



OpenLab EZChrom A.04.10

## Release Notes

# Notices

© Agilent Technologies, Inc. 2010-2020

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.

## Manual Part Number

M8201-90208  
Rev. B

## Edition

09/2020

Agilent Technologies, Inc.

## Software Revision

This guide is valid for Agilent OpenLab EZChrom version A.04.10

Microsoft ® is a U.S. registered trademark of Microsoft Corporation.

## 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

If software is for use in the performance of a U.S. Government prime contract or subcontract, Software is delivered and licensed as "Commercial computer software" as defined in DFAR 252.227-7014 (June 1995), or as a "commercial item" as defined in FAR 2.101(a) or as "Restricted computer software" as defined in FAR 52.227-19 (June 1987) or any equivalent agency regulation or contract clause. Use, duplication or disclosure of Software is subject to

Agilent Technologies' standard commercial license terms, and non-DOD Departments and Agencies of the U.S. Government will receive no greater than Restricted Rights as defined in FAR 52.227-19(c)(1-2) (June 1987). U.S. Government users will receive no greater than Limited Rights as defined in FAR 52.227-14 (June 1987) or DFAR 252.227-7015 (b)(2) (November 1995), as applicable in any technical data.

## 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

<b>Introduction</b>	<b>7</b>
For our Regulated Customers	7
<b>Version A.04.10</b>	<b>8</b>
Infrastructure Support Changes	8
Agilent LC Updates	8
Agilent GC updates	8
New for OpenLab EZChrom	8
<b>Version A.04.09</b>	<b>11</b>
Infrastructure Support Changes	11
Agilent LC Updates	11
Agilent GC updates	11
New for OpenLab CDS EZChrom	12
<b>Version A.04.08</b>	<b>13</b>
Infrastructure Support Changes	13
Agilent LC Updates	13
Agilent GC updates	13
New for OpenLab CDS EZChrom	13
<b>Version A.04.07 SR2</b>	<b>14</b>
Infrastructure Support Changes	14
Agilent LC Updates	14
Agilent GC updates	14

Third Party Instrument control drivers	14
New For OpenLab CDS EZChrom	15
<b>Version A.04.07 SR1</b>	<b>16</b>
<b>Version A.04.07</b>	<b>17</b>
Updated Agilent LC/CE RC.NET Driver, Version A.02.10, support of independent driver release A.02.11	17
Updated Agilent GC RC.NET Driver install package A.02.05	17
New Graphical Sample Entry (LC only)	17
New OpenLab CDS VL Product Option	18
Database Support	18
Operating System Support	18
OpenLab Control Panel	18
Classic Agilent 6890 GC	19
<b>Version A.04.06</b>	<b>20</b>
Licensing	20
Changes to Operating System support	20
M8620AA OpenLab Data Store Software A.02.01	20
OpenLab CDS Shared Services	21
M8370AA OpenLab Data Analysis A.01.02	21
Intelligent Reporting Enhancements	21
Rules and Alerts	22
Updated Agilent 78xx GC Driver, Version 5.03	22
Updated Agilent LC Driver (RC.NET Drivers A.02.09)	23
Updated Agilent 68xx GC Driver, Version 6.23	23

68xx Classic driver is available but no longer supported	23
3rd Party Instrument Driver Support	23
New Software Verification Tool	24
OpenLab CDS Upgrade	24
Product and instrument driver versions available in the Help “About” box	24
<b>Version A.04.05</b>	<b>25</b>
OpenLab Control Panel Improvements	25
EZChrom Platform Improvements	25
Agilent Parts Finder	26
Intelligent Report Enhancements	26
Display or Printing of chromatograms	26
Previewing Intelligent Reports	27
OpenLab Data Store A.01.02	27
Sample Submission	27
Lab Journal	28
Updated Agilent LC Drivers	28
Updated Agilent 78xx GC Driver (Version 5.01)	28
Barcode Scanning (7890B)	29
Resource Conservation (7890B)	29
Early Maintenance Feedback EMF (7890B)	29
5890 GC	30
Updated Hitachi drivers (LaChrom and Chromaster)	30
New M8370AA OpenLab Data Analysis A.01.01	30
<b>Version A.04.04</b>	<b>32</b>

OpenLab Control Panel/Shared Services Updates	32
The Data Store Storage Configuration	32
Master Installer Updates	33
Instrument-Project locking	33
Project Navigation Option	33
EP Signal/Noise Calculations	34
Intelligent Report Updates	34
ASCII Sequence Update	34
Spectral Fluorescence Data	34
Updated Instrument Control Drivers	35
Updated Agilent LC Control Driver	35
Updated Agilent GC Control Driver	35
Hitachi LC Control Driver	35
Updated 490 micro GC Control Driver	36
NEW html-page for accessing OpenLab Documentation and Manuals	36
<b>Version A.04.03</b>	<b>37</b>
Master Installer Updates	37
Reprocessing status while analyzing	37
Intelligent Report Update	37
Uploading Result Flag	37
New Product Introductions	38
OpenLab CDS EZChrom Value-Line	38
OpenLab CDS EZChrom Edition Compact	38
New Instrument Control Drivers	38

Updated Agilent LC Driver	38
New Agilent CTC RC.NET Driver	39
Updated GC Driver	39
<b>Version A.04.02</b>	<b>40</b>
Agilent OpenLab ECM connectivity	40
Application Virtualization	40
OpenLab Control Panel/Shared Services Updates	41
Master Installer	41
New Instrument Controls	41
iControl Panel Tool	42
<b>Version A.04.01</b>	<b>43</b>
Intelligent Reporting	43
Additional Workflows	43
New Result mode	43
Result Package Mode	44
Master Method mode	44
Instrument improvements	44
OpenLab Control Panel / OpenLab Shared Services	45
Instrument Management:	45
User Management:	45
License Management:	45
Project Management:	46
New Licensing	46
New Communication Layer	46

Printing	47
<b>Software Status and Release Bulletins</b>	<b>48</b>



# Introduction

This document provides a listing of the major feature modifications made in each release of the Agilent OpenLab EZChrom product. The following includes items modified in the Shared Services core product.

References to product documentation regarding known issues and workarounds are also provided.

## 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.

## Version A.04.10

### Infrastructure Support Changes

- Support for OpenLab Server 3.4
- Support for OpenLab ECM 3.6 (ECM supports an English system only)
- A.04.10 supports 7.15 [LTSR] version
- Windows Server 2019

### Agilent LC Updates

- Support for Agilent LC A.02.19 SR3 driver

### Agilent GC updates

- Support for Agilent GC 3.3.65 (DSA GC 7.3 [65]) driver

### New for OpenLab EZChrom

- **IR to CSV export:** A new workflow has been established with this update. EZChrom provides an option to specify the Result Set folder as the Export Path. This change allows a user to enter a tag <ResultSet> in the **Export Path** column of **Method > Advanced Reports**. The same can be achieved by keeping the **Export Path** field empty while defining the Method exports. EZChrom automatically fills the <ResultSet> tag when saving the method. The tag is not case sensitive.

This feature works for acquisition and result set processing. The report will not be exported to a result set folder when performing Simple Analysis (Analysis > Analyze).

The export file name format in the result name will be:  
<Data File Name Without Extension>\_<Line #>\_MAR\_<Template Name Without Extension>.<Export File Extension>

**Note:** Intelligent Reports can only be exported in Microsoft Excel (.xls) and CSV (.csv) file formats. Advanced Reports can only be exported in a Text (.txt) file format.

The table below lists the expected behavior of the Method Advanced Report export:

Report Type	Export Path		Export File Name/Extension	Expected behaviour
	<ResultSet>	External Path		
Intelligent Report (.rdl)	Yes	<Not Specified>	Blank (Note: Default extension is .csv)	Report is exported in Result Set folder by name format: <Data File Name Without Extension>_<Line #>_MAR_<Template Name Without Extension>.csv
Intelligent Report (.rdl)	Yes	<Not Specified>	File Name with extension specified (e.g. ExportFile.xls or .xls)	Report is exported in Result Set folder by name format: <Data File Name Without Extension>_<Line #>_MAR_<Template Name Without Extension>.xls  In case of export to Result Set, the file names have predefined file name format. File name specified by user is ignored, but extension is used.
Intelligent Report (.rdl)	<Not Specified>	Yes	This field cannot be blank. (Note: Default extension is .xls)	Report is exported to specified path with the file name extension given in Export File Name field. The same report is also exported to result set, name is predefined as mentioned in previous rows and extension is taken from Export File Name field.
Advanced Report (.tpl)	Yes	<Not Specified>	Blank (Note: Default extension is .txt)	Report is exported in Result Set folder by Name: <Data File Name Without Extension>_<Line #>_MAR_<Template Name Without Extension>.txt

Report Type	Export Path		Export File Name/Extension	Expected behaviour
	<ResultSet>	External Path		
Advanced Report (.tpl)	Yes	<Not Specified>	File Name with extension specified (e.g. ExportFile.txt)	Report is exported in Result Set folder by Name: <Data File Name Without Extension>_<Line #>_MAR_<Template Name Without Extension>.txt
Advanced Report (.tpl)	<Not Specified>	Yes	This field cannot be blank. (Note: Default extension is .txt)	Report is exported to specified path with the file name extension given in Export File Name field. The same report is also exported to result set, name is predefined as mentioned in previous rows and extension is taken from Export File Name field.

## Version A.04.09

### Infrastructure Support Changes

- Support for OpenLab Server 2.3
- Support for OpenLab ECM 3.5 (ECM supports an English system only)
- A.04.09 supports only Citrix 7.8 and 7.15 [LTSR] versions
- Windows Server 2016

### Agilent LC Updates

- Support for Agilent LC A.02.19 driver

### Agilent GC updates

- Support for Agilent GC B.01.04 driver

## New for OpenLab CDS EZChrom

- An “enable/disable multiple signature signing from the same level” toggle option in the Edit Signatures window in the Project Settings. With this feature, all new OpenLab CDS EZChrom projects have the option of multiple users with the same level to perform multiple signatures at the same level.
  - Result sets will be locked after final e-signature approval is done.
  - After e-signature is applied, no further change is allowed, especially changes to the raw data set due to the new technical implementation of the “marker file.” This was fixed in A.04.07 SR2 Update 6 and included in A.04.09 as a roll up.
- New snippet “ResultSetVersion” is available for Intelligent Report under Special Objects > Live System Values. “Result set version” can be reported for systems with ECM 3.x and ECM XT. “Result set version” for File System will be reported as “N/A”.

**Note:** The versioning format reported in Standard Report/Method Report is different than the format in Intelligent Report, where [Sequence version][modified state] is reported for ECM 3.x and [Sequence version][Result Set Version] is reported for ECM XT (Data Store/Content Management) in Standard Report.

## Version A.04.08

### Infrastructure Support Changes

- Support for Mixed Mode which allows OpenLab CDS EZChrom and OpenLab CDS 2.x to be run together on the same OpenLab Server
- Support for Windows 10
- Support for OpenLab Server 2.1
- Support for OpenLab ECM 3.5

### Agilent LC Updates

- Support for Agilent LC A.02.16 driver

### Agilent GC updates

- Support for Agilent GC B.01.02 driver

### New for OpenLab CDS EZChrom

- Option to enable global audit trail enforcement. With this feature all new OpenLab CDS EZChrom projects will have the audit trail automatically enabled. This feature was introduced in OpenLab CDS EZChrom A.04.07 SR2 Hotfix 2

## Version A.04.07 SR2

### Infrastructure Support Changes

- Support of OpenLab CDS A.02.02 SR2 and OpenLab CDS 2.0 sharing the same OpenLab Data Store server
- Support of OpenLab Data Store 2.0 and ECM 3.4.1 SP2

### Agilent LC Updates

- Includes Agilent LC/CE RC.NET Driver Version A.02.13

### Agilent GC updates

- Includes Agilent GC RC.Net Driver Version A.03.02
- 7890B supports Agilent 8355 Sulfur (G3488A) and 8255 Nitrogen (G3489A) Chemiluminescence Detectors.
- GC Method Editor provides a Navigation Tree of the available instrument devices such as inlets, detectors, oven, and signals for 78xx only.
- 7820A supports FPD, FPD+, COC and the TCD timed events.

### Third Party Instrument control drivers

- Third party drivers consolidated on disk 5



## New For OpenLab CDS EZChrom

- Support for Citrix 7.6 (client only)
- ATK refreshed with updated documentation and sample code.

## Version A.04.07 SR1

New workflow for editing a running sequence and editing a method in a running sequence.

## Version A.04.07

### **Updated Agilent LC/CE RC.NET Driver, Version A.02.10, support of independent driver release A.02.11**

- Support for the new Agilent 1290 Infinity II Multisampler (G7167A/B)
- With A.02.11: support of new Agilent Infinity II modules

### **Updated Agilent GC RC.NET Driver install package A.02.05**

- Agilent 78xx RC.Net GC driver, version 5.05
- Agilent 68xx RC.Net GC driver, version 6.25

### **New Graphical Sample Entry (LC only)**

Introduces a graphical representation of autosamplers and sample containers (tray and plate) to simplify the task of creating sequences or sample lists.

## New OpenLab CDS VL Product Option

An Agilent LC Core system driver is bundled with the OpenLab CDS VL core license. This driver can run a full 1260 Infinity LC Instrument with selected modules (for details, see the “Supported Instruments and Firmware” guide).

## Database Support

- Microsoft SQL Server 2012 SP2 Standard or Enterprise Ed.
- PostgreSQL 9.2
- Oracle 12c R1
- Microsoft SQL Server 2008 R2 SP2 (Standard, Enterprise, or Express Edition) supported for upgrades only
- Oracle 11g R2 supported for upgrades only

## Operating System Support

- Windows 7 SP1 professional or Enterprise (32-bit / 64-bit)
- Windows 8.1 professional or Enterprise (32-bit / 64-bit)
- Microsoft Windows Server 2012 R2

## OpenLab Control Panel

Lab-at-a-glance view now has a column called “Current Sample” which displays the name of the running sample.

## Classic Agilent 6890 GC

Beginning in OpenLab CDS EZChrom Edition A.04.06, the Classic 6890 GC is no longer supported and cannot be used to acquire data. The 68xx classic driver is only available to reprocess data previously created with a Classic 6890 GC.

There will be no new updates or hotfixes for the Classic 6890 GC.

## Version A.04.06

### Licensing

With OpenLab CDS A.02.01, Agilent is increasing the license version. With OpenLab CDS EZChrom Edition A.04.06, the license version will increment to 4.6. The license must be upgraded before installing the new OpenLab CDS EZChrom Edition A.04.06. For customers with a valid SMA, an upgrade button is shown in SubscribeNet, which allows upgrading the license to license version 4.6.

### Changes to Operating System support

- Windows 8.1 professional or Enterprise (32-bit / 64-bit)
- Windows 7 SP1 professional or Enterprise (32-bit / 64-bit)
- Microsoft Windows XP no longer supported

### M8620AA OpenLab Data Store Software A.02.01

- OpenLab Data Store has its own Installer in a separate DVD
- Support for PostgreSQL DB
- Support for up to 30 instruments. See Data Store hardware and software requirements guide for details.
- Support of OpenLab CDS EZChrom Edition and ICP-MS Workstations on the same server
- Lab Applications is not supported with this release

## OpenLab CDS Shared Services

- Updates to OpenLab Shared Services Login, Failover and Local Configuration Screens
- Support for PostgreSQL Database
- Microsoft SQL Express edition supported for upgrades only

## M8370AA OpenLab Data Analysis A.01.02

OpenLab Data Analysis is a product for chromatographic data from OpenLab CDS ChemStation and EZChrom Editions. OpenLab Data Analysis A.01.02 adds productivity, calibration and quantification enhancements, user interface improvements for data review and reprocessing and more. Refer to the Release Notes on Disk 7 for a detail list of enhancements.

## Intelligent Reporting Enhancements

- Enhanced printing of calibration curves
- Enhanced scaling and new coloring options for chromatograms
- Change of font properties for multiple report items at the same time
- Direct access to properties of a table column
- Expression support for chart axis scaling
- Simple peak filter for chart control (like table/matrix)
- Change of paper size and orientation at any time
- Improved alignment tools
- Date/time filter for the report template audit trail viewer

## Rules and Alerts

- Configurable pre-injection error handling (missing vial for a tray; plunger error for injector)
  - Skip the vial
  - Stop
- Configurable action after pressing GC stop button
  - Pushing the stop key at the GC keyboard will stop the current run and generate a report then continue to the next line in the sequence.
  - Abort

## Updated Agilent 78xx GC Driver, Version 5.03

- Hydrogen sensor support for calibration, reporting and hydrogen shutdown when hydrogen leak detected
- Improved operation following short LAN interruptions
- NPD adjust setup now available in the Maintenance dialog
- 7697A headspace to 7890B GC communication shown in status
- Improved direct communications between 7890B GC & 7697A headspace for EMF and Sleep/Wake and Vent methods



## Updated Agilent LC Driver (RC.NET Drivers A.02.09)

- ISET 3 (Intelligent System Emulation Technology) - using 1290 Quaternary Pump [G4204A]
- Support of the HDR feature for DAD detectors (G4212A/B)
- Support of the Valve Head 5067-4214 (2ps/4pt-4pt, 1200 bar) with the UVD G1170A and the Flexible Cube G4227A

## Updated Agilent 68xx GC Driver, Version 6.23

Ability to configure autosampler error handling available from the Configuration tab/ALS in the Method Editor.

## 68xx Classic driver is available but no longer supported

The 68xx classic driver is only available to reprocess data previously created with a 68xx GC. There will be no new updates or hotfixes for the 68xx Classic driver.

## 3rd Party Instrument Driver Support

- Integrated Shimadzu RC.NET LC Driver
- Integrated Hitachi Primaide Driver
- Integrated Hitachi ChromasterUltra Rs

## **New Software Verification Tool**

The IQT Report and Installation Qualification Report name have been updated to Software Verification and Software Verification Report. There is no change to functionality.

## **OpenLab CDS Upgrade**

OpenLab CDS A.02.01 allows for lab operation during an upgrade phase to minimize lab down time. This mode supports running different versions of OpenLab CDS on the same Networked or Distributed environment (for details, see “OpenLab CDS Administration” guide).

## **Product and instrument driver versions available in the Help “About” box**

Detailed product and driver versions are now available.

## Version A.04.05

### OpenLab Control Panel Improvements

- Lab-at-a-glance status now tracks sequence progress with repetitions
- Improved user editing in Administration
- Improved instrument status display defaults

### EZChrom Platform Improvements

- In “Result Review Mode”, when a user opens a method, sample prep, sequence, or data file, the last opened method and sequence file, if available, will be kept open rather than being set to un-assigned while closing the “Result Set”
- Update signal displayed in the online plot automatically whenever a method is downloaded to the instrument
- Instrument aux trace curves now available in Intelligent Reporter
- API support for automation of LC driver: Set instrument parameters from a 3<sup>rd</sup> party application (e.g., flow rate, solvent composition) for the Agilent LC driver
- Option to select Local Time or UTC in exported reports
- Display and positioning of windows for instrument setup values and instrument curves is now persisted
- Auto periodic integration during acquisition now functional
- Auto baseline subtract channel definition from sequence run type
- Easy Sequence files upload to ECM
- Cut, copy, paste functions added to method advanced
- Sequence description field display truncation extended

## Agilent Parts Finder

The new Agilent Parts Finder Tool is tightly integrated with OpenLab CDS EZChrom Edition. It can be called from the Instrument menu and allows users to quickly locate the part number and add it to the parts list or favorites and then print the parts list or save it to a file.

The parts finder supports the 7693A and 7650A ALS, 7890A and 7890B GC.

## Intelligent Report Enhancements

Table columns can now be formatted using conditional formatting based on expression results. Expressions may include custom variables and parameters.

## Display or Printing of chromatograms

- A new printing option optimizes the printout of chromatograms for black and white printers. All signals are drawn with black color.
- The drawing of baselines has been improved.
- For a series of overlaid signals, the user may now choose to limit the peak annotation to a single signal or annotate all.
- Users may turn off bitmap data compression for high detector sampling rates. This option does not affect the acquired raw data. Switching off data compression only improves the display and printing of the signal.
- The new reporting option "fraction delay" allows correction for the delay time between fraction collector and detector when marking the fractions in the chromatogram printout.

- Printing performance has been improved by now reading audit trail information from EZChrom register files instead of the ACAML file. The audit trail still gets reported as expected.

## Previewing Intelligent Reports

Interactively updated report previews can now be printed as a hard copy or saved as PDF.

## OpenLab Data Store A.01.02

With OpenLab Data Store A.01.02, the Lab Applications have been introduced to accommodate common laboratory workflows and includes two applications:

### Sample Submission

- Request an analysis - Prepare a chromatography sequence and assign it to the analyst through Data Store.
- Accept and complete the analysis - The analyst receives the sequence, runs it on an instrument using an Agilent CDS system, and stores the results into Data Store.
- Review results - Import the results, review, approve or reject them in Data Store.

## Lab Journal

- Manage lab assets, such as instruments and columns.
- Capture important lab events such as instrument service, replacements, calibrations, and etc.
- Review the operations that occurred on a specific instrument or column.
- Display simple usage statistics for instruments or columns.

In addition, the Data Store user interface has been redesigned to support the new applications.

## Updated Agilent LC Drivers

Support for Agilent 1260 Preparative LC Systems including:

- 1260 Isocratic Preparative Pump (G1361A)
- Cluster of up to four 1260 Isocratic Preparative Pumps (G1361A)
- Fraction Collector (G1364A)
- Cluster of up to three Fraction Collectors (G1364A/B/C or G5664A) and additionally one as Recovery Fraction Collector (G1364A/B/C or G5664A)
- 1260 Dual Loop Autosampler (G2258A)
- 1260 Preparative Autosampler (G2260A)

## Updated Agilent 78xx GC Driver (Version 5.01)

- Support for new Agilent 7890B GC instrument
- Improved handling of Autosampler (ALS) errors such as a missing vial; ALS abort/retry option provided in GC status user interface
- New Resource Conservation (Sleep/Wake) with 7890B
- New Early Maintenance Feedback (EMF) with 7890B

- New Barcode Scanning with Auto-Input with 7890B
- Updated Column Configuration, including the ability to lock the column
- New Syringe and Liner Configuration
- Support for the 7667A Mini Thermal Desorber (TD)
- GC Calculators integrated into the method editor
- Updated Columns Database
- New Syringe and Liner Databases
- New Keypad Lock Options

## Barcode Scanning (7890B)

Barcode scanning enables automatic transfer of column, liner and syringe information into the method.

## Resource Conservation (7890B)

Users may now reduce gas and power consumption with the new Resource Conservation functionality. The instrument can be put to sleep manually or by scheduling the time in the Wake and Sleep methods.

## Early Maintenance Feedback EMF (7890B)

User-defined EMF counters can now be set in OpenLab CDS based on the configuration of the GC. Warnings and Service Due notifications are provided based on the configured thresholds. EMF exceptions can be printed on Intelligent Reports.

## 5890 GC

With OpenLab CDS EZChrom Edition A.04.04, the 5890 GC was removed as an instrument selection in Agilent OpenLab Control Panel. It was no longer possible to configure a 5890 GC. The ability to configure a 5890 GC is now re-introduced in OpenLab CDS EZChrom Edition A.04.05 for backwards compatibility only.

The 5890 driver is only available to reprocess data previously created with a 5890 GC. The 5890 driver is not tested and no longer supported. There will be no new updates or hotfixes for the 5890 GC driver.

## Updated Hitachi drivers (LaChrom and Chromaster)

## New M8370AA OpenLab Data Analysis A.01.01

With OpenLab Data Analysis, Agilent introduces a new data analysis package to create sample reports, sequence summary reports, and cross-sequence summary reports. It is fully integrated and compatible with data from EZChrom and ChemStation Edition.

The first release is designed for data analysis in Chemical and Petrochemical laboratories and Hydrocarbon Processing Industry. Version A.01.01 features:

- Very fast reprocessing (more than 10 times faster than OpenLab CDS).
- Intuitive operation with "flat" user interface.
- Microsoft-style function ribbons and improved data selection tree allow fast access to main functions and data. Select data from multiple folders, load complete result sets, or select single samples.



- One-click peak integration tool for fast review.
- Easy sample review: New data viewing concept allows overlaying and comparing hundreds of signals. Work with both LC and GC instruments at the same time and use multiple methods and data sets in parallel.
- Automatically scale to a specific peak, ignore main peaks, or scale to the baseline. Supports customization of screen layout based on four predefined configurable layouts.

## Version A.04.04

### OpenLab Control Panel/Shared Services Updates

- Enhancements to the Licensing User Interface
- Generic projects
- Authentication for Data Store
- Data Store Synchronization

### The Data Store Storage Configuration

OpenLab CDS EZChrom Edition and Shared Services now support the use of a data store as the storage location for your scientific data. Use of data store storage is supported in networked and distributed configurations of OpenLab CDS and provides several benefits for your system:

- Central Storage with security
- Handles localized Chinese/Japanese content
- Web based data access for search, share, and review
- Result package version control
- 21 CFR 11 Compliant: e-signatures and audit trailing

## Master Installer Updates

Master Installer now supports installation on top of an existing A.04.03 system resulting in an automatic upgrade. Automatic upgrade from A.04.01 is not supported and requires uninstall.

Software installation repair is now supported using the repair utility found in the maintenance section of the master installer.

A new additional software installation tool is now available in the maintenance section of the master installer. This currently only supports the addition of data store components to an AIC or client upgraded from a previous revision. This function will support other add-ons in the future.

Scripted installation of clients and AICs is now supported. An XML file may be exported at the end of installer parameter definition which can then be used to install an identical configuration on other machines.

Master Installer now supports the installation of Data Store storage system components including servers, AICs, and clients.

## Instrument-Project locking

Projects now have a feature enabling the administrator to set a default project for instruments. This default project may also be set as the mandatory project for launching the instrument for all users.

## Project Navigation Option

Projects now have an option enabling the administrator to lock the users of the project within the defined project folders. This prevents users from opening or saving files to/from locations outside of the pre-defined project structure.

## EP Signal/Noise Calculations

The built in signal to noise custom parameter has been updated to reflect new European Pharmacopeia calculation requirements regarding baseline drift corrected noise calculation.

## Intelligent Report Updates

Reporting engine was updated to support v1.4 ACAML format. Method information from LC instrument parameters is now available in intelligent reports.

## ASCII Sequence Update

ASCII sequences now support the coding of injection volumes in the sequence to use the volume from the method. For the volume parameter insert the text "UseMethod" without quotes as the volume value in the standard ASCII sequence format.

## Spectral Fluorescence Data

OpenLab CDS EZChrom Edition now supports the capture and viewing of 3D data for Agilent LC fluorescence detectors (FLDs).

## Updated Instrument Control Drivers

### Updated Agilent LC Control Driver

- Improved status dashboard for enhanced control and interaction
- 1290 Quat Pump (G4204A)
- 1220 Infinity LC DAD (G4294B)
- Flexible Cube (G4227A)(Standalone Driver)
- UIB II (G1390B)

### Updated Agilent GC Control Driver

- Classic and Enhanced driver updates to fix minor issues
- New menu item to extend run for GC
- GC Tray user interface enhanced (RC.NET only)
- Method Audit Trail and Method Resolution Audit Trail

### Hitachi LC Control Driver

- Now built-in to be installed by default
- Chromaster LC Support
- LaChrom Elite and Ultra support on XP operating systems

## Updated 490 micro GC Control Driver

- Now enabled for use in Value-Line versions of EZChrom Edition

## NEW html-page for accessing OpenLab Documentation and Manuals

New html-page for OpenLab CDS Documentation can accessed from the START menu > All Programs > Agilent Technologies > OpenLab CDS Documentation.

## Version A.04.03

### Master Installer Updates

Master Installer now supports stack installation on top of an existing A.04.02 system resulting in an automatic upgrade. Automatic upgrade from A.04.01 is not supported.

### Reprocessing status while analyzing

Allows automation interfaces to get coefficient values of data collected after reprocessing and used data for calculations and/or integration with other applications such as LIMS. See automation documentation for details.

### Intelligent Report Update

Intelligent report engine has been updated to include the latest version (A.01.02). This provides improvements to report previews, report formats, and chromatogram displays. Engine was also updated to support v1.1, 1.2, and 1.3 ACAML formats. Newly generated ACAML information is now in the v.1.3 format.

### Uploading Result Flag

A new option is now available to upload result sets automatically to ECM as a result of reprocessing. The option can be selected or deselected.

## New Product Introductions

### OpenLab CDS EZChrom Value-Line

Provides lower cost option for workstation control of only compact instrumentation. See firmware support documentation for details on supported chromatography hardware.

### OpenLab CDS EZChrom Edition Compact

Provides lower cost option for workstation control of compact instruments and/or Micro 490GC instruments with a maximum of 2 simultaneous instrument sessions. See firmware support documentation for details on supported chromatography hardware

## New Instrument Control Drivers

### Updated Agilent LC Driver

- Now supports fraction collection modules
- Fraction Annotations
- Active Fraction Shading
- Insert a Fraction Report
- Running a Sequence with Fraction Collector
- Configuration changed notification
- Starting Location
- Delay Time



## New Agilent CTC RC.NET Driver

- Full control of Agilent CTC hardware
- Support CombiPAL, GC-PAL, LC-PAL, HTS-PAL, and HTS-PAL models
- Overlapping sample support
- Improved Configuration
- Custom cycle
- Improved Error handling
- Quick validation

## Updated GC Driver

- Variety of improvements based on reported issues
- Re-introduced Classic 6890 GC Driver  
Allows use of classic 6890 based methods from previous versions of EZChrom Software

## Version A.04.02

### Agilent OpenLab ECM connectivity

With the A.04.02 release of OpenLab CDS EZChrom Edition, ECM is now supported as a data storage location. With this feature you get the following enhancements for your data system:

Store and retrieve data directly in the Agilent OpenLab Enterprise Content Management (ECM) system from the CDS with ECM as a valid storage location and authentication provider. This provides the following benefits:

- Keyword searches for data, methods and sequences
- Web based data access for sharing and review
- Automated data archival
- Result package version control
- Enhanced security
- Consolidate CDS data with all laboratory content
- Thin client configuration support (ChemStation Edition)

### Application Virtualization

Support for OpenLab CDS clients being distributed via the following thin client solutions:

- Citrix XenApp 5.0 and 6.0
- Windows Terminal Services with Windows Server 2008 R2

## OpenLab Control Panel/Shared Services Updates

Support for injection count (injection X of Y) and remaining runtime column in the lab-at-a-glance view of the OpenLab Control Panel

Support for Oracle 11g as a database server for Shared Services

## Master Installer

Automated uninstallation of all OpenLab CDS components is now available from the master installer.

## New Instrument Controls

The following new or updated instrument control drivers are available for this release:

- Agilent 7890A with LTM support (Low Thermal Mass)
- Agilent 490/4900 Micro GC (formerly Varian)
- Varian 200X MicroGC
- Support for new Agilent 7697A Headspace Sampler
- Varian 3800/3900 GC
- SS420X A/D
- Perkin Elmer Nelson A/D
- Perkin Elmer Autosystem XL GC
- Perkin Elmer Series 200 LC

## iControl Panel Tool

This user contributed application allows for viewing of the OpenLab Control Panel information remotely. This allows users to see Instrument status as well as logbooks remotely from an iPhone, iPad or other smartphone. This application is provided on the support disc that is delivered with your OpenLab CDS order. See the installation guide provided in the Tools\OpenLab CDS iControlPanel folder on Disc 6 of the OpenLab CDS installation media for more information.

## Version A.04.01

### **Intelligent Reporting**

OpenLab CDS provides a new intelligent reporting feature. The new intelligent reporting is provided in addition to the method and advanced reporting. The new intelligent reporting allows you to create state-of-the-art reports in an industry standard reporting format. You can easily create and modify report templates with the new Report Template Editor built into the system that has an easy to use WYSIWYG interface. Report templates are saved in the standardized Report Definition Language (RDL) format that is also used by Microsoft Business Intelligence Studio.

### **Additional Workflows**

The workflows in the OpenLab CDS EZChrom Edition have been adapted to be more flexible and intuitive.

### **New Result mode**

OpenLab CDS EZChrom Edition now names a sequence of runs a result set after data has been collected. The new menu is under the File/Open dialog and will open up the entire sequence of runs in the navigation table for easy review.

## Result Package Mode

OpenLab CDS EZChrom Edition now has a mechanism to create a result package. This package will create a subdirectory and copy all necessary items for reprocessing the runs at any time in the future. Acquisition method, report templates, reports and sample prep templates can all be selected to be put into the package. In this mode, when reprocessing data any changes to the method are executed on the method contained in the result package, not the method in the methods folder. You can add additional runs to the result package at any time.

## Master Method mode

The result package mode also enables the ability to work with a “Master Method”. This is the method contained in the project/methods folder where the method that exists with the data in the result package becomes the “working” method. Within the system there are now new permissions that can restrict access to the Master Method while still allowing a user to make necessary changes to the working method.

## Instrument improvements

You can now have a single instrument configured with multiple injection sources and select the source in the method. This allows easier use of instruments such as a GC with an automatic liquid sampler, Headspace, and or valves to all be configured at the same time. The GC instrument configuration is also now part of the method.

## OpenLab Control Panel / OpenLab Shared Services

OpenLab Shared Services allow the management of users, instruments and licenses using a common interface for ChemStation and EZChrom editions of OpenLab CDS. OpenLab Shared Services is managed using the OpenLab Control Panel.

### Instrument Management:

OpenLab Control Panel allows you to set up and configure instruments individually or as groups for easier management. Instruments are launched from the OpenLab Control Panel. In addition, the Control Panel provides a lab-at-a-glance overview of the status of the instruments on a workstation or networked workstation.

### User Management:

OpenLab Control Panel allows you to manage users, groups, roles, and privileges. If you manage your users within an ECM system or a Windows domain, you can map those existing users into OpenLab Shared Services.

### License Management:

This service includes the administration of licenses that are required your instrument modules and Add-Ons. You can add or remove licenses and view the status of all licenses.

## Project Management:

OpenLab Control panel allows the creation of projects, their paths and their configuration settings for the system. You can easily group projects together and apply settings to the entire group.

## New Licensing

With OpenLab CDS EZChrom Edition, a new licensing strategy is introduced that helps you use your licenses more effectively. All instrument control, driver and Add-on licenses are floating licenses. Any instrument that starts up requests licenses from License Management, and when the instrument is closed, it returns the licenses. The new licensing is based on Flexera Software™ SubScribeNet® technology. It simplifies your software license management and license tracking and allows you to receive new or urgent software and patch download information immediately after release.

## New Communication Layer

OpenLab CDS EZChrom Edition now uses modern industry standard tools provided by Microsoft® for communication (WPF/WCF). This technology uses TCP/IP as its communication protocol and removes the necessity for the outdated DCOM protocol in systems configured with newer instruments. This allows normal firewall and port settings to be used without affecting the system allowing it to fit in a standard and safe IT environment.



## Printing

With OpenLab CDS EZChrom Edition A.04.01 a new mechanism has been implemented for generating printed results using a PDF printer and monitored folder. PDF copies of the results are always created and can be optionally printed using this new feature. This allows local and network printers to better be utilized and not have to be setup across the different components of the system to be used.

## Software Status and Release Bulletins

For known issues and workarounds in the OpenLab EZChrom software at the time of release see the Software Status Bulletins, which can be found online at:

<http://www.agilent.com/en-us/support/m82xxreva-04xssb>

You may also visit <http://www.agilent.com> for up to date issue information for all Agilent software products.

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

© Agilent Technologies, Inc. 2020

Edition 09/2020



M82 01-90208

