

Release Notes for Agilent Drivers for Thermo Scientific Chromeleon 7 (Rev. 1.2)

This document provides information on the major features and supported modules of the Agilent Drivers for Chromeleon 7, Revision 1.2.

Contents

Introduction	2
Agilent Drivers for Chromeleon 1.2 – What’s new?	3
Agilent Drivers for Chromeleon 1.1 – Features	4
Compatibility	5
Installation	6
Support Information/User Documentation	7
Online-Help	8
User Guides	8
Obtaining Technical Support	9
Supported Agilent LC Modules and Firmware	10



Introduction

The Agilent Drivers for Thermo Scientific Chromeleon 7 is an instrument control application for the Agilent LC instrument portfolio running in the Thermo Scientific Chromeleon 7.2 SR5 or higher environment.

References to product documentation for installation and usage are provided, as well as references to documentation regarding known issues and workarounds.

Table 1 Terms and Abbreviations used in this document

Terms	Description
Agilent Drivers	Agilent Drivers for Thermo Chromeleon 7
Chromeleon 7	Thermo Scientific Chromeleon 7 Chromatography Data System (CDS)
Chromeleon	Thermo Scientific and Chromeleon are registered trademarks of Thermo Fisher Scientific.
Thermo	Thermo Fisher Scientific

For our Regulated Customers

When any change is made to Agilent software, the validation status of the software needs to be re-established by the user. Whenever software is changed, a validation analysis should be conducted not just for the validation of an individual change, but also to determine the extent and impact of that change on the entire software system.

Agilent Drivers for Chromeleon 1.2 – What's new?

- *New LC instrumentation*
 - G4782A 1260 Infinity II SFC Binary Pump
 - G4767A 1260 Infinity II SFC Multisampler
 - G7161A 1260 Infinity II Preparative Binary Pump
 - G7161B 1290 Infinity II Preparative Binary Pump
 - G7170B 1290 Infinity II MS Flow Modulator
 - G7129C 1260 Infinity II Vialsampler
 - G7104C 1260 Infinity II Flexible Pump
 - G4260B 1260 Infinity II Evaporative Light Scattering Detector with LAN connection
 - Prep 6-column selector valve head 5067-4267
 - G4232E (5067-4283) 2pos/10port Valve head 800 bar
 - G4234D (5067-4284) 6 column selector 800 bar with different port layout
 - G4237D (5067-4279) 4-column selector 800 bar
 - G4231D (5067-4282) 2pos/6port Valve Head 800 bar
 - new characterizations for ISET 4.2 (support for G7104C)
(For more details, see [“Supported Agilent LC Modules and Firmware”](#) on page 10)
- *Shutdown Method*

Enables the user to shut down the pump and detector lamp after a run.
- *Direct Actions*

Enables the user (expert permissions) to execute direct actions via the command tree or to place them on an ePanel.
- *Improvements timetable handling*

Invalid timetable script changes are notified in more detail at method check.

Agilent Drivers for Chromeleon 1.1 – Features

The Agilent Drivers for Chromeleon 7 supports the following features:

- Support for the following Agilent LC Systems and Modules:
 - 1220/1260/1290 Agilent Infinity II LC
 - 1220/1260/1290 Agilent Infinity LC
 - 1100/1200 Series and 1120
- Chromeleon specific user experience to control Agilent LC equipment, as these drivers offer a new look and feel similar to the previous look and feel of the Chromeleon Native Driver for Agilent LC instrumentation:
 - *ePanels for each module class*

In addition to the Home ePanel offering the LC Instrument Control Dashboard module ePanels are now available. There is one ePanel available for each module class, which can be customized to fit the configuration in use.
 - *Independent module access in instrument method tree*

The instrument method offers all modules as a separate unit enabling direct access and allowing method creation via the guided method wizard.
 - *Chromeleon specific command handling*
 - The Command Tree offers Agilent method parameters.
 - The Chromeleon method script includes the Agilent instrument method parameters now in scripted format.
 - Timetable parameters listed in method script’s “Run stage” in order of their execution.
 - Enabling full feature set for custom variables handling in method script.
 - *Elimination of duplicate entries*
 - One entry for overall run time/stop time.
 - Injection volume defined by sequence only (manual overwrite with script possible).
 - *Alignment of status information*
 - Agilent LC Status Dashboard runtime / Chromeleon runtime information
 - Agilent LC Status Dashboard colored status / Chromeleon colored status information
 - *Audit trail captures changes falling within and outside a run on module level*
 - *Data Audit Trail for Instrument Method*

Using the enabled versioning and data audit trails is now possible to perform a method comparison of different versions via the scripted method and no longer in an ICF specific view.
 - *Enabling development of shutdown method to switch off lamps or pumps at end of the sequence*

- *Sequence handling*
 - Acts as master for injection volume (manual overwrite via method script possible).
 - Improved overlapped injection
- *Performance improvements*

The performance of the graphical user interface components, e.g. close/open method windows has been improved.
- *Partially available Method Migration*
 - User guidance migrating previously created methods using Instrument Control Framework (ICF) is given, the resulting method continues to work with the scripted method (minor manual interaction still required).
 - No complete method migration possible due to the revised command naming structure. However, all incompatible commands are marked red and can easily be transferred manually.

The Agilent Drivers for Chromeleon 7 does not support the following features:

- *Blend Assist*

This pump feature is offered by the LC driver, but it is not supported in this release.

Compatibility

The compatibility matrix provides information about installation and execution prerequisites with respect to hardware, firmware, and the operating system.

Supported Chromatographic Data System (CDS)

The following CDS versions support the Agilent Drivers for Chromeleon 7 (Rev. 1.2):

- Thermo Scientific Chromeleon 7.2 SR 5
- Thermo Scientific Chromeleon 7.2 SR 5 MUX
- Thermo Scientific Chromeleon 7.2.6
- Thermo Scientific Chromeleon 7.2.7
- Thermo Scientific Chromeleon 7.2.8
- Thermo Scientific Chromeleon 7.2.9

Supported Agilent Components

The following Agilent components are included in Agilent Drivers for Chromeleon 7 (Rev. 1.2):

- Agilent Instrument Control Framework A.02.05
- Agilent LC Drivers A.02.18
- Agilent ELSD Drivers 1.7
- Agilent LC Modules/Firmware (see [“Supported Agilent LC Modules and Firmware”](#) on page 10)

Supported Operating Systems

The supported operating system in use is determined by the hosting CDS.

The following operating systems have been tested by Agilent:

- Microsoft Windows 7 and 7 SP1, 32 bit and 64 bit
- Microsoft Windows 8.1, 64 bit
- Microsoft Windows 10, 64 bit
- Microsoft Windows 2012 Server R2, 64 bit

The LC Drivers have been optimized for the Windows default font size (100 %). Larger font sizes may require increasing the window size or may cause truncations.

Supported Language Settings

The Agilent Drivers for Chromeleon 7 are localized in English only.

Installation

The Agilent Drivers for Chromeleon 7 are delivered on the Chromeleon installation medium available at time of release. Installation prerequisites are outlined by Chromeleon (e.g. CPU, memory and hard drive space).

The preferred installation is the *automatic installation* using the advanced option of the Chromeleon installer. Please refer to the Chromeleon Installer documentation for installation, updates and uninstallation.

Manual installation is described in the Agilent User Guide (Users Guide Agilent Drivers for Chromeleon.pdf), available on the Chromeleon DVD in folder \Packages\Agilent Chromeleon Drivers\Documentation.

Installation Verification

Agilent offers a Software Verification Tool (SVT) to verify the correct installation of the software components. The tool is delivered along with the Agilent Drivers for Chromeleon 7.

- Using the Chromeleon installation routine, the tool is installed along with the Agilent Drivers.
- Manual installation of the Agilent Drivers for Chromeleon requires the Software Verification Tool (SVT) to be installed prior to the Agilent Drivers installation.

After the installation of the Software Verification Tool (SVT) execute the verification with the following steps:

- 1 Open **Start > Programs > Agilent Technologies** and select **Software Verification Tool**.
- 2 In the upcoming dialog box, select the required report type and the components of interest and click **Qualify**. The corresponding browser opens and shows the resulting files and passed or failed status of the installation.

NOTE

- The Chromeleon IQ does not start the Agilent SVT, the Agilent Software Verification Tool requires manual execution.
- Silent execution of the installation verification is possible, please refer to the User Guide for the Agilent Drivers for Chromeleon 7.
- The Agilent Drivers for Chromeleon 7 are not listed among the third party components in the Chromeleon IQ report.

NOTE

If the Agilent Drivers for Chromeleon are already installed on the system, the tool does not have to be upgraded or installed again.

Support Information/User Documentation

The Agilent Drivers for Chromeleon 7 include the following components:

- Agilent Driver for Thermo Chromeleon 7 (integration adapter)
- Agilent Instrument Control Framework (ICF)
- Agilent Instrument Control Framework - LC Driver

The following components documents are present in the folder structure \Packages\Agilent Chromeleon Drivers\Documentation of the Chromeleon DVD:

- *Release Notes*

The release notes document new and changed feature sets, important information on the required operating environment, supported modules and firmware, impact analysis etc.

- *Software Status Bulletin (SSB)*

The Software Status Bulletin lists known limitations and incompatibilities and information about available fixes or workarounds for this and previous versions

- The *LC Driver SSB* is updated regularly. Please visit our Website for the latest version:

http://www.chem.agilent.com/Library/Support/Patches/SSBs/LC_RC_Net.html

- The *ICF Driver SSB* is updated regularly. Please visit our Website for the latest version:

[http://www.agilent.com/cs/library/support/Patches/SSBs/Agilent_Instrument_Control_Framework_\(ICF\).html](http://www.agilent.com/cs/library/support/Patches/SSBs/Agilent_Instrument_Control_Framework_(ICF).html)

- The *Agilent Driver for Chromeleon 7 Driver SSB* is updated regularly. Please visit our Website for the latest version:

[http://www.agilent.com/en-us/support/liquid-chromatography/agilent-drivers-fo
r-chromeleon-7](http://www.agilent.com/en-us/support/liquid-chromatography/agilent-drivers-for-chromeleon-7)

- *Software Release Bulletin (SRB)*

The Software Release Bulletin is an excerpt from the SSB listing issues which have been fixed with this revision.

- *Validation Certificate/Declaration of Software Quality*

This document provides the assurance that the Agilent software product listed was developed and tested using Agilent's product development and lifecycle processes.

Online-Help

LC Drivers

Online help is available either via the help button present on the window screen or using the F1 button. F1 brings up online help even if there is no help button present. LC Driver Help explains the parameters present on the current window along with the possible parameter ranges, variables and allowed formats, which may be entered.

Agilent Drivers for Chromeleon

Online help for the Agilent Drivers for Chromeleon can only be accessed via the Help button below the LC dashboard. Help provides information on special handling required to run Agilent LC modules in a Chromeleon environment.

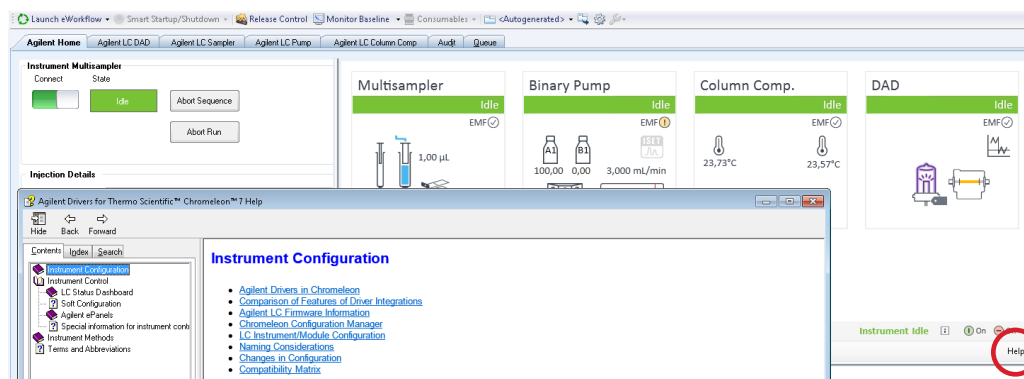


Figure 1 How to access Agilent Drivers for Chromeleon Help

User Guides

A User Guide is included in the installation for the Agilent Drivers for Chromeleon 7. Please navigate on the Chromeleon disk to the folder \Packages\Agilent Chromeleon Drivers\Documentation and open the Users Guide Agilent Drivers for Chromeleon.pdf.

The User Guide offers information on

- how to install the Agilent Drivers
- how to configure the instrument
- how to run injections
- the method handling (Agilent Method user interface/Chromeleon Script Editor)
- how to migrate methods based on native drivers and ICF based drivers
- how to troubleshoot

Obtaining Technical Support

For all technical support enquiries regarding Thermo Scientific Chromeleon 7 Chromatography Data System (CDS) software please contact your Thermo Fisher Scientific customer support organization as your first point of contact regarding any data system and instrument control enquiries.

If your question or problem is related directly to your Agilent LC instrument, please contact your local Agilent Sales & Support organization for assistance.

In any communication with the Thermo Fisher Scientific or Agilent Technologies support teams regarding a problem, please clearly state the following:

- Your name, address, e-mail address and telephone number.
- Your Chromeleon version number together with installed Chromeleon updates.
- Your instrument driver information is listed in the Agilent Software Verification Tool report. Please run this tool by navigating to **Start > AgilentTechnologies > Software Verification Tool** from your Windows operating system.
- Your instrument information can be found in Chromeleon 7 by accessing the LC Instrument Status dashboard and clicking **i**.
- A description of the problem including any errors that were displayed in the Instrument Audit Trail, what you were trying to do when the problem occurred and the frequency of the problem.

Supported Agilent LC Modules and Firmware

Pumps

Module No.	Module Name	Min. Firmware	Min. version of Agilent Drivers for Chromeleon 7
G1310A	1200 Isocratic Pump	A.06.10	1.1
G1310B	1260 Infinity Isocratic Pump	A.06.32	1.1
G1311A	1200 Series Quaternary Pump ¹	A.06.10	1.1
G1311B	1260 Infinity Quaternary Pump ¹	A.06.32	1.1
G1311C	1260 Infinity Quaternary Pump VL ¹	A.06.32	1.1
G1312A	1200 Series Binary Pump ¹	A.06.10	1.1
G1312B	1260 Infinity Binary Pump ¹	A.06.10	1.1
G1312C	1260 Infinity Binary Pump VL ¹	A.06.32	1.1
G1361A	1260 Infinity Preparative Pump	A.06.50	1.1
G1376A	1260 Infinity Capillary Pump	A.06.10	1.1
G2226A	1260 Infinity Nanoflow Pump	A.06.10	1.1
G4204A	1290 Infinity Quaternary Pump ¹	B.06.50	1.1
G4220A	1290 Infinity Binary Pump ¹	B.06.23	1.1
G4220B	1290 Infinity Binary Pump VL ¹	B.06.43	1.1
G4302A	1260 Infinity SFC Binary Pump ¹	A.06.32	1.1
G4782A	1260 Infinity II SFC Binary Pump ¹	D.07.13	1.2
G5611A	1260 Infinity Bio-inert Quaternary Pump ¹	A.06.32	1.1
G5654A	1260 Infinity II Bio-Inert Quaternary Pump ¹	D.07.01	1.1
G7104A	1290 Infinity II Flexible Pump ¹	B.06.71	1.1
G7104C	1290 Infinity II Flexible Pump ¹	D.07.20	1.2
G7110B	1260 Infinity II Isocratic Pump	D.07.01	1.1
G7111A	1260 Infinity II Quaternary Pump VL ¹	D.07.01	1.1
G7111B	1260 Infinity II Quaternary Pump VL ¹	D.07.01	1.1
G7112B	1260 Infinity II Binary Pump ¹	D.07.01	1.1
G7120A	1290 Infinity II High Speed Pump ¹	B.06.71	1.1
G7161A	1260 Infinity II Preparative Binary Pump		1.2
G7161B	1290 Infinity II Preparative Binary Pump		1.2
<i>Cluster</i>			
N/A	Pumps marked with ¹ are able to build a pump valve cluster with up to two valves of type G1160A and/or G1170A with 5067-4159 or 5067-4147	See modules	1.1
N/A	1260 Infinity Preparative Pump Cluster with up to four G1361As	A.06.50	1.1

Sampling Systems

Module No.	Module Name	Min. Firmware	Min. version of Agilent Drivers for Chromeleon 7
G1328A/B	Manual Injector	N/A	No
G1330A/B	Thermostat for Agilent Sampler	N/A	1.1
G1313A	1100 Series Autosampler	A.06.10	1.1
G1329A	1200 Series Standard Autosampler	A.06.10	1.1
G1329B	1260 Infinity Standard Autosampler	A.06.10	1.1
G1367A	1100 Series Well-plate Autosampler	A.06.31	1.1
G1367B	1200 Series High Performance Autosampler	A.06.31	1.1
G1367C	1200 Series High Performance Autosampler SL	A.06.31	1.1
G1367D	1200 Series High Performance Autosampler SL+	A.06.31	1.1
G1367E	1260 Infinity High Performance Autosampler	A.06.32	1.1
G1377A	1260 Infinity High Performance Micro Autosampler	A.06.12	1.1
G1389A	1100 Series Micro Thermostatted Autosampler	A.06.10	1.1
G2258A	1260 Infinity Dual-Loop Autosampler	A.06.50	1.1
G2260A	1260 Infinity Preparative Autosampler (High flow)	A.06.50	1.1
G4226A	1290 Infinity Autosampler	A.06.31	1.1
G4303A	1260 Infinity SFC Standard Autosampler	A.06.54	1.1
G4767A	1260 Infinity II SFC Multisampler	D.07.13	1.2
G5667A	1260 Infinity Bio-inert High Performance Autosampler	A.06.32	1.1
G5668A	1260 Infinity II Bio-inert Multisampler	D.07.01	1.1
G7129A	1260 Infinity II Vialsampler	D.06.60	1.1
G7129B	1290 Infinity II Vialsampler	D.06.60	1.1
G7129C	1260 Infinity II Vialsampler	D.07.20	1.2
G7157A	1260 Infinity II Preparative Autosampler	D.07.01	1.2
G7167A	1260 Infinity II Multisampler	D.06.60	1.1
G7167B	1290 Infinity II Multisampler	D.06.60	1.1

Detectors

Module No.	Module Name	Min. Firmware	Min. version of Agilent Drivers for Chromeleon 7
G1314A	1100 Series Variable Wavelength Detector	A.06.10	1.1
G1314B	1260 Infinity Variable Wavelength Detector VL	A.06.10	1.1
G1314C	1260 Infinity Variable Wavelength Detector VL+	A.06.10	1.1
G1314D	1200 Series Variable Wavelength Detector	B.06.32	1.1
G1314E	1290 Infinity Variable Wavelength Detector	B.06.32	1.1
G1314F	1260 Infinity Variable Wavelength Detector	B.06.32	1.1
G1315A	1100 Series Diode Array Detector	A.06.10	1.1
G1315B	1200 Series Diode Array Detector	A.06.10	1.1
G1315C	1260 Infinity Diode Array Detector VL+	B.06.30	1.1
G1315D	1260 Infinity Diode Array Detector VL	B.06.30	1.1
G1365A	1100 Series Multiple Wavelength Detector	A.06.10	1.1
G1365B	1200 Series Multiple Wavelength Detector	A.06.10	1.1
G1365C	1260 Infinity Multiple Wavelength Detector	B.06.30	1.1
G1365D	1260 Infinity Multiple Wavelength Detector VL	B.06.30	1.1
G1321A	1200 Series Fluorescence Detector (FLD)	A.06.10	1.1
G1321B	1260 Infinity Fluorescence Detector Spectra	A.06.32	1.1
G1321C	1260 Infinity Fluorescence Detector	A.06.54	1.1
G1362A	1260 Infinity Refractive Index Detector	A.06.10	1.1
G4212A	1290 Infinity Diode Array Detector	B.06.30	1.1
G4212B	1260 Infinity Diode Array Detector	B.06.30	1.1
G7114A	1260 Infinity II Variable Wavelength Detector	D.07.01	1.1
G7114B	1290 Infinity II Variable Wavelength Detector	D.06.70	1.1
G7115A	1260 Infinity II Diode Array Detector WR	D.07.01	1.1
G7117A	1290 Infinity II Diode Array Detector FS	D.06.70	1.1
G7117B	1290 Infinity II Diode Array Detector	D.06.70	1.1
G7117C	1260 Infinity II Diode Array Detector HS	D.07.01	1.1
G7121A	1260 Infinity II Fluorescence Detector	D.07.01	1.1
G7121B	1260 Infinity II Fluorescence Detector Spectra	D.07.01	1.1
G7165A	1260 Infinity II Multi Wavelength Detector	D.07.01	1.1
G4218A	1260 Infinity Evaporative Light Scattering Detector	1.3	No
G4260A	380-ELSD	25.00	1.1
G4261A	385-ELSD	25.00	1.1
G4260B	1260 Infinity II Evaporative Light Scattering Detector	30.35	1.1
G4261B	1290 Infinity Evaporative Light Scattering Detector	30.35	1.1
G7102A	1290 Infinity II Evaporative Light Scattering Detector	30.42	1.1
G7162A	1260 Infinity II Refractive Index Detector	D.06.76	1.1
G7162B	1290 Infinity II Refractive Index Detector	D.06.76	1.1

Module No.	Module Name	Min. Firmware	Min. version of Agilent Drivers for Chromeleon 7
<i>Cluster</i>			
HDR-DAD Cluster	2 × G4212A or 2 × G4212B or a combination of 1 × G4212A and 1 × G4212B	B.06.57	1.1
HDR-DAD Cluster	2 × G7117A or 2 × G7117B or a combination of 1 × G7117A and 1 × G7117BB	B.06.70	1.1

Column Compartments

Module No.	Module Name	Min. Firmware	Min. version of Agilent Drivers for Chromeleon 7
G1316A	1260 Infinity Thermostatted Column Compartment	A.06.10	1.1
G1316B	1200 Series Thermostatted Column Compartment SL	A.06.10	1.1
G1316C	1200 Series Thermostatted Column Compartment SL	A.06.14	1.1
G7116A	1260 Infinity II Multicolumn Thermostat	D.07.01	1.1
G7116B	1290 Infinity II Multicolumn Thermostat (Host with firmware B.06.75/D.06.75 required)	C.06.70	1.1
G7130A	Integrated Column Compartment ICC (option to G7129A/B)	C.06.76	1.1
<i>Cluster</i>			
N/A	Cluster with up to three G1316C with integrated 8-pos/9-port valves (product G4230A/B) or a minimum of two G13161C TCCs; the third TCC can be a G1316A, B or C.	See module	1.1

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.

Module	Min. Module Firmware	Min. Host Module Firmware	Min. version of Agilent Drivers for Chromeleon 7
G7116B	C.06.75	B.06.75/D.06.75	1.1
G1170A	C.06.75	B.06.75/D.06.75	1.1
G7130A (option of G7129A/B)	D.06.76	N/A	1.1
G1316C	A.06.55	N/A	1.1
G1316A/B	A.06.10	N/A	1.1

Quick Change Valves for Agilent LC Modules

Module No.	Module Name	Min. Firmware	Min. version of Agilent Drivers for Chromeleon 7
G1157A	1200 Series 2-Position/10-Port Valve	A.06.02	1.1
G1158A	1200 Series 2-Position/6-Port Valve	A.06.02	1.1
G1158B	1200 Series 2-Position/6-Port Valve (600bar)	A.06.02	1.1
G1159A	1200 Series 6-Position Selection Valve	A.06.02	1.1
G1160A	1100 Series Multiple Purpose Switching Valve (12-Position/13-Port)	A.06.02	1.1
G1162A	1200 Series 2-Position/6-Port Micro Valve	A.06.02	1.1
G1163A	1200 Series 2-Position/10-Port Micro Valve	A.06.02	1.1
G1170A	1290 Infinity Valve Drive (Host required with firmware B.06.40/D.06.060)	C.06.40	1.1

Other Modules

Module No.	Module Name	Min. Firmware	Min. version of Agilent Drivers for Chromeleon 7
G1390A	1100 Series Universal Interface Box (UIB)	A.06.02	1.1
G1390B	1200 Infinity Series Universal Interface Box II (Host required with B.06.53 firmware)	C.06.53	1.1
G4227A	1290 Infinity Flexible Cube (Host required with B.06.52 firmware)	C.06.52	1.1
G1364A	1100 Series Automatic Fraction Collector	A.06.53	No
G1364B	1260 Infinity Fraction Collector (preparative-scale)	A.06.53	No
G1364C	1260 Infinity Fraction Collector (analytical-scale)	A.06.53	No
G1364D	1100 Series Micro Fraction Collector	A.06.53	No
G5664A	1260 Infinity Bio-inert fraction collector AS	A.06.53	No
G4240A	1260 Infinity Chip Cube MS Interface	A.06.36	No
G4301A	1260 Infinity Analytical SFC System	A.03.07	1.1
G7170B	1290 Infinity II MS Flow Modulator		1.2
<i>Cluster</i>			
N/A	Any combination of G1364A/B/C or G5664A plus a fourth G1364A/B/C or G5664A for recovery can be clustered. Multiple single Fraction Collectors are not supported	See module	No

Agilent LC Systems (1120 Compact, 1220 Infinity LC System, 1220 Infinity II LC System)

Module No.	Module Name	Min. Firmware	Min. version of Agilent Drivers for Chromeleon 7
G4286A	1120 LC Isocratic	B.06.50	1.1
G4286B	1220 LC System Isocratic, Man. Inj., VWD, 600 bar	B.06.50	No
G4287A	1120 LC Isocratic with Oven and ALS	B.06.50	1.1
G4287B	1220 LC Isocratic, ALS, TCC, VWD, 600 bar	B.06.50	1.1
G4288A	1120 LC Gradient	B.06.50	1.1
G4288B	1220 LC Gradient, Man. Inj., VWD, 600 bar	B.06.50	No
G4289A	1120 LC Gradient with Oven	B.06.50	1.1
G4289B	1220 LC Gradient, ALS, TCC, VWD, 600 bar	B.06.50	1.1
G4290A	1120 LC Gradient with oven and ALS	B.06.50	1.1
G4290B	1220 LC Gradient, ALS, Man. Inj., TCC, VWD, 600 bar	B.06.50	1.1
G4291B	1220 LC Isocratic, Man. Inj., TCC, VWD, 600 bar	B.06.50	No
G4292B	1220 LC Isocratic, ALS, VWD, 600 bar	B.06.50	1.1
G4293B	1220 LC Gradient, ALS, VWD, 600 bar	B.06.50	1.1
G4294B	1220 LC Gradient, ALS, TCC, DAD, 600 bar	B.06.50	1.1
G4288C	1220 LC System VL Gradient, Man. Inj. VWD, 400 bar	B.06.50	No
G4289C	1220 LC System VL Gradient, Man. Inj. VWD, 400 bar	B.06.50	No
G4290C	1220 LC System VL Gradient, ALS, TCC, VWD, 400 bar	B.06.50	1.1
G4293C	1220 LC System VL Gradient, ALS, VWD, 400 bar	B.06.50	1.1

Capillary Electrophoresis

Module No.	Module Name	Min. Firmware	Agilent Drivers for Chromeleon 7
G7150A	G7100 Capillary Electrophoresis II	B.06.25	No
G7151A	Diode Array Detector for CE	B.06.25	No

Driver Add-Ons and Special Driver Features

Module No.	Module Name	Min. Firmware	Min. version of Agilent Drivers for Chromeleon 7
Additional Driver Features	External Contacts Board G1351A		1.1
Additional Driver Features	Blend Assist		Not supported
Additional Driver Features	ISET G2197AA I		1.1
	ISET G2197AA II		1.1
	ISET G2197AA III		1.1
	ISET G2197AA IV		1.1

Module No.	Module Name	Min. Firmware	Min. version of Agilent Drivers for Chromeleon 7
Special Solutions	Buffer Advisor (G5617AA)	No	1.1 (Import Buffer Files)
	2DLC (G2198AA)	No	No
	Method Scouting Wizard (G2196AA)	No	No
	Automated Purification Software (M8368/M8369AA)	No	No

