Mass spectrometer	Pump	Time (days)	1 year
	Filament 1	Time (days)	1 year
	Filament 2	Time (days)	1 year
	Source (time since last cleaning)	Time (days)	1 year
	EMV at last tune	V (volts)	2600

Viewing Maintenance Counters

To view the maintenance counters:

1. Select the **Maintenance** tab.
2. Select the desired component type. The Status column lists the counter for the corresponding component.
3. Scroll to view additional components, as applicable.

To Enable, Reset, or Change a Limit for an EMF Counter

When using the GC without a data system, enable or change the limit for a counter as follows:

1. Locate the counter you wish to change. See [Viewing Maintenance Counters](#).
2. Select the **Details** listing for the counter you want to change. The dialog includes a trend plot for past reset events for this counter.
The trend plot legend includes the following controls:
 - **User Reset**: Select to toggle manual counter resets.
 - **Maintenance Reset**: Select to toggle counter resets performed automatically when using an automated maintenance procedure.
 - **Service Due**: Select to toggle the Service Due limit on the plot.
 - **Service Warning**: Select to toggle the Service Warning limit on the plot.
3. To modify a threshold:
 - a. Select the threshold entry. A data entry dialog box appears.
 - b. Enter the desired value. See [Default Thresholds](#).
4. To enable or disable a warning, select or deselect **Enable** for the corresponding counter.
5. Select **Apply**. The dialog box closes. The entered value is displayed in the corresponding field.
6. To reset the counter:
 - a. Select **Reset Counter**. A confirmation dialog box appears.
 - b. Select **Yes**. The confirmation dialog box closes.
7. Select **Apply**.

The GC logs each counter reset. Access this data from the maintenance log.

Note that any change to a **Service Due** or **Service Warning** limit does not appear on the plot until the next counter reset.

EMF Counters for Columns

Because GC columns are sometimes transferred from GC to GC or from front to back on the same GC, the EMF counters for GC columns are associated not with the GC, but with the column's serial number. This serial number is entered during column configuration or is found in the columns database if using an Agilent data system.

- If the column is reconfigured on the same GC, for example, changed from column #1 to column #2, the GC will transfer the EMF data to the new column position, and will continue to cumulatively track EMF data.
- If a column is completely removed from the GC (the current GC configuration no longer includes that column serial number), the GC will track the EMF data for any new column using the new column's data.
- When applicable, the GC stores column EMF data in the columns database. The data in the columns database do not get reset by the GC.

Column EMF data will not transfer between GCs unless using an Agilent data system with a columns database. If using an Agilent data system, column EMF data may not transfer between data system installations (each data system usually has its own columns database).

EMF Counters for Autosamplers

The GC provides access to the counters for the autosampler. The functionality for ALS counters depends on the ALS model and firmware version. In all cases, the GC shows EMF counter status and allows you to enable, disable, and reset the counters using the Browser Interface.

Counters for 7693A and 7650 ALS with EMF-enabled firmware

If using an Agilent 7693 injector with firmware version G4513A.10.8 (or higher), or a 7650 injector with firmware version G4567A.10.2 (or higher), each injector independently tracks its EMF counters.

- The injector counters will increment as long as the injector is used on an 8860B Series GC. You can change positions on the same GC or install the injector on a different GC without losing the current ALS counter data.
- The ALS will report an exceeded limit only when mounted on an 8860B Series GC.

Counters for ALS with earlier firmware

If using a 7693 or 7650 injector with earlier firmware, or if using another injector model, the GC tracks the counters for that injector. The GC uses the injector serial number to distinguish between installed injectors, but only maintains up to two sets of counters—one for the front injector, and one for the back injector.

- The GC will track injector counters regardless of installed position (front or back inlet). Because the GC tracks the injector serial number, you can change the injector position without losing the counters as long as the injector remains installed on the GC.
- Each time the GC detects a new injector (different model or different serial number), the GC resets the ALS counters in the new injector's position.
- In this case, ALS EMF counters are not shared between GCs.

EMF Counters for MS Instruments

For GC-MS, GC-HS, and GC-MS-HS systems, all EMF counters are available at the Browser Interface. In addition, most counters can be reset at the Browser Interface. Some counter types, for example, a counter which requires a calibration to be performed on a headspace sampler, cannot be reset at the Browser Interface, but can still be viewed.

9

Logs

Logs View

Maintenance logs

Run log

System log

This section describes the log features available on the Agilent GC.

Logs View

The Logs view provides listings of GC events including maintenance events, run events, system events, sequence events, run history, and logs for certain connected instruments (such as an 8697 headspace sampler), sorted by date/time.

Selecting one of the buttons on the Logs view brings up the corresponding log page. .

For Maintenance and System log items, items are sorted by date and time. For Run log items, relative time (from the start of the run) is used.

Use the scroll buttons to scroll through the log entries.

Select **Cancel** to return to the Logs view.

Maintenance logs

The maintenance log contains entries made by the system when:

- A system event occurs (for example, a detector shutdown)
- Any component counter reaches a monitored limit

The log entry contains a description of the maintenance event and the date/time that the event occurred. In addition, each user task related to the counter is recorded in the log, including resetting, enabling, or disabling monitoring, and changing limits or units (cycles or duration).

Run log

The run log is cleared at the start of each new run. During the run, any deviations from the planned method (including touchscreen, control button, or Browser Interface intervention) are listed in the run log table.

System log

The system log records significant events that occur during GC operation. Some of the events also appear in the run log if they occur during a run.

Sequence log

The Sequence Log record events associated with browser interface sequences (not data system sequences).

Run history log

The Run History Log records general run history, including runs started from connected data systems.

10

Settings

About Settings

Service Mode

About

Calibration

Maintaining EPC calibration—inlets, detectors, PCM, PSD, and AUX

To zero a specific flow or pressure sensor

System Settings

Configuring the IP address for the GC

To change the system locale

To access stored run data

To control Browser Interface access

Tools (Browser Interface)

Performing a column compensation run

About Settings

The Settings view provides access to configuration and system settings for the GC.

- Select **Scheduler** to access the GC Instrument Schedule settings. See [Resource Conservation](#).
- Select **Service Mode** to access the service mode settings for the GC. See [Service Mode](#).
- Select **About** to get information about this GC. See [About](#).
- Select **Calibration** to access the calibration functions. See [Calibration](#).
- Select **System Settings** to access the system settings for the GC, including setting the network address, system date and time, touchscreen settings, system setup information, and so on. See [System Settings](#).
- Select **Tools** to access the Tools page. See [Tools \(Browser Interface\)](#).

Service Mode

The Service Mode feature allows you to view the installed GC system component details. This includes serial numbers, firmware versions, voltages, currents, temperatures, and so on.

To view the specifications for the various components of your instrument:

1. Select the desired component type. The Service Mode page for the selected component appears.
2. Use the page selection buttons on the left side of the page to display related functional information.

About

The About feature allows you to view details about the GC.

The About screen lists the GC manufacturing date, serial number, firmware revision, and help and information revision.

Select **Close** on the About page to return to the Settings view.

Calibration

Calibration allows you to adjust the following items (when available):

- ALS
- Inlets
- Oven
- Detectors
- EPC modules
- Hydrogen sensor (optional). [Hydrogen Sensor](#)

To change the calibration settings:

1. Use the page selection buttons on the left side of the page to display related functional information.
2. Make changes to calibration settings as desired. See [Maintaining EPC calibration—inlets, detectors, PCM, and AUX, To zero a specific flow or pressure sensor](#) for more information.
3. Select **Apply**. The entered changes are saved to the GC.

Maintaining EPC calibration—inlets, detectors, PCM, and AUX

The EPC gas control modules contain flow and/or pressure sensors that are calibrated at the factory. Sensitivity (slope of the curve) is quite stable, but zero offset requires periodic updating.

Flow sensors

All inlet modules use flow sensors, as do PSDs, and channel 1 of a PCM. If the **Auto flow zero** feature is selected, they are zeroed automatically after each run. This is the recommended setting. They can also be zeroed manually—see [To zero a specific flow or pressure sensor](#).

Pressure sensors

All EPC control modules use pressure sensors. They must be zeroed individually. There is no automatic zero for pressure sensors.

Auto Zero flow

A useful calibration option is **Auto Zero flow**. When selected, after the end of a run, the GC shuts down the flow of gases to an inlet, waits for the flow to drop to zero, measures and stores the flow sensor output, and turns the gas back on. This takes about two seconds. The zero offset is used to correct future flow measurements.

Auto zero septum purge

This is similar to **Auto Zero flow**, but is for the septum purge flow.

Zero conditions

Flow sensors are zeroed with the carrier gas connected and flowing.

Pressure sensors are zeroed with the supply gas line disconnected from the gas control module.

Zero intervals

Table 14. Flow and Pressure Sensor Zero Intervals

Sensor type	Module type	Zero interval
Flow	All	Use Auto flow zero and/or Auto zero septum purge
Pressure	Inlets	
	Small capillary columns (id $\leq$ 0.32 mm)	Every 12 months
	Large capillary columns (id $>$ 0.32 mm)	At 3 months, at 6 months, then every 12 months
	Auxiliary channels	Every 12 months
	Detector gases	Every 12 months

To zero a specific flow or pressure sensor

1. Select **Settings > Calibration > Detectors**, and select the desired detector.
2. Select **On** next to the desired sensor to zero it.
3. For **Flow sensors**. Verify that the gas is connected and flowing (turned on).
4. For **Pressure sensors**. Disconnect the gas supply line at the back of the GC. Turning it off is not adequate; the valve may leak.
5. Reconnect any gas line disconnected in the previous step and restore operating flows.

System Settings

System settings include setting the network address, system date and time (GC clock), locale setting, and system setup information.

To change the GC clock (date and time), run the system setup wizard. In the Browser Interface, select **System Setup > Start System Setup**. Continue through the wizard until it reaches the date and time setup. Set the GC time and date, then continue through the wizard until it completes.

Select **Apply** to save any changes made to the GC.

Configuring the IP address for the GC

For network (LAN) operation, the GC needs an IP address. It can get this from a DHCP server, or it can be entered directly from the touchscreen. In either case, see your LAN administrator for appropriate settings.

To use a DHCP server

1. From the System Settings page, select the **Network** page selection button. The Network Configuration page appears.
2. For DHCP, select **Enabled**.
3. Select **Apply**.
4. If prompted, restart the GC.

To set the LAN address at the touchscreen

1. Select the System Settings page.
2. If **Enabled DHCP** is selected:
 - a. Deselect **Enabled DHCP**.
 - b. When prompted, restart the GC.
 - c. Return to the System Settings page
3. Enter the **IP Address** in its corresponding field
4. Enter the **Gateway** number in the corresponding field.
5. Enter the subnet mask in the **Net Mask** field.
6. Select **Apply**.
7. If prompted, restart the GC.

To change the system locale

1. From the **Settings** icon, select **MISC** tab to select your locale. See [Figure 19](#).

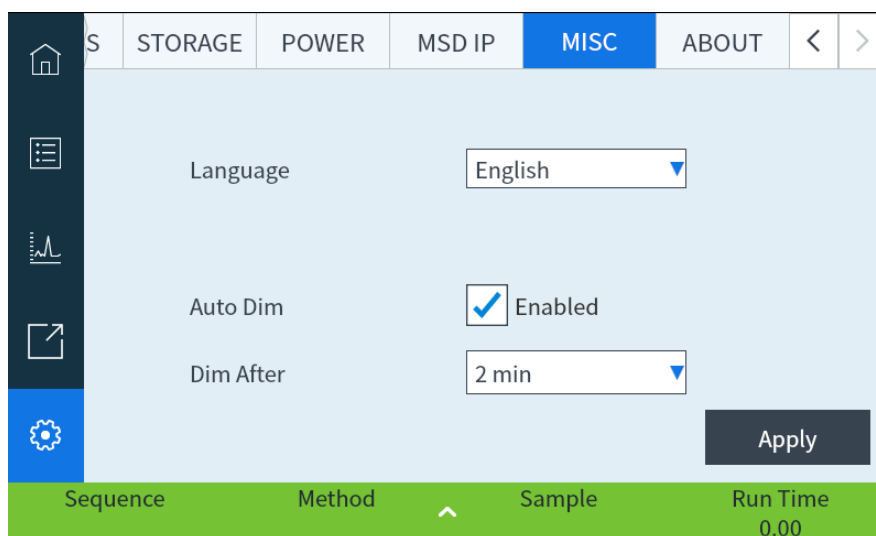


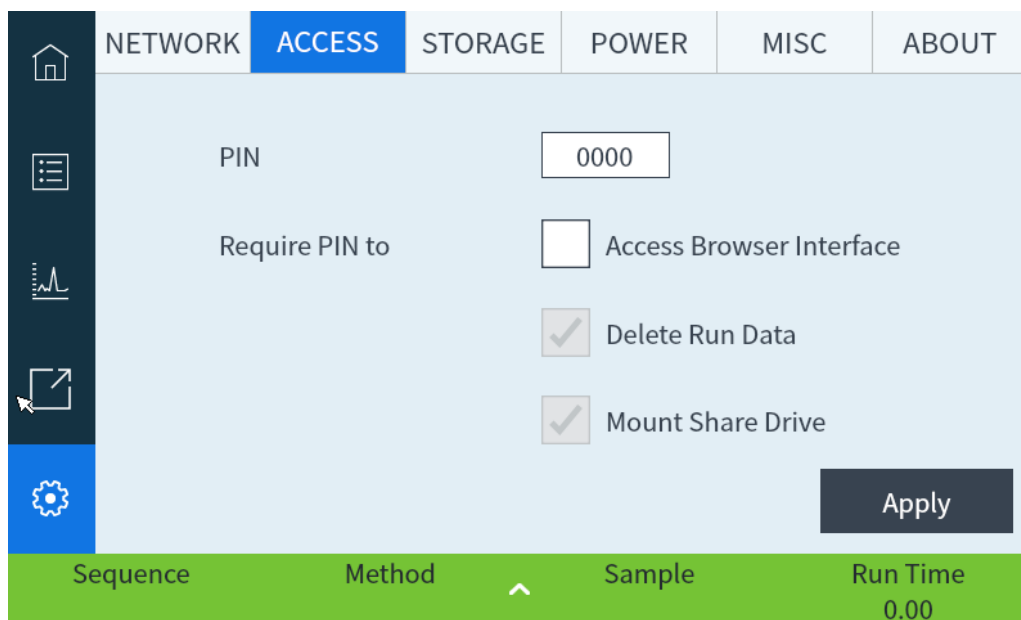
Figure 19. Settings

2. Choose the desired **Language** from the corresponding drop-down list box.
3. Select **Apply**. The GC saves the change made. The system is changed to the selected locale. This may take a few moments.

To access stored run data

If using the Browser Interface to perform runs and collect data, the GC stores the result data internally. To access that data:

1. From the **Settings** page, select the **Access** tab then enter the appropriate PIN.



2. Select the **Storage** tab. Note the path to the GC share.

Settings

System Settings

- On your PC, map a network drive to the GC share. When prompted, connect using the credentials:
user: results
password: the PIN (default: 0000).

To control Browser Interface access

The GC is set so that a four-digit PIN must be used to perform the following actions for your GC:

- Delete run data.
- Mount a share drive.

By default, the PIN is set to 0000. Additionally, you may choose to require the PIN to access the Browser Interface. To set a new PIN:

- From the **Settings** page, select the **Access** tab.
- Select the PIN field and input a new PIN.
- If desired, select the checkbox next to Access Browser Interface to require the PIN for Browser Interface access.

The screenshot shows the 'Settings' application with the 'ACCESS' tab selected. The interface includes a sidebar with icons for home, list, graph, share, and settings. The main content area has tabs for NETWORK, ACCESS (selected), STORAGE, POWER, MISC, and ABOUT. Under the ACCESS tab, there is a 'PIN' field with the value '0000'. Below this is a 'Require PIN to' section with three checkboxes: 'Access Browser Interface' (unchecked), 'Delete Run Data' (checked), and 'Mount Share Drive' (checked). An 'Apply' button is located at the bottom right of the settings area. At the bottom of the screen, a green status bar displays 'Sequence', 'Method', 'Sample', and 'Run Time 0.00'.

Tools (Browser Interface)

The Tools page in the browser interface allows you to perform column compensation runs for the installed columns on the GC.

In temperature programmed analysis, bleed from the column increases as the oven temperature rises. This causes a rising baseline which makes peak detection and integration more difficult. Column compensation corrects for this baseline rise.

A column compensation run is made with no sample injected. The GC collects an array of data points from any installed detectors. If a detector is not installed or is turned off, that part of the array is filled with zeros.

Each array defines a set of curves, one for each detector, that can be subtracted from the real run to produce a flat baseline.

When a connected data system is used, the raw signal, and the column compensation data, is output to the data system so that a compensated, and uncompensated, signal are available for analysis.

Performing a column compensation run

All conditions must be identical in the column compensation run and the real run. The same detector and column must be used, operating under the same temperature and gas flow conditions.

Up to four column compensation runs can be made. The GC retains the results of these runs for later use.

Any column compensation run can then be used to compensate a rising baseline during a run.

1. Go to **Settings > Tools**. With the Tools page displayed, select the desired **Column Compensation** in the **Start Specified Run** column. The GC performs the column compensation run. No injection occurs as a part of this run.
2. Use the connected data system to edit the method. Set the detector to **Subtract from Signal: Column compensation Curve #x** (where **x** is the number of the column compensation run).
3. Run the method. The results use the column compensation run data to compensate for baseline changes in the column.

About Configuration

Configuration Changes

Valve Configuration

Inlet Configuration

Columns

- To configure a single column

- To configure a packed column

- Additional notes on column configuration

Oven

Detector Configuration

Analog Output Settings

MSD Configuration

- MSD configuration

- To enable or disable MS communications

- To use the GC when the MS is shut down

Headspace Sampler (8697)

Valve Box

PCMs

About Configuration

A device's configuration properties are constant for an instrument hardware setup, unlike method settings which can change from sample run to sample run. Two example configuration settings are the gas type flowing through a pneumatic device and the operating temperature limit of a device. The GC will try to set any configuration properties necessary for any device that it finds installed. For example, if the GC knows it is connected to an MSD, it will automatically configure the MSD heated transfer line. The only settings that need to be configured manually are items that can be changed depending on your analysis, such as installed columns and carrier gas type, or an item which may be installed in a non-standard location. For example, an auxiliary EPC module installed in an inlet EPC module position may require manual configuration.

Ensure all devices have been configured correctly prior to running any samples.

Configuration Changes

To change the configuration properties for a device:

1. Select **Method**.
2. Select the desired device type from the list on the left side of the screen. The properties for the selected device type appear on the right side of the screen.
3. Scroll to the device setting and change the property. This can involve making a selection from a list, or entering a numeric value.
4. When all required changes have been made, select the **Save** icon .

Valve Configuration

Valve configuration provides the ability to specify valve types, loop volumes, step times, and BCD inversion settings. BCD inversion allows for changing the BCD input (1's become 0's and 0's become 1's). This feature accommodates coding convention differences between valve manufacturers.

Note that the Valves page appears regardless of whether any valves are currently installed in the GC.

To configure valves

1. Select **Method > Valves**.
2. Use the **Valve Type** drop-down list box to select the corresponding valve type. The possible valve types are:
 - **Gas Sampling** Two-position (load and inject) valve. In load position, an external sample stream flows through an attached (gas sampling) or internal (liquid sampling) loop and out to waste. In inject position, the filled sampling loop is inserted into the carrier gas stream. When the valve switches from Load to Inject, a run starts if one is not already in progress.
 - **Switching** Two-position valve with four, six, or more ports. These are general-purpose valves used for such tasks as column selection, column isolation, and many others.
 - **Other** Something else.
 - **Not installed** Self-explanatory.
3. To enter the loop volume and step time, select **Method > Configuration > Miscellaneous**, and set the values as desired.
4. Select **Apply**. The entered changes are saved to the GC.
5. When all required changes have been made, select the Save icon .

Inlet Configuration

1. Select **Method > Inlets >** for your desired inlet. The properties for the selected device type appear on the right side of the screen.
2. Select the desired gas type from the **Carrier Gas Type** drop-down list box.
3. Select **Readiness** from the list of devices on the left, and then click the checkbox next to your inlet to enable its readiness. If Readiness is disabled, the GC becomes ready even if this device has not reached or cannot reach its setpoints.
4. Adjust the settings for the second inlet as necessary.
5. When all required changes have been made, select the Save icon .

Columns

Length

The length, in meters, of a capillary column. Enter 0 for a packed column or if the length is not known.

Diameter

The inside diameter, in millimeters, of a capillary column. Enter 0 for a packed column.

Film thickness

The thickness, in microns, of the stationary phase for capillary columns.

Inlet

Identifies the source of gas for the column.

Outlet

Identifies the device into which the column effluent flows.

Thermal zone

Identifies the device that controls the temperature of the column.

To configure a single column

You define a capillary column by entering its length, diameter, and film thickness. You then enter the device controlling the pressure at the Inlet (end of the column), the device controlling the pressure at the column Outlet, and the Thermal zone that controls its temperature.

With this information, the instrument can calculate the flow through the column. This has great advantages when using capillary columns because it becomes possible to:

- Enter split ratios directly and have the instrument calculate and set the appropriate flow rates.
- Enter flow rate or head pressure or average linear velocity. The instrument calculates the pressure needed to achieve the flow rate or velocity, sets that, and reports all three values. (Split/splitless inlet only.)
- Perform splitless injections with no need to measure gas flows.
- Choose any column mode. Except for the simplest configurations, such as a column connected to a specific inlet and detector, we recommend that you begin by making a sketch of how the column will be connected.

To configure a capillary column:

1. Select **Method > Configuration > Columns**.
2. Double click the column you want to configure.
3. Under Column Type, select **Capillary**.

4. Input the Length, Diameter, and Film Thickness into the appropriate fields.
If you do not know the column dimensions—they are usually supplied with the column—or if you do not wish to use the GC calculating features, enter **0** for either **Length** or **Diameter**. The column will be not defined.

NOTE

Agilent recommends always defining capillary columns.

5. Under Inlet connection, select the desired inlet from the dropdown menu. Selections include the installed GC inlets and installed PCM channels.
6. Under Outlet connection, select the desired outlet from the dropdown menu.
 - When a detector is selected, the outlet end of the column is controlled at 0 psig for the FID, TCD, FPD+, NPD, and ECD or vacuum for the MSD.
 - If the column exhausts into a nonstandard detector or environment (neither ambient pressure nor complete vacuum), select Other and enter the outlet pressure.
7. Under Heated By, select the desired thermal zone from the dropdown menu.
8. Set the Min Temp, Max Temp, and Max Program Temp for your column.
9. Set the column's
 - Manufacturer
 - Serial Number
10. Click **OK** to confirm your changes.

To configure a packed column

To configure a packed column:

1. Select **Method > Configuration > Columns**.
2. Double click the column you want to configure.
3. Under Column Type, select **Packed**.
4. Set the Min Temp, Max Temp, and Max Program Temp for your column.
5. Set the column's
 - Manufacturer
 - Serial Number
6. Under Inlet connection, select the desired inlet from the dropdown menu. Selections include the installed GC inlets and installed PCM channels.
7. Under Outlet connection, select the desired outlet from the dropdown menu.
 - When a detector is selected, the outlet end of the column is controlled at 0 psig for the FID, TCD, FPD+, NPD, and ECD or vacuum for the MSD.
 - If the column exhausts into a nonstandard detector or environment (neither ambient pressure nor complete vacuum), select **Other** and enter the outlet pressure.
8. Under **Heated By**, select the desired thermal zone from the dropdown menu.
9. Select **Apply**. This saves the changes to the GC.

Additional notes on column configuration

You should check configurations for all columns to verify that they specify the correct pressure control device at each end. The GC uses this information to determine the flow path of the carrier

gas. Only configure columns that are in current use in your GC's carrier gas flow path. Unused columns configured with the same pressure control device as a column in the current flow path cause incorrect flow results.

It is possible, and sometimes appropriate, to configure both installed columns to the same inlet.

Some pneumatic setpoints change with oven temperature because of changes in gas viscosity. This may confuse users who observe pneumatics setpoints changing when their oven temperature changes. However, the flow condition in the column remains as specified by the column mode (constant flow or pressure, ramped flow or pressure) and the initial setpoint values.

Oven

Readiness

When enabled, the GC will not become ready until the oven reaches its setpoints.

Oven Standby

When enabled, the oven temperature changes to the specified Temperature after the GC has been idle for the entered Time. A usage example would be to set the standby temperature to a relatively high value to keep the system clean between runs (reduce carryover) when the GC may sit idle for long periods.

Maximum temperature

Sets an upper limit to the oven temperature. Used to prevent accidental damage to columns. The range is 70 to 425 °C . See the column manufacturer's recommendations.

To configure the oven

1. Select **Method > Oven**.
2. Use the oven configuration window to set the **Maximum Oven Temperature, Equilibration Time**, and oven Temperatures.

NOTE

If the GC uses cryogenic cooling, it automatically configures the coolant type.
3. When all required changes have been made, select the Save icon .

Detector Configuration

If you are operating with column defined, you have a choice of two makeup gas modes. To set the makeup gas for a detector:

1. Select **Method > Detectors**, then select the detector you wish to configure.
2. If available, select the makeup gas type that is plumbed to the detector. Depending on your detector type and the carrier gas selection, the gas type may be already set for you.
3. Make or enter settings for the detector, as applicable:
 - **Lit Offset.** See **Detectors**.
 - **Readiness.** Enable if using this detector. When disabled, the GC becomes ready even if this device has not reached or cannot reach its setpoints.
 - **Target Offset.** NPD only. Enter the default target offset. The GC will slowly try to reach this target value after the detector flows are read, the temperature is stable, and the dry bead time has elapsed.
4. When all required changes have been made, select the Save icon .

Analog Output Settings

When connecting the GC to an integrator or recorder, the device must be fast enough to process data coming from the GC. If it cannot keep up with the GC, the data may be incorrect or corrupted. This usually shows up as broadened peaks and loss of resolution.

Speed is measured in terms of bandwidth. The recorder or integrator should have a bandwidth twice that of the signal being measured.

The GC can operate at two speeds. The faster speed allows minimum peak widths of 0.004 minutes (8 Hz bandwidth), while the standard speed allows minimum peak widths of 0.01 minutes (1.6 Hz bandwidth).

If using fast peaks feature, the integrator should operate at approximately 15 Hz.

NOTE

Agilent does not recommend using Fast peaks with a TCD detector. Since the gas streams switch at 5 Hz, the gain in peak width is offset by increased noise.

To change the analog output settings:

1. Select **Method > Analog Out**, and then select either **Analog Out 1** or **Analog Out 2**.
2. To use fast peaks for either analog output channel, select the corresponding **Fast Peaks** check box.

MSD Configuration

MSD configuration

The method for configuring a connected MSD varies based on the model of MSD being used.

MSD

To change the MSD settings:

1. From the touchscreen, select the settings page, and then the MSD config tab. See [Figure 20](#).

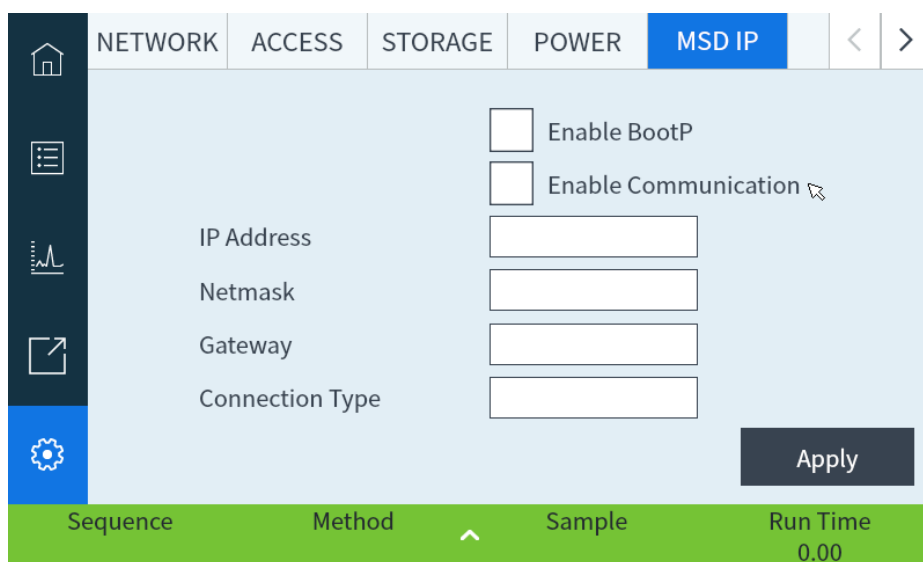


Figure 20. Detector MSD settings page

2. Select **LCOMM** from the dropdown menu under **Connection Type**.
3. Enter details to control the MSD.
4. Select **Apply**. The entered changes are saved to the GC.

To enable or disable MS communications

To temporarily disable GC-MS communications:

1. From the Browser Interface, select **Method > Detectors**, and select your MS. Alternatively, from the touchscreen, select the settings page, and then the **MSD IP** tab.
2. Scroll until the **Enable Communication** checkbox is visible.
3. Select or deselect the checkbox next to **Enable Communication** to enable or disable communication with the MS.

To use the GC when the MS is shut down

To use the GC while an MS is being repaired or maintained, do the following:

Be careful to avoid settings that send carrier gas into the MS, or which raise the temperature of parts which might cause burns when working on the MS.

Configuration**MSD Configuration**

1. Disable MS communications. See [To enable or disable MS communications](#).
2. If needed, completely uninstall the MS from the GC.

Headspace Sampler (8697)

The GC considers the Agilent 8697 headspace sampler to be an integrated part. Once installed, no further setup is needed for communication. To configure the headspace sampler, go to **Method > Configuration > Headspace**. See the 8697 documentation for details.

Readiness

The states of the various hardware components are among the factors that determine whether the GC is Ready for analysis.

Under some circumstances, you may not wish to have a specific component's readiness considered in the GC readiness determination. This parameter lets you make that choice. The following components allow readiness to be ignored: inlets, detectors, the oven, and PCM.

For example, suppose an inlet heater is defective but you don't plan to use that inlet today. By deselecting the **Enable** checkbox under **Readiness** for that inlet, you can use the rest of the GC. After the heater is repaired, reselect the checkbox or the run could start before that inlet's conditions are ready.

To ignore a component's readiness, click **Method > Readiness**. Click the checkbox next to the desired component to remove its checkmark.

To re-enable a component's readiness, click **Method > Readiness**. Click the checkbox next to the desired component to restore its checkmark.

Valve Box

The valve box mounts on top of the column oven. It may contain up to three valves mounted on heated blocks.

Valve positions on the blocks are numbered. We suggest that valves be installed in the blocks in numeric order.

All heated valves in a valve box are controlled by the same temperature setpoint.

To Configure a Valve:

1. Select **Method > Configuration > Miscellaneous**.
2. For each valve installed, select the valve type, then set the loop volume, step time, and BCD inversion as applicable to the valve type.
3. When all required changes have been made, select the Save icon .

PCMs

To configure a PCM:

1. Select **Settings > Configuration > PCMs**.
2. Select the PCMs Aux Mode:
 - Forward Pressure: Pressure is being sensed downstream from the flow proportional valve.
 - Back pressure: Pressure is being sensed upstream from the flow proportional valve.
3. Channel 1 of the PCM provides forward pressure or forward flow control. It can be used to provide column flow or pressure for valved systems, sample preparation devices, or advanced flow devices like splitters or switches.
4. Channel 2, or the auxiliary channel, can only provide forward pressure regulation when plumbed normally, or back pressure regulation when reversed. Thus, channel 2 (reversed) can be used as a controlled leak. If the input pressure drops below setpoint, the regulator can close down to restore it. If the input pressure rises above setpoint, the regulator can open to restore it.

Aux EPCs

An auxiliary pressure controller provides three channels of forward-pressure regulation. Three modules can be installed for a total of nine channels.

The numbering of the channels depends on where the controller is installed. Within a single module, channels are numbered and labeled.

To configure an Aux EPC:

1. Select **Settings > Configuration > Aux EPCs**.
2. For each channel, select the **Gas Type**.



12 Resource Conservation

Resource Conservation

Sleep methods

Wake and condition methods

To Set the GC to Conserve Resources

This section describes the resource-saving features of the GC.

Resource Conservation

The GC provides an instrument schedule to conserve resources such as electricity and gases. Using the instrument schedule, you can assign sleep, wake, and conditioning methods that allow you to program resource usage. A Sleep method sets low flows and temperatures. A Wake method sets new flows and temperatures, typically to restore operating conditions. A Condition method sets flows and temperatures for a specific run time, typically high enough to clean out contamination if present.

Load the sleep method at a specified time during the day to reduce flows and temperatures. Load the wake or condition method to restore analytical settings before operating the GC again. For example, load the sleep method at the end of each day or work week, then load the wake or condition method an hour or so before arriving to work the next day.

Sleep methods

Use a connected data system to create a sleep method to reduce gas and power usage during times of reduced activity.

When creating a sleep method, consider the following:

- The detector. While you can reduce temperatures and gas usage, consider the stabilization time required to prepare the detector for use. The power savings is minimal.
- Connected devices. If connected to an external device such as a mass spectrometer, set compatible flows and temperatures.
- The inlets. Maintain sufficient flow to prevent contamination.

See [Table 15](#) for general recommendations.

Table 15. Sleep method recommendations

GC Component	Comment
Columns	• Maintain some carrier gas flow to protect the columns.
Oven	• Reduce temperature to save power. • Turn off to save the most power.
Inlets	For all inlets: • Reduce temperatures. Reduce temperatures to 40 °C or Off to save the most power.
Split/splitless	• Use split mode to prevent diffusion of contamination from the vent line. Use reduced split ratio. • Reduce pressure. Consider using current Gas Saver levels, if used.
PPIP/PCI	• Use split mode to prevent diffusion of contamination from the vent line. Use reduced split ratio. • Reduce pressure. Consider using current Gas Saver levels, if used. • Reduce temperature.
Detectors	
FID	• Turn off the flame. (This turns off hydrogen and air flows.) • Reduce temperatures. (Keep at or above 150 °C to reduce contamination) • Turn off makeup flow.

GC Component	Comment
FPD+	• Turn off the flame. (This turns off hydrogen and air flows.) • Reduce temperatures. • Leave emission block at 125-175 °C. • Reduce transfer line to 150 °C. • Turn off makeup flow.
ECD	• Reduce makeup flow. Try using 15–20 mL/min and test results. • Maintain temperature to avoid long recovery/stabilization times.
NPD	• Maintain flows and temperatures. Sleep not recommended due to recovery times and also thermal cycling can reduce bead life.
TCD	• Leave filament on. • Leave block temperature on. • Reduce reference and makeup flows.
Other devices	
Valve box	• Reduce temperature. (Keep valve box temperature high enough to prevent sample condensation, if applicable.)
Aux thermal zones	• Reduce or turn off. Also refer to the manuals for any connected device (for example, a connected MSD).
Aux pressures or flows	• Reduce or turn off as appropriate for connected columns, transfer lines, and so forth. Always refer to the manual for any connected device or instrument (for example, a connected MSD), to maintain at least the minimum recommended flows or pressures.

Wake and condition methods

The GC can be programmed to wake in one of several ways:

- By loading the last Active Method used before going to sleep
- By loading the **Wake** method
- By running a method called **Condition**, then loading the last Active Method
- By running a method called **Condition**, then loading the **Wake** method

NOTE

The GC can also store wake, sleep, and condition methods which have been created by a connected data system. Although these methods are not visually displayed on the GC, once downloaded to the GC from the data system, they can be used by the GC scheduler functionality.

These choices provide flexibility in how you prepare the GC after a sleep cycle.

A Wake method sets a temperatures and flows. The oven temperature program is isothermal since the GC does not start a run. When the GC loads a **Wake** method, it remains at those settings until you load another method using the touchscreen, data system, or by starting a sequence.

A **Wake** method can include any settings, however it typically will do the following:

- Restore inlet, detector, column, and transfer line flows.
- Restore temperatures.
- Ignite the FID or FPD+ flame.
- Restore inlet modes.

A Condition method sets flows and temperatures for the duration of the method's oven program. When the program ends, the GC loads either the **Wake** method or the last Active Method before sleep, as specified in the instrument schedule (or when manually exiting the sleep state).

One possible use for a condition method is to set higher than normal temperatures and flows to bake out any possible contamination that may have collected in the GC during sleep.

To Set the GC to Conserve Resources

To set the GC to conserve resources by creating and using an Instrument Schedule:

1. Select the **Settings > Scheduler**.
2. If needed, select the **Sleep/Wake** tab.
3. Create the **Instrument Schedule**. You do not have to program events for every day. For example, you can program the GC to sleep on Friday evening, then wake on Monday morning, keeping it continuously at operating conditions during weekdays.
4. Choose **Set Wake Method** for each desired day, as applicable. This will cause the Wake method to be run when the GC wakes on the selected days.
5. Enter a **Wake Time** for each desired day, based on a 24-hour clock (for example, 9:00 AM is 09:00, 9:00 PM is 21:00).
6. Choose **Set Sleep Method** for each desired day, as applicable. This will cause the Sleep method to be run prior to the GC going to sleep on the selected days.
7. Enter a **Sleep Time** for each desired day, based on a 24-hour clock.
8. If desired, edit the sleep, wake or conditioning methods. Select **Edit Wake Method**, **Edit Conditioning Method**, or **Edit Sleep Method** as needed.
9. Select **Scheduler Options**.
10. Decide how to restore flows. Choose the desired options:
 - **Wake to last active method before sleep**: At the specified time, the GC will restore the last active method used before it went to sleep.
 - **Perform a conditioning run before waking**: At the specified time, the GC will load the conditioning method. This method runs once.
11. Select **Apply**. The settings are saved to the GC.



13 Programming

Clock Time Programming

- Using clock time events

- Adding events to the clock table

- Deleting clock time events

Clock Time Programming

Clock time programming allows certain setpoints to change automatically at a specified time during a 24-hour day. Thus, an event programmed to occur at 14:35 hours will occur at 2:35 in the afternoon. A running analysis or sequence has precedence over any clock table events occurring during this time. When this happens, such events are not executed.

Possible clock time events include:

- Valve control
- Method and sequence loading
- Starting sequences
- Initiating blank and prep runs
- Column compensation changes
- Adjustments of the detector offset

Using clock time events

The Clock Table function allows you to program events to occur during a day based on the 24-hour clock. Clock table events that would occur during a run or sequence are ignored.

For example, the clock table could be used to make a blank run before you even get to work in the morning.

Adding events to the clock table

1. Select **Settings > Scheduler > Clock Table**.
2. Select **Add**.
3. Choose your **Clock Type** and **Frequency** from their respective dropdown menus.
4. Set the **Time** you want this event to occur at.
5. Select **Add** to add this entry to the clock table.
6. Repeat this process until all entries are added.

Deleting clock time events

1. Select **Settings > Scheduler > Clock Table**.
2. Touch or click **Select**.
3. Touch or click the event to delete, then select **Delete**.
4. Select **Done** to accept the change.
5. Select **Apply**.

14

Flow and Pressure Modules

About Flow and Pressure Control

Maximum Operating Pressure

Auxiliary Pressure Controllers

Restrictors

Selecting a frit

Example: Using the PCM Channels

PIDs

About Flow and Pressure Control

The GC uses four types of electronic flow or pressure controllers: inlet modules, detector modules, pressure control modules (PCMs), and auxiliary pressure controllers (Aux EPCs).

All of these modules mount in the slots at the top rear of the GC. The slots are identified by numbers, as shown here.

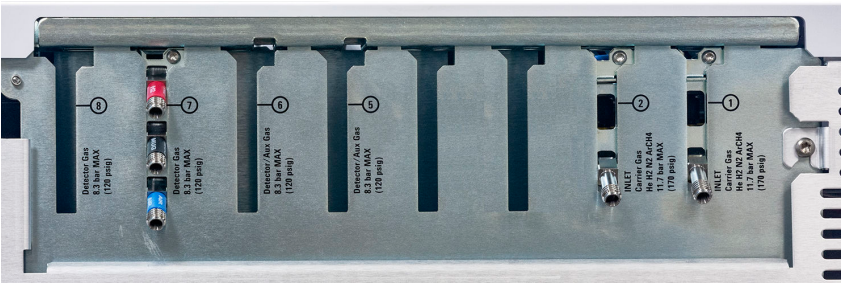


Figure 21. 8860 EPC module slots

Table 16. EPC Modules

Number	Usage
1	Inlet module
2	Inlet module
3	blank
4	blank
5	Aux EPC
6	Detector module/AuxEPC
7	Detector module
8	Detector module

Maximum Operating Pressure

We recommend a maximum continuous operating pressure of 170 psi to avoid excessive wear and leaks.

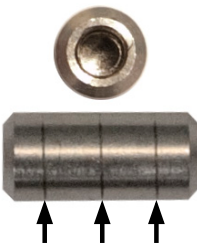
Auxiliary Pressure Controllers

The Auxiliary Pressure Controller (Aux epc) is also a general purpose device. It has three independent forward-pressure regulated channels. Channels are designated by numbers 1 through 3.

Restrictors

The Aux EPC and aux PCM channel use frit-type restrictors to allow accurate flow control. To work properly, there must be adequate flow resistance downstream of the pressure sensor. Each channel provides a frit-type restrictor. Four frits are available.

Table 17. Auxiliary channel frits

Frit marking		Flow resistance	Flow characteristic	Often used with
Three rings Blue		High	3.33 ± 0.3 SCCM @ 15 PSIG	NPD Hydrogen
Two rings Red		Medium	30 ± 1.5 SCCM H ₂ @ 15 PSIG	FID Hydrogen
One ring Brown		Low	400 ± 30 SCCM AIR @ 40 PSIG	FID Air, FPD+ Air, QuickSwap, Splitter, Deans Switch
None (brass tube)		Zero	No restriction	Headspace vial pressurization

The one ring frit (low resistance, high flow) is installed in all channels in the AUX EPC when the instrument (or accessory) is shipped. No frit ships in the PCM Aux channel.

When installing or replacing a frit, always use a new O-ring (5180-4181, 12/pk).

Selecting a frit

The frits change the control range of the channels. The objective is to find a frit that allows the required range of flows at reasonable source pressures.

- For an auxiliary channel ordered as an option (part of the GC order), use the frit supplied by the factory.

- For an auxiliary channel ordered as an accessory (separate from the GC order), see the instruction information supplied with the accessory.
- For a non-Agilent instrument, you must experiment to find the appropriate frit.

When you change a frit, you change the physical characteristics of the channel. It may be desirable (or necessary) to change the PID constants for that channel. See **PIDs**.

Example: Using the PCM Channels

The two channels in a PCM are different. Channel 1 is used to supply a pressure. Channel 2 may be used in the same way, but can also be used to maintain a pressure by reversing the input and output connections.

Channel 1: Forward-pressure

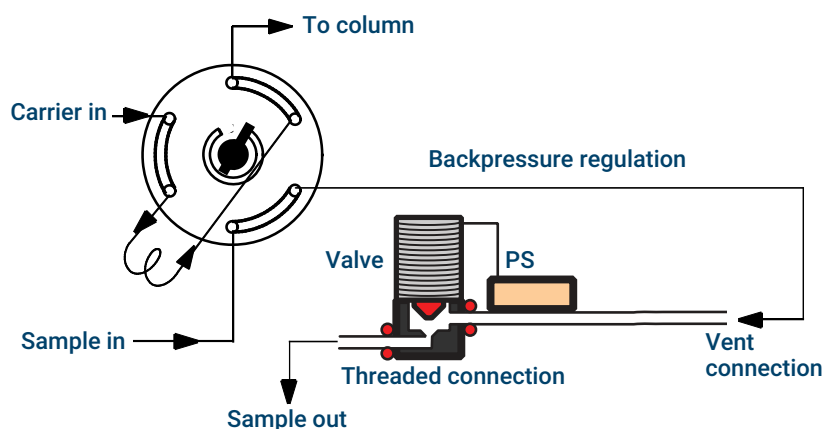
This is identical to the carrier gas channel for the packed column inlet.

Channel 2: Two-way channel

If gas is supplied at the threaded connection and delivered by the tubing, this operates as a forward pressure regulator. But the connections can be reversed—requiring some fittings—so that it will maintain the gas supplied to it at a fixed pressure. In this mode it behaves as a controlled leak.

This can be used to maintain the gas in a gas sampling valve at a fixed pressure, even if the supply varies somewhat. The result is improved reproducibility in sample amount.

The figure shows the connections with the valve in the Load position.



PIDs

The behavior of a pressure control module is governed by a set of three constants, called **P** (proportional), **I** (integral), and **D** (differential).

Certain gases or special applications (such as headspace vial pressurization, flow splitter, and backflush applications) require different PIDs than those provided at the factory.

The table summarizes custom PID values required for selected applications. Note that if you are updating an AUX EPC module, you will need to change the frit for the channel used. See also **"Restrictors."**

Table 18. PIDs and frits

Application	Module	AUX frit	Select Available PID Values
Headspace vial pressurization	AUX EPC	No color or rings	AUX_EPC_Headspace
Headspace sampling loop	PCM in backpressure control		PCM_Headspace

15 Inlets

Inlet Overview

About the Split/Splitless Inlet

Selecting the correct S/SL inlet liner

About the Purged Packed Column Inlet

About the Packed Column Inlet

About the Cool On-Column Inlet

Setup modes of the COC inlet

Retention gaps

Inlet Overview

Table 19. Comparing inlets

Inlet	Column	Mode	Sample concentration	Comments	Sample to column
Split/splitless	Capillary	Split	High		Very little
		Pulsed split ¹	High		Very little
		Splitless	Low		All
		Pulsed splitless ¹	Low		All
Purge packed column	Packed	n/a	Any	OK if resolution not critical	All
	Large capillary	n/a	Any		All
Packed column	Packed	n/a	Any		All
Cool on-column	Capillary	n/a	Low or labile	Minimal discrimination and decomposition	All

¹Pulsed split and pulsed splitless modes are not available for split/splitless inlets on EPR (electronic pneumatics regulation) equipped GCs.

About the Split/Splitless Inlet

This inlet is used for split, splitless, pulsed splitless, or pulsed split analyses. You can choose the operating mode from the inlet parameter list. The split mode is generally used for major component analyses, while the splitless mode is used for trace analyses. The pulsed splitless and pulsed split modes are used for the same type of analyses as split or splitless, but allow you to inject larger samples. Sleep mode is used for creating sleep methods, or if all gas needs to flow onto the column.

NOTE

Pulsed split and pulsed splitless modes are not supported on EPR (electronic pneumatics regulation) equipped GCs.

This inlet is available in both EPC (electronic pneumatics control) and EPR (electronic pneumatics regulation) equipped GCs.

Selecting the correct S/SL inlet liner

Split liner

A good liner for split mode operation will offer very little restriction to the split flow path between the bottom of the liner and the inlet gold seal and between the outside of the liner and the inside of the injection port body. The preferred Agilent split liner, part number 5183-4647, incorporates a glass positioning bead on the bottom to facilitate this. It also has glass wool inside the liner that provides for complete sample vaporization across the boiling point range of the sample. Select an appropriate liner from ???.

Table 20. Split mode liners

Liner	Description	Volume (μL)	Mode	Deactivated	Part Number
	Low Pressure Drop – Positioning Bead	870	Split – Fast Injection	Yes	5183-4647
	4 mm id, Glass Wool	990	Split – Fast Injection	No	19251-60540
	Low Pressure Drop, Glass Wool	870	Split	Yes	5190-2295

Splitless liner

NOTE

Pulsed split and pulsed splitless modes are not supported on EPR (electronic pneumatics regulation) equipped GCs.

The liner volume must contain the solvent vapor. The liner should be deactivated to minimize sample breakdown during the splitless time. Solvent vapor volume can be reduced by using Pulsed Splitless mode. Use the Vapor Volume Calculator to determine vapor volume requirements.

Vapor volume < 300 μL

Use 2 mm liner (250 μL volume), 5181-8818 or similar.

Vapor volume 225 – 300 μL

Consider pulsed splitless mode to reduce vapor volume.

Vapor volume > 300 μL

Use 4 mm liner, 5062-3587 or similar.

Vapor volume > 800 μL

Consider pulsed splitless mode to reduce vapor volume.

For thermally labile or reactive samples, use G1544-80700 (open top) or G1544-80730 (top taper) liners.

Table 21. Splitless mode liners

Liner	Description	Volume (uL)	Mode	Deactivated	Part Number
	Single Taper Glass Wool	900	Splitless	Yes	5062-3587
	Single Taper	900	Splitless	Yes	5181-3316
	Dual Taper	800	Splitless	Yes	5181-3315
	2 mm Quartz	250	Splitless	No	18740-80220
	2 mm Quartz	250	Splitless	Yes	5181-8818
	1.5 mm	140	Direct Inject, Purge and Trap, Headspace	No	18740-80200
	Single Taper Glass Wool	900	Splitless	Yes	5062-3587
	Single Taper	900	Splitless	Yes	5181-3316
	4 mm Single Taper	Direct column connect		Yes	G1544-80730
	4 mm Dual Taper	Direct column connect		Yes	G1544-80700
	Single Taper	900	Splitless	Yes	5190-2293

About the Purged Packed Column Inlet

This inlet is used with packed columns when high-efficiency separations are not required. It can also be used with wide-bore capillary columns, if flows greater than 10 mL/min are acceptable.

The inlet is flow controlled, whether using an undefined column (packed column) or a defined capillary column.

About the Packed Column Inlet

This inlet is used with packed columns when high-efficiency separations are not required.

The inlet is flow controlled regardless of whether the column is defined or not.

This inlet is available in EPR (electronic pneumatics regulation) equipped GCs only.

About the Cool On-Column Inlet

This inlet introduces liquid sample directly onto a capillary column. To do this, both the inlet and the oven must be cool at injection, either at or below the boiling point of the solvent.

Because the sample does not vaporize immediately in the inlet, problems with sample discrimination and sample alteration are minimized. If done properly, cool-on column injection also provides accurate and precise results.

You can operate the inlet in track oven mode, where the inlet temperature follows the column oven, or you can program up to three temperature ramps.

Setup modes of the COC inlet

The COC inlet hardware must be set up for one of three usages, depending on the type of injection and column size.

- 0.25 mm or 0.32 mm automated on-column.
- 0.53 mm automatic on-column or retention gap
- 0.2 mm manual

Retention gaps

Because the sample is injected directly onto the column, it is strongly suggested that a retention gap—or guard column—be used to protect your column. A retention gap is a deactivated column that is connected between the inlet and the analytical column. If you choose to use one, it is suggested that at least 1 m of retention gap be installed per 1 mL of sample injected. Information on ordering retention gaps can be found in the Agilent catalog for consumables and supplies.



16 Columns and Oven

Hydrogen Sensor

Instrument logs

Calibration

Status information

Operation with an Agilent data system

Hydrogen Sensor

The optional G6598A hydrogen sensor module checks for uncombusted free hydrogen in the GC column oven. During normal operation with hydrogen as a carrier gas, leaks from the inlets or detectors could possibly put hydrogen gas directly into the oven. Hydrogen–air mixtures are potentially explosive in concentrations of 4–74.2% hydrogen by volume. The sensor monitors the free hydrogen level in the oven and will trigger a shutdown of all hydrogen gas flows before the hydrogen level in the oven reaches 1%.

In the event of a hydrogen shutdown, the GC records the event in its system log. See the Troubleshooting manual for more information about GC shutdown events and how to clear them.

The GC can only shut down hydrogen gas flows that have been configured properly. Always configure the gas types used for the inlets, detectors, and so forth. The hydrogen sensor has possible cross-sensitivity to helium in rare cases.

NOTE

Accuracy of the hydrogen sensor may be degraded above 1829 m (6,000 ft) elevation.

Usage

Intended Use: The G6598A hydrogen sensor is intended for indication only. The hydrogen sensor is not intended as a safety device.

At startup, the GC automatically purges the oven (opens the flaps and runs the fan). After a few minutes, the GC automatically zeroes the hydrogen sensor and begins to monitor the hydrogen content in the GC oven.

Audible Alarm

If the GC is configured to sound audible alerts, the GC will beep whenever the hydrogen sensor triggers a shutdown. To clear the shutdown and stop the audible alert (beeping), power cycle the GC. Upon startup, the GC opens the oven flaps and runs the fan to purge the oven.

Instrument logs

The GC will log the following hydrogen sensor events in its event logs:

- Hydrogen shutdowns initiated by the hydrogen sensor
- Calibrations
- Hydrogen sensor tests

Calibration

The G6598A Hydrogen Sensor requires periodic calibration after it is installed. Agilent recommends calibrating the hydrogen sensor every 6 months. Use the GC's EMF feature to set up a scheduled reminder to perform calibration.

1. Go to **Maintenance > Hydrogen Sensor** and select the **On time since calibration** parameter.

2. Select **Details**.
3. Enable **Service Due**. As the value, enter the desired number of days between calibrations (default is 6 months) and select **Apply**. If desired, add a **Service Warning** for a shorter number of days, to receive an additional prompt.

As with other EMFs, when the **Service Warning** or **Service Due** counter expires, the GC raises a Condition to alert you that it is time to repeat the calibration. See [Calibrate the sensor](#).

When the reminder appears, use the GC's automated calibration procedure. If a calibration fails for any reason (for example, a lack of calibration gas), the sensor continues to use its existing calibration data.

Calibrate the sensor

Calibration requires several minutes, and includes steps to cool the GC oven, install the calibration gas, and set the calibration gas flow rate. The GC will walk you through the needed steps.

To calibrate the sensor:

Parts needed:

- Hydrogen sensor calibration gas cylinder (5190-6890)
- Flow meter capable of reading flows of 30 mL/min

You may also need:

- Regulator (G3440-80153)
- Copper tubing and 1/8-inch Swagelok nut and ferrules
- Open-end wrenches to connect the calibration gas to the GC

1. Select **Settings > Calibration > H2 Sensor**.
2. Select **Start Calibration**.
3. Follow the prompts. The GC will cool and purge the oven, guide you through connecting the calibration gas to the sensor and setting the calibration gas flow rate, and then perform the calibration.

When calibration completes, you may shut off and disconnect the calibration gas. If calibration fails, check the calibration gas supply.

Status information

If a calibration fails, a notification will appear on the diagnostics tab.

Operation with an Agilent data system

Using the hydrogen sensor with an Agilent data system provides additional features. Use the data system to:

- Print calibration reports. The report includes a plot of all calibration data stored in the GC.
- Store lot number and expiration date information for the calibration gas cylinder.

- View the hydrogen sensor status information in the GC status user interface. The status shows the current percent hydrogen level and any messages related to the hydrogen sensor.
- Plot the measured hydrogen level as a diagnostic signal, if desired.
- View and print all logged entries for calibrations, cylinder information, and shutdowns.
- Start a hydrogen sensor calibration.

Chromatographic Checkout

About Chromatographic Checkout

To Prepare for Chromatographic Checkout

To check FID performance with a purged packed, split splitless or cool-on column inlet

To check FID performance with a packed column inlet (PCI)

To check TCD performance with a purged packed, split splitless or cool-on column inlet

To check TCD performance with a packed column inlet (PCI)

To Check NPD Performance

To Check ECD Performance

To Check FPD+ Performance (Sample 5188-5953)

Preparation

Phosphorus performance

Sulfur performance

To Check FPD+ Performance (Sample 5188-5245, Japan)

Preparation

Phosphorus performance

Sulfur performance

This section describes the general procedure for verifying performance against the original factory standards. The checkout procedures described here assume a GC that has been in use for some period of time. Therefore the procedures ask that you perform bakeouts, replace consumable hardware, install the checkout column, and so forth.

About Chromatographic Checkout

The tests described in this section provide basic confirmation that the GC and detector can perform comparably to factory condition. However, as detectors and the other parts of the GC age, detector performance can change. The results presented here represent typical outputs for typical operating conditions and are not specifications.

The tests assume the following:

- Use of an automatic liquid sampler. If not available, use a suitable manual syringe instead of the syringe listed.
- Use of a 10 µL syringe in most cases. However, a 5 µL syringe is an acceptable substitute.
- Use of the septa and other hardware (liners, jets, adapters, and so forth) described. If you substitute other hardware, performance can vary.

Checkout can also be performed as a diagnostic test if your GC uses a simple flow path (inlet to column to detector, for example). You will need to prepare the GC as described in the following sections, including preparation of the checkout samples, gas types, and consumable parts. The automated procedure will walk you through the checkout process and create the checkout method for you. If your system is more complex, you can always manually perform a checkout as described in the following sections. To run the automatic procedure, go to **Diagnostics > Diagnostic Tests > System Checkout**.

To Prepare for Chromatographic Checkout

Because of the differences in chromatographic performance associated with different consumables, Agilent strongly recommends using the parts listed here for all checkout tests. Agilent also recommends installing new consumable parts whenever the quality of the installed ones is not known. For example, installing a new liner and septum ensures that they will not contribute any contamination to the results.

When the GC is delivered from the factory, these consumable parts are new and do not need replacement.

NOTE

For a new GC, check the installed inlet liner. The liner shipped in the inlet may not be the liner recommended for checkout.

1. Check the indicators/dates on any gas supply traps. Replace/recondition expended traps.
2. Install new consumable parts for the inlet and prepare the correct injector syringe (and needle, as needed).

Table 22. Recommended parts for checkout by inlet type

Recommended part for checkout	Part number
Split splitless inlet	
Syringe, 10- μ L	5181-1267
O-ring	5188-5365
Septum	5183-4757
Liner	5190-2295
Purged packed column inlet	
Syringe, 10- μ L	5181-1267
O-ring	5080-8898
Septum	5183-4757
Packed column inlet	
Syringe, 10- μ L	5181-1267
O-ring	5080-8898
Septum	5183-4757
Cool on-column inlet	
Septum	5183-4758
Septum nut	19245-80521
Syringe, 5- μ L on-column	5182-0836
0.32-mm needle for 5- μ L syringe	5182-0831
7693A ALS: Needle support insert, COC	G4513-40529
Insert, fused silica, 0.32-mm id	19245-20525

Table 23. Checkout standards

Standard	Part number	Number of ampoules
FID checkout	5188-5372	3
TCD checkout	18710-60170	3

Chromatographic Checkout

To Prepare for Chromatographic Checkout

Standard	Part number	Number of ampoules
ECD checkout	18713-60040	3
NPD checkout	18789-60060	3
FPD+ checkout (methyl parathion)	5188-5953	3
OQ/PV Headspace checkout	5182-9733	1

To check FID performance with a purged packed, split splitless or cool-on column inlet

FID performance is checked differently depending on the inlet being used. For GCs equipped with a packed column inlet (PCI), use **To check FID performance with a packed column inlet (PCI)**.

- Gather the following:
 - Evaluation column, HP-5 30 m × 0.32 mm × 0.25 µm (19091J-413)
 - FID performance evaluation (checkout) sample (5188-5372)
 - Chromatographic-grade isooctane
 - 4-mL solvent and waste bottles or equivalent for autoinjector
 - 2-mL sample vials or equivalent for sample
 - Inlet and injector hardware (See **"To Prepare for Chromatographic Checkout."**)
- Verify the following:
 - Capillary column jet installed. If not, select and install a capillary column jet.
 - Capillary column adapter installed. If not, install it.
 - Chromatographic-grade gases plumbed and configured: helium as carrier gas, nitrogen, hydrogen, and air.
 - Empty waste vials loaded in sample turret.
 - 4-mL solvent vial with diffusion cap filled with isooctane and inserted in Solvent A injector position.
- Replace consumable parts (liner, septum, traps, syringe, and so forth) as needed for the checkout. See **"To Prepare for Chromatographic Checkout."**
- Install the evaluation column.
 - Bake out the evaluation column for at least 30 min at 180 °C.
 - Be sure to configure the column.
- Check the FID baseline output. The output should be between 5 pA and 20 pA and relatively stable. (If using a gas generator or ultra pure gas, the signal may stabilize below 5 pA.) If the output is outside this range or unstable, resolve this problem before continuing.
- If the output is too low:
 - Check that the electrometer is on.
 - Check that the flame is lit.
 - Check that the signal is set to the correct detector.
- Create or load a method with the parameter values listed in **Table 24**.

Table 24. FID Checkout Conditions

Column and sample	
Type	HP-5, 30 m × 0.32 mm × 0.25 µm (19091J-413)
Sample	FID checkout 5188-5372
Column flow	6.5 mL/min
Column mode	Constant flow for EPC equipped GCs. Constant pressure mode (30 psi) for EPR equipped GCs.

Chromatographic Checkout

To check FID performance with a purged packed, split splitless or cool-on column inlet

Split/splitless inlet

Temperature	250 °C
Mode	Splitless
Split flow	40 mL/min
Purge time	0.5 min
Gas saver	ON
Gas saver flow	20 mL/min
Gas saver time	3 min

Purged packed column inlet

Temperature	250 °C
Septum purge	3 mL/min

Cool on-column inlet

Temperature	Track oven
Septum purge	15 mL/min

Detector

Temperature	300 °C
H2 flow	30 mL/min
Air flow	400 mL/min
Makeup flow (N2)	25 mL/min
Lit offset	Typically 2 pA

Oven

Initial temp	75 °C
Initial time	0.5 min
Rate 1	20 °C/min
Final temp	190 °C
Final time	0 min

ALS settings (if installed)

Sample washes	2
Sample pumps	6
Sample wash volume	8 (maximum)
Injection volume	1 µL
Syringe size	10 µL
Solvent A pre washes	2
Solvent A post washes	2
Solvent A wash volume	8
Solvent B pre washes	0
Solvent B post washes	0
Solvent B wash volume	0
Injection mode (7693A)	Normal
Airgap Volume (7693A)	0.20
Viscosity delay	0
Inject Dispense Speed (7693A)	6000
PreInjection dwell	0
PostInjection dwell	0

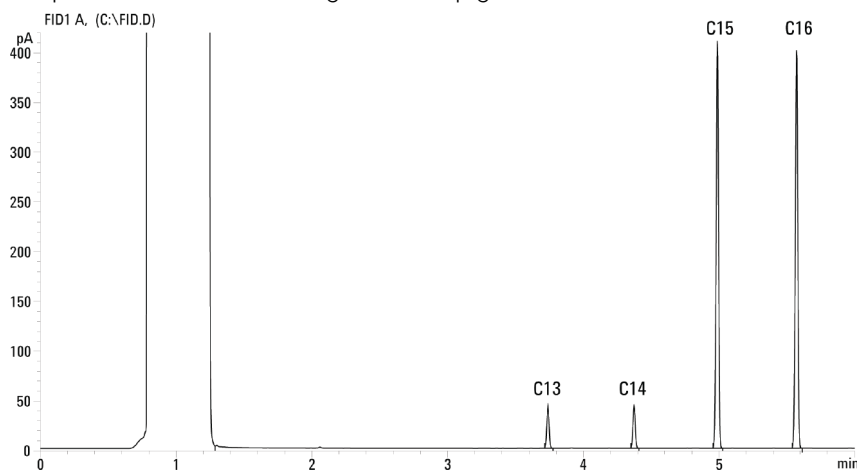
Manual injection

Chromatographic Checkout

To check FID performance with a purged packed, split splitless or cool-on column inlet

Injection volume	1 μ L
Data system	
Data rate	5 Hz

8. If using a data system, prepare the data system to perform one run using the loaded checkout method. Make sure that the data system will output a chromatogram.
If not using a data system, create a one sample sequence using the GC Browser Interface.
9. Start the run.
If performing an injection using an autosampler, start the run using the data system, or press the "Start" button.
If performing a manual injection (with or without a data system):
 - a. Press the "Prep Run" button to prepare the inlet for splitless injection.
 - b. When the GC becomes ready, inject 1 μ L of the checkout sample and press "Start" on the GC.
 - c. The following chromatogram shows typical results for a new detector with new consumable parts installed and nitrogen makeup gas.



To check FID performance with a packed column inlet (PCI)

FID performance is checked differently depending on the inlet being used. This procedure applies to GCs equipped with a packed column inlet (PCI). For all other inlet types, use [To check FID performance with a purged packed, split splitless or cool-on column inlet](#).

1. Gather the following:
 - Evaluation column, 10% OV-101, 5 ft, OD 1/8, ID 2 mm (G3591-81093)
 - FID MDL sample (5188-5372)
 - Chromatographic-grade isooctane
 - 4-mL solvent and waste bottles or equivalent for autoinjector
 - 2-mL sample vials or equivalent for sample
 - Inlet and injector hardware (See [“To Prepare for Chromatographic Checkout.”](#))
2. Verify the following:
 - Packed column jet installed. If not, select and install a packed column jet.
 - Packed column adapter installed. If not, install it.
 - Chromatographic-grade gases plumbed and configured: helium as carrier gas, nitrogen, hydrogen, and air.
 - Empty waste vials loaded in sample turret.
 - 4-mL solvent vial with diffusion cap filled with isooctane and inserted in Solvent A injector position.
3. Replace consumable parts (liner, septum, traps, syringe, and so forth) as needed for the checkout. See [To Prepare for Chromatographic Checkout](#).
4. Install the evaluation column. (See the procedure for the PCI in the Maintenance manual.)
 - Bake out the evaluation column for at least 30 min at 180 °C. (See the procedure for the PCI in the Maintenance manual.)
 - Be sure to configure the column.
5. Check the FID baseline output. The output should be between 5 pA and 20 pA and relatively stable. (If using a gas generator or ultra pure gas, the signal may stabilize below 5 pA.) If the output is outside this range or unstable, resolve this problem before continuing.
6. If the output is too low:
 - Check that the electrometer is on.
 - Check that the flame is lit.
 - Check that the signal is set to the correct detector.
7. Create or load a method with the parameter values listed in [Table 25](#).

Table 25. FID Checkout Conditions- Packed Column Inlet

Column and sample	
Type	10% OV-101, 5 ft, OD 1/8, ID 2 mm (G3591-81093)
Sample	FID MDL sample (5188-5372)

Chromatographic Checkout

To check FID performance with a packed column inlet (PCI)

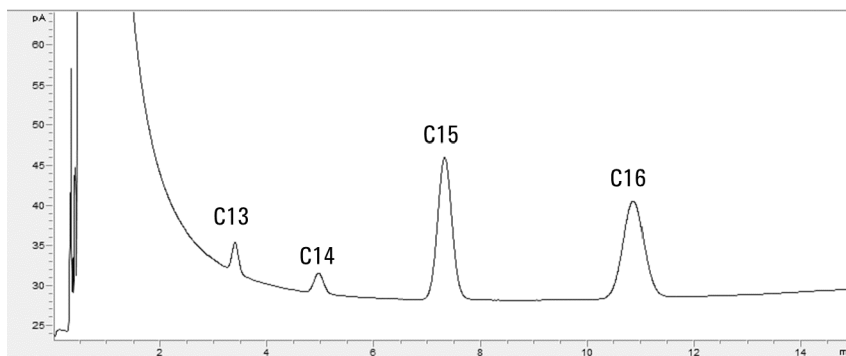
Column flow	20 mL/min
Column mode	Flow mode
Packed column inlet	
Temperature	250 °C
Detector	
Temperature	300 °C
H2 flow	30 mL/min
Air flow	400 mL/min
Makeup flow (N ₂)	OFF
Mode	Constant makeup flow OFF
Flame	On
Lit offset	Typically 2 pA
Oven	
Constant temperature	180 °C
Time	15 min
ALS settings (if installed)	
Sample washes	2
Sample pumps	6
Sample wash volume	8 µL
Injection volume	1 µL
Syringe size	10 µL
Solvent A pre washes	2
Solvent A post washes	2
Solvent A wash volume	8 µL
Injection mode (7693A)	Normal
Airgap Volume (7693A)	0.20
Viscosity delay	0
Inject Dispense Speed (7693A)	6000
PreInjection dwell	0
PostInjection dwell	0
Manual injection	
Injection volume	1 µL
Data system	
Data rate	5 Hz

8. If using a data system, prepare the data system to perform one run using the loaded checkout method. Make sure that the data system will output a chromatogram.
If not using a data system, create a one sample sequence using the Browser Interface.
9. Start the run.
If performing an injection using an autosampler, start the run using the data system, or press  on the Browser Interface.
If performing a manual injection (with or without a data system):
 - a. Press the "Prep Run" button to prepare the inlet for splitless injection.
 - b. When the GC becomes ready, inject 1 µL of the checkout sample and press the "Start" button.

Chromatographic Checkout

To check FID performance with a packed column inlet (PCI)

- c. The following chromatogram shows typical results for a new detector with new consumable parts installed and nitrogen makeup gas.



To check TCD performance with a purged packed, split splitless or cool-on column inlet

TCD performance is checked differently depending on the inlet being used. For GCs equipped with a packed column inlet (PCI), use **To check TCD performance with a packed column inlet (PCI)**.

- Gather the following:
 - Evaluation column, HP-5 30 m × 0.32 mm × 0.25 µm (19091J-413)
 - FID/TCD performance evaluation (checkout) sample (18710-60170)
 - 4-mL solvent and waste bottles or equivalent for autoinjector
 - Chromatographic-grade hexane
 - 2-mL sample vials or equivalent for sample
 - Chromatographic-grade helium as carrier, makeup, and reference gas
 - Inlet and injector hardware See **"To Prepare for Chromatographic Checkout."**
- Verify the following:
 - Chromatographic-grade gases plumbed and configured: helium as carrier gas and reference gas.
 - Empty waste vials loaded in sample turret.
 - 4-mL solvent vial with diffusion cap filled with hexane and inserted in Solvent A injector position.
- Replace consumable parts (liner, septum, traps, syringe, and so forth) as needed for the checkout. See **"To Prepare for Chromatographic Checkout."**
- Install the evaluation column.
 - Bake out the evaluation column for at least 30 min at 180 °C.
 - Configure the column
- Create or load a method with the parameter values listed in **Table 26**.

Table 26. TCD Checkout Conditions

Column and sample	
Type	HP-5, 30 m × 0.32 mm × 0.25 µm (19091J-413)
Sample	FID/TCD checkout 18710-60170
Column flow	6.5 mL/min
Column mode	Constant flow for EPC equipped GCs. Constant pressure mode (30 psi) for EPR equipped GCs.
Split/splitless inlet	
Temperature	250 °C
Mode	Splitless
Purge flow	60 mL/min
Purge time	0.75 min
Gas saver	ON
Gas saver flow	20 mL/min

Chromatographic Checkout

To check TCD performance with a purged packed, split splitless or cool-on column inlet

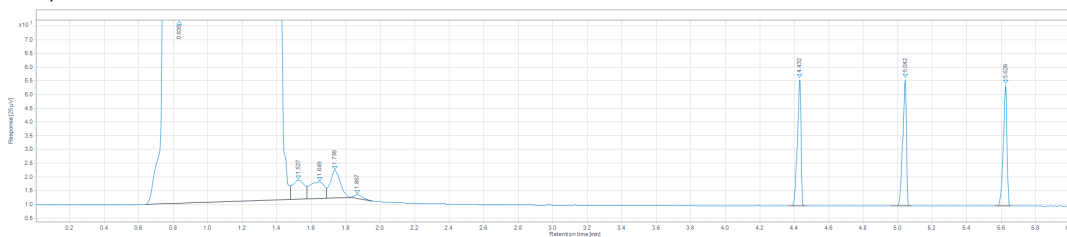
Gas saver time	3 min
Purged packed column inlet	
Temperature	250 °C
Cool on-column inlet	
Temperature	Oven track
Septum purge	15 mL/min
Detector	
Temperature	300 °C
Reference flow (He)	30 mL/min
Makeup flow (He)	2 mL/min
Baseline output	< 30 display counts on Agilent OpenLab CDS ChemStation Edition (< 750 μ V)
Oven	
Initial temp	75 °C
Initial time	0.5 min
Rate 1	20 °C/min
Final temp	190 °C
Final time	0 min
ALS settings (if installed)	
Sample washes	2
Sample pumps	6
Sample wash volume	8 (maximum)
Injection volume	1 μ L
Syringe size	10 μ L
Solvent A pre washes	2
Solvent A post washes	2
Solvent A wash volume	8
Solvent B pre washes	0
Solvent B post washes	0
Solvent B wash volume	0
Injection mode (7693A)	Normal
Airgap Volume (7693A)	0.20
Viscosity delay	0
Inject Dispense Speed (7693A)	6000
PreInjection dwell	0
PostInjection dwell	0
Manual injection	
Injection volume	1 μ L
Data system	
Data rate	5 Hz

- Display the signal output. A stable output at any value between 12.5 and 750 μ V (inclusive) is acceptable.
 - If the baseline output is < 0.5 display units (< 12.5 μ V), verify that the detector filament is on. If the offset is still < 0.5 display units (< 12.5 μ V), your detector requires service.

Chromatographic Checkout

To check TCD performance with a purged packed, split splitless or cool-on column inlet

- If baseline output is > 30 display units (> 750 μV), there may be chemical contamination contributing to the signal. Bakeout the TCD. If repeated cleanings do not give an acceptable signal, check gas purity. Use higher purity gases and/or install traps.
7. If using a data system, prepare the data system to perform one run using the loaded checkout method. Make sure that the data system will output a chromatogram.
 8. Start the run.
If performing an injection using an autosampler, start the run using the data system or press the "Start" button.
If performing a manual injection (with or without a data system):
 - a. Press the "Prep Run" button to prepare the inlet for splitless injection.
 - b. When the GC becomes ready, inject 1 μL of the checkout sample and press the "Start" button.
 - c. The following chromatogram shows typical results for a new detector with new consumable parts installed.



To check TCD performance with a packed column inlet (PCI)

TCD performance is checked differently depending on the inlet being used. This procedure applies to GCs equipped with a packed column inlet (PCI). For all other inlet types, use [To check TCD performance with a purged packed, split splitless or cool-on column inlet](#).

- Gather the following:
 - Evaluation column, 10% OV-101, 5 ft, OD 1/8, ID 2 mm (G3591-81093)
 - FID/TCD performance evaluation (checkout) sample (18710-60170)
 - 4-mL solvent and waste bottles or equivalent for autoinjector
 - Chromatographic-grade hexane
 - 2-mL sample vials or equivalent for sample
 - Chromatographic-grade helium as carrier, makeup, and reference gas
 - Inlet and injector hardware (See [“To Prepare for Chromatographic Checkout.”](#))
- Verify the following:
 - Chromatographic-grade gases plumbed and configured: helium as carrier gas and reference gas.
 - Empty waste vials loaded in sample turret.
 - 4-mL solvent vial with diffusion cap filled with hexane and inserted in Solvent A injector position.
- Replace consumable parts (liner, septum, traps, syringe, and so forth) as needed for the checkout. See [“To Prepare for Chromatographic Checkout.”](#)
- Install the evaluation column. (See the procedure for the PCI in the Maintenance manual.)
 - Bake out the evaluation column for at least 30 min at 180 °C. (See the procedure for the PCI in the Maintenance manual.)
 - Configure the column
- Create or load a method with the parameter values listed in [Table 27](#).

Table 27. TCD Checkout Conditions- Packed Column Inlet

Column and sample	
Type	10% OV-101, 5 ft, OD 1/8, ID 2 mm (G3591-81093)
Sample	FID/TCD checkout 18710-60170
Column flow	20 mL/min
Column mode	Flow mode
Packed column inlet	
Temperature	250 °C
Detector	
Temperature	300 °C
Reference flow (He)	20 mL/min
Makeup flow (He)	OFF

Chromatographic Checkout

To check TCD performance with a packed column inlet (PCI)

Baseline output	< 30 display counts on Agilent OpenLAB CDS ChemStation Edition (< 750 μ V)
Oven	
Constant temp	180 °C
Time	15 min
ALS settings (if installed)	
Sample washes	2
Sample pumps	6
Sample wash volume	8 μ L
Injection volume	1 μ L
Syringe size	10 μ L
Solvent A pre washes	2
Solvent A post washes	2
Solvent A wash volume	8 μ L
Injection mode (7693A)	Normal
Airgap Volume (7693A)	0.20
Viscosity delay	0
Inject Dispense Speed (7693A)	6000
PreInjection dwell	0
PostInjection dwell	0
Manual injection	
Injection volume	2 μ L
Data system	
Data rate	5 Hz

6. Display the signal output. A stable output at any value between 12.5 and 750 μ V (inclusive) is acceptable.
 - If the baseline output is < 0.5 display units (< 12.5 μ V), verify that the detector filament is on. If the offset is still < 0.5 display units (< 12.5 μ V), your detector requires service.
 - If baseline output is > 30 display units (> 750 μ V), there may be chemical contamination contributing to the signal. Bakeout the TCD. If repeated cleanings do not give an acceptable signal, check gas purity. Use higher purity gases and/or install traps.
7. If using a data system, prepare the data system to perform one run using the loaded checkout method. Make sure that the data system will output a chromatogram.
8. Start the run.

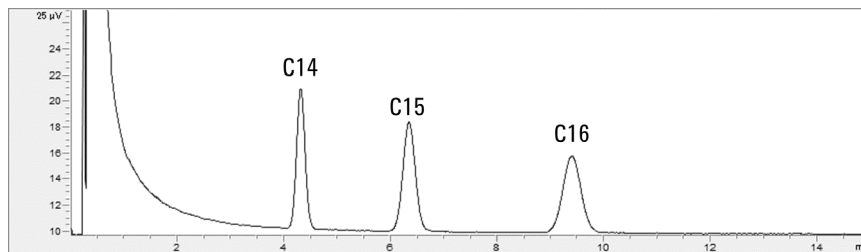
If performing an injection using an autosampler, start the run using the data system or press the "Start" button on the GC.

If performing a manual injection (with or without a data system):

 - a. Press the "Prep Run" button to prepare the inlet for splitless injection.
 - b. When the GC becomes ready, inject 1 μ L of the checkout sample and press the "Start" button on the GC.
 - c. The following chromatogram shows typical results for a new detector with new consumable parts installed.

Chromatographic Checkout

To check TCD performance with a packed column inlet (PCI)



To Check NPD Performance

- Gather the following:
 - Evaluation column, HP-5 30 m × 0.32 mm × 0.25 µm (19091J-413)
 - NPD performance evaluation (checkout) sample (18789-60060)
 - 4-mL solvent and waste bottles or equivalent for autoinjector.
 - Chromatographic-grade isooctane
 - 2-mL sample vials or equivalent for sample.
 - Inlet and injector hardware See [To Prepare for Chromatographic Checkout](#).
- Verify the following:
 - Capillary column jet installed. If not, select and install a capillary column jet.
 - Chromatographic-grade gases plumbed and configured: helium as carrier gas, nitrogen, hydrogen, and air.
 - Empty waste vials loaded in sample turret.
 - 4-mL vial with diffusion cap filled with isooctane and inserted in Solvent A injector position.
- Replace consumable parts (liner, septum, traps, syringe, and so forth) as needed for the checkout. See [To Prepare for Chromatographic Checkout](#).
- If present, remove any protective caps from the inlet manifold vents.
- Install the evaluation column.
 - Bake out the evaluation column for at least 30 min at 180 °C.
 - Be sure to configure the column
- Create or load a method with the parameter values listed in [Table 28](#).

Table 28. NPD Checkout Conditions

Column and sample	
Type	HP-5, 30 m × 0.32 mm × 0.25 µm (19091J-413)
Sample	NPD checkout 18789-60060
Column mode	Constant flow
Column flow	6.5 mL/min (helium)
Split/splitless inlet	
Temperature	200 °C
Mode	Splitless
Purge flow	60 mL/min
Purge time	0.75 min
Gas saver	ON
Gas saver flow	20 mL/min
Gas saver time	3 min
Purged packed column inlet	
Temperature	200 °C
Cool on-column inlet	
Temperature	Track oven

Chromatographic Checkout

To Check NPD Performance

Septum purge	15 mL/min
Detector	
Temperature	300 °C
H2 flow	3 mL/min
Air flow	60 mL/min
Makeup flow (N2)	3 mL/min
Carrier Gas Flow Correction	None (constant makeup and fuel flow)
Output	20 display units (20 pA)
Oven	
Initial temp	60 °C
Initial time	0 min
Rate 1	20 °C/min
Final temp	200 °C
Final time	3 min
ALS settings (if installed)	
Sample washes	2
Sample pumps	6
Sample wash volume	8 (maximum)
Injection volume	1 µL
Syringe size	10 µL
Solvent A pre washes	2
Solvent A post washes	2
Solvent A wash volume	8
Solvent B pre washes	0
Solvent B post washes	0
Solvent B wash volume	0
Injection mode (7693A)	Normal
Airgap Volume (7693A)	0.20
Viscosity delay	0
Inject Dispense Speed (7693A)	6000
PreInjection dwell	0
PostInjection dwell	0
Manual injection	
Injection volume	1 µL
Data system	
Data rate	5 Hz

7. If using a data system, prepare the data system to perform one run using the loaded checkout method. Make sure that the data system will output a chromatogram.
8. Start the run.

If performing an injection using an autosampler, start the run using the data system, or creating a one sample sequence and pressing  on the Browser Interface.

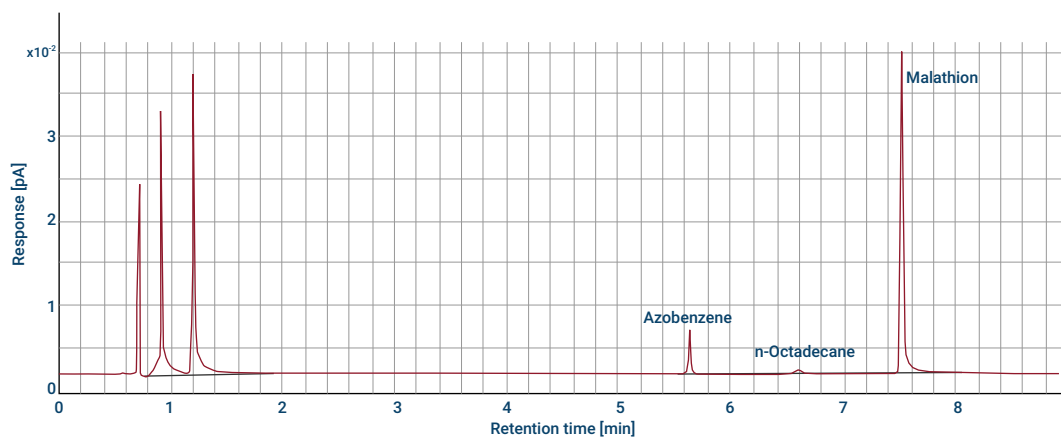
If performing a manual injection (with or without a data system):

 - a. Press the "Prep Run" button to prepare the inlet for splitless injection.
 - b. When the GC becomes ready, inject 1 µL of the checkout sample and press the "Start" button.

Chromatographic Checkout

To Check NPD Performance

- c. The following chromatogram shows typical results for a new detector with new consumable parts installed.



To Check ECD Performance

- Gather the following:
 - Evaluation column, HP-5 30 m × 0.32 mm × 0.25 μm (19091J-413)
 - ECD performance evaluation (checkout) sample (18713-60040, Japan: 5183-0379)
 - 4-mL solvent and waste bottles or equivalent for autoinjector.
 - Chromatographic-grade isooctane
 - 2-mL sample vials or equivalent for sample.
 - Inlet and injector hardware See **To Prepare for Chromatographic Checkout**.
- Verify the following:
 - Clean fused silica indented mixing liner installed. If not, install it.
 - Chromatographic-grade gases plumbed and configured: helium for carrier gas, nitrogen for makeup.
 - Empty waste vials loaded in sample turret.
 - 4-mL vial with diffusion cap filled with hexane and inserted in Solvent A injector position.
- Replace consumable parts (liner, septum, traps, syringe, and so forth) as needed for the checkout. See **To Prepare for Chromatographic Checkout**.
- Install the evaluation column.
 - Bake out the evaluation column for at least 30 minutes at 180 °C.
 - Be sure to configure the column.
- Display the signal output to determine baseline output. A stable baseline output at any value between 0.5 and 1000 Hz (OpenLab CDS ChemStation Edition display units) (inclusive) is acceptable.
 - If the baseline output is < 0.5 Hz, verify that the electrometer is on. If the offset is still < 0.5 Hz, your detector requires service.
 - If the baseline output is > 1000 Hz, there may be chemical contamination contributing to the signal. Bakeout the ECD. If repeated cleanings do not give an acceptable signal, check gas purity. Use higher purity gases and/or install traps.
- Create or load a method with the parameter values listed in **Table 29**.

Table 29. ECD Checkout Conditions

Column and sample	
Type	HP-5, 30 m × 0.32 mm × 0.25 μm (19091J-413)
Sample	ECD checkout (18713-60040 or Japan: 5183-0379)
Column mode	Constant flow
Column flow	6.5 mL/min (helium)
Split/splitless inlet	
Temperature	250 °C
Mode	Splitless
Purge flow	60 mL/min
Purge time	0.75 min

Chromatographic Checkout

To Check ECD Performance

Gas saver	ON
Gas saver flow	20 mL/min
Gas saver time	3 min
Purged packed column inlet	
Temperature	250 °C
Cool on-column inlet	
Temperature	Track oven
Septum purge	15 mL/min
Detector	
Temperature	300 °C
Makeup flow (N2)	25 mL/min (constant + makeup)
Baseline output	Should be < 1000 display counts. In Agilent OpenLab CDS ChemStation Edition (< 1000 Hz)
Oven	
Initial temp	80 °C
Initial time	0 min
Rate 1	15 °C/min
Final temp	180 °C
Final time	10 min
ALS settings (if installed)	
Sample washes	2
Sample pumps	6
Sample wash volume	8 (maximum)
Injection volume	1 µL
Syringe size	10 µL
Solvent A pre washes	2
Solvent A post washes	2
Solvent A wash volume	8
Solvent B pre washes	0
Solvent B post washes	0
Solvent B wash volume	0
Injection mode (7693A)	Normal
Airgap Volume (7693A)	0.20
Viscosity delay	0
Inject Dispense Speed (7693A)	6000
PreInjection dwell	0
PostInjection dwell	0
Manual injection	
Injection volume	1 µL
Data system	
Data rate	5 Hz

7. If using a data system, prepare the data system to perform one run using the loaded checkout method. Make sure that the data system will output a chromatogram.
8. Start the run.

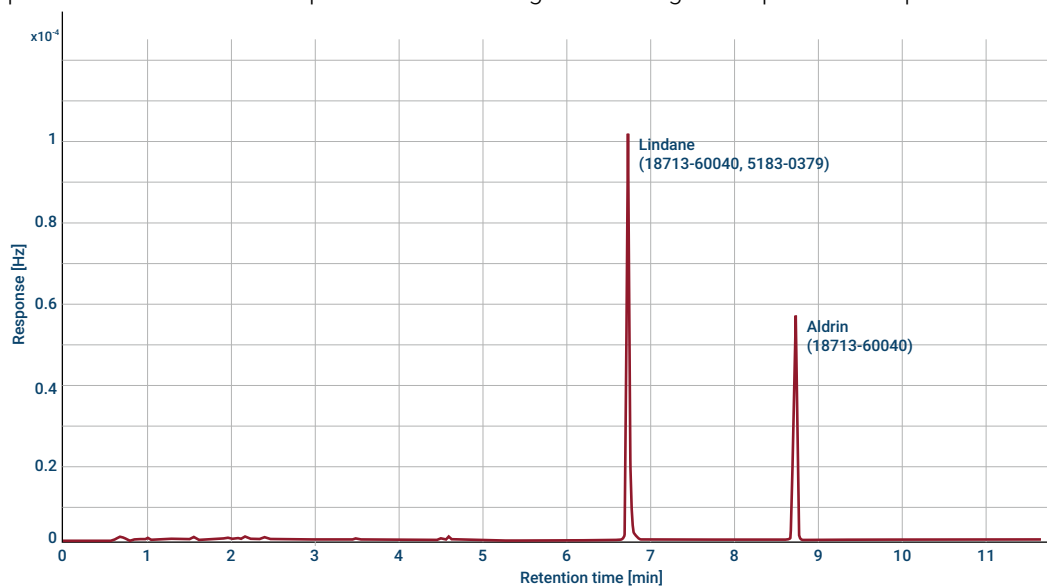
Chromatographic Checkout

To Check ECD Performance

If performing an injection using an autosampler, start the run using the data system or press the "Start" button.

If performing a manual injection (with or without a data system):

- a. Press the "Prep Run" button to prepare the inlet for splitless injection.
 - b. When the GC becomes ready, inject 1 μL of the checkout sample and press the "Start" button.
9. The following chromatogram shows typical results for a new detector with new consumable parts installed. The Aldrin peak will be missing when using the Japanese sample 5183-0379.



To Check FPD+ Performance (Sample 5188-5953)

To check FPD+ performance, first check the phosphorus performance, then the sulfur performance.

Preparation

- Gather the following:
 - Evaluation column, HP-5 30 m × 0.32 mm × 0.25 µm (19091J-413)
 - FPD+ performance evaluation (checkout) sample (5188-5953), 2.5 mg/L (± 0.5%) methylparathion in isooctane
 - Phosphorus filter
 - Sulfur filter and filter spacer
 - 4-mL solvent and waste bottles or equivalent for autoinjector.
 - 2-mL sample vials or equivalent for sample.
 - Chromatographic-grade isooctane for syringe wash solvent.
 - Inlet and injector hardware. See [To Prepare for Chromatographic Checkout](#).
- Verify the following:
 - Chromatographic-grade gases plumbed and configured: helium as carrier gas, nitrogen, hydrogen, and air.
 - Empty waste vials loaded in sample turret.
 - 4-mL vial with diffusion cap filled with isooctane and inserted in Solvent A injector position.
- Replace consumable parts (liner, septum, traps, syringe, and so forth) as needed for the checkout. See [To Prepare for Chromatographic Checkout](#).
- Verify that the Lit Offset is set appropriately. Typically, it should be about 2.0 pA for the checkout method.
- Install the evaluation column.
- Set the oven, inlet, and detector to 250 °C and bake out for at least 15 minutes.
- Be sure to configure the column.

Phosphorus performance

- If it is not already installed, install the phosphorus filter.
- Create or load a method with the parameter values listed in [Table 30](#).

Table 30. FPD+ Checkout Conditions (P)

Column and sample	
Type	HP-5, 30 m × 0.32 mm × 0.25 µm (19091J-413)
Sample	FPD checkout (5188-5953)
Column mode	Constant flow
Column flow	6.5 mL/min
Split/splitless inlet	

Chromatographic Checkout

To Check FPD+ Performance (Sample 5188-5953)

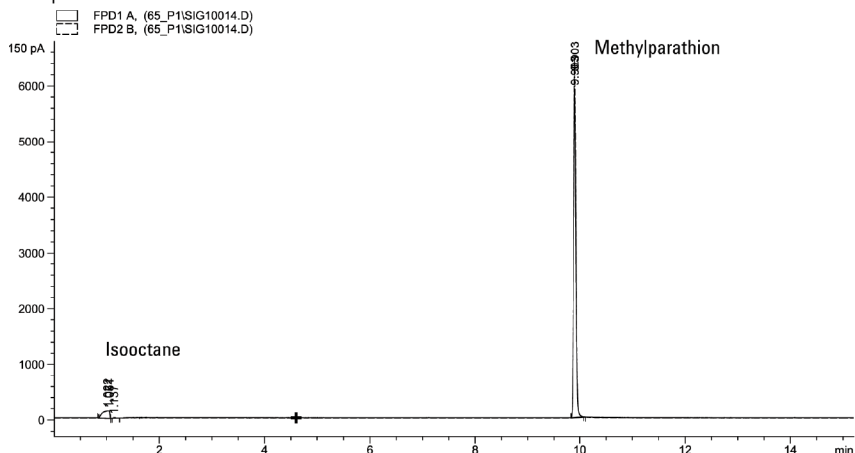
Temperature	180 °C Split/splitless
Mode	Splitless
Purge flow	60 mL/min
Purge time	0.75 min
Gas saver	ON
Gas saver flow	20 mL/min
Gas saver time	3 min
Purged packed column inlet	
Temperature	180 °C
Cool on-column inlet	
Temperature	Track oven
Septum purge	15 mL/min
Detector	
Heater and Aux Transfer Line Temperature	200 °C (On)
Emission Block Temperature	125 °C (On)
Hydrogen flow	60 mL/min (On)
Air (Oxidizer) flow	60 mL/min (On)
Carrier Gas Flow Correction	None (constant makeup and fuel flow)
Makeup flow	60 mL/min (On)
Makeup gas type	Nitrogen
Flame	On
Lit offset	Typically 2 pA
PMT voltage	On
Emission Block	125 °C
Oven	
Initial temp	70 °C
Initial time	0 min
Rate 1	25 °C/min
Final temp 1	150 °C
Final time 1	0 min
Rate 2	5 °C/min
Final temp 2	190 °C
Final time 2	7 min
ALS settings (if installed)	
Sample washes	2
Sample pumps	6
Sample wash volume	8 (maximum)
Injection volume	1 µL
Syringe size	10 µL
Solvent A pre washes	2
Solvent A post washes	2
Solvent A wash volume	8
Solvent B pre washes	0
Solvent B post washes	0
Solvent B wash volume	0

Chromatographic Checkout

To Check FPD+ Performance (Sample 5188-5953)

Injection mode (7693A)	Normal
Airgap Volume (7693A)	0.20
Viscosity delay	0
Inject Dispense Speed (7693A)	6000
PreInjection dwell	0
PostInjection dwell	0
Manual injection	
Injection volume	1 µL
Data system	
Data rate	5 Hz

- Ignite the FPD+ flame, if not lit.
- Display the signal output and monitor. This output typically runs around 10. Wait for the output to stabilize. This takes approximately 1 hour.
If the baseline output is too high:
 - Check column installation. If installed too high, the stationary phase burns in the flame and increases measured output.
 - Check for leaks.
 - Bake out the detector and column at 250 °C.
 - Wrong flows set for installed filter.
 If the baseline output is zero, verify the electrometer is on and the flame is lit.
- If using a data system, prepare the data system to perform one run using the loaded checkout method. Make sure that the data system will output a chromatogram.
- Start the run.
If performing an injection using an autosampler, start the run using the data system or press the "Start" button.
If performing a manual injection (with or without a data system):
 - Press the "Prep Run" button to prepare the inlet for splitless injection.
 - When the GC becomes ready, inject 1 µL of the checkout sample and press the "Start" button.
 - The following chromatogram shows typical results for a new detector with new consumable parts installed.



Sulfur performance

1. Install the sulfur filter and filter spacer.
2. Ignite the FPD+ flame if not lit.
3. Display the signal output and monitor. This output typically runs between 50 and 60 but can be as high as 70. Wait for the output to stabilize. This takes approximately 1 hour.

If the baseline output is too high:

- Check column installation. If installed too high, the stationary phase burns in the flame and increases measured output.
- Check for leaks.
- Bake out the detector and column at 250 °C.
- Wrong flows set for installed filter.

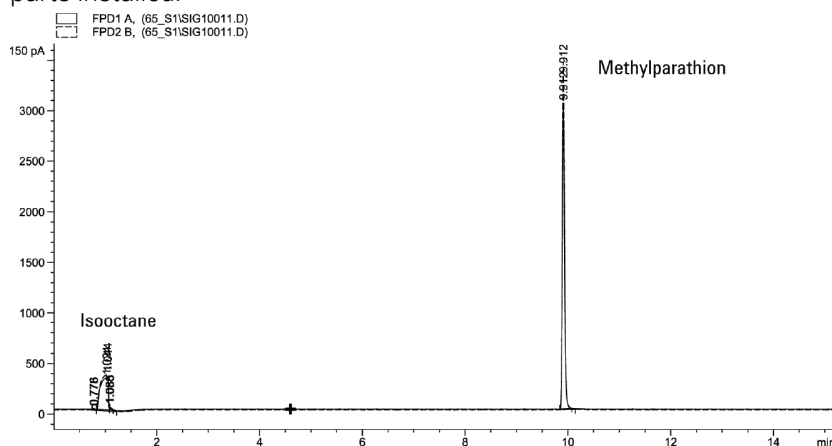
If the baseline output is zero, verify the electrometer is on and the flame is lit.

4. If using a data system, prepare the data system to perform one run using the loaded checkout method. Make sure that the data system will output a chromatogram.
5. Start the run.
If performing an injection using an autosampler, start the run using the data system or press the "Start" button.

If performing a manual injection (with or without a data system):

- a. Press the "Prep Run" button to prepare the inlet for splitless injection.
- b. When the GC becomes ready, inject 1 µL of the checkout sample and press the "Start" button.

6. The following chromatogram shows typical results for a new detector with new consumable parts installed.



To Check FPD+ Performance (Sample 5188-5245, Japan)

To verify FPD+ performance, first check the phosphorus performance, then the sulfur performance.

Preparation

- Gather the following:
 - Evaluation column, DB5 15 m × 0.32 mm × 1.0 μm (123-5513)
 - FPD performance evaluation (checkout) sample (5188-5245, Japan), composition: n-Dodecane 7499 mg/L (± 5%), Dodecanethiol 2.0 mg/L (± 5%), Tributyl Phosphate 2.0 mg/L (± 5%), tert-Butyldisulfide 1.0 mg/L (± 5%), in isooctane as solvent
 - Phosphorus filter
 - Sulfur filter and filter spacer
 - 4-mL solvent and waste bottles or equivalent for autoinjector.
 - 2-mL sample vials or equivalent for sample.
 - Chromatographic-grade isooctane for syringe wash solvent.
 - Inlet and injector hardware See [To Prepare for Chromatographic Checkout](#).
- Verify the following:
 - Chromatographic-grade gases plumbed and configured: helium as carrier gas, nitrogen, hydrogen, and air.
 - Empty waste vials loaded in sample turret.
 - 4-mL vial with diffusion cap filled with isooctane and inserted in Solvent A injector position.
- Replace consumable parts (liner, septum, traps, syringe, and so forth) as needed for the checkout. See [To Prepare for Chromatographic Checkout](#).
- Verify the lit offset is set appropriately. Typically, it should be about 2.0 pA for the checkout method.
- Install the evaluation column.
- Set the oven, inlet, and detector to 250 °C and bake out for at least 15 minutes.
- Configure the column.

Phosphorus performance

- If it is not already installed, install the phosphorus filter.
- Create or load a method with the parameter values listed in [Table 31](#).

Table 31. FPD+ Phosphorus Checkout Conditions

Column and sample	
Type	DB-5MS, 15 m × 0.32 mm × 1.0 μm (123-5513)
Sample	FPD checkout (5188-5245)

Chromatographic Checkout

To Check FPD+ Performance (Sample 5188-5245, Japan)

Column mode	Constant flow
Column flow	7.5 mL/min
Split/splitless inlet	
Temperature	250 °C
Mode	Splitless
Total purge flow	69.5 mL/min
Purge flow	60 mL/min
Purge time	0.75 min
Gas saver	ON
Gas saver flow	20 mL/min
Gas saver time	3 min
Purged packed column inlet	
Temperature	250 °C
Cool on-column inlet	
Temperature	Oven track
Septum purge	15 mL/min
Detector	
Heater and Aux Transfer Line Temperature	200 °C (On)
Emission Block Temperature	125 °C (On)
Hydrogen flow	60 mL/min (On)
Air (Oxidizer) flow	60 mL/min (On)
Carrier Gas Flow Correction	None (constant makeup and fuel flow)
Makeup flow	60 mL/min (On)
Makeup gas type	Nitrogen
Flame	On
Lit offset	Typically 2 pA
PMT voltage	On
Emission Block	125 °C
Oven	
Initial temp	70 °C
Initial time	0 min
Rate 1	10 °C/min
Final temp	105 °C
Final time	0 min
Rate 2	20 °C/min
Final temp 2	190 °C
Final time 2	7.25 min for sulfur
	12.25 min for phosphorus
ALS settings (if installed)	
Sample washes	2
Sample pumps	6
Sample wash volume	8 (maximum)
Injection volume	1 µL
Syringe size	10 µL

Chromatographic Checkout

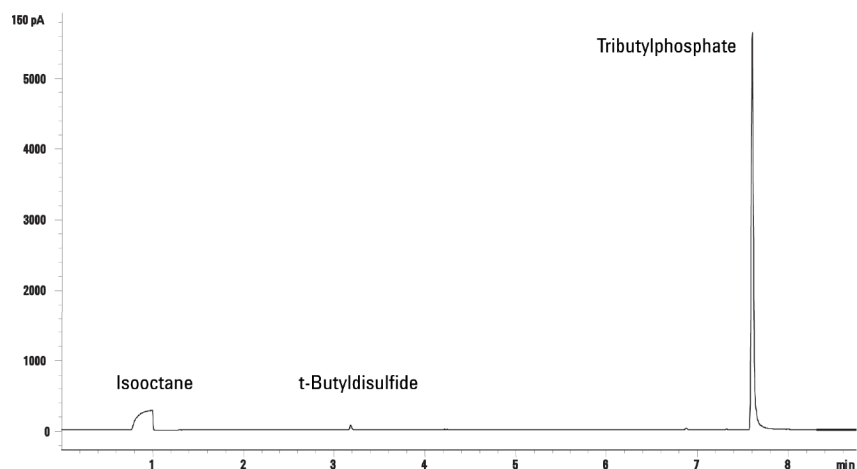
To Check FPD+ Performance (Sample 5188-5245, Japan)

Solvent A pre washes	2
Solvent A post washes	2
Solvent A wash volume	8
Solvent B pre washes	0
Solvent B post washes	0
Solvent B wash volume	0
Injection mode (7693A)	Normal
Airgap Volume (7693A)	0.20
Viscosity delay	0
Inject Dispense Speed (7693A)	6000
PreInjection dwell	0
PostInjection dwell	0
Manual injection	
Injection volume	1 µL
Data System	
Data rate	5 Hz

- Ignite the FPD+ flame, if not lit.
- Display the signal output and monitor. This output typically runs around 10. Wait for the output to stabilize. This takes approximately 1 hour.
If the baseline output is too high:
 - Check column installation. If installed too high, the stationary phase burns in the flame and increases measured output.
 - Check for leaks.
 - Bake out the detector and column at 250 °C.
 - Wrong flows set for installed filter.
 If the baseline output is zero, verify the electrometer is on and the flame is lit.
- If using a data system, prepare the data system to perform one run using the loaded checkout method. Make sure that the data system will output a chromatogram.
- Start the run.
If performing an injection using an autosampler, start the run using the data system or press the "Start" button.
If performing a manual injection (with or without a data system):
 - Press the "Prep Run" button to prepare the inlet for splitless injection.
 - When the GC becomes ready, inject 1 µL of the checkout sample and press the "Start" button.
- The following chromatogram shows typical results for a new detector with new consumable parts installed.

Chromatographic Checkout

To Check FPD+ Performance (Sample 5188-5245, Japan)

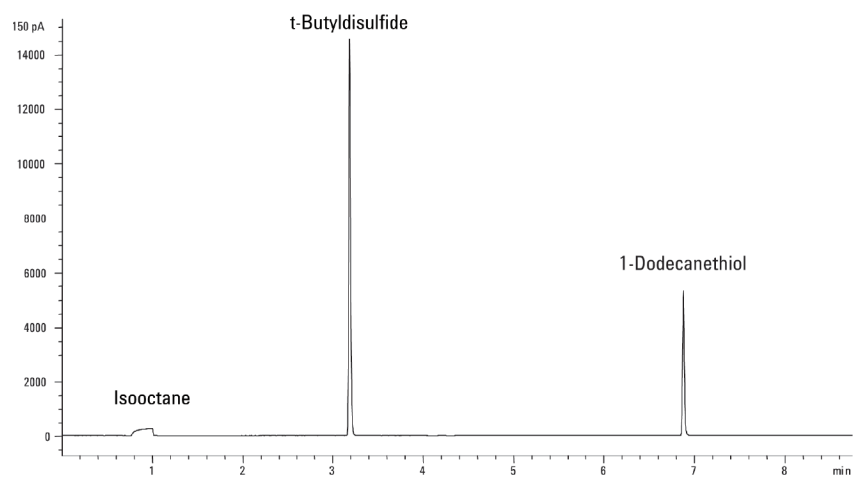


Sulfur performance

1. Install the sulfur filter.
2. Ignite the FPD+ flame, if not lit.
3. Display the signal output and monitor. This output typically runs between 50 and 60 but can be as high as 70. Wait for the output to stabilize. This takes approximately 2 hours.
If the baseline output is too high:
 - Check column installation. If installed too high, the stationary phase burns in the flame and increases measured output.
 - Check for leaks.
 - Bake out the detector and column at 250 °C.
 - Wrong flows set for installed filter
 If the baseline output is zero, verify the electrometer is on and the flame is lit.
4. If using a data system, prepare the data system to perform one run using the loaded checkout method. Make sure the data system will output a chromatogram.
5. Start the run.
If performing an injection using an autosampler, start the run using the data system or press the "Start" button.
If performing a manual injection (with or without a data system):
 - a. Press the "Prep Run" button to prepare the inlet for splitless injection.
 - b. When the GC becomes ready, inject 1 μ L of the checkout sample and press the "Start" button.
6. The following chromatogram shows typical results for a new detector with new consumable parts installed.

Chromatographic Checkout

To Check FPD+ Performance (Sample 5188-5245, Japan)



FPD+ and ECD Unit Conversion Factors

Conversion factors for the FPD+

Conversion factor for the ECD

Using the Conversion Factors

References

The GC complies with the following company standard: Q31/0115000033C005-2016-02.

China Metrology testing of the 8860B GC is performed in accordance with company standard Q31/0115000033C005-2016-02. This chapter provides information and techniques to correctly identify noise and drift when testing an FPD+ or ECD.

FPD+ and ECD Unit Conversion Factors

At the time of publication, China Metrology testing requires noise and drift metrics as shown below:

Detector	Reporting units
FID	A
TCD	mV
NPD	A
FPD+	A
ECD	mV or Hz

However, data collection is required to come through the digital output available through the GC and data system. For the FID, NPD, and TCD, the data system provides data in the desired reporting units. However, for the ECD and FPD+, Agilent reports output to its data systems in “display units,” (DU). This section describes how to accurately convert/scale the FPD+ and ECD digital results to make them consistent with Chinese Metrology requirements.

The conversion factors for the FPD+ and ECD take the display unit output from the Agilent data system digital path to an absolute value for current or voltage. Agilent developed the conversion factors empirically, based on measurements from a single system that simultaneously output both the digital and analog data. The conversion factors also incorporate:

- The scaling applied to analog versus digital signals
- An analog signal range setting of 5 (2⁵) at the GC
- The unique filtering applied by the 35900 ADC
- The differences in bandwidth (BW) associated with the GC digital channel (5 Hz) and the 35900 ADC analog path (3 Hz)

The differences in channel bandwidth between the analog and digital signal paths can be taken into consideration as follows:

$$BW = 35900 \text{ ADC path} / \text{GC digital path} = \sqrt{(3 \text{ Hz} / 5 \text{ Hz})} = 0.7$$

Conversion factors for the FPD+

For the FPD+, the conversion factor is the same whether the phosphorus or sulfur filter is used:

$$\text{FPD+: } 1 \text{ DU} = 150 \text{ pA}$$

Conversion factor for the ECD

$$1 \text{ DU} = 1 \text{ Hz}$$

Using the Conversion Factors

To use the conversion factors, multiply the ASTM noise reported from the Agilent data system for the GC digital signal path by the appropriate conversion factor.

For example, consider applying the FPD+ and ECD conversion factors to a statistical sampling of digital noise performances measured for both detectors at Agilent:

Average FPD+ ASTM noise, DU^{1,2}: 1.54

Average ECD ASTM noise, DU³: 0.16

Applying the conversion factors:

FPD+: $1.54 \text{ DU} \times 150 \text{ pA/1DU} = 231 \text{ pA}$

ECD: $0.16 \text{ DU} \times 1 \text{ Hz/1 DU} = 0.16 \text{ Hz}$

¹Agilent data for FPD+ noise in this example represents sulfur mode only.

²Data collected for purposes of comparison should be acquired with a nominal FPD+ offset of < 100 DU in sulfur mode and < 20 DU in phosphorus mode and at a data rate of 5 Hz.

³Data collected for purposes of comparison should be acquired with a nominal ECD baseline at or below 150 DU and at a data rate of 5 Hz.

References

"Calculation of Performance Factors for the HP 6890 Gas Chromatograph Using Different Data Handling Devices" Agilent Technologies publication 5964-0282E.

"Calculation of Performance Factors for the HP 6890 Gas Chromatograph Using Different Data Handling Devices" Agilent Technologies publication 5091-9207E.

"Calculation of Performance Factors for the HP 6890 Gas Chromatograph Using Different Data Handling Devices" Agilent Technologies publication 5965 8901E.

The terms in **Table 32** are used in discussing this product. They are collected here for your convenience.

Table 32. Terms

Term	Definition
ADC	Analog to digital converter
ALS	Automatic liquid sampler
AS	Autosampler
BCD	Binary coded decimal
COC	Cool on column inlet
DHCP	Dynamic host creation protocol
ECD	Electron capture detector
ELVDS	Port for external communications with Agilent MSDs
EMF	Early maintenance feedback
EPC	Electronic pneumatics control
EPR	Electronic pneumatics regulation
FID	Flame ionization detector
FPD+	Flame photometric detector plus
GC	Gas chromatograph
HS	Headspace sampler
LAN	Local area network
LUI	Local User Interface
LVDS	Low-voltage differential signaling
MS	Mass spectrometer
MSD	Mass selective detector
NCD	Nitrogen chemiluminescence detector
NPD	Nitrogen-phosphorus detector
NTP	Normal temperature and pressure (25 °C and 1 atmosphere)
PCI	Packed column inlet
PCM	Pressure Control Module
PID	Proportional integral and differential
PP	Purged packed inlet
PTFE	Polytetrafluoroethylene
SCD	Sulfur chemiluminescence detector
SSL	Split/splitless inlet
TCD	Thermal conductivity detector

www.agilent.com

© Agilent Technologies, Inc. 2026

G2790-90022



May 2026
Printed in USA

