

OpenLab EZChrom A.04.10

Supported Instruments and Firmware Guide

Notices

Manual Part Number

M8201-90038 Rev. C
September 2020

Copyright

© Agilent Technologies, Inc. 2020
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This guide is valid for the A.04.10 revision or higher of Agilent (Product Name), until superseded.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.

Contents

OpenLab CDS Driver Information	5
Minimum required Firmware	5
Operating Systems	5
Driver types	5
Instrument Communication Information	6
Instrument LAN Interface Card	6
Agilent HPLC/LC Instrument and Firmware Compatibility	7
General Firmware Support Statement	7
Support statement for HPLC/LC Modules	8
Agilent LC Sampling Systems	8
Agilent LC Pumps	9
Agilent LC Column Compartments	10
Agilent LC Detectors	11
Agilent LC Fraction Collector (Some features may not be available in OpenLab EZChrom)	12
Agilent LC – Valves, Valve Drives and Clusters	13
Agilent LC Valve Thermostat Cluster (VTC)	13
Agilent LC – Other Module types	14
Agilent 218/SD1 Prep LC A.01.02	14
Agilent 1120 Compact LC and 1220 Infinity LC Systems	15
Lab Advisor Software Support for LC	15
Agilent Analog-Digital Converter Information	16
Agilent A/D Converter	16
Agilent GC Instrument Compatibility	17
General Support Statement	17
Supported Agilent GC hardware	17
GC Sampler, Autosampler and Tray Compatibility	18
7650 ALS hardware	18
7683B ALS hardware	18
7683A GC ALS hardware	18

7693 ALS hardware	19
Headspace Support with GC	19
Control of PAL Autosampler with GC	19
Agilent GC software support	20
OpenLab CDS MatchCompare	20
Hardware not supported with the current OpenLab EZChrom edition	21
Supported Non-Agilent GC Devices	22
Supported Non-Agilent LC Devices	23
OpenLab CDS Compact and VL Licenses	24
OpenLab CDS VL	24
OpenLab EZChrom VL rev. A.04.10	24
OpenLab EZChrom Compact	24
Supported Configurations for OpenLab EZChrom VL	25
A. Control of one Agilent GC, using the respective GC instrument type	25
B. Control of one Agilent 1260 Infinity LC instrument, configure modules using instrument type "Agilent LC Core System".	25
C. Control of one 1120 Compact LC, or one 1220 Infinity LC system using Instrument type "Agilent Compact LC".	26
Agilent OpenLab EZChrom Compact Support Information	28
OpenLab EZChrom Compact Compatibility with GC.	28
OpenLab EZChrom Compact Compatibility using instrument type "Agilent Compact LC"	28
In This Book	30

OpenLab CDS Driver Information

Minimum required Firmware

Agilent recommends always using the most recent firmware revisions. This will provide the highest level of system capability and ensure latest firmware features and improvements are present.

Drivers are forward compatible with respect to firmware, i.e. the firmware can be updated without the need of updating the driver or CDS.

In the following sections this guide lists the instruments and modules compatible with the current version of OpenLab EZChrom. Whenever you use new hardware you have to update your firmware to the current revision.

Even though Agilent recommends using the most recent firmware revisions, older instruments can be operated with lower firmware revisions. Note, that in this case you need to ensure compatibility to the current CDS and between the firmware versions of individual components of your system. Please see the following sections in this guide for more detail.

The minimum instrument driver and firmware versions listed have been verified to work with OpenLab EZChrom version A.04.10. Please note, that minimum required firmware lower than the most recent firmware might not be apt to support the latest features.

Whenever you use new hardware you have to update your firmware to the current revision.

Operating Systems

Unless otherwise noted all instrument drivers are supported on

- Windows 10 (64-bit)
- Windows Server 2016
- Windows Server 2019

For more information on support of server operating systems please see the *OpenLab EZChrom Requirements* guide.

Driver types

Current Agilent OpenLab CDS drivers are based on the RC.NET architecture. Information in this guide refers to this driver type, unless otherwise noted.

Instrument Communication Information

Instrument LAN Interface Card

Product	Hardware	Minimum required Firmware Revision	Notes
G1369C	LAN Interface card	B.06.40	CAN support for hosted modules/CAN slave (e.g. G1170A Universal Valve Drive) SW 7 and SW 8 must be in the OFF position 10/100BaseT
G1369B	LAN Interface card	A.01.01	SW 7 and SW 8 must be in the OFF position 10/100BaseT
G1369A	LAN Interface card	A.01.10	2 nd version – Board revision C.03.00 – firmware A.01.10 – Supported with 35900E and GC 1 st version – firmware A.01.05 – Not Supported with GC and 35900E

J4100A Jet Direct Card and 82357B USB /GBIB interface are no longer compatible with the current version of OpenLab EZChrom.

Agilent HPLC/LC Instrument and Firmware Compatibility

General Firmware Support Statement

Agilent recommends using current LC firmware sets with OpenLab EZChrom A.04.10. The latest LC firmware sets are 7.01, 6.50, and 6.30.

Firmware Set 7.01 can be downloaded from <http://www.agilent.com/en-us/firmwareDownload?whid=99818>.

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.

Modules combined into one LC instrument always need to have firmware from the same set. Firmware of one set is fully compatible with your CDS that supports this firmware set. However, if one module uses a firmware revision lower than the specified Minimum Firmware Revision some functionality might not be supported.

Do not mix firmware revisions from one set with older or newer sets. Firmware is not tested across set borders!

You need to upgrade ALL existing modules to the latest version only if

- you add a new LC module to the existing system (recommended).
- one of the existing modules requires an upgrade due to a bug fix solved with the latest release.

For more information on downloading the current LC firmware, please refer to the Agilent website [LC Firmware News and Downloads](#) and [Firmware Update Tools & Procedures](#).

EZChrom A.04.10 will support LC driver release A.02.19. If you plan to use a more recent driver revision, please refer to the respective *Release Note for Agilent LC and CE drivers* delivered with the drivers for more information.

Support statement for HPLC/LC Modules

Some of the modules are also supported with Firmware revisions lower than the most recent one. The firmware revision listed below or higher can be used. For more detail on firmware revisions required, please refer to "Release Note for Agilent LC and CE Drivers Revision A.02.19".

Agilent LC Sampling Systems

Product Number	Module Name	Minimum Firmware Rev. (RC.NET Driver A.02.19)	Comments
G1313A	1100 Standard Autosampler	A.06.10	<i>Exception:</i> no support for G1313A#21 multi-draw kit
G1329A	1200 Series Standard Autosampler	A.06.10	
G1329B	1260 Infinity Standard Autosampler	A.06.10	
G1367A	1100 Series Well-plate Autosampler	A.06.31	
G1367B	1200 Series High Performance Autosampler	A.06.31	
G1367C	1200 Series High Performance Autosampler SL	A.06.31	
G1367D	1200 Series High Performance Autosampler SL+	A.06.31	
G1367E	1260 Infinity High Performance Autosampler	A.06.32	
G1377A	1260 Infinity High Performance Micro Autosampler	A.06.12	
G1389A	1100 Series Micro Thermostatted Autosampler	A.06.10	
G2258A	1260 Infinity Dual-Loop Autosampler	A.06.50	
G2260A	1260 Infinity Preparative Autosampler (High flow)	A.06.50	
G4226A	1290 Infinity Autosampler	A.06.31	Based on G1367D
G4303A	1260 Infinity SFC standard autosampler	A.06.54	
G5667A	1260 Infinity Bio-inert High Performance Autosampler	A.06.32	
G5668A	1260 Infinity II Bio-inert Multisampler	D.07.13	
G7167A	1260 Infinity II Multisampler	D.07.13	
G7167B	1290 Infinity II Multisampler	D.07.13	
G7129A	1260 Infinity II Vialsampler	D.06.76	
G7129B	1290 Infinity II Vialsampler	D.06.76	
G7129C	1260 Infinity II Vialsampler	D.07.20	

Product Number	Module Name	Minimum Firmware Rev. (RC.NET Driver A.02.19)	Comments
G7157A	1260 Infinity II Preparative Autosampler	D.07.01	
G4767A	1260 Infinity II SFC Multisampler	D.07.13	

Agilent LC Pumps

Product Number	Module Name	Minimum Firmware Rev. (RC.NET Driver A.02.19)	Comments
G1310A	1200 Series Isocratic Pump	A.06.10	
G1310B	1260 Infinity Isocratic Pump	A.06.32	
G1311A	1200 Series Quaternary Pump*	A.06.10	G1354A (bundle)
G1311B	1260 Infinity Quaternary Pump*	A.06.32	
G1311C	1260 Infinity Quaternary Pump VL*	A.06.32	
G1312A	1200 Series Binary Pump*	A.06.10	
G1312B	1260 Infinity Binary Pump*	A.06.10	
G1312C	1260 Infinity Binary Pump VL*	A.06.32	
G1361A	1260 Infinity Preparative Pump	A.06.50	G1391A (gradient extension)
G1361A cluster	1260 Infinity Preparative Pump Cluster	A.06.50	Cluster with up to 4 pumps
G1376A	1260 Infinity Capillary Pump	A.06.10	G1382A (includes degasser)
G2226A	1260 Infinity Nanoflow Pump	A.06.10	
G4204A	1290 Infinity Quaternary Pump*	B.06.50	
G4220A	1290 Infinity Binary Pump*	B.06.23	
G4220A cluster	1290 Infinity Binary Pump* Cluster	B.06.23	Pump-Valve-Cluster: up to two G1160A valves can be clustered together with a pump
G4220B	1290 Infinity Binary Pump VL*	B.06.43	
G4302A	1260 Infinity SFC Binary Pump*	A.06.32	
G4782A	1260 Infinity II SFC Binary Pump*	D.07.13	
G5611A	1260 Infinity Bio-inert Quaternary Pump*	A.06.32	
G5654A	1260 Infinity II Bio-inert Quaternary Pump*	D.07.01	
G7104A	1290 Infinity II Flexible Pump*	B06.71	

Product Number	Module Name	Minimum Firmware Rev. (RC.NET Driver A.02.19)	Comments
G7104C	1290 Infinity II Flexible Pump*	D.07.20	
G7110B	1260 Infinity II Isocratic Pump	D.07.01	
G7111A	1260 Infinity II Quaternary Pump VL*	D.07.01	
G7111B	1260 Infinity II Quaternary Pump*	D.07.01	
G7112B	1260 Infinity II Binary Pump*	D.07.01	
G7120A	1290 Infinity II High Speed Pump*	B.06.71	
G7161A	1260 Infinity II Preparative Binary Pump	D.07.20	
G7161B	1290 Infinity II Preparative Binary Pump	D.07.20	

*Pump valve clusters are possible for marked pumps with up to 2 valves of type G1160A and/or G1170A with 5067-4159 or 5067-4147.

Agilent LC Column Compartments

Product Number	Module Name	Minimum Firmware Rev. (RC.NET Driver A.02.19)	Comments
G1316A	1260 Infinity Thermostatted Column Compartment	A.06.10	no support for G1316A#56 microvalve control.
G1316B	1200 Series Thermostatted Column Compartment SL	A.06.10	no support for G1316B#56 microvalve control.
G1316C	1200 Series Thermostatted Column Compartment SL*	A.06.14	no support for new 2/10 and 2/6 valves with Classic.
G1316C cluster	1200 Series Thermostatted Column Compartment SL* cluster	D.06.14	no support for new 2pos/6port or 2pos/10port valves
G7116A	1260 Infinity II Multicolumn Thermostat (firmware for host module in brackets)	C.07.01 (B.07.01/D.07.01)	
G7116B	1290 Infinity II Multicolumn Thermostat (firmware for host module in brackets)	C.06.75 (B.06.75/D.06.75)	
G7130A	Integrated Column Compartment ICC	D.06.76	

*Cluster with up to three G1316C with integrated 8pos/9port valves (products G4230A/B). Minimum two G1316C TCCs, the third TCC can be a G1316A, B or C.

Agilent LC Detectors

Product Number	Module Name	Minimum Firmware Rev. (RC.NET Driver A.02.19)
G1314A	1100 Series Variable Wavelength Detector	A.06.10
G1314B	1260 infinity Variable Wavelength Detector VL	A.06.10
G1314C	1260 Infinity Variable Wavelength Detector VL+	A.06.10
G1314D	1200 Series Variable Wavelength Detector	B.06.32
G1314E	1290 Infinity Variable Wavelength Detector	B.06.32
G1314F	1260 Infinity Variable Wavelength Detector	B.06.32
G1315A	1100 Series Diode Array Detector	A.06.10
G1315B	1200 Series Diode Array Detector	A.06.10
G1315C	1260 Infinity Diode Array Detector VL+	B.06.30
G1315D	1260 Infinity Diode Array Detector VL	B.06.30
G1365A	1100 Series Multiple Wavelength Detector	A.06.10
G1365B	1200 Series Multiple Wavelength Detector	A.06.10
G1365C	1260 Infinity Multiple Wavelength Detector	B.06.30
G1365D	1260 Infinity Multiple Wavelength Detector VL	B.06.30
G1321A	1200 Series Fluorescence Detector (FLD)	A.06.10
G1321B	1260 Infinity Fluorescence Detector Spectra	A.06.32
G1321C	1260 Infinity Fluorescence Detector	A.06.54
G1362A	1260 Infinity Refractive Index Detector	A.06.10
G4212A	1290 Infinity Diode Array Detector	B.06.30
G4212B	1260 Infinity Diode Array Detector	B.06.30
G4212A/B HDR-DAD Cluster	2x G4212A or 2x G4212B or a combination of 1x G4212A and 1x G4212B	B.06.57
G7121A	1260 Infinity II Fluorescence Detector	D.07.01
G7121B	1260 Infinity II Fluorescence Detector Spectra	D.07.01
G7165A	1260 Infinity II Multiple Wavelength Detector	D.07.01
G4218A	1260 Infinity Evaporative Light Scattering Detector	1.3
G4260A	Agilent 380-ELSD	25.00
G4260B	1260 Infinity II Evaporative Light Scattering Detector	32.06

Product Number	Module Name	Minimum Firmware Rev. (RC.NET Driver A.02.19)
G4261A	Agilent 385-ELSD	25.00
G4261B	Agilent 1290 Infinity ELSD	32.06
G7102A	1290 Infinity II Evaporative Light Scattering Detector	32.06
G7162A	1260 Infinity II Refractive Index Detector	D.06.76
G7162B	1290 Infinity II Refractive Index Detector	D.06.76

Agilent LC Fraction Collector (Some features may not be available in OpenLab EZChrom)

Product Number	Module Name	Minimum Firmware Rev. (RC.NET Driver A.02.19)
G1364A	1100 Series Automatic Fraction Collector	A.06.53
G1364A cluster*	1100 Series Automatic Fraction Collector Cluster of up to 3	A.06.53
G1364B	1260 Infinity Fraction Collector (preparative-scale)	A.06.53
G1364B cluster*	1260 Infinity Fraction Collector (preparative-scale) Cluster of up to 3	A.06.53
G1364C	1260 Infinity Fraction Collector (analytical-scale)	A.06.53
G1364C cluster*	1260 Infinity Fraction Collector (analytical-scale) Cluster of up to 3	A.06.53
G1364D	1100 Series Micro Fraction Collector	A.06.53
G1364E	1260 Infinity II Preparative Fraction Collector**	D.07.23
G1364F	1260 Infinity II Analytical Fraction Collector**	D.07.23
G5664A	1260 Infinity Bio-inert fraction collector AS	A.06.53
G7159B	1290 Infinity II Preparative Open-Bed Fraction Collector***	D.07.23
G7166A	1260 Infinity II Preparative Valve-Based Collector (firmware for host module in brackets)	C.07.10 (B.07.23/D.07.23)

Note: *Any combination of G1364A/B/C or G5664A plus a fourth G1364A/B/C or G5664A for recovery can be clustered. Multiple individual Fraction Collectors are not supported.

**Can be clustered with a G7166A or the same module type for recovery collection.

***Can be clustered with a G7166A for Recovery.

Agilent LC – Valves, Valve Drives and Clusters

Product Number	Module Name	Minimum Firmware Rev. (RC.NET Driver A.02.19)
G1156A	1200 Series 6 Position / 7 Port Valve (400 bar)	A.06.02
G1157A	1200 Series 2 Position / 10 Port Valve	A.06.02
G1158A	1200 Series 2 Position / 6 Port Valve	A.06.02
G1158B	1200 Series 2 Position / 6 Port Valve (600bar)	A.06.02
G1159A	1200 Series 6 Position Selection Valve	A.06.02
G1160A	1100 Series Multiple Purpose Switching Valve (12 Position / 13 Port)	A.06.02
G1162A	1200 Series 2 Position/ 6 Port Micro Valve	A.06.02
G1163A	G1163A 2Pos/10Port Valve (micro)	A.06.02
G1170A	1290 Infinity Valve Drive (firmware for host module in brackets)	C.06.40 (B.06.40/D.06.60)
G9322A	1260 Infinity II Clustering Valve	D.07.23
VTC	Combinations of G7116B, G1170A and G1316C (valve or column hosts) and G1361A/B and G7130A (column hosts)	see table below
FCC	Combinations of preparative Fraction Collectors (G7159B, G1364E, G7116A) and a Clustering Valve G9322A or analytical Fraction Collectors (G1364F, G5664B) and a Quick Change Valve (G1170A with 5067-4159 or 5067-4194)	

Agilent LC Valve Thermostat Cluster (VTC)

Product Number	Minimum module FW	Minimum host module FW
G7116B	C.06.75	B.06.75/D.06.75
G1170A	C.06.75	B.06.75/D.06.75
G7130A (within G7129A/B)	D.06.76	n/a
G1316C	A.06.55	n/a
G1316A/B	A.06.10	n/a

Agilent LC – Other Module types

Product Number	Module Name	Minimum Firmware Rev. (RC.NET Driver A.02.16)	Comments
G1390A*	1100 Series Universal Interface Box (UIB)	A.06.02	Not supported on Chinese EZChrom
G1390B*	1200 Infinity Series Universal Interface Box II (firmware for host module in brackets)	C.06.53 (B.06.53/D.06.60)	
G4227A	1290 Infinity Flexible Cube (firmware for host module in brackets)	C.06.52 (B.06.52/D.06.60)	
G4240A	Chip Cube	A.06.36	
G4301A	1260 Infinity Analytical SFC System	A.03.09	
G7170B	1290 Infinity II MS Flow Modulator (firmware for host module in brackets)	C.07.20 (B.07.20/D.07.20)	

Note: *The UIB is not compatible with CE modules.

Agilent 218/SD1 Prep LC A.01.02

Product Number	Module	Minimum Firmware Rev. (RC.NET Driver A.02.16)	Communications
Agilent 210	Pump	1.73 rev C	RS232
Agilent 218	Pump	1.1 rev B	RS232
Agilent SD1	Pump	2.2	RS232
Agilent 325	UV/VIS Detector	2.07	LAN
Agilent 410	Autosampler	2.07	RS232
Agilent 440	Fraction Collector	1.31	RS232

Agilent 1120 Compact LC and 1220 Infinity LC Systems

NOTE

Agilent 1120 and 1220 instruments are configurable using the instrument type "Agilent Compact LC" only.

Product Number	1120 System	Minimum Firmware Rev. (RC.NET Driver A.02.19)
G4286A	1120 Compact LC, Isocratic pump	B.06.50
G4287A	1120 compact LC. Isocratic with Oven and ALS	B.06.50
G4288A	1120 Compact LC, Gradient	B.06.50
G4289A	1120 Compact LC, Gradient with oven	B.06.50
G4290A	1120 Compact LC, Gradient with oven and ALS	B.06.50

Figure 1.

Product Number	1220 System	Minimum Firmware Rev. (RC.NET Driver A.02.19)
G4286B	1220 Infinity LC System Isocratic, Man. Inj., VWD, 600 bar	B.06.50
G4287B	1220 Infinity LC Isocratic, ALS, TCC, VWD, 600 bar	B.06.50
G4288B	1220 Infinity LC Gradient, Man. Inj., VWD, 600 bar	B.06.50
G4289B	1220 Infinity LC Gradient, ALS, TCC, VWD, 600 bar	B.06.50
G4290B	1220 Infinity LC Gradient, ALS, Man. Inj., TCC, VWD, 600 bar	B.06.50
G4291B	1220 Infinity LC Isocratic, Man. Inj., TCC, VWD, 600 bar	B.06.50
G4292B	1220 Infinity LC Isocratic, ALS, VWD, 600 bar	B.06.50
G4293B	1220 Infinity LC Gradient, ALS, VWD, 600 bar	B.06.50
G4294B	1220 Infinity LC Gradient, ALS, TCC, DAD, 600 bar	B.06.50
G4288C	1220 Infinity LC System VL, Gradient, Man. Inj. VWD, 400 bar	B.06.50
G4289C	1220 Infinity LC System VL, Gradient, Man. Inj. VWD, 400 bar	B.06.50
G4290C	1220 Infinity LC System VL, Gradient, ALS, TCC, VWD, 400 bar	B.06.50
G4293C	1220 Infinity LC System VL, Gradient, ALS, VWD, 400 bar	B.06.50

Lab Advisor Software Support for LC

Model	Description	Lab Advisor
M8550A or M8555A	Co-existence of Lab Advisor and EZChrom on the same PC is supported.	B.02.01 and later

Agilent Analog-Digital Converter Information

Agilent A/D Converter

Instrument	Supported Firmware Version (minimum)	Comment
Agilent 35900E Dual Channel Interface	E.01.02	RC.NET TCP/IP, LAN only
35900E Series II	E.02.04.40	RC.net driver
SS420X	A.01.01	

3rd and 4th Signal, supported configurations:

Instrument	35900E Classic Driver	35900E RC.NET Driver
7890A 7820A	Not supported	yes
68xx RC.net	Not supported	Yes
68xx classic	Not supported	Not supported
LC classic	Not supported	Not supported
LC RC.net	Yes (LAN only)	Yes

Agilent 35900E RC.Net driver supports separating the channels between 2 instruments on an AIC distributed system

Agilent GC Instrument Compatibility

General Support Statement

Agilent releases GC firmware updates independently of software releases. All Agilent GC instrument driver revisions have been designed to be backward compatible to the installed instrument base.

Agilent recommends always using the latest module firmware revision to provide the highest level of system capability. A firmware upgrade should be done only if you face problems or want to add system capability to your GC.

Only RC.Net drivers version B.01.04 are supported for Agilent GC instruments in A.04.10. Classic Agilent GC drivers are no longer supported and cannot be used to acquire data. The Classic Agilent 6890 GC driver is available to reprocess data previously created with a 6890 GC only.

NOTE

Communication is handled via LAN, unless stated otherwise

Supported Agilent GC hardware¹

Product Number	Module Name	Comments
G3950A, G3952A, G3953A	Agilent Intuvo 9000 GC	7683 ALS family not supported
G3450A, G3542A, G3543A, G3545A	8890A	
G2790A	8860A	
G3440A G3445A	7890A	RTL, Backflush wizard, method transfer, and Easy SamplePrep are not supported
G3440B G3445B	7890B	RTL, Backflush wizard, method transfer, and Easy SamplePrep are not supported
G4350A	7820A	RTL, Backflush wizard, method transfer, and Easy SamplePrep are not supported
G1530N G1540N	6890N	RC.NET Driver support only
G1530A G1540A	6890A 6890Plus	RC.NET Driver support only
G2629A	6850 Handheld Controller	

¹ For more detail please check current Agilent GC Drivers Release Note

G2630A	6850 serial #<= US00003200 6850 serial # >= US10243001	RC.NET Driver support only
G3581A	490 Micro GC	TCP-IP
Product Number	Module Name	Comments
5890	5890 Series II	
35900E	35900E	Thoroughly tested
35900E Series II	35900E	Thoroughly tested

GC Sampler, Autosampler and Tray Compatibility

7650 ALS hardware

Model	Description	Support Statement
G4567A	Injector	Thoroughly tested

7683B ALS hardware

Model	Description	Support Statement
G2913A	Injector	Thoroughly tested
G2614A	Tray /Controller	Thoroughly tested
G2615A	BCR/Mixer	Thoroughly tested, Mixing only

7683A GC ALS hardware

Model	Description	Support Statement
G2613A	Injector	Thoroughly tested
G2614A	Tray/Controller	Thoroughly tested
G2615A	BCR/Mixer	Thoroughly tested, Mixing Only

7693 ALS hardware

Model	Description	Support Statement
G4513A	Injector	Thoroughly tested
G4514A	Tray	Thoroughly tested
G4515A	BCR/Mixer	Thoroughly tested, Heater, Mixing Only
G4516A	External Controller	Thoroughly tested
G4517A	6890Plus ALS card upgrade	Thoroughly tested
G4521A	LVI Syringe Carriage	Thoroughly tested
G4522A	Cooling Accessory	Thoroughly tested
G4520A	Tray with BCR/Mixer	Thoroughly tested, Heater, Mixing Only

Headspace Support with GC

Model	Description	Support Statement
7697A	Headspace	Thoroughly tested
G1888A	Headspace	Thoroughly tested

Control of PAL Autosampler with GC

Model	Description	FW Revision	SW Driver	Support Statement
G6501B G6502B G6509B	Agilent PAL-xt CTC Sampler	Agilent 4.3.0	G3382AA B.01.08	Thoroughly tested.
G6501B, G6502B, G6509B	Agilent GC Sampler, Injector	Agilent 4.3.0	G3382AA B.01.08	Thoroughly tested PAL-xt method parameters are available to Intelligent Reporting beginning in PAL Software rev. B.01.07. Note: License number no longer required for installation. The software registration packet will still include a license number.

Agilent GC software support

Product Number	Module Name	Revision	Comments
G3382AA	PAL Control Software	B.01.08	Thoroughly tested
G7388AA	PAL Control Software	A.01.03	Thoroughly tested

OpenLab CDS MatchCompare

Application	Revision	Comments
OpenLab CDS Match Compare (M8350AA)	A.01.03	Supported with Workstation Standalone

Hardware not supported with the current OpenLab EZChrom edition

List may not be complete!

Model	Support Statement
5890 Series II	Not supported, but in code.
19405A/B Sampler Event Control Module (SECM)	Not supported
G2403A LAN/RS232 converter Dudley box	Not supported
18596C Tray (7673C)	Not supported
G1512A Controller (7673C)	Not supported
7673A family (18593A, 18594A, 18596A)	Not supported
7673B family (18593B, 18594B, 18596B)	Not supported
7673C family (G1512A, G1513A, 18596C)	Not supported
G1926A	Not supported
5890A	Not Supported
G1176A or G1180A (6820 GC)	Not Supported
3000 micro GC, M200, M400	Not Supported
35900C/D	Not Supported
G1289A, G1290A, G1883A (7694E), 7695A Headspace	Not Supported; Run stand-alone only. Headspace A and E models are not supported
7673A family (18593A, 18594A, 18596A)	Not Supported
7673B family (18593B, 18594B, 18596B)	Not Supported
G1926A	Not Supported
G7360A Archon Purge and Trap Autosampler for Water Only	Not Supported
G7361A The Archon Purge and Trap Autosampler for Needle Sparge	Not Supported

Supported Non-Agilent GC Devices

Please refer to the A.04.10_Non_Agilent_Compatibility_Matrix.

Supported Non-Agilent LC Devices

Please refer to the A.04.10_Non_Agilent_Compatibility_Matrix.

OpenLab CDS Compact and VL Licenses

OpenLab CDS VL

OpenLab EZChrom offers a VL version. The VL versions provide the full functionality of the *Workstation* configuration including the new OpenLab Data Analysis, while restricting instrument control.

The VL license allows to control one LC or GC instrument only and restricts the use of the instruments to standard modules. The supported modules and firmware options for VL Systems are listed in more detail on the following pages of this document. LC-MS configurations are not supported.

OpenLab EZChrom VL rev. A.04.10

Allows control of one of the below:

- **One** 1260 Isocratic Infinity LC system, or one 1260 Quaternary Infinity LC system with selected 1260 Infinity LC modules (use instrument type “*Agilent LC Core System*”).
- **One** Agilent 1220 Infinity LC system or one 1120 Compact LC system (use Instrument type “*Agilent Compact LC*”). Control of supplemental Infinity LC modules is supported for all LC modules with the exception of modular pumps.
- **One** 7820A GC, 490 Micro GC, 490-GC Micro-GC, or one 4900 Micro GC.
- **One** 8860 GC or 990 Micro GC

OpenLab EZChrom Compact

OpenLab EZChrom offers a Compact version. The license is a fixed configuration with a limited feature set that includes instrument control and data analysis. The supported modules for Compact are shown in the following sections of this document.

It provides total control of a maximum of **two** instruments of :

- Agilent 1120 and 1220 LC
- 7820A GC and 8860 GC
- Agilent 490 Micro GC, 490-GC Micro-GC, CP-4900 Micro-GC, and 990 Micro GC

Supported Configurations for OpenLab EZChrom VL

A VL license allows to control one of the configurations described below: A, B or C.

A. Control of one Agilent GC, using the respective GC instrument type

The appropriate instrument drivers are bundled with the instrument hardware.

Table 2 GC instruments supported with OpenLab VL EZChrom

Model	Description	Min. FW Revision	SW Driver	Support Statement
G4350B	7820A	A.01.15.012	RC.NET	Thoroughly tested; No support with Software Add-ons
G3581A	490 Micro GC	3.32	1.9	Supported
	CP-4900 Micro GC	1.40 build 45		Supported
	490-GC Micro GC	1.40 build 45		Supported
G2790A	8860A	3.3	RC.NET	Supported
	990-GC Micro GC	1.40 build 45		Supported

B. Control of one Agilent 1260 Infinity LC instrument, configure modules using instrument type “Agilent LC Core System”.

The Agilent LC modules listed below are supported with Value Line software license and may be part of the 1260 Infinity LC instrument.

For compatible driver and support information, please refer to section 3 of this guide.

Product Number	Module Name
G1329B	Autosampler
G1310B	1260 Series Isocratic Pump
G1311B	1260 Infinity Quaternary Pump
G1311C	1260 Infinity Quaternary Pump VL
G1314B	1260 Infinity Variable Wavelength Detector VL
G1314C	1260 Infinity Variable Wavelength Detector VL+
G1314F	1260 Infinity Variable Wavelength Detector
G1315C	1260 Infinity Diode Array Detector VL+
G1315D	1260 Infinity Diode Array Detector VL
G4212B	1260 Infinity Diode Array Detector

Product Number	Module Name
G1365C	1260 Infinity Refractive Index Detector
G1365D	1260 Infinity Refractive Index Detector
G1321B	1260 Infinity Fluorescence Detector Spectra
G1321C	1260 Infinity Fluorescence Detector
G1362A	1260 Infinity Refractive Index Detector
G1390B	1200 Infinity Series Universal Interface Box II

C. Control of one 1120 Compact LC, or one 1220 Infinity LC system using Instrument type “Agilent Compact LC”.

The appropriate instrument drivers are bundled with the instrument hardware.

Systems	
G4286A	1120 Compact LC, Isocratic
G4286B	1220 Infinity LC System Isocratic, Man. Inj., VWD, 600 bar
G4287A	1120 Compact LC, Isocratic with Oven and ALS
G4287B	1220 Infinity LC Isocratic, ALS, TCC, VWD, 600 bar
G4288A	1120 Compact LC, Gradient
G4288B	1220 Infinity LC Gradient, Man. Inj., VWD, 600 bar
G4288C	1220 Infinity LC System VL, Gradient, Man. Inj. VWD, 400 bar
G4289A	1120 Compact LC, Gradient with Oven
G4289B	1220 Infinity LC Gradient, ALS, TCC, VWD, 600 bar
G4289C	1220 Infinity LC System VL, Gradient, Man. Inj. VWD, 400 bar
G4290A	1120 Compact LC, Gradient with oven and ALS
G4290B	1220 Infinity LC Gradient, ALS, Man. Inj., TCC, VWD, 600 bar
G4290C	1220 Infinity LC System VL, Gradient, ALS, TCC, VWD, 400 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

G4293C	1220 Infinity LC System VL, Gradient, ALS, VWD, 400 bar
G4294B	1220 Infinity LC Gradient, ALS, TCC, DAD, 600 bar
G1321A	FLD
G1362A	Refractive Index Detector
G4218A	ELSD
G4260A	Agilent 380 ELSD
G4261A	Agilent 385 ELSD
G4260B	Agilent 1260 Infinity ELSD
G4261B	Agilent 1290 Infinity ELSD

NOTE

Using the "Agilent Compact LC" instrument type you will be able to configure selected Infinity LC modules along with the 1220 instrument. Please note that the number of hardware configurations is limited due to functionality incongruity. Please contact support before adding an 1100/1200/1260 Infinity module to an Agilent 1220 Infinity System to ensure compatibility.

Agilent OpenLab EZChrom Compact Support Information

OpenLab EZChrom Compact Compatibility with GC.

Table 6 EZChrom Compact compatibility with GC Hardware

Model	Description	Min. FW Revision	SW Driver	Support Statement
G4350B	7820A	A.01.15.012	RC.NET	Thoroughly tested; No support with Software Add-ons
G3581A	490 Micro GC	3.32	1.9	Supported
	CP-4900 Micro GC	1.40 build 45		Supported
	490-GC Micro GC	1.40 build 45		Supported
G2790A	8860A	3.3	RC.NET	Supported
	990-GC Micro GC	1.40 build 45		Supported

OpenLab EZChrom Compact Compatibility using instrument type “Agilent Compact LC”

Table 7 EZChrom Compact compatibility with Agilent 1120 Compact LC or Agilent 1220 Infinity LC Systems

Product Number	Module 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
G4286B	1220 Infinity LC System 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

<i>Product Number</i>	<i>Module Name</i>
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
G1321A	FLD
G1362A	Refractive Index Detector
G4218A	ELSD
G4260A	Agilent 380 ELSD
G4261A	Agilent 385 ELSD
G4260B	Agilent 1260 Infinity ELSD
G4261B	Agilent 1290 Infinity ELSD

NOTE

Using the “Agilent Compact LC” instrument type you will be able to configure selected Infinity LC modules along with the 1220 instrument. Please note that the number of hardware configurations is limited due to functionality incongruity. Please contact support before adding an 1100/1200/1260 Infinity module to an Agilent 1220 Infinity System to ensure compatibility.

In This Book

This guide summarizes the instruments supported for operation with the Agilent OpenLab EZChrom A.04.10.

