

Agilent OpenLAB Chromatography Data System (CDS)

Workstation Installation Guide



Agilent Technologies

Notices

© Agilent Technologies, Inc. 2010-2012, 2013

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

M8301-90005

Edition

01/2013

Printed in Germany

Agilent Technologies
Hewlett-Packard-Strasse 8
76337 Waldbronn

This product may be used as a component of an in vitro diagnostic system if the system is registered with the appropriate authorities and complies with the relevant regulations. Otherwise, it is intended only for general laboratory use.

Software Revision

This guide is valid for revision A.01.05 of Agilent OpenLAB CDS.

Microsoft ® and Windows Server ® are U.S. registered trademarks 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.

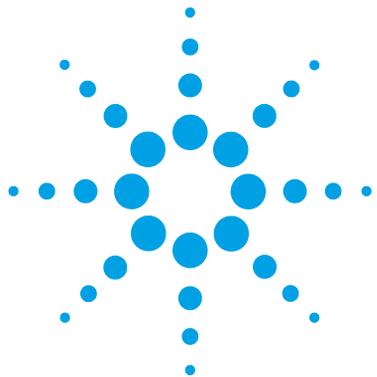
WARNING

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

1	Introduction	5
	Who Should Read This Guide?	6
	The Agilent OpenLAB CDS Software	7
	Before You Begin	8
2	Install the Software	11
	Step 1: Run the Application	12
	Step 2: Copy DVDs to a Centralized Folder for Installation (Optional)	13
	Step 3: Run the System Configuration Checker	14
	Step 4: Prepare Third Party Tools	16
	Step 5: Run the OpenLAB CDS Installation Wizard	18
	Post-Installation Tasks	23
	Optional: Install OpenLAB Data Analysis as Add-on or Standalone Application	24
	What to do Next	26
3	Uninstall the Software	27
	About Uninstallation	28
	Run the Master Installer	29
	Run the OpenLAB CDS Uninstallation Wizard	30
4	Scripted Installation	31
	About Scripted Installation	32
	Export as XML	33
	Parameters and Return Codes	34
	Installation, Upgrade, or Repair	37
	Uninstallation	38
	Logging and Tracing	39
5	Optional Procedures	41
	Run the IQ After Software Installation	42
	Upgrade OpenLAB CDS Workstation to Latest Version	43

Contents



1 Introduction

Who Should Read This Guide? 6

The Agilent OpenLAB CDS Software 7

Before You Begin 8

This chapter gives you an overview of the OpenLAB CDS software. It also includes certain requirements before starting the installation.



1 Introduction

Who Should Read This Guide?

Who Should Read This Guide?

This installation guide is designed to help system administrators and other users install the Agilent OpenLAB Chromatography Data System (CDS) to individual workstations quickly and correctly.



Figure 1 Sample workstation hardware configuration

See your Agilent representative for other system configuration installation guides.

The Agilent OpenLAB CDS Software

The Agilent OpenLAB CDS, along with OpenLAB Enterprise Content Manager (ECM), and OpenLAB Data Store With Lab Applications plus several add-ons make up the Agilent OpenLAB software suite.

These integrated solutions form a system designed to grow with customers' needs, from a small standalone analytical lab to a global network with hundreds of locations. Using the OpenLAB software, instruments from multiple vendors can connect to any size network – from individual workstations and workgroups to global enterprises. With full instrument control, centralized administration and ease-of-support, users can achieve lower operational costs and better instrument use. See your Agilent representative to learn how other products in the OpenLAB suite can help you.

The Agilent OpenLAB CDS software is provided on several edition-specific disks. Table 1 lists the disks you will receive for your software edition.

Table 1 OpenLAB CDS Components

Disk Name	ChemStation Edition	EZChrom Edition
OpenLAB CDS Shared Services Software	X	X
OpenLAB CDS ChemStation Edition Core Software	X	
OpenLAB CDS ChemStation Edition Drivers	X	
OpenLAB CDS EZChrom Edition Core Software		X
OpenLAB CDS EZChrom Edition Drivers		X
OpenLAB CDS Support	X	X
OpenLAB CDS Data Store Software	X	X
OpenLAB Data Analysis	X	X

Before You Begin

To simplify installation of the software, it is helpful to decide on some configuration options before you begin the actual software installation.

- 1 For installing OpenLAB CDS, you need to have administrator privileges for all servers and clients. Power user privileges are not sufficient (the installation does not start).
- 2 Decide on a directory location to store all files related to the data system software, including data, methods, sequences, and configurations. The directory must always be accessible to the PC running the software.
- 3 If you will be using OpenLAB ECM with your system, obtain the ECM server name.

NOTE

Make sure you have administrator privileges for both ECM and OpenLAB.

- 4 Decide on the software delivery approach you want to use:
 - *Install directly from the DVDs to your computer* – Load the disks as required directly to the workstation disk drive.
 - *Copy DVDs to a centralized location for installation* – You can use the utility to copy DVD content to a network share folder or USB drive, and run an uninterrupted installation from that location. This approach will save time; when you install from the folder, you won't have to reload the disks.
- 5 Remember that when you launch the application, you will be able to review the following PDFs in the installer **Planning** and **Resources** menus before you install the software.

Under the **Installer Planning** menu, you will find:

- *OpenLAB CDS Hardware and Software Requirements* – Use this PDF to determine whether your hardware and software will support the system.
- *Workstation Installation Guide* – An electronic copy of this installation guide is provided in PDF format for your convenience.

Under the **Installer Resources** menu, you will find:

- *OpenLAB CDS Network Requirements* –Use this PDF to check that your settings comply with the network requirements.

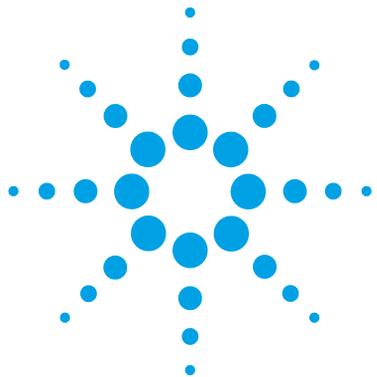
6 Set up your workstation environment:

If you are installing OpenLAB CDS to Windows 7 go to the **Windows Control Panel** and switch off **User Account Control**. After installation is complete, you can switch the controls back on.

7 Install all required hardware, including any A/D boards, cables, GPIB boards, instrument detectors, and communication cables.

1 Introduction

Before You Begin



2 Install the Software

Step 1: Run the Application	12
Step 2: Copy DVDs to a Centralized Folder for Installation (Optional)	13
Step 3: Run the System Configuration Checker	14
Step 4: Prepare Third Party Tools	16
Install Adobe PDF Reader	16
Install .NET Framework 4.0	17
Step 5: Run the OpenLAB CDS Installation Wizard	18
License Agreement Screen	18
Installation Folder Screen	19
Installation Type Screens	19
Summary Screen	21
Post-Installation Tasks	23
Optional: Install OpenLAB Data Analysis as Add-on or Standalone Application	24
What to do Next	26

The installation is automated for you by the included OpenLAB CDS Master Installer. This tool installs the various components of OpenLAB CDS. OpenLAB Data Analysis can also be installed as add-on version or standalone application.



2 Install the Software

Step 1: Run the Application

Step 1: Run the Application

To run the application for the first time on your workstation:

- 1 For the direct DVD approach, insert the OpenLAB CDS Installation disk. Autorun.inf will automatically run Agilent. OpenLABCDSSetupFromDVD.exe and display the installer **Planning** screen.
- 2 For the portable data storage device approach, insert the device in a computer USB port, navigate to \Disk1\Setup.bat and run the application to proceed to the **Planning** screen.
- 3 For the share file approach, copy all DVDs to a centralized folder as described in Step 2 below.

Step 2: Copy DVDs to a Centralized Folder for Installation (Optional)

Completing this step will enable you to run an uninterrupted installation directly from a folder—i.e., you will not have to load individual disks during actual installation.

- 1 From the Master Installer **Planning** screen, select **Installation** from the sidebar menu.
- 2 Select **Preparation of an Installation from Network Share**.
- 3 At the **Network Share** screen, browse to a directory and create a destination folder as follows:

NOTE

Installations into the root of a drive may cause problems during operations and are not supported.

- a Select the button with the three dots.
 - b Navigate to the directory where you want to create the folder (you can navigate to a USB drive port to create the folder on a USB drive).
 - c Select **Make New Folder**.
 - d Type in the folder name.
 - e Select **OK**. The system will return you to the **Network Share** screen, with the path displayed.
 - f Select the disks you want to copy to the folder: **All disks**, **ChemStation disks**, **EZChrom disks**, or **OpenLAB Data Analysis disks**.
 - g Select **Start**.
 - h Follow the prompts to insert new disks as they appear.
- 4 When processing is complete, copy the files to the local drive or map the location to a network drive.
 - 5 Close the application and navigate to the directory and folder you created. Open the folder.
 - 6 Select the Disk 1 folder, then execute Setup.bat to run the application. The system will display the installer **Planning** screen.

Step 3: Run the System Configuration Checker

- 1 At the **Planning** screen, select **System Configuration Checker**. If .NET 3.5 SP1 or higher is not on your system, the **Agilent OpenLAB CDS Prerequisites Setup Screen** appears.
 - a Read the licensing agreement.
 - b Select **Accept** to accept the agreement and begin .NET 3.5 SP1 installation.

NOTE

If you select **Don't Accept** at this screen, the install program will exit. You will have to install .NET 3.5 SP1 or higher manually to proceed with OpenLAB CDS installation.

- 2 The **Site Preparation Tool** opens. Select the correct edition and setup of the OpenLAB CDS software you are installing from the drop-down list:
 - **OpenLAB CDS ChemStation Edition C.01.XX**
 - **OpenLAB CDS EZChrom Edition A.04.XX**
 - **OpenLAB Data Analysis A.01.XX**
- 3 Select **OK**.
- 4 Complete page 1 of the **Contact Information—System details** by typing in the fields provided.
 - System Location fields
 - System Information fields
 - Configuration fields

NOTE

The regional settings must be configured to use the dot ('.') as a decimal separator.

- 5 Review the system details and make any necessary entries. The system will follow the paths specified.
- 6 Select the green check-mark icon in the top left corner of the screen to begin the software check. A summary report is displayed showing the results for each check category. Results are expressed as **Pass**, **Warning**, **Critical Warning**, or **Fail**.

Step 3: Run the System Configuration Checker

Fail results must be corrected before continuing with the installation. Agilent recommends investigating and correcting any **Critical Warnings** and **Warnings** whenever possible before proceeding.

- 7 To view details of the report, select the appropriate link: **System Hardware Details**, **Operating System and Software Details**, or **Manual Verification Required**.
- 8 To save the report, select the **Save** icon at the top left of the screen.
- 9 E-mail the saved report to your Agilent Service Representative for evaluation, and for validation of your personal computer for Agilent Software Systems Installs.

Step 4: Prepare Third Party Tools

The **OpenLAB CDS Master Installer** offers a list of tools that can be installed directly from the **Installation** screen.

Install Adobe PDF Reader

In order to view site prep or administrative reports (such as system reports), or to use the new **Report Viewer** feature (ChemStation Edition only) you need *version 11* of Adobe PDF Reader. If Adobe PDF Reader is not installed on your system, or if an older version is installed:

- 1 Select **Third Party Tools** and then **Adobe PDF Reader**.
 - a The Adobe Reader setup screen appears. Click **Install** to continue.
 - b If Adobe Reader was successfully installed, click **Finish** to exit the setup screen.

Alternatively, you can install Adobe Reader version 11 from the OpenLAB CDS installation DVDs. It is available on Disk1/Tools/Adobe Reader in 4 language versions (English, Japanese, Chinese and Portuguese).

Run the self-extracting exe-file of your choice and follow the instructions of the Adobe Reader X Setup wizard.

NOTE

If you install Adobe PDF Reader directly from the DVDs: When OpenLAB CDS users open a PDF file for the first time, they will be asked to confirm the Adobe Reader license agreement. This dialog will appear for each newly configured instrument.

Install .NET Framework 4.0

If you want to install OpenLAB Data Analysis, .NET Framework 4.0 must be installed on your system.

- 1 Select **Third Party Tools** and then **.NET Framework 4.0**.
- 2 The **.NET Framework 4 Setup** window opens. Accept the license terms to continue.
Click **Install** to continue.
- 3 If .NET 4.0 was successfully installed, click **Finish** to return to the **Installation** screen.

Step 5: Run the OpenLAB CDS Installation Wizard

At this point, you'll either be running your installation using the DVDs directly, or from a centralized location (USB drive or share folder). The installation steps are identical, until you start the installation wizard: *you won't be prompted to load the DVDs if you're installing from a centralized location.*

License Agreement Screen

- 1 From the OpenLAB CDS Master Installer screen, select **Installation**.
- 2 If you want to install OpenLAB Data Analysis, .NET Framework 4.0 must be installed on your system (see [“Install .NET Framework 4.0”](#) on page 17).
- 3 Select **OpenLAB CDS**.
- 4 If .NET 3.5 SP1 or higher is not on your system, the **Agilent OpenLAB CDS Prerequisites Setup Screen** appears.
 - a Read the licensing agreement.
 - b Select **Accept** to accept the agreement and begin .NET 3.5 SP1 installation.

NOTE

If you select **Don't Accept** at this screen, the install program will exit. You will have to install .NET 3.5 SP1 or higher at a later time to proceed with OpenLAB CDS installation.

- 5 The **OpenLAB CDS Installation Wizard** opens. Read the terms of the **License Agreement**. Master Installer provides a printable PDF of the license agreement under the **Resources** option of the main menu.
- 6 Select **I agree with the terms and conditions**. You cannot proceed with installation unless you agree to these terms.
- 7 Select **Next** to proceed to the **Installation Folder** screen.

Installation Folder Screen

- 1 Type the folder name or browse to the directory where you want to store the application components.

NOTE

Installations into the root of a drive may cause problems during operation and are not supported.

- 2 To run an installation qualification as part of this installation, select **Run Installation Qualification**. The Installation Qualification (IQ) provides documentary evidence that your system has been built and installed correctly, and that all design specifications have been met. You can run the IQ at a later time if you prefer (see [“Run the IQ After Software Installation”](#) on page 42).
- 3 Select **Next** to proceed to the **Installation type** screen.

Installation Type Screens

- 1 Under **Installation Type**, select **Standalone Workstation**.
- 2 Select **Next** to proceed to the CDS Edition screen. Select the edition of your data system.
- 3 Select **Next** to proceed to the installation screens specific to your edition.
- 4 Choose the CDS edition you wish to install.

If you chose **OpenLAB ChemStation Edition** as your edition:

- Under **Installation folder**, type the folder name or browse to the directory where you want to store the application components. Folder names must be entered without spaces.

NOTE

ChemStation Edition may not be installed into the same directory as the *OpenLAB CDS Shared Server Core*.

- Select the **Install OpenLAB Data Analysis** check box if you want to install this application in addition to your CDS edition. OpenLAB Data Analysis can be installed as add-on version at a later time if you prefer (see [“Optional: Install OpenLAB Data Analysis as Add-on or Standalone Application”](#) on page 24).

2 Install the Software

Step 5: Run the OpenLAB CDS Installation Wizard

NOTE

OpenLAB Data Analysis is not supported for installations with central data storage.

- Select the **Install IO Library** check box if you want to install the Agilent IO library suite. The suite can be installed manually at a later time if you prefer.

If you chose *EZChrom* as your edition:

- Select the **Install OpenLAB Data Analysis** check box if you want to install this application in addition to your CDS edition. OpenLAB Data Analysis can be installed as add-on version at a later time if you prefer (see [“Optional: Install OpenLAB Data Analysis as Add-on or Standalone Application”](#) on page 24).

NOTE

OpenLAB Data Analysis is not supported for installations with central data storage.

- Confirm the installation of the print server.

5 Select **Next** to proceed to the **Additional items** screen.

NOTE

If you have selected the **Install OpenLAB Data Analysis** check box, you will proceed directly to the **Summary** screen.

6 If you want to use *OpenLAB ECM* with your data system:

- a Check the **EM Server** option button. Type in a server name and press the **Test Connection...** button.
- b The system will perform a connectivity check to verify access to a functional OpenLAB ECM server. If the connectivity check is successful, the message **Connection succeeded** appears. Click **OK** to continue. If the connectivity test fails, you will be returned to the **Additional items** screen. From here you can select **Next** to run the test again. If the test is still unsuccessful:
 - Enter a new OpenLAB ECM server and try another test.
 - Call internal support for assistance if you cannot connect to an OpenLAB ECM server.

- You can uncheck the box and run the installation without OpenLAB ECM at this time. You will be able to add it to your data system at a later time, when a server is determined.

NOTE

The storage type **Data Store Server** is only available for Networked Workstation or higher.

- 7 Select **Next** to proceed to the **Summary** screen.

Summary Screen

- 1 Review the installation settings that you have selected in the preceding steps. Select **Back** as necessary to change installation settings, or **Cancel** to cancel the installation.

NOTE

Before starting or canceling the installation, you can save an XML file with your installation settings. This XML can then be used for a scripted installation (see [“About Scripted Installation”](#) on page 32).

- 2 Select **Start** to begin installation.

The system performs an automated system check before it proceeds with the listed activities.

If a *system check passed* message appears, installation continues.

If a *system check failed* message appears, you can either:

- Decline to view the system report, and continue installation.
- View the system report, and decide to continue installation.
- View the system report and postpone installation until the problem is fixed.

NOTE

To view the system report as PDF file, Adobe PDF Reader must be installed (see [“Install Adobe PDF Reader”](#) on page 16).

- Decline to view the system report and postpone.

2 Install the Software

Step 5: Run the OpenLAB CDS Installation Wizard

- 3 Your installation settings determine which disks need to be copied during installation. If you are installing directly from the DVDs, follow the **Insert Disk** prompts as they appear.

If you are installing from a local drive, you won't insert disks or browse to disk folders – the entire installation will run automatically, based on your installation settings.

If the **OpenLAB CDS Installation Wizard** cannot locate a necessary file (whether on a disk or in a folder), a **Disk or File not found** prompt will appear.

For disks, follow the prompts to:

- **Retry** using the same disk.
- Insert the correct disk (if you made an error), and select **OK**.
- Select **Cancel** to cancel the installation. The system will uninstall any components installed so far.

If a file in a centralized folder cannot be found, you may have to recreate the folder. See “[Step 2: Copy DVDs to a Centralized Folder for Installation \(Optional\)](#)” on page 13 above.

If the installation of any vital application or component fails, the installation will roll back (uninstall). If a failed component can be installed at a later time without affecting basic installation (e.g. ECM API, Driver) the installation will proceed. In this case you will be notified of any failure(s), and that further steps are necessary to fully complete installation.

- 4 If an IQ was completed as part of this installation, review the *Installation Qualification Report*. If the report indicates failure, verify the computer requirements and reinstall the data system. Do not use the system until an Installation Qualification Report gives a ‘pass’ result.
- 5 Click **Finish** to close the installation wizard.

Post-Installation Tasks

To make sure that all the net.tcp components are properly initialized, non-http activation must be enabled for your server and workstations.

- 1 Select **Start > Control Panel > Programs > Programs and Features** and click on **Turn Windows features on or off**.

The **Windows Features** window opens.

- 2 Expand the **Microsoft .NET Framework 3.5.1** node and select the **Windows Communication Foundation Non-HTTP Activation** check box.
- 3 Then click **OK**.

Optional: Install OpenLAB Data Analysis as Add-on or Standalone Application

If you have not installed OpenLAB Data Analysis together with OpenLAB CDS installation, you can install OpenLAB Data Analysis as an add-on to OpenLAB CDS or as a standalone application.

- 1 Run the Master Installer.
- 2 From the **Master Installer** screen, select **Installation**.
- 3 OpenLAB Data Analysis requires .NET Framework 4.0 on your system (see “[Install .NET Framework 4.0](#)” on page 17). If this application is already installed on your PC, you will find **Microsoft.NET Framework 4** components under **Start > Control Panel > Programs > Programs and Features**.
- 4 Select **OpenLAB Data Analysis**.
- 5 The **OpenLAB Data Analysis Installation Wizard** opens. Read the terms of the **License Agreement**. The Master Installer provides a printable PDF of the license agreement under the **Resources** option of the main menu.
- 6 Select **I agree with the terms and conditions**. You cannot proceed with installation unless you agree to these terms.
- 7 Click **Next** to proceed to the **Installation Folder** screen.
- 8 If you have installed OpenLAB CDS on your PC, OpenLAB Data Analysis will be installed as an add-on.
 - a To run an installation qualification as part of this installation, select **Run Installation Qualification**. The Installation Qualification (IQ) provides documentary evidence that your system has been built and installed correctly. You can run the IQ at a later time if you prefer.
 - b Click **Next** to proceed to the **Summary** screen.
 - c Continue with step 10.

Optional: Install OpenLAB Data Analysis as Add-on or Standalone Application

- 9 If you have not installed OpenLAB CDS on your PC, OpenLAB Data Analysis will be installed as a standalone application.
 - a You can use the default folder, type a new folder name or browse to the directory where you want to install OpenLAB Data Analysis.
 - b To run an installation qualification as part of this installation, select **Run Installation Qualification**. The Installation Qualification (IQ) provides documentary evidence that your system has been built and installed correctly. You can run the IQ at a later time if you prefer.
 - c Click **Next** to proceed to the **Installation Type** screen.
 - d Under **Data Analysis Installation Type**, choose **Standalone Workstation**.
 - e Click **Next** to proceed to the **Summary** screen.
- 10 Click **Start** to begin installation.
- 11 When the installation is finished, click **Next** to proceed to the **Installed Features** screen.
- 12 Click **Finish** to close the installation wizard.

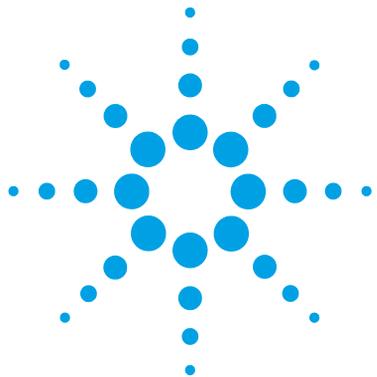
What to do Next

The basic installation of the data system software is complete.

There is a *60-day Startup License* for this system, and the expiration period starts with your first launch of an application.

To request and download your *final software license*, and add the *license file* to your system, see the *Software License Installation Guide*.

After you have acquired and installed your *final software license*, you will continue to prepare your data system for operation by end users by configuring projects, users, and instruments. This is accomplished through the *OpenLAB CDS Control Panel*, which features an *online help* resource. For more information, please contact your Agilent representative.



3 Uninstall the Software

About Uninstallation 28

Run the Master Installer 29

Run the OpenLAB CDS Uninstallation Wizard 30

This chapter contains information on the uninstallation by using the OpenLAB CDS Uninstallation Wizard.



About Uninstallation

Like the installation, the uninstallation of OpenLAB CDS is automated by the OpenLAB CDS Master Installer.

For your convenience, Master Installer uses the same user interfaces for the software uninstallation of all OpenLAB CDS configurations (standalone or networked workstation). The **OpenLAB CDS Uninstallation Wizard** under the **Maintenance** section of the Master Installer guides you through the uninstallation steps.

OpenLAB Data Analysis can be uninstalled separately (see *OpenLAB Data Analysis - Getting Started*).

Run the Master Installer

To run the Master Installer on your workstation choose one of the following approaches:

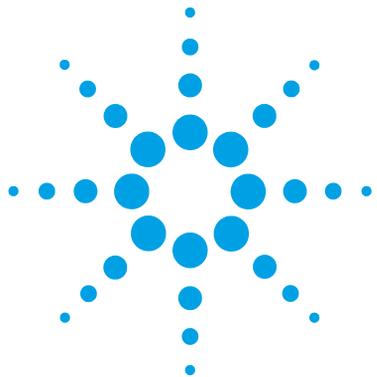
- To run the uninstallation from disk, insert the OpenLAB CDS Installation disk (Disk 1).
Autostart will automatically run Agilent.OpenLABCDSSetupFromDVD.exe and display the installer **Planning** screen.
- For the *portable data storage device* approach, insert the device in a computer USB port, navigate to \Disk1 and run Setup.bat to proceed to the **Planning** screen.
- For the *share file* approach, copy the contents to the local drive and navigate to the respective share file folder. Then navigate to \Disk1 and run Setup.bat to proceed to the **Planning** screen.

3 Uninstall the Software

Run the OpenLAB CDS Uninstallation Wizard

Run the OpenLAB CDS Uninstallation Wizard

- 1 From the Master Installer **Planning** screen, select **Maintenance** from the side bar menu.
- 2 Select **OpenLAB CDS Uninstallation**.
The **OpenLAB CDS Uninstallation Wizard** opens.
- 3 In the **Shared Components** screen, check the **Uninstall Installation Qualification** check box.
Note: IQ needs to be uninstalled if you wish to re-install OpenLAB CDS at a later time
- 4 In the **Summary** screen under **Uninstallation of OpenLAB CDS Components**, there is a list of the components you want to uninstall.
- 5 Select **Start** to start the uninstallation.
If you want to abort the uninstallation, select **Cancel**. If you want to change any settings, select **Back**.
All listed components are automatically uninstalled, one after another.
- 6 When the uninstallation has finished, click **Finish** to close the uninstallation wizard.



4 Scripted Installation

About Scripted Installation	32
Export as XML	33
Parameters and Return Codes	34
Installation, Upgrade, or Repair	37
Uninstallation	38
Logging and Tracing	39

This chapter describes the syntax and parameters for an installation or uninstallation in command line mode.



About Scripted Installation

The OpenLAB CDS Master Installer supports a command line mode for installation, the so-called *scripted installation*. This mode supports installation, upgrade, repair, and uninstallation. You can execute scripted installations either manually or as part of software management systems such as LANDesk or HP CM. With the corresponding parameter (-q), the scripted installation is completely UI-less.

Export as XML

The Master Installer supports a feature to export the installation parameters into an XML file which you can then use for the scripted installation.

This feature is also supported for upgrade and repair. However, for these cases the exported installation XML file is not appropriate. For scripted repair and upgrade, you must prepare specific XML files using the respective Master Installer wizards.

- 1 Launch the OpenLab CDS Installation Wizard.
- 2 Follow the instructions as described under *Install the Software* in this manual.
- 3 When you have reached the **Summary** screen, click the icon  on the top right corner to export the installation parameters to XML. Save the file on a physical drive.

NOTE

Installation file and XML file must not be in the same file path.

You can now use the XML file for the scripted installation.

Parameters and Return Codes

Parameters

You can call Agilent.OpenLAB.CDSInstaller.exe in command line mode with the following parameters:

- *-i*
Installation, upgrade or repair
- *-u*
Uninstallation
- *ConfigurationXML="<ConfigurationXMLFilePath>"*
The XML file contains all required inputs of the Master Installer to install, upgrade, or repair a certain topology (see “Export as XML” on page 33). Replace <ConfigurationXMLFilePath> with the correct file path and XML file name.

NOTE

Do not enter a blank before or after the equals (=) sign. The scripted installation and uninstallation mode will not work as expected.

- *KeepComponents*
Optional parameter for the uninstallation process, which can contain one or more shared components that should stay on your system. Without this parameter, all OpenLAB CDS components will be removed from your system. To keep certain shared components, list the according IDs from the table below in double quotes and separated by comma.

Component Name	Id
Installation Qualification	IQT
Microsoft SQL Server	SQLServer
IO Library	IOLibraries

- *-q*
Silent mode – no installation or uninstallation wizard will be shown.

Return Codes

After installation, uninstallation, upgrade, or repair in the command line mode, the system will return a number code which is explained below.

Table 2 Return codes

Error/Return Code	Return value
Unknown (default)	-1
Success	0
CoreComponentFailure	1
NonCoreComponentFailure	2
TestConnectivityFailure	3
ExpectedWindowsInstallerNotInstalled (WI 4.5 missing)	4
ParameterMismatchError	5
CannotProceedWithFreshInstallation	6
CannotProceedWithUpgrade	7
CannotProceedWithUninstallation	8
CannotProceedWithRepair	9
CannotProceedWithReRegistration	10
ReRegistrationNotSupported	11
IncompleteTopologyFound	12
InvalidUNCPath	13
MissingInstallable	14
NotAStrongPassword	15
DowngradeNotSupported	16
RestartRequired	17
RegistryCleanupError	18
InvalidInputXML	19
InvalidMode	20

4 Scripted Installation

Parameters and Return Codes

Table 2 Return codes

Error/Return Code	Return value
SitePrepFailure	21
DatabaseConnectionFailed	22
DotNetFramework4NotInstalled	23
OLSSConnectionFailed	24
PDFReaderNotInstalled	25
AllComponentsInstallationFailed	26
SomeComponentsInstallationFailed	27
Failed	28
AddOnListEmpty	29
EULANotAccepted	30
ScriptedNotSupported	31

Installation, Upgrade, or Repair

In installation mode, the OpenLAB CDS Master Installer checks if .Net Framework is present on your system. If not, it will automatically be installed. Select **Accept** to agree with the license agreement.

The Master Installer evaluates the products already installed on your system. Depending on the installed components, the Master Installer will offer one of the following options:

- Start a fresh installation
- Upgrade
- Repair

If a required installable is missing, the Master Installer will create an entry in a log file, and, depending on the component type, will continue or rollback the installation. A corresponding error code will be returned in such scenarios.

Preparations

You must have copied all disks to a centralized folder (see [“Step 2: Copy DVDs to a Centralized Folder for Installation \(Optional\)”](#) on page 13). This step is mandatory for scripted installation.

- 1 Right-click the executable of the command prompt or Power shell prompt, and run it as administrator.

You will get a return code for the scripted installation only if you start it as administrator.

- 2 Navigate to the drive where you have saved the disks.

For example: C:\CDS_DVD

- 3 To start the installation, call `Agilent.OpenLab.CDSInstaller.exe` with the following syntax:

```
Agilent.OpenLab.CDSInstaller.exe -i ConfigurationXML="<path to xml file>" -q
```

For example:

```
Agilent.OpenLab.CDSInstaller.exe -i ConfigurationXML="c:\settings\  
ConfigurationXML.xml" -q
```

With this command, you start the installation wizard without a user interface.

Uninstallation

- 1 Right-click the executable of the command prompt or Power shell prompt, and run it as administrator.

You will get a return code for the scripted uninstallation only if you start it as administrator.

- 2 Navigate to the drive where you have saved the disks.

For example: C:\CDS_DVD

- 3 To start the uninstallation, call `Agilent.OpenLab.CDSInstaller.exe` with the following syntax:

Agilent.OpenLab.CDSInstaller.exe -u KeepComponents="<list of components>" -q

For Example:

`Agilent.OpenLab.CDSInstaller.exe -u KeepComponents="IQT,IOLibraries" -q`

With the `KeepComponents` parameter, you can specify a list of shared components that you want to keep on the system (see “[Parameters](#)” on page 34). With the command given in the example, the OpenLAB CDS components Installation Qualification (IQT) and IO Library (IOLibraries) will be kept.

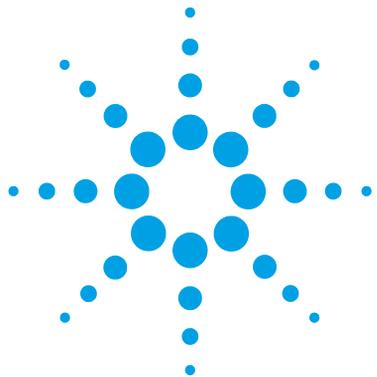
Logging and Tracing

All exceptions, errors and information messages are logged in the following locations:

- During installation, upgrade, or repair: under <BaseInstallDirectory>\Logs
- During uninstallation: under <User's Temp>\<Company Name>\Logs\<Log folder>\<Wizard Name>.txt

4 **Scripted Installation**

Logging and Tracing



5 Optional Procedures

Run the IQ After Software Installation [42](#)

Upgrade OpenLAB CDS Workstation to Latest Version [43](#)

This chapter includes information on how to use the Installation Qualification Tool, and how to upgrade your OpenLAB CDS Workstation to a higher version.



Run the IQ After Software Installation

The installation qualification (IQ) provides documentary evidence that your system has been built and installed correctly, and that all design specifications have been met.

1 Using your Windows operating system, go to **Start > All Programs > Agilent Technologies > Installation Qualification Tool**.

2 Select **Qualify**.

The system will run the application and generate an Installation Qualification Report.

3 If the report indicates failure, verify the computer requirements and reinstall the data system.

Do not use the system until an Installation Qualification Report gives a 'pass' result.

Upgrade OpenLAB CDS Workstation to Latest Version

Using the Master Installer, you can easily upgrade your OpenLAB CDS Workstation to a higher version.

To upgrade your workstation to OpenLAB CDS A.01.05, perform the following steps:

- 1 From the Master Installer **Planning** screen, switch to the **Installation** screen.
- 2 Select **OpenLAB CDS Installation**.
If OpenLAB CDS is already installed, this automatically opens the **OpenLAB CDS Upgrade Wizard**.
In the **Summary** screen of the Upgrade Wizard, the components for the upgrade are listed.
- 3 Select **Upgrading** to start the upgrade.
If an error occurs during the upgrade, an error message appears.
- 4 After the upgrade is finished, select **Finish** to close the **OpenLAB CDS Upgrade Wizard**.

www.agilent.com

In This Book

This installation guide is designed to help system administrators install the Agilent OpenLAB Chromatography Data System (CDS) to individual workstations quickly and correctly.

© Agilent Technologies 2010-2012, 2013

Printed in Germany
01/2013



M8301-90005



Agilent Technologies