



OpenLab ChemStation

## Instrument support



# Notices

## Manual Part Number

D0013761 Rev. B  
Edition 09/2022

## Copyright

© Agilent Technologies, Inc. 2010-2020, 2022

No part of this manual may be reproduced in any form or by any means (including electronic storage and retrieval or translation into a foreign language) without prior agreement and written consent from Agilent Technologies, Inc. as governed by United States and international copyright laws.

Agilent Technologies, Inc.  
5301 Stevens Creek Blvd.  
Santa Clara, CA 95051  
USA

## Software Revision

This guide is valid for the LTS 01.11 revision or higher of Agilent OpenLab ChemStation, until superseded.

## 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.

## Table of Contents

<b>1</b>	<b>General Driver Information</b>	<b>5</b>
	Operating Systems	5
	Instrument Driver types	5
	Instrument Driver Revisions	6
	Firmware Revisions	6
	Instrument Communication Information	7
<b>2</b>	<b>Agilent HPLC and LC Instrument Compatibility</b>	<b>8</b>
	Recommended Firmware for Agilent HPLC/LC Modules	8
	Supported Agilent HPLC/LC Modules	10
	Agilent LC Sampling Systems	10
	CTC PAL Autosamplers (Agilent and CTC) with LC	12
	Agilent LC – Column Compartments and Valve Thermostat Cluster	12
	Agilent LC – Pumps	13
	Agilent LC – Detectors	14
	Agilent LC – Valves	16
	Agilent LC – Fraction Collectors	17
	Agilent LC – Other Module types	18
	Agilent 218/SD1 Prep LC driver (formerly Varian Prep LC)	18
	Agilent 1120 Compact LC Systems	19
	Agilent 1220 Infinity LC Systems	19
<b>3</b>	<b>Agilent LC/MS Module Compatibility</b>	<b>20</b>
<b>4</b>	<b>Agilent CE Module Compatibility</b>	<b>22</b>
	Agilent Capillary Electrophoresis	22
<b>5</b>	<b>Agilent Analog/Digital Converter Information</b>	<b>23</b>
	Agilent A/D Converter	23
	Other	23

<b>6</b>	<b>Agilent GC Instrument Compatibility</b>	<b>24</b>
	Agilent GC support	24
	Agilent Micro GC support	25
	Other, unsupported instruments	25
	GC Sampler, Autosampler and Tray Compatibility	26
	7650 ALS hardware:	26
	7693 ALS hardware	26
	7683B ALS hardware	26
	7683A ALS hardware:	27
	Other, unsupported instruments	27
	Headspace support with GC	28
	Control of PAL sampler with GC	28
<b>7</b>	<b>Compatibility with Waters instruments</b>	<b>29</b>
	M8505AA Waters LC Control	29
	Compatibility with Waters ACQUITY and ACQUITY H-Class	29
	Compatibility with Waters Alliance instruments:	30
<b>8</b>	<b>OpenLab ChemStation Workstation VL</b>	<b>31</b>
	Overview	31
	Supported Instruments with OpenLab ChemStation Workstation VL	32
	GC instrument	32
	Agilent 1260 Infinity II LC instrument	32
	Agilent 1220 Infinity II LC system	34
	Agilent 1120 Compact LC	35
<b>9</b>	<b>Classic Driver Support Information</b>	<b>36</b>

# 1

## General Driver Information

### Operating Systems

All instrument drivers are supported on the following OS unless otherwise noted

- Windows 11 21H2 Enterprise and Professional (64-bit)
- Windows 10 21H2 Enterprise and Professional (64-bit)

For more detail and support information on server operating systems, please see the [OpenLab ChemStation Requirements guide](#).

### Instrument Driver types

Agilent OpenLab drivers are based on the RC.Net architecture. Information in this guide refers to this driver type, unless otherwise noted. RC.Net 3.6.1 is integral part of OpenLab CDS ChemStation LTS.

#### NOTE

Instrument driver versions must always match in a network system

For single quad Mass Spectrometer, only “Classic Plus” drivers are available with OpenLab ChemStation. The Classic Plus driver types can be combined with RC.Net drivers via their CIC ports.

# Instrument Driver Revisions

The following Agilent instrument drivers are part of the OpenLab ChemStation LTS 01.11 software package and are installed with the software:

**Table 1** Drivers provided with OpenLab ChemStation LTS 01.11

RC.Net Instrument Driver	Driver Software Revision
Agilent LC	3.5
Agilent ELSD	1.8
Agilent GC	3.7
Agilent Micro GC	2.3
Agilent 35900E A/D Converter	2.3

If you are looking for information on the build number of the included drivers, please refer to *Driver Revisions*, accessible from the **Resource Platform > Administration > Configuration**.

## Firmware Revisions

Agilent recommends using the most recent firmware revisions to provide the highest level of system capability.

Agilent and other vendors may release RC.net drivers independent of the OpenLab ChemStation releases. Agilent driver software is forward compatible regarding firmware, i.e., the firmware can be updated without the need of updating the driver or the Chromatography Data System.

To check the firmware of your instrument, do the following:

In the ChemStation **Method and Run Control** view,

### LC Instruments:

In the menu, go to **Instrument > Revisions & Serial #s**

### GC Instruments:

- 1 In the menu, go to **Instrument > Instrument Configuration**.
- 2 Select the GC instrument under **Selected modules**,
- 3 Click the **configure** button, then select the configuration tab.

## Instrument Communication Information

### NOTE

On Agilent's *Analytical Instrument Controllers* (AICs), OpenLab ChemStation exclusively supports instruments and modules connected via LAN.

GPIB, RS-232, USB, or any converters are not supported on AICs.

---

The G2403A LAN/RS232 converter Dudley box is not supported with OpenLab ChemStation LTS

## 2

# Agilent HPLC and LC Instrument Compatibility

OpenLab ChemStation can control all Agilent LC modules in support using LC & CE driver release 3.5 or higher.

Starting with the LC & CE Drivers release 3.5 you may restrict the access to menu items that affect the instrument configuration to users with the ChemStation permission **Modify instrument configuration**.

The tables in this chapter list the supported LC modules. Please note that some of the driver features such as valve-thermostat clusters or new temperature control modes require current firmware versions.

## Recommended Firmware for Agilent HPLC/LC Modules

Agilent recommends using current LC firmware sets with your OpenLab ChemStation. The latest LC firmware sets are 7.01, 6.50, and 6.30.

Some of the modules are supported with firmware revisions lower than the ones available in the current firmware set. Please see the respective Release Note for Agilent LC and CE drivers for more information.

**Table 2** Recommended firmware per LC instrument series

Device	Recommended Firmware
Agilent 1100 Series, 1200 Series and 1200 Infinity	A.07.01 or later
<b>NOTE:</b> 1100 Series Modules are supported on "best effort" basis.	
Agilent 1200 Series, 1200 Infinity, and 1120 Compact LC	B.07.36 or later
Agilent 1200 Infinity, hosted modules	C.07.30 or later
Agilent 1260/1290 Infinity II	D.07.36 or later

## Agilent HPLC and LC Instrument Compatibility

Please visit the Agilent site [LC Firmware News and Downloads](#) to download current Agilent LC firmware.

Firmware Set 7.xx can be downloaded from <http://www.agilent.com/en-us/firmwareDownload?whid=99818>.

For more information, please see the respective LC driver release notes. The Release Notes for Agilent LC and CE drivers (ReleaseNotes\_Driver\_LC-CE.pdf) is available on your ChemStation software media.

### NOTE

LC and CE firmware revisions are grouped into sets for each module or system. Firmware sets include just the latest firmware of each module.

A firmware update within set A/B/C/D.07.01 is required for all modules in that stack, not only new modules, as for example the fraction collector uses new detector features.

Please note that the LC Driver releases are independent from ChemStation Releases. As a rule, OpenLab ChemStation is forward compatible with new driver software.

### Supported Agilent HPLC/LC Modules

The listed Agilent LC modules are supported to be controlled with OpenLab ChemStation unless stated otherwise.

The Agilent LC-drivers are backwards compatible.

Modules with identical product numbers are supported, even if the tables below list only the name of the current model version. 1100 Series models are supported on a best effort basis only.

For more detail, please see the driver Release Notes of the driver revision you are using.

#### Agilent LC Sampling Systems

Product Number	Module Name	Comment
G1313A	1100 Series Autosampler	Best effort
G1327A	Thermostatted Autosampler	G1329A with G1330A/B (Thermostat for ALS)
G1329A	1200 Series Standard Autosampler	Supported
G1329B	1260 Infinity Standard Autosampler	Supported
G1367A	1100 Series Well-Plate Sampler	Best effort
G1367B	1200 Series High Performance Autosampler	Supported
G1367C	1200 Series High Performance Autosampler SL	Supported
G1367D	1200 Series High Performance Autosampler SL+	Supported
G1367E	1260 Infinity High Performance Autosampler	Supported
G1368A	Thermostatted Well plate sampler	G1367A with G1330A/B
G1373A	Purge Kit <sup>1</sup>	Supported
G1377A	1260 Infinity High Performance Micro Autosampler	Supported

<sup>1</sup> Includes G1156A 6Pos/7Port valve, cabinet top kit with rail, accessories and ChemStation SW, for use with 1200 Wellplate sampler

## Agilent LC Sampling Systems

Product Number	Module Name	Comment
G1378A	Thermostatted micro well-plate sampler	Supported G1377A with G1330A/B
G1389A	1100 Micro-Scale Autosampler	Best effort
G2258A	1260 Infinity Dual-Loop Autosampler	Supported
G2260A	1260 Infinity Preparative Autosampler (High flow)	Supported
G2261A	Thermostatted High Flow ALS	Supported G2260A with G1330A/B
G4226A	1290 Infinity Autosampler	Supported
G4303A	1260 Infinity SFC Standard Autosampler	Supported
G4767A	1260 Infinity II SFC Multisampler	Supported
G5667A	1260 Infinity Bio-Inert High-Performance Autosampler	Supported
G5668A	1260 Infinity II Bio-Inert Multisampler	Supported
G7129A	1260 Infinity II Vialsampler	Supported
G7129B	1290 Infinity II Vialsampler	Supported
G7129C	1260 Infinity II Vialsampler	Supported
G7157A	1260 Infinity II Preparative Autosampler	Supported
G7158B	1290 Infinity II Open-bed Sampler / Collector	Supported Instrument configuration shows two modules: G7169B (injector) and G7159B (fraction collector)
G7167A	1260 Infinity II Multisampler	Supported
G7167B	1290 Infinity II Multisampler	Supported
G7169B	Agilent 1290 Infinity II Open-bed Sampler / Fraction Collector - Sampler Driver	Supported

## CTC PAL Autosamplers (Agilent and CTC) with LC

Product Number	Module Name	Minimum required Revision	
		CTC Firmware	Software
G4277A	Agilent 1290 Infinity LC Injector HTS	4.3.0	B.01.08
G4278A	Agilent 1290 Infinity LC Injector HTC	4.3.0	B.01.08
G4270-CTC	CTC Analytics HTC PAL Injection System1	2.5.2 (for legacy mainboard) 2.6.8 (for legacy mainboard)	B.01.08 Cycle Parameters in English
G4271-CTC	CTC Analytics HTS PAL Injection System1	4.3.0 (for new XT mainboard)	

## Agilent LC – Column Compartments and Valve Thermostat Cluster (VTC)

Product Number	Module Name	Comment
G1316A	1260 Infinity Thermostatted Column Compartment	Supported
G1316B	1200 Thermostatted Column Compartment SL	Supported
G1316C	1290 Infinity Thermostatted Column Compartment	Supported
G7116A	1260 Infinity II Multicolumn Thermostat (MCT)	Supported
G7116B	1290 Infinity II Multicolumn Thermostat (MCT)	Thermal Equilibration Device
G7130A	Integrated Column Compartment ICC (option for G7129 A/B)	Supported
TCC cluster	Up to three G1316C with integrated 8pos/9port valves (products G4230A/B) can be clustered; minimum two G1316C TCCs; the third TCC can be a G1316A or B	Supported Exception: No support for new 2pos/6port or 2pos/10port valves
VTC	The Valve Thermostat Cluster is a combination of G7116B, G1170A and G1316C as valve or column hosts and G1316A/B and G7130A as column hosts.	Supported

## Agilent LC – Pumps

Product Number	Module Name	Comment
G1310A	1200 Series Isocratic Pump	Supported
G1310B	1260 Infinity Isocratic Pump	Supported
G1311A	1200 Series Quaternary Pump <sup>2</sup>	Supported
G1311B	1260 Infinity Quaternary Pump <sup>1</sup>	Supported
G1311C	1260 Infinity Quaternary Pump VL <sup>1</sup>	Supported
G1312A	1260 Infinity Binary Pump <sup>1</sup>	Supported
G1312B	1260 Infinity Binary Pump SL <sup>1</sup>	Supported
G1312C	1260 Infinity Binary Pump VL <sup>1</sup>	Supported
G1361A	1260 Infinity Preparative Pump <sup>3</sup>	Supported
G1376A	1260 Infinity Capillary Pump	Supported
G2226A	1260 Infinity Nanoflow Pump	Supported
G4204A	1290 Quaternary Pump <sup>1</sup>	Supported
G4220A	1290 Infinity Binary Pump <sup>1</sup>	Supported
G4220B	1290 Infinity Binary Pump VL <sup>1</sup>	Supported
G4302A	1260 Infinity SFC Binary Pump <sup>1</sup>	Supported
G4782A	1260 Infinity II SFC Binary Pump	Supported
G5611A	1260 Infinity Bio-inert Quaternary Pump <sup>1</sup>	Supported
G5654A	1260 Infinity II Bio-inert Pump <sup>1</sup>	Supported
G7104A	1290 Infinity II Flexible Pump	Supported
G7104C	1260 Infinity II Flexible Pump	Supported
G7110B	1260 Infinity II Isocratic Pump	Supported
G7111A	1260 Infinity II Quaternary Pump VL <sup>1</sup>	Supported

<sup>2</sup> 1200 Pump Valve Clusters are supported with G1311, G1312, G4220, G4204, G4302 or (G5611 + 5067-4147). Pump valve clusters are possible with up to 2 valves of type G1160A and/or G1170A,

<sup>3</sup> Prep Pump cluster supported with up to four G1361A per cluster

## Agilent HPLC and LC Instrument Compatibility

Product Number	Module Name	Comment
G7111B	1260 Infinity II Quaternary Pump <sup>1</sup>	Supported
G7112B	1260 Infinity II Binary Pump <sup>1</sup>	Supported
G7120A	1290 Infinity II High Speed Pump	Supported
G7161A	1260 Infinity II Preparative Binary Pump	Supported
G7161B	1290 Infinity II Preparative Binary Pump	Supported

### Agilent LC – Detectors

Product Number	Module Name	Comment
<b>VWD</b>		Supported
G1314A	1100 Series Variable Wavelength Detector	Supported
G1314B	1260 Infinity Variable Wavelength Detector VL	Supported
G1314C	1260 Infinity Variable Wavelength Detector VL+	Supported
G1314D	1200 Series Variable Wavelength Detector	Supported
G1314E	1290 Infinity Variable Wavelength Detector	Supported
G1314F	1260 Infinity Variable Wavelength Detector	Supported
G7114A	1260 Infinity II Variable Wavelength Detector	Supported
G7114B	1290 Infinity II Variable Wavelength Detector	Supported
<b>MWD</b>		
G1365A	1100 Series Multiple Wavelength Detector	Best effort
G1365B	1200 Series Multiple Wavelength Detector	Supported
G1365C	1260 Infinity Multiple Wavelength Detector	Supported
G1365D	1260 Infinity Multiple Wavelength Detector VL	Supported
G7165A	1260 Infinity II Multiple Wavelength Detector	Supported

## Agilent HPLC and LC Instrument Compatibility

Product Number	Module Name	Comment
<b>FLD</b>		
G1321A	1100 / 1200 Fluorescence Detector	Supported
G1321B	1260 Infinity Fluorescence Detector Spectra	Supported
G1321C	1260 Infinity Fluorescence Detector	Supported
G7121A	1260 Infinity II Fluorescence Detector	Supported
G7121B	1260 Infinity II Fluorescence Detector Spectra	Supported
<b>DAD</b>		
G1315A	1100 Series Diode Array Detector	Supported
G1315B	1200 Series Diode Array Detector	Supported
G1315C	1200 Series Diode Array Detector VL+	Supported
G1315D	1200 Series Diode Array Detector VL	Supported
G4212A	1290 Infinity Diode Array Detector	Supported
G4212B	1260 Infinity Diode Array Detector	Supported
G4212A/B HDR-DAD Cluster	2x G4212A or 2x G4212B or a combination of 1x G4212A and 1x G4212B	No add-on software required. HDR function included in RC.net drivers
G7115A	1260 Infinity II Diode Array Detector WR	Supported
G7117A	1290 Infinity II Diode Array Detector FS	Supported
G7117B	1290 Infinity II Diode Array Detector	Supported
G7117C	1260 Infinity II Diode Array Detector HS	Supported
G7117A/B HDR-DAD Cluster	2x G7117A or 2x G7117B or a combination of 1x G7117A and 1x G7117B	No add-on software required. HDR function included in RC.net drivers
<b>RID</b>		
G1362A	1100/1200 Infinity Refractive Index Detector	Supported
G7162A	1260 Infinity II Refractive Index Detector	Supported
G7162B	1290 Infinity II Refractive Index Detector	Supported

## Agilent HPLC and LC Instrument Compatibility

Product Number	Module Name	Comment
<b>ELSD</b>		
G4218A	1260 Infinity Evaporative Light Scattering Detector	Not supported
G4260A	1260 Infinity 380 Evaporative Light Scattering Detector	Supported English only NOT localized for Japanese and Chinese ChemStation
G4261A	1290 Infinity 385 Evaporative Light Scattering Detector	
G4260B	1260 Infinity II Evaporative Light Scattering Detector	
G4261B	1290 Infinity Evaporative Light Scattering Detector	RS232, not supported on AIC.
G7102A	1290 Infinity II Evaporative Light Scattering Detector	Supported

### Agilent LC – Valves

#### NOTE

The below table lists just selected valves, valve drives and clusters. Please refer to the LC&CE Driver Release Note of your driver version for a complete list of supported Agilent valves.

Product Number	Module Name	Comment
G1156A	1200 6-Position / 7 Port Valve (400 bar)	Not configurable in ChemStation, requires an ALS
G1157A	1200 2-Position / 10 Port Valve	Supported
G1158A	1200 2-Position / 6 Port Valve	Supported
G1158B	1200 2-Position / 6 Port Valve (600bar)	Supported
G1159A	1200 6-Position Selection Valve	Supported
G1160A	1200 12-Position/13-Port Valve	Supported
G1162A	1200 2-Position/ 6 Port Micro Valve	Supported
G1163A	1200 2-Position/ 10 Port Micro Valve	Supported
G1170A	1290 Infinity II Valve Drive	Supported
G1323B	Local User Interface	Supported

## Agilent HPLC and LC Instrument Compatibility

Product Number	Module Name	Comment
G2229A	Nano Flow LC System <sup>4</sup>	Supported
G4208A	Instant System Pilot	Supported
G4734A	Hosted valve: 6-Column-selector	requires valve host G1170A or G7116A/B
G9322A	1260 Infinity II Clustering Valve	Supported
G4230A/B	Valve Kits	Supported
G4234A		Supported
G4234C		Supported
G4240A	Chip Cube interface	Not supported
Valve Thermostat Cluster	The Valve Thermostat Cluster is a combination of G7116B, G1170A and G1316C as valve or column hosts and G1316A/B and G7130A as column hosts.	Supported

### Agilent LC – Fraction Collectors

Product Number	Module Name	Comment
G1364A	1100 Series Automatic Fraction Collector	Best effort
G1364B	1260 Infinity Preparative-Scale Fraction Collector	Supported
G1364C	1260 Infinity Analytical-Scale Fraction Collector	Supported
G1364D	1260 Infinity Micro-Scale Fraction Collector/Spotter	Supported
G1364E	1260 Infinity II Preparative-Scale Fraction Collector	Supported
G1364F	1260 Infinity II Analytical-Scale Fraction Collector	Supported
G5664A	1260 Infinity Bio-Inert Fraction Collector	Supported
G5664B	1260 II Infinity Bio-Inert Fraction Collector	Supported
G7159B	1290 Infinity II Preparative Open-Bed Fraction Collector	Supported
G7166A	1260 Infinity II Preparative Valve-Based Fraction Collector	Supported

<sup>4</sup>: Includes nanoflow pump, micro degasser, micro well-plate sampler, handheld controller (G1377A + G2226A)

## Agilent HPLC and LC Instrument Compatibility

Product Number	Module Name	Comment
Fraction Collector Cluster	Any combination of up to three G1364A/B/C or G5664A plus a fourth G1364A/B/C or G5664A for recovery can be clustered. controlling multiple un-clustered Fraction Collectors is not supported. G1364 E /F can be clustered with a G7166A or the same module type for recovery collection. G7159 can be clustered with a G7166A for recovery	

### Agilent LC – Other Module types

Product Number	Module Name	Comment
G1390A	1100 Series Universal Interface Box (UIB)	Supported
G1390B	InfinityLab Universal Interface Box	Supported
G2254A	1260 Automation Interface	Supported
G4227A	1290 Infinity Flexible Cube	Supported
G4240A	1260 Chip Cube MS Interface	Not supported
G4301A	1260 Infinity Analytical SFC System	Supported
G7170B	1290 Infinity II MS Flow Modulator	Supported

### Agilent 218/SD1 Prep LC driver (formerly Varian Prep LC)

Product Number	Module Name	Communication	Comment
Agilent 210	Pump	RS232	Not supported
Agilent 218	Pump	RS232	Not supported
Agilent SD-1	Pump	RS232	Not supported
Agilent 325	UV/VIS Detector	LAN	Not supported
Agilent 410	Autosampler	RS232	Not supported
Agilent 440	Fraction Collector	RS232	Not supported

## Agilent 1120 Compact LC Systems

Product Number	Module Name	Comment
G4286A	1120 Compact LC, Isocratic pump	Supported
G4287A	1120 compact LC. Isocratic with Oven and ALS	Supported
G4288A	1120 Compact LC, Gradient	Supported
G4289A	1120 Compact LC, Gradient with oven	Supported
G4290A	1120 Compact LC, Gradient with oven and ALS	Supported

## Agilent 1220 Infinity LC Systems

Product Number	Module Name	Comment
G4286B	1220 Infinity LC System Isocratic, Man. Inj., VWD, 600 bar	Supported
G4287B	1220 Infinity LC Isocratic, ALS, TCC, VWD, 600 bar	Supported
G4288B	1220 Infinity LC Gradient, Man. Inj., VWD, 600 bar	Supported
G4289B	1220 Infinity LC Gradient, ALS, TCC, VWD, 600 bar	Supported
G4290B	1220 Infinity LC Gradient, ALS, Man. Inj., TCC, VWD, 600 bar	Supported
G4291B	1220 Infinity LC Isocratic, Man. Inj., TCC, VWD, 600 bar	Supported
G4292B	1220 Infinity LC Isocratic, ALS, VWD, 600 bar	Supported
G4293B	1220 Infinity LC Gradient, ALS, VWD, 600 bar	Supported
G4294B	1220 Infinity LC Gradient, ALS, TCC, DAD, 600 bar	Supported
G4288C	1220 Infinity LC System VL, Gradient, Man. Inj. VWD, 400 bar	Supported
G4289C	1220 Infinity LC System VL, Gradient, TCC, Man. Inj. VWD, 400 bar	Supported
G4290C	1220 Infinity LC System VL, Gradient, ALS, TCC, VWD, 400 bar	Supported
G4293C	1220 Infinity LC System VL, Gradient, ALS, VWD, 400 bar	Supported
G4294B	1220 LC Infinity Gradient System	Supported

## 3 Agilent LC/MS Module Compatibility

Single Quadrupole Modules listed in the table below are supported, unless stated otherwise.

Requires Keysight I/O libraries.

Product Number	Module Name	Comment
G6110A G6120A G6130A G6140A	6100 Series Single Quadrupole LC/MS System	Not supported
G6120B	6100 Series Single Quadrupole LC/MS System	Supported
G6125B	LC/MSD	Supported
G6130B	6100 Series Single Quadrupole LC/MS System	Supported
G6135B	LC/MSD XT	Supported
G6150B	6100 Series Single Quadrupole LC/MS System	Supported
G6120C G6130C G6125C G6135C	6100 Series Single Quadrupole LC/MS System LC/MSD LC/MSD XT	Not supported need OpenLab CDS v2.x
G6160A	LC/MSD iQ Single Quadrupole LC/MS System	Not supported needs OpenLab CDS v2.3 or higher

### LC Single-Quad Sources

G1947A G1971A G1948A	APCI LC SQ Source APPI LC SQ Source ESI LC SQ Source	Not supported
G1947B G1971B G1948B	APCI LC SQ Source APPI LC SQ Source ESI LC SQ Source	Supported

## Agilent LC/MS Module Compatibility

Product Number	Module Name	Comment
G1978B	Multimode Source	Supported
G4240A	Chip Cube Interface	Not supported
G1968C	Active Splitter Kit	Supported on English ChemStation only
	CTC Autosampler	Classic Driver only. Releases B.01.01 or later are NOT compatible with the Cycle Composer software from CTC. Cannot co-reside. Not supported on Japanese and Chinese ChemStation
G1979A	SIDEBox	Not supported
G1958B	Agilent Jet Stream for Single Quad	Supported
G1951A	Analog Output Accessory	Supported

## 4 Agilent CE Module Compatibility

### NOTE

When OpenLab ChemStation is used to control an Agilent 3D Capillary Electrophoresis system, no additional instruments are supported on the PC system.

When OpenLab ChemStation is used to control the Agilent LC/MS, in conjunction with the Agilent 3D Capillary Electrophoresis system, no additional instruments are supported on the PC system (the one PC controls only the CE and MS system).

CE/MS systems are supported with OpenLab ChemStation when configured as CE/MS system, workstation, and networked workstation only. Co-installation is possible, but no Co-execution.

### Agilent Capillary Electrophoresis

Product Number	Module Name	Comment
G1600 CE G1600 CE/MS	CE and CE/MS families	Not supported
G7150A G7151A	G7100 Capillary Electrophoresis II Diode Array Detector for CE II	Configure using Instrument Type "Agilent 7100 CE System."
35900E	A/D converter	RC.net only, LAN only!
G1607B	CE-ESI-MS Sprayer kit AJS compatible	Only when configured as CE/MS system, workstation, and networked workstation.

## 5 Agilent Analog/Digital Converter Information

### Agilent A/D Converter

Module Name	Comment
35900C/D	Not supported
35900E 35900E Series II	Supported

### Other

Module Name	Comment
SS420x A/D interface	Not supported

## 6 Agilent GC Instrument Compatibility

Agilent recommends using the latest firmware revision to provide the highest level of system capability. See current Hardware Service Notes for latest firmware available.

### NOTE

Communication is handled via LAN. GCs using GPIB must update to LAN.

### Agilent GC support

Product Number	Module Name	Comment
G3950A G3952A G3953A	Intuvo 9000 GC system	Supported (7683 ALS family not supported)
G3540A G3542A G3543A G3545A	8890 Series GC System	Supported (7683 ALS family not supported)
G2790A	8860 Series GC System	Supported (7683 ALS family not supported)
G3440A G3445A	7890A	Supported
G3440B G3442B G3443B G3445B	7890B	Supported
G4350A G4350B	7820A	Supported (Backflush and Easy SamplePrep are not supported with 7820)
G1530N G1540N	6890N	Supported (Manual inlets and manual detectors are not supported)

Product Number	Module Name	Comment
G1530A G1540A	6890A 6890Plus	Supported (Manual inlets and manual detectors are not supported)
G2629A	6850 Handheld Controller	Not Supported Out of Support
G2630A G2630B	6850	Supported

### Agilent Micro GC support

Product Number	Module Name	Comment
G3588A	990 Micro GC	Supported
G3581A	490 Micro GC	Supported TCP/IP. For legacy mainboard: minimum required FW is 1.40
G3582A	490 Micro GC Analyzer	Supported
M200 M400	3000 micro GC	Not Supported

### Other, unsupported instruments

Product Number	Module Name	Comment
G1176A G1180A	6820 GC	Not Supported
5890A 5890 Series II 4890D	GC systems	Not Supported

## GC Sampler, Autosampler and Tray Compatibility

**NOTE**

Communication for GC samplers and tray is handled through the GC.

### 7650 ALS hardware:

Product Number	Model Name	Comment
G4567A	Injector	Supported

### 7693 ALS hardware

Product Number	Model Name	Comment
G4513A	7693 Injector	Supported
G4514A	7693 Tray	Supported
G4515A	BCR/Mixer/Heater	Supported
G4516A	ALS Controller Board for 6890 Plus GC	Supported
G4517A	ALS Controller for 6890A GC upgrade	Supported

### 7683B ALS hardware

Product Number	Model Name	Comment
G2912A	7683B ALS Controller for 6890	Supported
G2913A	7683B Injector	Supported
G2614A	Tray	Supported
G2615A	BCR/Mixer	Supported

**7683A ALS hardware:**

Product Number	Model Name	Comment
G2612A	7683B ALS Controller for 6890	Supported
G2613A	7683A Injector	Supported
G2614A	Tray	Supported
G2615A	BCR/Mixer	Supported
G2880A	Injector	Supported

**Other, unsupported instruments**

Product Number	Model Name	Comment
19405A	Sampler Event Control Module (SECM)	Not supported
G1926A	Barcode Reader/Mixer	Not supported
G8130A	WorkBench Tray	Not supported
G7360A	Archon Purge and Trap Auto sampler for Water only	Not supported
G7361A	Archon Purge and Trap Auto sampler for Needle Sparge	Not supported

## Headspace support with GC

**NOTE**

License number no longer required for installation. The software registration packet will still include a license number.

**NOTE**

Communication is handled via LAN

Product Number	Model Name	Comment
G4511A	8697 Headspace	Communication is handled through the GC
G4556A, G4557A	7697A Headspace	Supported
G1888A	G1888 Headspace	Supported
G1289B, G1290B, G1883A	7694A 7694B  7694E Headspace	Not supported 7694B headspace was last supported with version C.01.05 with Headspace software rev B.01.04.

## Control of PAL sampler with GC

Product Number	Model Name	Comment
G6501B, G6502B, G6509B	Agilent PAL-xt CTC Sampler	Supported PAL-xt method parameters are available to Intelligent Reporting beginning in PAL Software Rev. B.01.07
G7366A, G7367A, G7368A, G7370A	Agilent PAL-3 CTC Autosampler	Supported

# 7

## Compatibility with Waters instruments

Supported driver for Waters instruments with OpenLab ChemStation:

### M8505AA Waters LC Control for OpenLab EZChrom and ChemStation

It includes the following drivers:

- Waters ACQUITY Driver rev. A.02.03
- Waters Alliance /eAlliance LC Driver rev. 6.3.1

#### NOTE

English language only

For further information please consult the respective driver user- and service documentation, your local Agilent sales representative or check into SubscribeNet.

### Compatibility with Waters ACQUITY and ACQUITY H-Class instruments

Instruments Element	Sub-element	Revision	Support Statement
UPLC	Sample Manager (SM)	1.65	Please check for driver release service note
	Sample Organizer*	1.60	
	Binary Solvent Manager (BSM)	1.72	
	Column Manager (CM)	1.40.74	*To be configured via the Sample Manager
	TUV Detector	1.70	
	PDA Detector	1.65	
	PDA eLambda Detector	1.65	
	FLR Detector	1.69	
ELS Detector	1.40		
I-Class	Sample Manager (SM-FL)	1.70	
	Sample Organizer*	1.60	
	Binary Solvent Manager (BSM)	1.65	
	Column Manager Active (CM-A)	1.69	
	TUV Detector	1.70	

## Compatibility with Waters instruments

	PDA Detector	1.65	
	PDA eLambda Detector	1.65	
	FLR Detector	1.69	
	ELS Detector	1.40	
<b>H-Class</b>	Sample Manager FTN	1.71	*To be configured via the Sample Manager
	Sample Organizer*	1.60	
	Quaternary Solvent Manager (QSM)	1.65	
	Column Manager Active (CM-A)	1.69	
	Column Manager Auxiliary (CM-AUX)	1.69	
	TUV Detector	1.70	
	PDA Detector	1.65	
	PDA eLambda Detector	1.65	
	FLR Detector	1.69	
	ELS Detector	1.40	
<b>H-Class Bio</b>	Sample Manager FTN Bio	1.71	
	Sample Organizer*	1.60	
	Quaternary Solvent Manager Bio	1.65	
	Column Manager Active (CM-A)	1.69	
	Column Manager Auxiliary (CM-AUX)	1.69	
	TUV Detector	1.70	
	PDA Detector	1.65	
	PDA eLambda Detector	1.65	
	FLR Detector	1.69	
	ELS Detector	1.40	

## Compatibility with Waters Alliance instruments:

Instruments Element	Sub-element	Revision	Support Statement
<b>Integrated LC (Autosampler + Pump)</b>	Waters Alliance 2690	1.20	
	Waters Alliance 2695	2.04	
	Waters Alliance e2695	3.00	
<b>UV Detector</b>	Waters 2487 UV	1.01 or 1.02	Please check for driver release service note
	Waters 2489 UV	1.65.2517	
<b>PDA Detector</b>	Waters 996 PDA	1.8	
	Waters 2996 PDA	R3.5	
	Waters 2998 PDA	1.65.6205	
<b>RI Detector</b>	Waters 2414 RI	2.10	
<b>FLR Detector</b>	Waters 2475 FLR	2.20	

## 8 OpenLab ChemStation Workstation VL

### Overview

OpenLab ChemStation offers a VL version for Workstations. It provides the full functionality of the Workstation configuration while restricting instrument control to one LC or GC instrument only.

In addition, it restricts the use of the instruments to standard modules. LC/MS configurations are not supported. The supported modules for VL Systems are listed in more detail in tables 4 to 7 below.

### **OpenLab ChemStation Workstation VL allows control of *one* of the below instruments:**

- *One* 1260 Isocratic Infinity (II) LC system, or one 1260 Quaternary Infinity (II) LC system with selected 1260 Infinity (II) LC modules (see table 5).

#### NOTE

1290 series LC modules are not compatible with VL.

- *One* Agilent 1220 Infinity LC System or one 1120 Compact LC. Control of supplemental Infinity LC modules is supported except for modular pumps.
- *One* 8860 or 7820 GC System, or *one* 490 or 990 Micro GC instrument.

## Supported Instruments with OpenLab ChemStation Workstation VL

### GC instrument

Use instrument type **Agilent GC Core System (VL)** to configure your GC instrument.

**Table 4** GC instruments supported with OpenLab ChemStation Workstation VL

Product Number	Model Name	Instrument Driver
G2790A	8860 GC System	G2790A GC VL driver
G4350B	7820A GC System	Instrument driver is bundled with the instrument hardware
G3581A	490 Micro GC System	M8238AA Agilent Micro GC Instrument Driver
G3588A	990 Micro GC System	

### Agilent 1260 Infinity II LC instrument

Use instrument type **Agilent LC Core System** to configure your modular 1260 LC instrument

The following table summarizes all 1260 Infinity LC modules that can be added to a 1260 Infinity LC system with a OpenLab ChemStation Workstation VL license.

**Table 5** Agilent LC modules that may be part of the 1260 LC instrument

Product Number	Model Name	
G1310B	1260 Isocratic Pump	
G1311B,	1260 Quaternary Pump	
G1311C	1260 Quaternary Pump VL	
G7110B G7111A G7111B	1260 Infinity II Pumps:	Isocratic Quaternary Quaternary
G1316A	1260 Column Compartment	
G1328C	1260 Manual Injector	

## OpenLab ChemStation VL – Support Information

Product Number	Model Name	
G7116A	1260 Infinity II Multicolumn Thermostat	
G1328C	1260 Manual Injector	
G1329B,	1260 Autosampler	may add G1330B Thermostat
G7129A G7129C	1260 Infinity II Vialsamplers	Includes ICC: G7130A
G1315C, G1315D, G4212B	1260 Diode Array Detectors	DAD
G1365C, G1365D	1260 Multiple Wavelength Detectors.	MWD
G1314B, G1314C	1260 Variable Wavelength Detectors.	VWD
G1362A, G7162A	1260 Refractive Index Detector	RID
G1321B, G1321C	1260 Fluorescence Detector	FLD
G7114A G7115A, G7121A, G7162A G7165A	1260 Infinity II detectors <sup>33</sup>	VWD, DAD, FLD, RID, MWD
G1390B	1200 UIB II	

## Agilent 1220 Infinity II LC system

To configure your 1220 LC System, use instrument type **Agilent 1220 LC System**.

The appropriate instrument drivers are bundled with the instrument hardware.

**Table 6** Agilent 1220 (II) Infinity LC Systems configurable with OpenLab ChemStation Workstation VL license

Product Number	Model Name
G4286B	1220 Infinity LC - Isocratic, Man. Inj., VWD, 600 bar
G4287B	1220 Infinity LC - Isocratic, ALS, TCC, VWD, 600 bar
G4288B	1220 Infinity LC - Gradient, Man. Inj., VWD, 600 bar
G4289B	1220 Infinity LC - Gradient, ALS, TCC, VWD, 600 bar
G4290B	1220 Infinity LC - Gradient, ALS, Man. Inj., TCC, VWD, 600 bar
G4291B	1220 Infinity LC - Isocratic, Man. Inj., TCC, VWD, 600 bar
G4292B	1220 Infinity LC - Isocratic, ALS, VWD, 600 bar
G4293B	1220 Infinity LC - Gradient, ALS, VWD, 600 bar
G4294B	1220 Infinity LC - Gradient, ALS, TCC, DAD, 600 bar
G4288C	1220 Infinity LC - System VL, Gradient, Man. Inj. VWD, 400 bar
G4289C	1220 Infinity LC - System VL, Gradient, Man. Inj. VWD, 400 bar
G4290C	1220 Infinity LC - System VL, Gradient, ALS, TCC, VWD, 400 bar
G4293C	1220 Infinity LC - System VL, Gradient, ALS, VWD, 400 bar

With exception of modular pumps all Agilent LC modules may be connected as supplemental module to your 1220 LC system.

## Agilent 1120 Compact LC

Use instrument type **Agilent 1220 LC System** to configure your 1120 Compact LC.

The appropriate instrument drivers are bundled with the instrument hardware.

**Table 7** Agilent 1120 Compact LCs configurable with ChemStation Workstation VL

Product Number	Model Name
G4286A	1120 Compact LC, Isocratic
G4287A	1120 Compact LC Isocratic with oven and ALS
G4288A	1120 Compact LC, gradient
G4289A	1120 Compact LC, gradient with oven
G4290A	1120 Compact LC gradient with oven and ALS
G1321A	1100/1200 FLD
G1362A	1100/1200 RID
G4260A	ELSD 380 Detector
G4261A	ELSD 385 Detector
G4260B	ELSD 1260 Infinity Detector
G4261B	ELSD 1290 Detector

## 9 Classic Driver Support Information

With exception of the MSD drivers, classic drivers are *not supported* with OpenLab ChemStation.

For information how to reconfigure instruments with classic drivers to use RC.net drivers please see the [OpenLab ChemStation Configuration Guide \(CDS\\_CS\\_Configure.pdf\)](#), Chapter 5, section "Reconfigure Instruments Using Classic Drivers After Upgrade to ChemStation C.01.08 or Higher".

## In This Book

The manual describes the compatibility of OpenLab ChemStation software with the following instruments and drivers:

- Agilent HPLC and LC instruments
- Agilent LC/MS Modules
- Agilent CE Modules
- Agilent Analog/Digital converters
- Agilent GC instruments
- Waters LC ACQUITY/ Alliance

It includes information on

- OpenLab ChemStation Workstation VL
- Classic Driver Support

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

© Agilent Technologies, Inc. 2010- 2020, 2022

Edition 09/2022

D0013761 Rev. B

