



OpenLab CDS Clients and Instrument Controllers

## Installation and Configuration

# Notices

## Document Information

Document No: D0007072 Rev. A  
Edition: 04/2021

## Copyright

© Agilent Technologies, Inc. 2015-2021

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

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

USA

## Software Revision

This guide is valid for revision 2.6 of Agilent OpenLab CDS.

## Warranty

The material contained in this document is provided "as is," and is subject to being changed, without notice, in future editions. Further, to the maximum extent permitted by applicable law, Agilent disclaims all warranties, either express or implied, with regard to this manual and any information contained herein, including but not limited to the implied warranties of merchantability and fitness for a particular purpose. Agilent shall not be liable for errors or for incidental or consequential damages in connection with the furnishing, use, or performance of this document or of any information contained herein. Should Agilent and the user have a separate written agreement with warranty terms covering the material in this document that conflict with these terms, the warranty terms in the separate agreement shall control.

## Technology Licenses

The hardware and/or software described in this document are furnished under a license and may be used or copied only in accordance with the terms of such license.

## Restricted Rights Legend

U.S. Government Restricted Rights. Software and technical data rights granted to the federal government include only those rights customarily provided to end user customers. Agilent provides this customary commercial license in Software and technical data pursuant to FAR 12.211 (Technical Data) and 12.212 (Computer Software) and, for the Department of Defense, DFARS 252.227-7015 (Technical Data - Commercial Items) and DFARS 227.7202-3 (Rights in Commercial Computer Software or Computer Software Documentation).

## Safety Notices

### CAUTION

A **CAUTION** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in damage to the product or loss of important data. Do not proceed beyond a **CAUTION** notice until the indicated conditions are fully understood and met.

### WARNING

A **WARNING** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in personal injury or death. Do not proceed beyond a **WARNING** notice until the indicated conditions are fully understood and met.

# In this Guide ...

This document provides instructions for installation, configuration, administration, and maintenance of OpenLab CDS Clients and Instrument Controllers. It includes information on the license generation with SubscribeNet and operating system configuration.

**Table 1      Terms and abbreviations used in this document**

Term	Description
Content Management	Component of OpenLab Server used to manage your analytical data, including a database. Always used in Client/Server systems, optional for Workstations.
AIC	Agilent's Analytical Instrument Controller
Control Panel	Control Panel for Agilent OpenLab software
Microsoft Control Panel	Part of the Microsoft Windows operating system
Shared Services	Set of administrative services that control, for example, the security policy and the central configuration of OpenLab CDS. Shared services are accessed via the Control Panel.

## 1    Prepare your PC

This chapter describes how to configure a non-Agilent-delivered PC. Agilent-delivered PC Bundle systems are delivered with the supported pre-installed Windows operating system and are configured for optimum performance. Non-Agilent PCs require some manual configuration changes in order to provide optimum performance.

## 2    Install OpenLab CDS

This chapter describes the installation of the OpenLab CDS software in a Client/Server environment.

## 3    Post Installation Tasks

This chapter describes tasks that are relevant after finishing the installation.

## **4 Optional Procedures**

This chapter describes the installation or upgrade of additional software. It also contains information on the installation of OpenLab Help and Learning only, and on performance improvement on offline machines.

## **5 Licensing**

This chapter provides basic information on OpenLab licensing and describes how you generate a license file with SubscribeNet.

## **6 Configure OpenLab CDS**

This chapter describes the initial configuration steps after installing the software. All configuration tasks are performed in the Control Panel. For more details, refer to the Control Panel section in OpenLab Help & Learning.

## **7 Set Up NIST Library Search**

This chapter describes the installation of NIST software components, and the configuration steps in OpenLab CDS that are required for using the NIST MS library search.

## **8 About the OpenLab CDS Software**

This chapter contains an overview of the basic software features.

## **9 System Setup and Maintenance**

This chapter contains information on the Control Panel and Shared Services Maintenance. In addition, it contains information on maintenance procedures.

## **10 Upgrade to the latest OpenLab CDS Version**

This chapter describes the upgrade of the OpenLab CDS software in a Client/Server environment.

## **11 Uninstall OpenLab CDS**

This chapter describes the uninstallation of the OpenLab CDS software.

# Contents

## 1 Prepare your PC 8

- Install and Update Windows 9
- Install the .NET 3.5 and 4.x Framework 10
- Before You Begin the Installation of OpenLab CDS 11
- Run the System Preparation Tool 12

## 2 Install OpenLab CDS 19

- Installation Workflow Overview 20
- Prepare for Installation 22
- Install an OpenLab Server 23
- Install OpenLab CDS (Client or Instrument Controller) 24
- Install CDS Client on Terminal Server 33
- Silent Installation 34
- Install or Upgrade Instrument Driver Software 38

## 3 Post Installation Tasks 39

- Run a Software Verification after Software Installation 40
- Create Account for Network Access and Printing 41
- Configure the Antivirus Program 44

## 4 Optional Procedures 47

- Install OpenLab Help and Learning Only 48
- Improve Performance on Offline Machines 49

## 5 Licensing 50

- About OpenLab CDS Licensing 51
- Get a License 52
- Install Your License 57

<b>6</b>	<b>Configure OpenLab CDS</b>	<b>58</b>
	Configure Authentication	59
	Configure Security Policy	60
	Configure users and roles	61
	Configure Initial Project	66
	Configure Initial Instrument	66
	Other settings in the Control Panel	67
<b>7</b>	<b>Set Up NIST Library Search</b>	<b>68</b>
	Install Software and Library Components	69
	Configure MS Library in OpenLab CDS	79
<b>8</b>	<b>About the OpenLab CDS Software</b>	<b>80</b>
	Software Architecture	81
	21 CFR Part 11 Compliance	85
	Customization	87
<b>9</b>	<b>System Setup and Maintenance</b>	<b>89</b>
	Control Panel	90
	Other Maintenance Procedures	95
<b>10</b>	<b>Upgrade to the latest OpenLab CDS Version</b>	<b>96</b>
	License Upgrade	97
	Upgrade Process	99
	Step 1: Upgrade the OpenLab Server	101
	Step 2: Upgrade an Analytical Instrument Controller	101
	Step 3: Upgrade an OpenLab CDS Client	103
	Step 4: Upgrade Drivers	105
	Register a Client or Instrument Controller on the Server	106
	Sample Scheduler for OpenLab	109
<b>11</b>	<b>Uninstall OpenLab CDS</b>	<b>110</b>
	Uninstall a Client or Analytical Instrument Controller (AIC)	111

Uninstall the OpenLab Server	111
Uninstall OpenLab Help and Learning Only	112

## **12 Appendix 113**

Privileges in the Control Panel	114
Sales and Support Assistance	123

# 1

## Prepare your PC

Install and Update Windows	9
Install the .NET 3.5 and 4.x Framework	10
Before You Begin the Installation of OpenLab CDS	11
Run the System Preparation Tool	12
Reference of SPT Checks	17

This chapter describes how to configure a non-Agilent-delivered PC. Agilent-delivered PC Bundle systems are delivered with the supported pre-installed Windows operating system and are configured for optimum performance. Non-Agilent PCs require some manual configuration changes in order to provide optimum performance.



## Install and Update Windows

- 1 Install the Windows operating system from the Microsoft installation media or qualified PC image media provided by your IT department. During the setup, provide the computer name, administrator password and network settings.

For the computer name, do not use underscores; installation is not possible if the computer name contains an underscore.

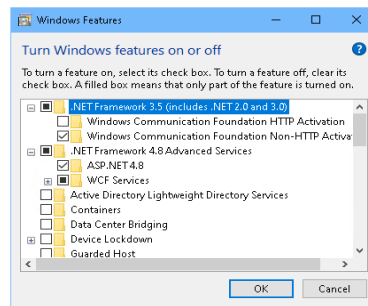
Both clients and instrument controllers must be in a Windows domain; systems in workgroup mode are not supported.

- 2 In the Microsoft Control Panel under **System > Windows activation**, click **Change product key**. Enter a valid value to activate Windows.
- 3 Update to the recommended version of Windows 10 (also known as a semi-annual Feature Update). See the FAQ section of the OpenLab CDS page at <https://www.agilent.com/>.
- 4 Check for Windows quality updates, and apply all critical security patches. Do not choose to install a newer version of Windows 10 (see previous step). Make sure the Windows quality updates have finished installing before proceeding to install OpenLab CDS.
- 5 Check that your computer meets all requirements. You can access the *Requirements and Supported Instruments for Instrument Controllers, Clients, or Workstations* guide (CDS\_Requirements.pdf) from the Setup\Docs folder on the installation medium.  
  
Use this PDF to check that your settings comply with the network requirements, and to determine whether your hardware and software will support the system.
- 6 To secure your system against viruses please install an antivirus program. Be sure to open the firewall ports listed in the Firewall Settings in the *Requirements and Supported Instruments Instrument Controllers, Clients, or Workstations* guide (CDS\_Requirements.pdf).
- 7 To configure remote settings: In the Microsoft Control Panel navigate to **System > Remote settings**. On the **Remote** tab. In the **Remote Desktop** section, select **Don't allow connections to this computer**.
- 8 In the Microsoft Control Panel under **Date and Time**: Choose the time zone of your regional location.

## Install the .NET 3.5 and 4.x Framework

If .NET 3.5 and .NET 4.x are not installed on your system, their installation will automatically be triggered by the installation wizard. However, this may require a system reboot. To avoid the system reboot during installation, install .NET in advance.

- 1 Go to the Microsoft Control Panel.  
In the Windows Start menu, enter "Control Panel" in the **Search programs and files** field (alternatively, press [Win+R] and enter "Control panel"). To view all items in the Control Panel view, select **Small icons** in the **View by** field.
- 2 Go to **Programs and Features**.
- 3 Go to **Turn Windows features on or off**.
- 4 Enable .NET 3.5 as follows:
  - a Expand the **.NET Framework 3.5 (includes .NET 2.0 and 3.0)** node.
  - b Select the **Windows Communication Foundation Non-HTTP Activation** check box.



**Figure 1** Enable .NET 3.5 (Win 10)

### NOTE

This requires an internet connection. If it does not work as expected, or the computer has no internet access, install .NET 3.5 from the Windows installation media. For Windows 10, see details 10 under <https://support.microsoft.com/en-us/kb/2734782>.

If you do not have installation media, create them as described under <https://www.microsoft.com/en-us/software-download/windows10>.

- 5 Select the **.NET Framework 4.8 Advanced Services** check box. Use the default values for sub items.

## Before You Begin the Installation of OpenLab CDS

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

**1** Decide on a computer name.

The computer name will be reflected in the license. To avoid effort, it is recommend to keep the computer name unchanged after installing OpenLab CDS.

### NOTE

To make sure that a DNS server can resolve the computer name, follow the internet standard for protocols (*RFC952*) and use only the following characters:

- Letters (a-z, A-Z)
- Digits (0-9)
- Hyphen (-)

Do not use an underscore. Installation is not possible if the computer name contains an underscore.

- 2** Install all required hardware, including any cables, instrument detectors, and communication cables. GPIB interfaces may be required for some non-Agilent instruments.
- 3** Prepare an account with administrative privileges to run the installation.
- 4** To allow the successful installation and activation of all components, make sure that the LAN-cable is connected and a network is available. A local network is sufficient. If you install the workstation without a network cable, the activation of the Content Management component will fail, and trying to run the database will return an error.
- 5** Make sure that the antivirus software is disabled during the installation.
- 6** Check the Agilent Service Notes for software updates for your installation package and software products. Service Notes are available from your Agilent support representative.
- 7** If you use Trend Micro<sup>TM</sup> as an antivirus software, turn off **Web Reputation** to allow the installation of all components.
- 8** If you plan to upgrade from a previous version of OpenLab CDS please refer to the *Upgrade OpenLab CDS* chapter.

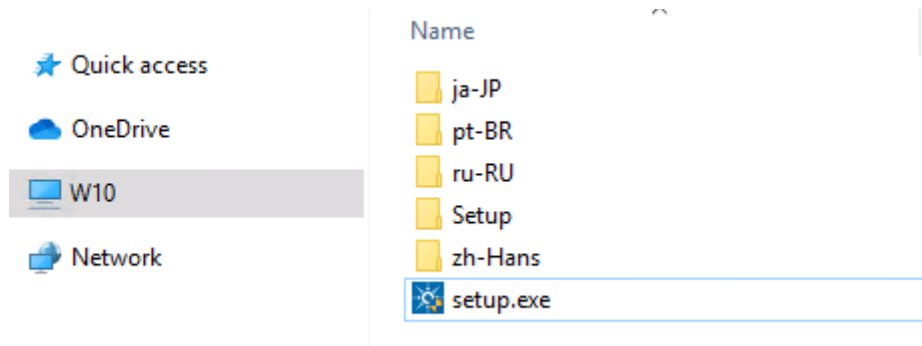
## Run the System Preparation Tool

The System Preparation Tool (SPT) checks and applies Windows settings on your machine. The settings are also applied automatically when you run the OpenLab installer. Running the SPT in advance helps you to shorten the installation process. For an overview of both mandatory and recommended settings, see [Table 2](#) on page 17.

- 1 If you want to keep your setting in the System Preparation Tool from one run to the next: Copy the entire content of the USB media to a local drive or centralized folder, then remove the USB media from the PC.
- 2 To open the installer, right-click the setup.exe file, and run it as administrator.

### NOTE

If User Account Control (UAC) is switched on, this step requires active confirmation to continue.

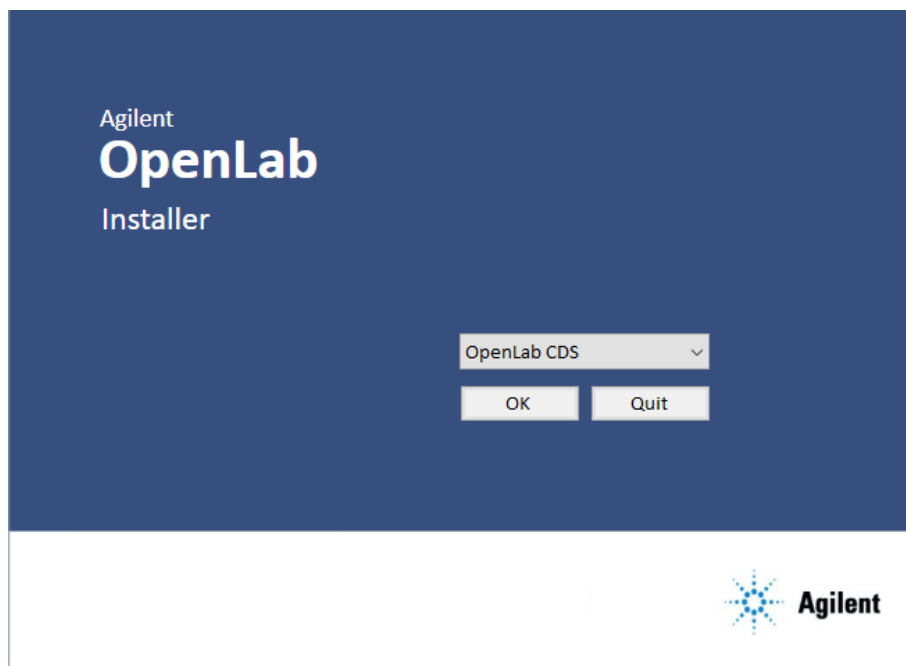


1

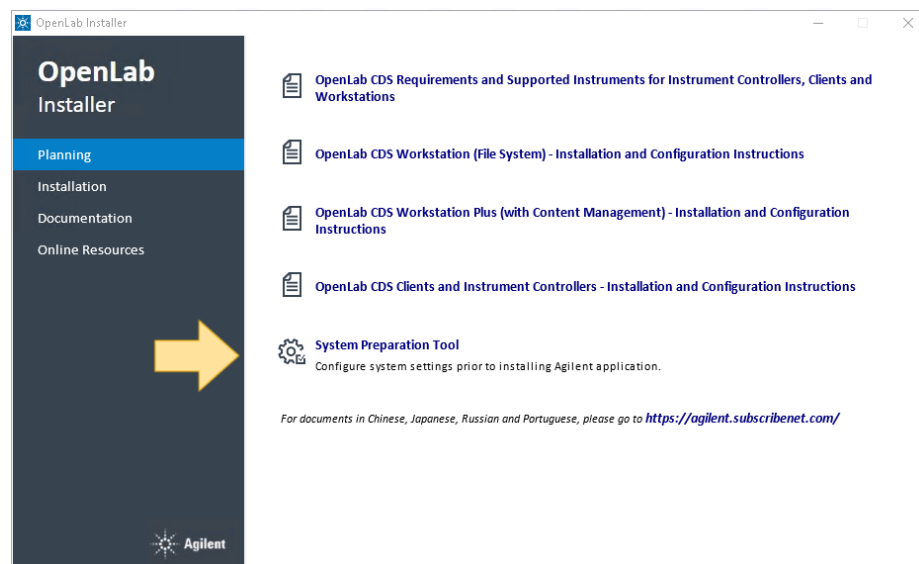
## Prepare your PC

### Run the System Preparation Tool

- 3 On the start screen, select **OpenLab CDS**, and click **OK**.



- 4 From the **Planning** tab, select **System Preparation Tool**.



The **System Preparation Tool** window opens.

## Prepare your PC

### Run the System Preparation Tool

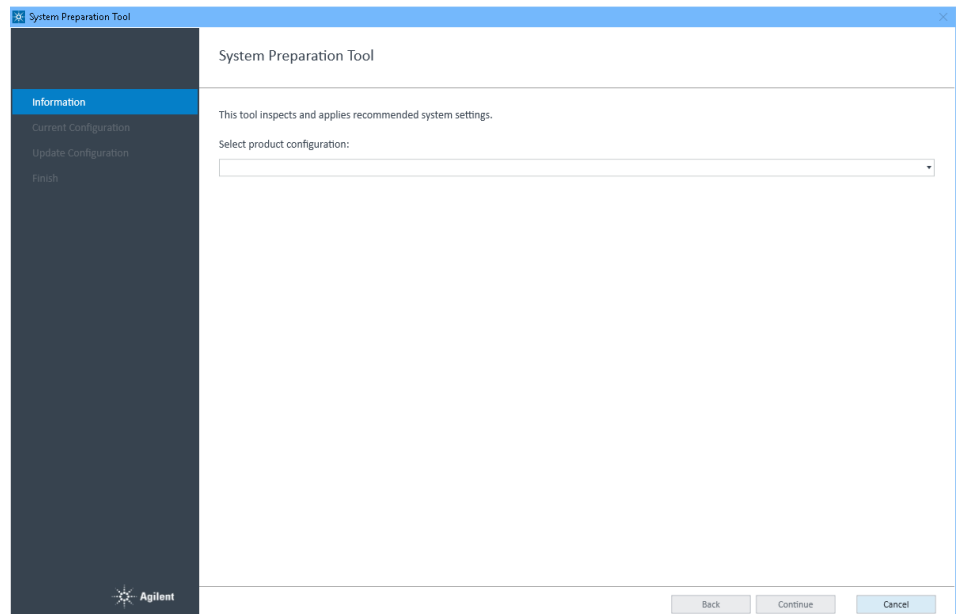
- 5 Select the product configuration corresponding to your system.

For AICs:

- **OpenLab CDS~2.6~AIC~Win10**
- **OpenLab CDS~2.6~AIC~Win2016**
- **OpenLab CDS~2.6~AIC~Win2019**

For clients:

- **OpenLab (CDS, ECMXT)~2.6~(Client, CMServices)~Win10**
- **OpenLab (CDS, ECMXT)~2.6~(Client, CMServices)~Win2016**
- **OpenLab (CDS, ECMXT)~2.6~(Client, CMServices)~Win2019**



Click **Continue**. The installer automatically applies all mandatory Windows settings to ensure proper installation.

- 6 Select which recommended settings to apply to the system.

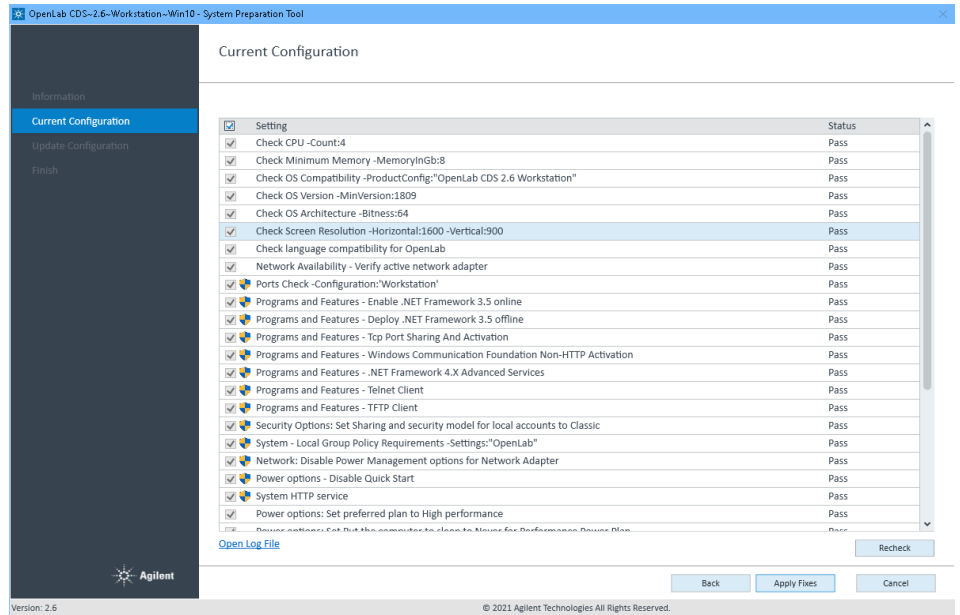
There are several settings that are recommended that can improve the performance and stability of your system, but do not need to be completed to deploy the application. The recommended settings are listed after the mandatory settings.

## Prepare your PC

### Run the System Preparation Tool

You can clear the check boxes for recommended settings. Mandatory settings cannot be cleared. Recommended actions are selected by default and will be applied unless they are cleared.

For more information on mandatory and recommended settings, see [Table 2](#) on page 17.



- 7 Click **Apply Fixes** to apply the correct settings.

The System Preparation Tool attempts to fix the selected settings and displays the new status on the **Update Configuration** page. All actions are saved to a log file.

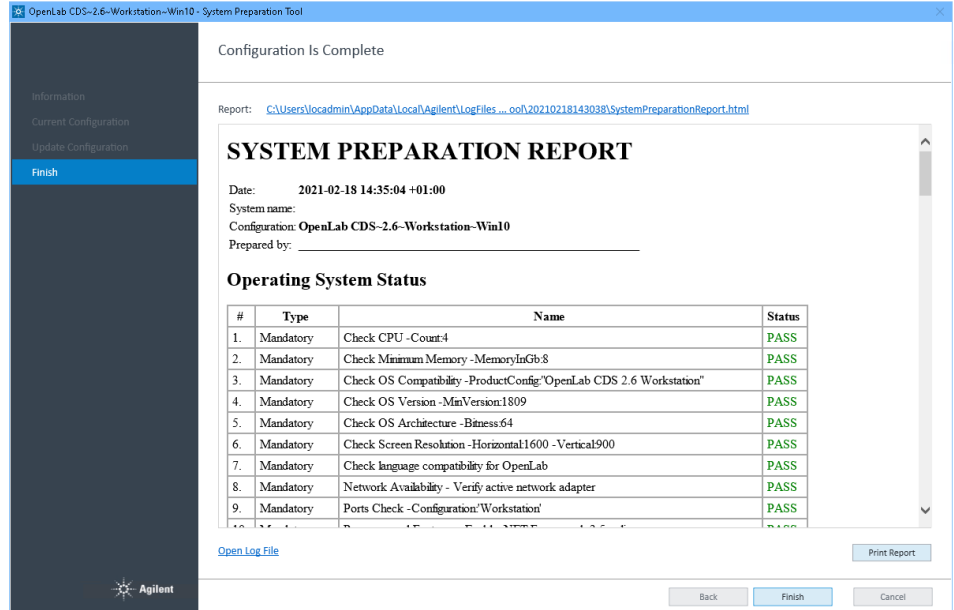
- 8 Click **Next** to proceed to the **System Preparation Report** page.

The System Preparation Report is displayed. It lists the new status for all selected settings.

The System Preparation Report for each run is saved separately to disk. Its location is shown at the top of the page.

## Prepare your PC

### Run the System Preparation Tool



9 Click **Print Report** to print the *System Preparation Report*.

You may print to a file, for example, using the *Adobe PDF* printer, and add comments.

10 The System Preparation Report lists any mandatory or recommended settings that are not automatically updated by the System Preparation Tool. Follow the instructions provided in the System Preparation report to manually update operating system settings.

11 Click **Finish**.

12 Reboot your system if requested to do so.



## Reference of SPT Checks

**Table 2** Checks and Setting updates by the System Preparation Tool

Name	Workstation Workstation Plus	Client CM Services	OpenLab Server ECM XT OLSS Server for ECM 3.x	AIC
<b>Mandatory</b>				
Check Minimum Memory	●	●	●	●
Check CPU	●	●	●	●
Check OS Compatibility	●	●	●	●
Check OS Version	●			
Check OS Architecture (64 bit)	●			
Check Screen Resolution	●	●	●	●
Check language compatibility for OpenLab	●	●	●	●
Network Availability - Verify active network adapter	●	●	●	●
Ports Check	●	●	●	●
Programs and Features - Enable/deploy .NET Framework 3.5	●	●	●	●
Programs and Features - Tcp Port Sharing And Activation	●	●	●	●
Programs and Features - Windows Communication Foundation Non-HTTP Activation	●	●	●	●
Programs and Features - .NET Framework 4.X Advanced Services	●			
Programs and Features - Telnet Client	●	●		●
Programs and Features - TFTP Client	●	●		●
System - Local Group Policy Requirements <sup>1</sup>	●	●	●	●
Power options - Disable Quick Start	●	●	●	●
System HTTP service	●	●	●	●

**Table 2** Checks and Setting updates by the System Preparation Tool

Name	Workstation Workstation Plus	Client CM Services	OpenLab Server ECM XT OLSS Server for ECM 3.x	AIC
<b>Mandatory, manual</b>				
Windows Update - Apply pending updates	●	●	●	●
Windows Activation	●	●	●	●
System domain membership	●	●	●	●
Region - Change system locale	●	●	●	●
Indexing options: Disable indexing Options for all drivers and locations			●	
<b>Recommended</b>				
System - Startup and Recovery Settings	●	●		●
System - Turn off system protection (restore points) for all drives	●			
Windows Update - Disable Windows Update service	●	●	●	●
Personalization - Disable advertising info	●	●		●
Personalization - Combine Taskbar Buttons	●	●		●
<b>Recommended, manual</b>				
File Explorer - Display Settings	●	●		●
Recycle Bin - Set Recycle Bin properties	●	●		●
System - Performance Settings	●	●		●
Privacy - Set privacy settings	●	●		●
Apps - Default browser	●	●		●
Personalization - Turn off show lockscreen background picture	●	●		●

<sup>1</sup> Local Group Policy Requirements for OpenLab: - Set "Hide entry points for Fast User Switching" to Enabled - Adds "Users" group to "Access this computer from the network" setting

## 2

# Install OpenLab CDS

Installation Workflow Overview	20
Prepare for Installation	22
Install an OpenLab Server	23
Install OpenLab CDS (Client or Instrument Controller)	24
Sample Scheduler for OpenLab	32
Install CDS Client on Terminal Server	33
Silent Installation	34
Export Properties File	34
Run Silent Installation	35
Parameters and Return Codes	36
Logging and Tracing	37
Install or Upgrade Instrument Driver Software	38

This chapter describes the installation of the OpenLab CDS software in a Client/Server environment.

## Installation Workflow Overview

### NOTE

Decide if you want to go for a cloud deployment. In this case, check with your Agilent representative or refer to the manual *OpenLab CDS Configuration in the Cloud* on the Agilent Software Service Desk:

<https://servicedesk.li.agilent.com/plugins/servlet/desk>.

### Install OpenLab Server

1. Set up the Windows Operating System
2. Install OpenLab Server including Content Management
3. Run Installation Verification

If you use the Content Management component provided by OpenLab Server or OpenLab ECM XT: Refer to the *Agilent OpenLab Server and OpenLab ECM XT* installation guide.

If you use an existing ECM 3.X system to store you data, refer to the *Configuring OpenLab CDS with ECM 3.x* guide.



### Prepare

- Check OpenLab CDS Requirements and Supported Instruments for Client/Server Systems
- Run System Preparation Tool
- Configure Windows (manual steps)

Running the System Preparation Tool before starting the installation is recommended. Running it helps to reduce or avoid reboots during the installation.

See “Run the System Preparation Tool” on page 12.

If you plan scripted installations, see “Silent Installation” on page 34.



#### Install Clients and AICs

1. Install OpenLab CDS  
on AICs and clients
2. Post Installation  
Software verification  
Set account to enable automatic printing
3. Optional  
Improve performance on offline machines  
Install Add-ons

It is recommended to install an AIC first before installing the clients.

- See "Install OpenLab CDS (Client or Instrument Controller)" on page 24
- See "Post Installation Tasks" on page 39
- See "Optional Procedures" on page 47



#### Get Licenses

1. Obtain licenses via SubscribeNet:
  - OpenLab CDS
  - Instrument products
  - Add-ons
2. Install your licenses

See "Licensing" on page 50



#### Configure

- Authentication
- Projects, incl. audit trail settings
- Instruments

See "Configure OpenLab CDS" on page 58. All configuration tasks are performed in the Control Panel. For more details, refer to the Control Panel section in OpenLab Help & Learning.

## Prepare for Installation

- 1 Make sure that the antivirus software is disabled during the installation.
- 2 Do not run the Windows Update Service during installation. Make sure that no Windows updates are performed during the installation.
- 3 Make sure that no system reboot is pending.  
Pending reboots are indicated in the System Preparation Tool (see [“Run the System Preparation Tool”](#) on page 12).
- 4 Ensure that all required ports can be used by OpenLab CDS. For details, see *Requirements and Supported Instruments for Instrument Controllers, Clients, or Workstations* (CDS\_Requirements.pdf).

## Install an OpenLab Server

If you use the Content Management component provided by OpenLab Server or OpenLab ECM XT: Refer to the *Agilent OpenLab Server and OpenLab ECM XT Installation* guide (ECM\_XT\_InstallationGuide.pdf).

If you use an existing ECM 3.X system to store you data, refer to the *Configuring OpenLab CDS with ECM 3.x* guide (CDS\_configure-with-ECM.pdf).

## Install OpenLab CDS (Client or Instrument Controller)

This procedure describes the installation of the OpenLab CDS software on a client or Analytical Instrument Controller (AIC). Carry out this procedure on each client PC and each AIC machine.

### Prerequisites

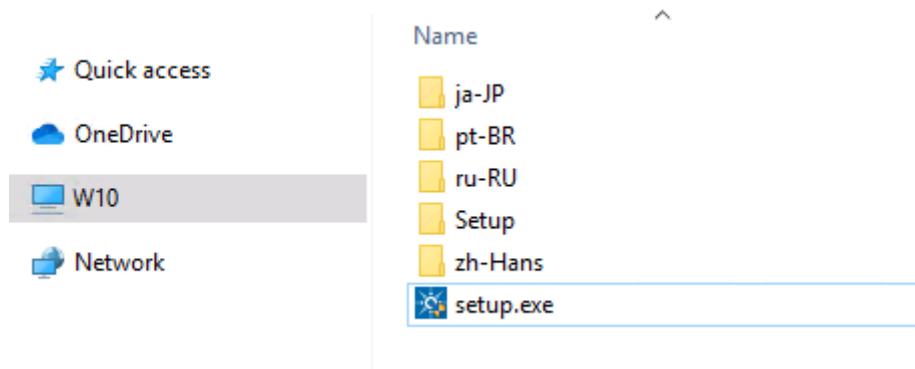
An OpenLab Server is already installed and available in your network.

OpenLab CDS versions on clients and AICs must not be higher than the OpenLab Server version.

### NOTE

If a client needs to be installed on a scalable system, connect it to virtual host (typically, the configured hostname of the load balancer) during the installation.

- 1 To open the installer, right-click the setup.exe file, and run it as administrator.

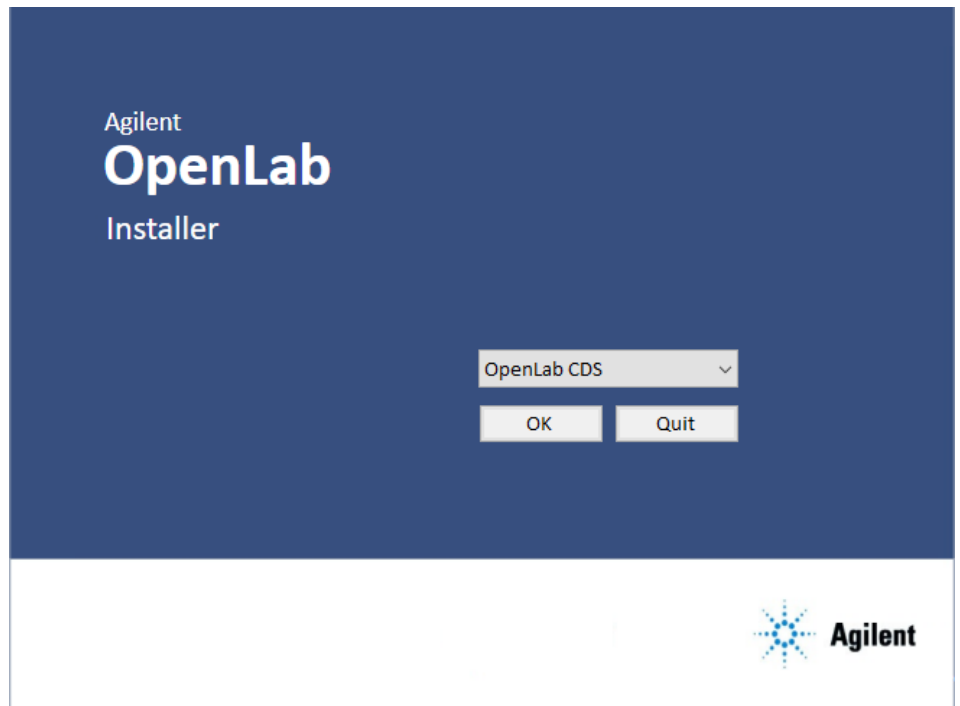




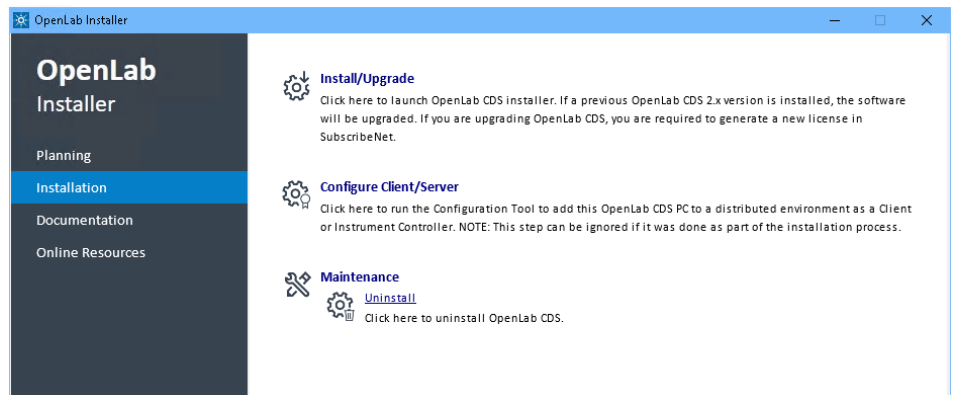
## Install OpenLab CDS

### Install OpenLab CDS (Client or Instrument Controller)

- 2 On the start screen, select **OpenLab CDS**, and click **OK**.



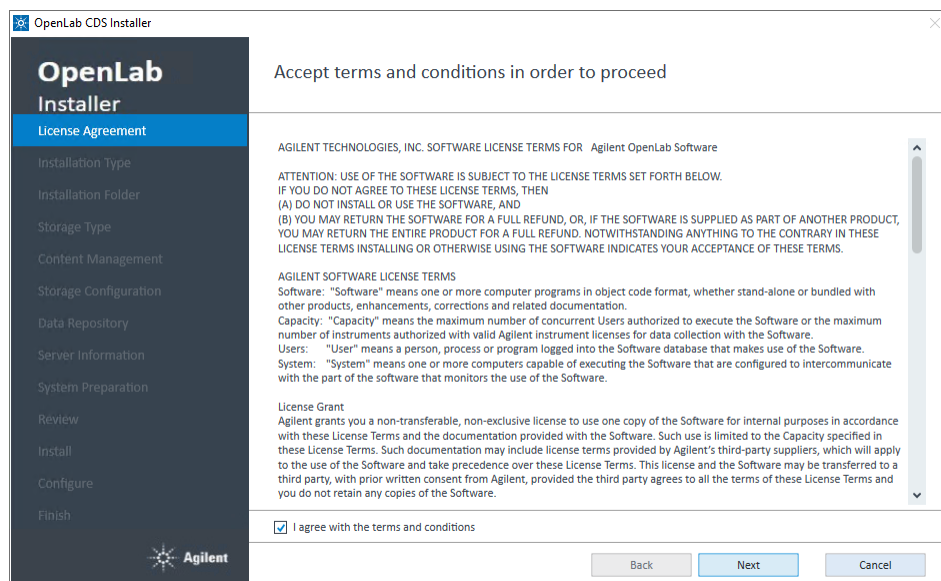
- 3 From the **Installation** tab, select **Install/Upgrade**.



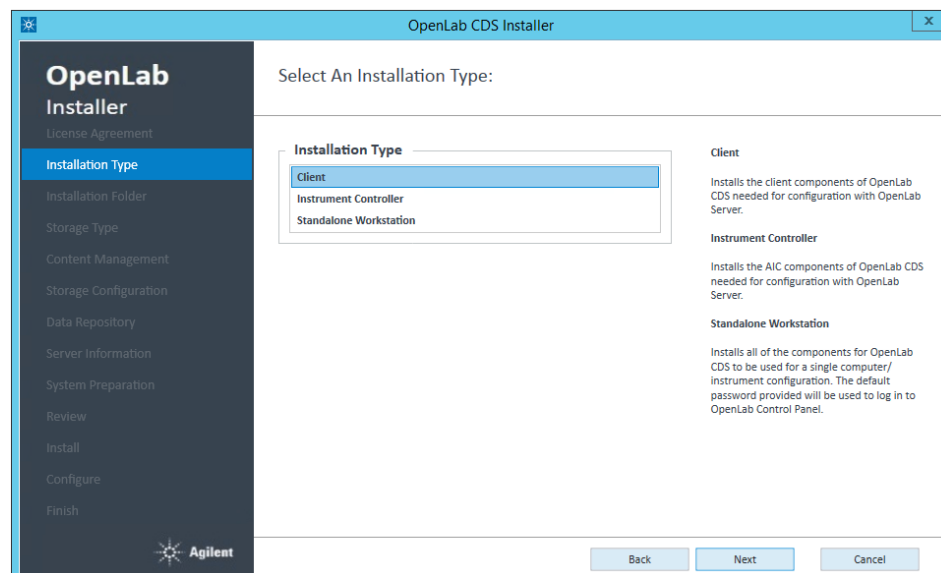
## Install OpenLab CDS

### Install OpenLab CDS (Client or Instrument Controller)

#### 4 License Agreement: Read and confirm Agilent terms and conditions.



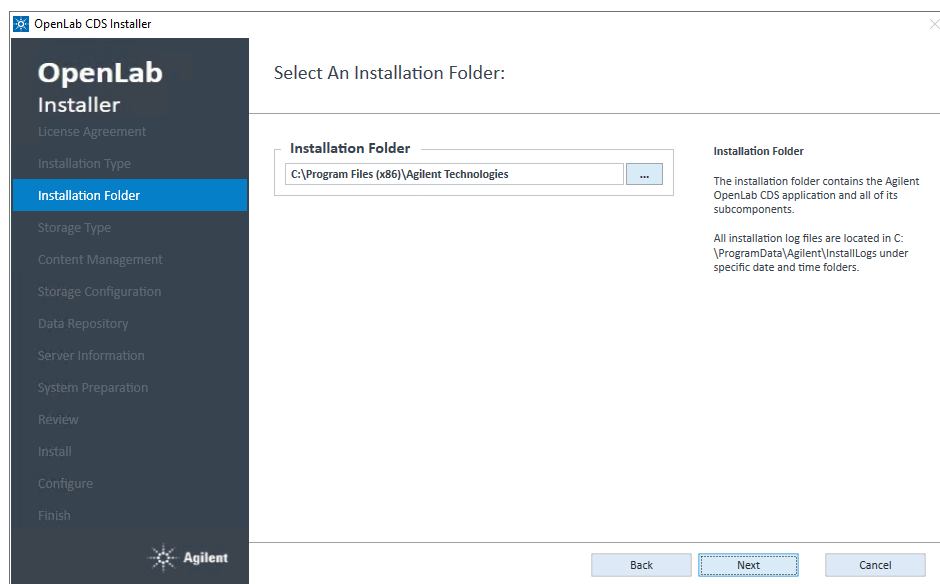
#### 5 Installation Type: Select **Client** or **Instrument Controller**, depending on what you are installing.



## Install OpenLab CDS

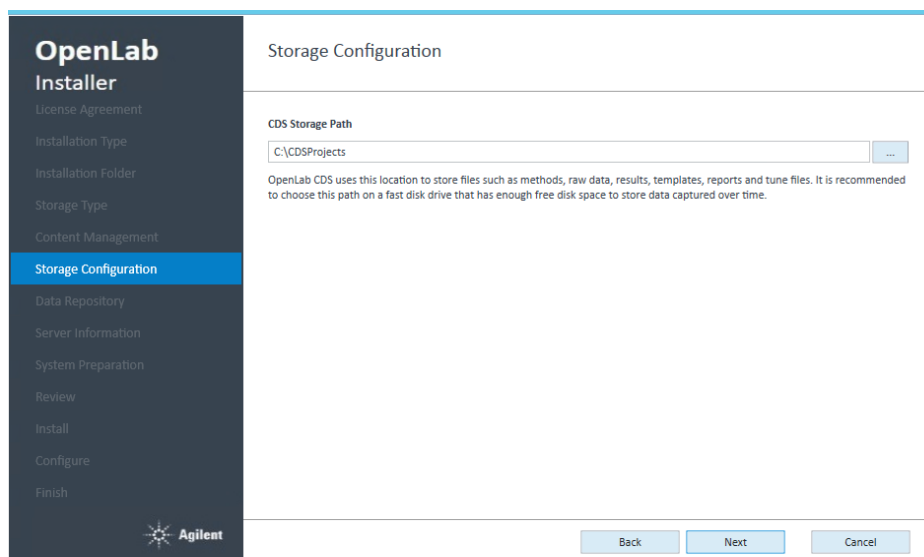
### Install OpenLab CDS (Client or Instrument Controller)

- 6 **Installation Folder:** Provide an installation folder for OpenLab CDS. Do not use the root folder of any drive.



The screenshot shows the 'OpenLab CDS Installer' window. On the left is a dark sidebar with a list of steps: License Agreement, Installation Type, **Installation Folder** (highlighted in blue), Storage Type, Content Management, Storage Configuration, Data Repository, Server Information, System Preparation, Review, Install, Configure, and Finish. The main area is titled 'Select An Installation Folder:'. It features a text box labeled 'Installation Folder' containing 'C:\Program Files (x86)\Agilent Technologies' with a browse button (...). To the right, there is explanatory text: 'The installation folder contains the Agilent OpenLab CDS application and all of its subcomponents.' and 'All installation log files are located in C:\ProgramData\Agilent\installLogs under specific date and time folders.' At the bottom right are three buttons: 'Back', 'Next' (which is highlighted with a dashed border), and 'Cancel'.

- 7 If you are installing an Instrument Controller: Provide a Projects Root Path that can be used to store projects data in a failover scenario.



The screenshot shows the 'OpenLab CDS Installer' window at the 'Storage Configuration' step. The sidebar on the left is the same as in the previous screen, but 'Storage Configuration' is now highlighted in blue. The main area is titled 'Storage Configuration'. It contains a section 'CDS Storage Path' with a text box showing 'C:\CDSProjects' and a browse button (...). Below this, a note states: 'OpenLab CDS uses this location to store files such as methods, raw data, results, templates, reports and tune files. It is recommended to choose this path on a fast disk drive that has enough free disk space to store data captured over time.' At the bottom right are three buttons: 'Back', 'Next', and 'Cancel'.

- 8 **Server Information:** Enter the hostname of the Shared Services server you would like to connect to, and click **Connect**. If you install an Instrument Controller, you must additionally provide the OpenLab Control Panel login credentials of an administrator for the specified server.

**OpenLab Installer**

- License Agreement
- Installation Type
- Installation Folder
- Storage Type
- Content Management
- Storage Configuration
- Data Repository
- Server Information**
- System Preparation
- Review
- Install
- Configure
- Finish

**Server Information**

Server Name

Enter the hostname of Shared Services server and click the Connect button.

**Connect**

**Administrator Authentication**

Enter administrator login information for the server.

Username

Password

Domain

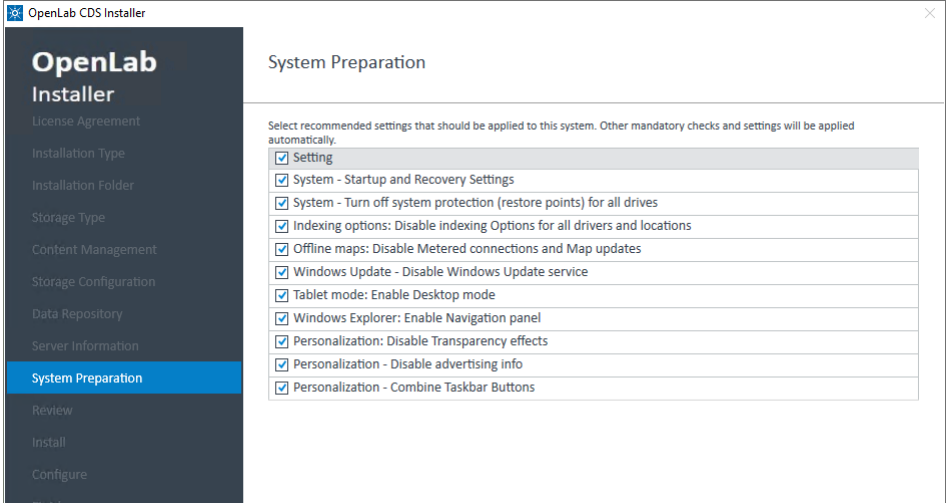
**Agilent**

**Back** **Next** **Cancel**

- 9 **System Preparation:** The installer shows the list of recommended settings for the system. You may unselect items that you do not want to apply on the system. Other mandatory settings will be applied automatically during installation. For an overview of both mandatory and recommended settings, see [Table 2](#) on page 17.

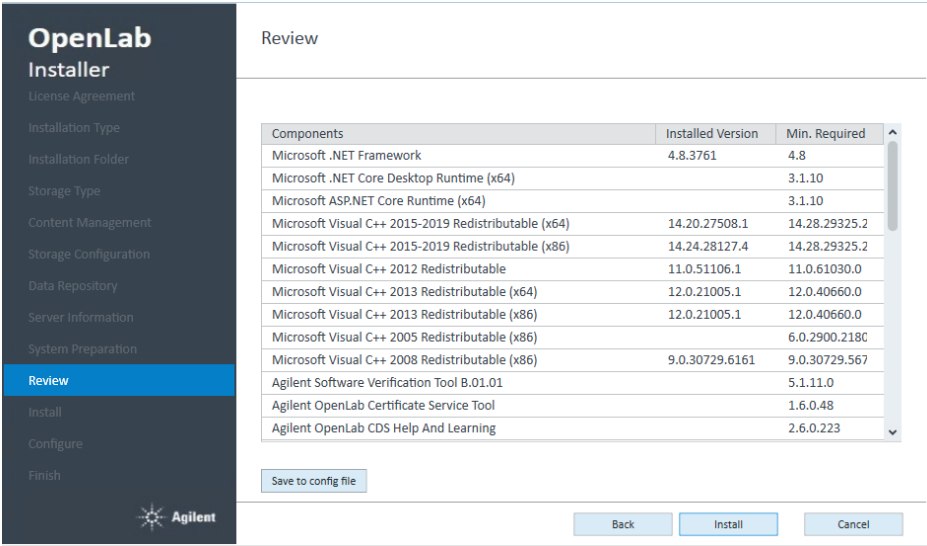
Click **Next** to proceed to the **Review** page. The settings will be applied as part of the installation.

Install OpenLab CDS  
Install OpenLab CDS (Client or Instrument Controller)



**10 Review:** All components (with their version numbers) that will be installed, and all settings that will be applied are listed here. The screenshot shows the review for an OpenLab CDS client.

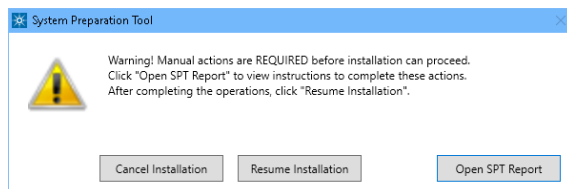
- To save a properties file for a future silent installation (see “[Silent Installation](#)” on page 34), click **Save to config File**.
- To start the installation, click **Install**.



## Install OpenLab CDS

### Install OpenLab CDS (Client or Instrument Controller)

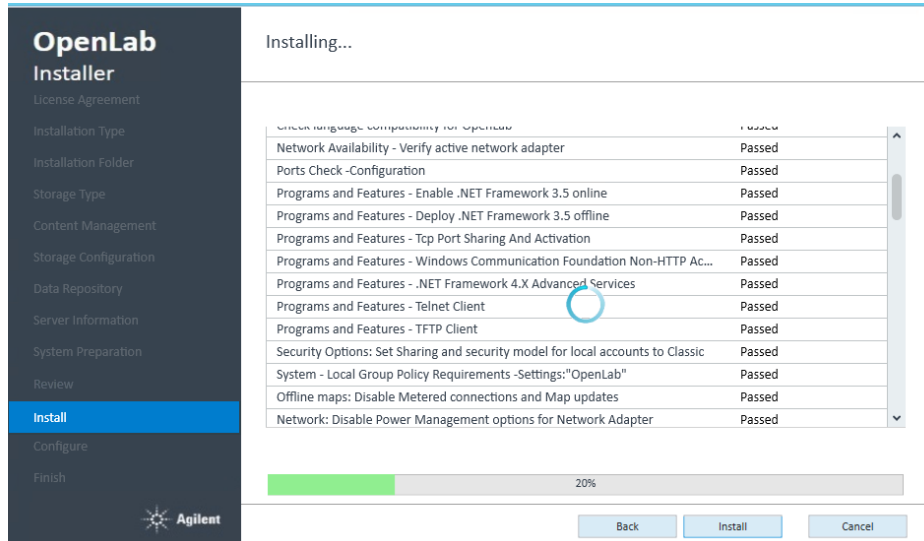
- 11 If there are still manual actions required before installation, a warning is shown.



Options to proceed

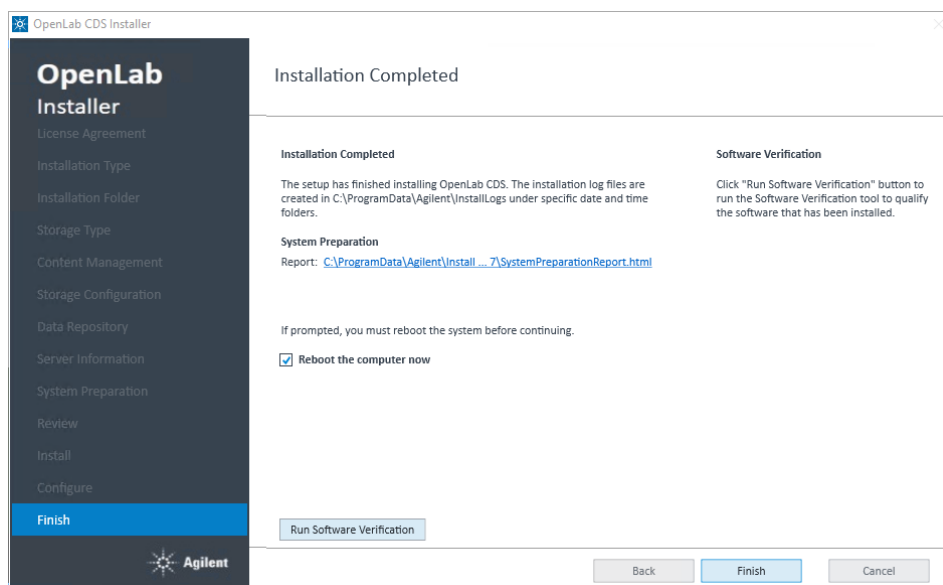
- *Recommended:* Click **Open SPT Report** to view instructions to complete these actions. After completing the operations, click **Resume Installation**.
- **Cancel Installation:** Installation is aborted. Make necessary updates and restart installation.
- **Resume Installation:** The dialog is closed, and installation continues with the installation of the CDS components even if a setting was not applied, or the Operating System does not meet requirements.

- 12 **Install:** After the installation has completed, click **Next**.



**13 Finish:**

- To view the report of the applied Windows settings, click the link under **System Preparation**. If you still need to apply settings manually, a warning will be displayed. You will find details in the linked report.
- To confirm everything has been installed correctly, click **Run Software Verification**<sup>1</sup>.
- To complete the installation, select the **Reboot the computer now** check box, and click **Finish**.



<sup>1</sup> To start the tool separately at a later point in time, select *Start > Agilent Technologies > Software Verification Tool*.

- 14** In case of errors during the installation: Check the installation log files under C:\ProgramData\Agilent\InstallLogs\<date and time>. Note that ProgramData is a hidden folder.

As of OpenLab CDS version 2.5, registration of the CDS application is no longer done at the end of the CDS Client installation. For regulated customers, we recommend to:

- Perform the Operational Qualification (OQ) of the CDS Clients after having installed at least *one Analytical Instrument Controller (AIC)*.
- or
- For configuration with no AIC to run the OpenLab Configuration tool after installing the first CDS Client (see “[Register a Client or Instrument Controller on the Server](#)” on page 106).

The installation includes a set of standard instrument connections. If you need other instrument driver software, install it in a separate step. See “[Install or Upgrade Instrument Driver Software](#)” on page 38.

## Sample Scheduler for OpenLab

As of OpenLab CDS version 2.6, Sample Scheduler for OpenLab is installed by default with an OpenLab CDS Client/Server installation. The following components are available in the Client/Server topology after upgrading to 2.6:

- Server: Sample Scheduler Database as part of the Data Repository, LIMS Agent, Web Server
- Instrument Controller: Sample Scheduler Agent
- Client: Sample Scheduler Client, Sample Scheduler Configuration

Sample Schedule services are disabled after installation. To try out Sample Scheduler for OpenLab select **Agilent Technologies > Agilent Sample Scheduler** from the Start menu. Click **Access free dashboard**, and follow the instructions.

If you decide to regularly use Sample Scheduler or want to evaluate more capabilities please activate it first. For details on activation see information page that is launched.



## Install CDS Client on Terminal Server

- 1 Set up a Terminal Server.
- 2 Install the CDS client as described under “[Install OpenLab CDS \(Client or Instrument Controller\)](#)” on page 24.
- 3 Publish the **Control Panel** RemoteApp program from the Terminal Server setup.
- 4 From a thin client, access the Terminal Server published link

## Silent Installation

OpenLab CDS supports a command-line mode for installation, also referred to as *silent installation*. This mode supports installation, upgrade, and uninstallation. You can execute silent installations either manually or as part of software management systems such as LANDesk or HP CM.

### Export Properties File

The OpenLab Installer supports a feature to export the installation parameters into a properties file which you can then use for the silent installation.

- 1 Launch the OpenLab Installer.
- 2 Follow the instructions of the wizard.
- 3 When you have reached the **Review** screen, click **Save to config file**.  
Save the file to a suitable location. The file will automatically be saved as a .properties file.

You can now use the properties file for the silent installation.

## Run Silent Installation

### Prerequisites

- You have prepared a properties file for silent installation. See [“Export Properties File”](#) on page 34.
- .Net Framework is present on your system.  
If it is not present, it will automatically be installed. You must then manually select **Accept** to agree with the license agreement.

### Install application on client or AIC

- 1 Copy the content of the USB media to a centralized folder.
- 2 Copy the properties file to the same directory as the CDSInstaller.exe.
- 3 Right-click the executable of the command prompt or Power shell prompt, and run it as administrator.
- 4 Navigate to the location where you have saved the installation files.  
For example: C:\CDS
- 5 To start the installation, call CDSInstaller.exe with the following syntax:  
*CDSInstaller.exe -s -c <PropertiesFile>*  
For example:  
*CDSInstaller.exe -s -c Silent.Properties*  
With this command, you start the OpenLab Installer without a user interface.
- 6 Wait about 5 minutes while the installation takes place. To check the process of installation, look at the log files under %ProgramData%\Agilent\InstallLogs.  
If a required installable is missing, the OpenLab Installer will create an entry in a log file, and, depending on the component type, will continue or roll back the installation. An error code will be returned in such scenarios.
- 7 After the installation has finished, reboot the PC.

### Register AIC on the server

- 1 Call registercds.exe with the required parameters.  
For example, register an AIC on a server named *server01* with the user *admin* and password *admin* in the domain *domain01*:  
*registercds.exe -registerapp -regaic -olssserver="server01" -olssuser="admin" -olsspw="admin" -olssdomain="domain01" -log="C:\temp\log.log"*

## Parameters and Return Codes

### Parameters

Use the following parameters when calling CDSInstaller.exe in command line mode:

- `-s`  
Silent mode - no user interface will be shown.
- `-c`  
Configuration file - a properties file contains all parameters.
- `<PropertiesFile>`  
The properties file contains all required inputs for the installer. Replace `<PropertiesFile>` with the correct file path and file name. The file must be located in the same directory as the CDSInstaller.exe.
- `-uninstall`  
Uninstall the currently installed CDS.

Use the following parameters when calling registercds.exe in command line mode:

- `-registerapp`  
Register the OpenLab CDS software on the server. This adds permissions, permission groups, and roles; it enables activity log, creates the RecoveryData folder structure and adds firewall exceptions.
- `-regaic`  
Register the machine as an AIC. This adds the name, hostname, port, and description of the machine to the list of AICs on the server. It also registers the installed drivers, registers the instrument icons, and adds support for multiple instruments.
- `-olssserver="nameOfServer"`  
Connect to the server specified and set default connection.
- `-olssuser="nameOfUser"`  
Username for connecting to the OLSS server specified by the `-olssserver` flag
- `-olsspw="password"`  
Password for connecting to the OLSS server specified by the `-olssserver` flag

- *-olssdomain="nameOfDomain"*  
Domain for the user connecting to the OLSS server specified by the *-olssserver* flag
- *-log="PathAndNameToLogfile"*  
Log the actions performed by the tool

## Return Codes

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

**Table 3**      **Return codes**

Error/return code	Return value
Success. You can see all of the information in the log file.	0
Reboot required. After reboot you need to execute installation command again.	3010 1641 350
Failure. Verify against the log file to see what failed.	any other number

## Logging and Tracing

All exceptions, errors and information messages are logged under C:\ProgramData\Agilent\InstallLogs\<date and time>. Note that ProgramData is a hidden folder.

## Install or Upgrade Instrument Driver Software

The following driver software packages are automatically installed and configured with OpenLab CDS. For details, see chapter *Supported Instruments* in the *OpenLab CDS Requirements and Supported Instruments for Client/Server Systems* guide (CDS\_ClientServer\_Requirements.pdf).

- Agilent GC & GC/MS
- Agilent LC & LC/MS
- Agilent 35900 A/D
- Agilent SS420x
- Agilent Data Player (Virtual Instruments)

Other instrument driver software, add-on software, or upgrades to existing driver software, must be installed and configured manually. Instrument driver software and add-on software can be found on the installation media under Setup\Packages\Add-Ons.

The latest Agilent drivers are available in SubscribeNet. In the Product List, select **OpenLab Software > OpenLab Agilent Instrument Drivers**.

### On the Analytical Instrument Controller (AIC)

- 1 Run the driver installer package, and follow the installation wizard.  
For details on the installation or upgrade procedures, refer to the respective driver documentation.
- 2 Register the new drivers on the server. The procedure is identical to registering an existing OpenLab PC as a Client or AIC on the server. See [“Register a Client or Instrument Controller on the Server”](#) on page 106.  
New instrument types are now available in the Control Panel.

#### NOTE

Both steps can alternatively be done via silent installation. For details, see [“Run Silent Installation”](#) on page 35.

### On the relevant clients

- 1 Run the installer package on each client from which you will access the instrument.



## 3 Post Installation Tasks

Run a Software Verification after Software Installation	40
Create Account for Network Access and Printing	41
Set up the domain user account	41
Enable automatic printing	43
Configure the Antivirus Program	44

This chapter describes tasks that are relevant after finishing the installation.

## Run a Software Verification after Software Installation

The Software Verification Tool (SVT) provides documentary evidence that your system has been built and installed correctly, and that all design specifications have been met. You do not need to run the software verification again if it has run successfully at the end of the installation.

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

**2** Select the components to qualify.

**3** Select **Qualify**.

The system will run the application and generate a Software Verification Report. Reports are automatically saved to C:\SVReports.

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

Do not use the system until the Software Verification Report gives a 'pass' result.



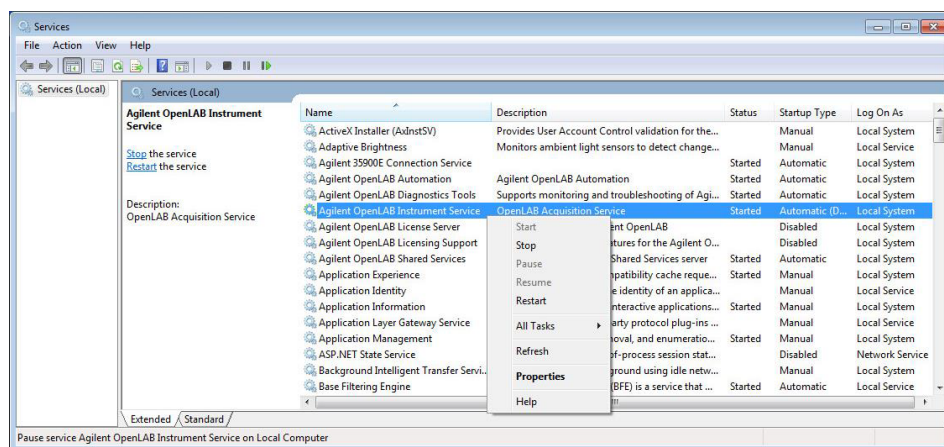
## Create Account for Network Access and Printing

OpenLab CDS allows you to automatically process your data during acquisition, without opening Data Analysis. During this processing specific domain user privileges to access a network share/network printer and log on as a service are required:

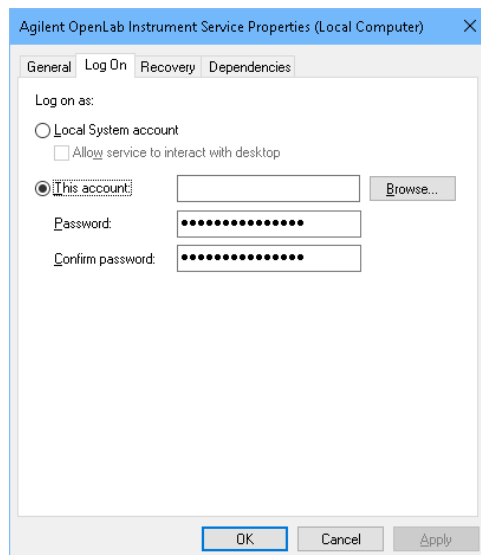
- You can generate reports and print them to printer or save them as files. Exporting the reports to a network share is a typical way how they are sent to an external system (for example, LIMS).
- You can export raw data or results during a run, and save the files to a network share.

## Set up the domain user account

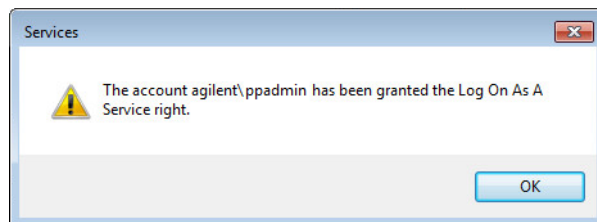
- 1 Log in as Windows domain user who has local administrative privileges.
- 2 Go to **Control Panel > All Control Panel Items > Administrative Tools** and double-click **Services**. Make sure the status of **Agilent OpenLAB Instrument Service** is **Started** or **Running**. If not, right-click and select **Start**.
- 3 Right-click **Agilent OpenLAB Instrument Service** and select **Properties**.



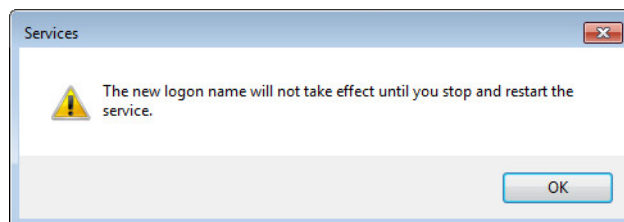
- 4 Click the **Log On** tab, select **This account** and enter the login credentials of a domain user who has network printing privileges. Click **OK**.



- 5 The **Services** window confirms that the account has been granted the **Log On As A Service** right. Click **OK**.



- 6 Click **OK** in the Services window to acknowledge that **The new logon name will not take effect until you stop and restart the service.**



- 7 Restart the PC. This will enable the *Log On* user specified in Instrument Service to automatically print reports in Data Acquisition.
- 8 After the PC reboots, log in using the same domain user account (for example, `agilent\ppadmin`) specified in the **Log On** tab of **Agilent OpenLab Instrument Service Properties** window.

## Enable automatic printing

If you plan to acquire a single sample or a sequence that specifies a processing method that has **Printer** report destination, add a default printer to the PC to enable automatic printing of reports.

### Prerequisites

You have set up a domain user account (see ["Set up the domain user account"](#) on page 41).

- 1 Go to **Control Panel > All Control Panel Items > Devices and Printers** and click **Add a printer**.
- 2 Select **Add a network, wireless or Bluetooth printer** and browse or type a shared printer name that you want to add. Once the printer is added, print a test page and set it as a default printer.

## Configure the Antivirus Program

- 1 Be sure to open the firewall ports listed in the Firewall Settings in the *OpenLab CDS Requirements and Supported Instruments (Client/Server)* guide (CDS\_ClientServer\_Requirements.pdf).
- 2 The following folders should be excluded from Antivirus scan. If you want to have these folders scanned, you should do this while the system is not acquiring or doing data Analysis as scanning may cause slowness and runs to be aborted due to concurrent access to the same file by the AV program and the CDS Application.
  - [C:\]DsData\DsArchive
  - [C:\]DsData\DsContent
  - [C:\]DsData\DsIndex
  - [C:\]Program Files (x86)\Agilent Technologies
  - [C:\]ProgramData\Agilent
  - [C:\]ProgramData\Agilent IPB Files
  - [C:\]ProgramData\Agilent Technologies
  - [C:\]ProgramData\ChromatographySystem
  - [C:\]ProgramData\Firebird
  - [C:\]ProgramData\IsolatedStorage

Refer to your specific antivirus software documentation on how to configure folder exclusions.

### Settings for Trend Micro™ antivirus software

OpenLab CDS can be used with other antivirus programs as well. If you use Trend Micro™, the following settings are recommended to optimize system performance.

- 1 If your version of Trend Micro has **Web Reputation**: Turn off to maximize performance.

The risk of turning off Web Reputation is that web traffic through browsing from the machine will not be checked.

Ensure that there is another URL/web scanner on the gateway level to protect the endpoint, or ensure that the endpoints have limited access to Internet. These production machines should not have access to Internet websites where most of the infections are coming from.

- 2 Real time scan:** Add exclusions, and modify scan direction from **Created/Modified/Retrieved** to **Created/Modified**.

Exclusions ensure that the working directory of Agilent Technologies will not be scanned, thus improving performance.

The risk is that only files that are created and changed on this machine are scanned. Files that are just accessed will be bypassed. Dormant Files that got infected without being noticed at the time they were created or written to the machine will not be scanned.

Increase scheduled scan to daily to ensure all files on the machine are being checked for infections that are dormant or not moving.

- 3 Behavior Monitoring:** Add below list of programs to **Approved programs**.

- OpenLab\Services\Distributed Transaction Coordinator Service\  
Agilent.OpenLab.DistributedTransactionCoordinator.Rest.exe
- OpenLab Backup Utility\Monitoring Service\  
Agilent.OpenLab.BackupRestore.BackupMonitoringService.exe
- OpenLab Backup Utility\Notification Service\  
Agilent.OpenLab.BackupRestore.NotificationService.exe
- OpenLab Backup Utility\Task Status Cache Service\  
Agilent.OpenLab.BackupRestore.TaskStatusCacheService.exe
- OpenLab\Services\Electronic Signature Service\  
Agilent.OpenLab.ESignature.Rest.exe
- OpenLab Acquisition\  
Agilent.OpenLab.Acquisition.AcqInstrumentService.exe
- OpenLab Acquisition\Agilent.OpenLab.AcquisitionClient.exe
- OpenLab Data Analysis\Bin\  
Agilent.Chromatography.DataAnalysis.Processing.ProcessingServer.exe
- OpenLab Data Analysis\Bin\  
Agilent.Chromatography.DataAnalysis.UI.CustomCalculationDesigner.exe
- OpenLab Data Analysis\Bin\Agilent.OpenLab.DataAnalysis.exe
- OpenLab Data Analysis\Bin\  
Agilent.OpenLab.DataAnalysis.Api.ApplicationService.exe
- OpenLab Data Analysis\Bin\Reporting\  
Agilent.OpenLab.Reporting.RdlDescriptor.exe
- OpenLab Data Analysis\Bin\Reporting\  
Agilent.OpenLab.Reporting.RdlDescriptorContextMenu.exe
- OpenLab Data Analysis\Bin\Reporting\  
IntelligentReporting.RenderServiceHost.exe

- OpenLab Data Analysis\Bin\Reporting\TemplateDocumentation.exe
- OpenLab Platform\Data Repository\Data Repository\Base\BaseService\Agilent.OpenLab.DR.BaseService.exe
- OpenLab Reverse Proxy Configuration Service\ConfigurationService\Agilent.OpenLab.ReverseProxy.ConfigurationService.exe
- OpenLab Services\Distributed Transaction Coordinator Service\RegistrationTool\Agilent.OpenLab.DataRepository.RegistrationTool.exe
- OpenLab Services\Automation\AutomationServerHost.exe
- OpenLab Services\Diagnostics\DiagnosticsToolsServiceHost.exe
- OpenLab Services\Licensing\Flexera\ladmin.exe
- OpenLab Services\Licensing\Licensing.Service.Host.exe
- OpenLab Services\Server\SharedServicesHost.exe
- OpenLab Services\UI\Agilent.OpenLab.ControlPanel.exe
- Test Services\Agilent.TestServices.WebService.exe
- Test Services\Central Management Service\Agilent.TestServices.Server.Main.exe

C:\Program Files (x86)\Agilent Technologies\...

The risk is that if any of the excluded files get infected, it will not be detected. For example, trigger a schedule on a daily basis to cover these files.

- 4 Realtime monitoring:** Add below folder to the exclusion list of Realtime Monitoring setting:

C:\Program Files (x86)\Agilent Technologies\



## 4 Optional Procedures

Install OpenLab Help and Learning Only 48

Improve Performance on Offline Machines 49

This chapter describes the installation or upgrade of additional software. It also contains information on the installation of OpenLab Help and Learning only, and on performance improvement on offline machines.

## Install OpenLab Help and Learning Only

Use this option to install OpenLab Help and Learning content without installing OpenLab CDS applications.

Do not use this option on a machine where OpenLab CDS is, or will be, installed.

- 1 Insert the USB media, right-click the setup.exe file, and run it as administrator.
- 2 On the start screen, select **OpenLab CDS**, and click **OK**.
- 3 In the OpenLab CDS Installer, click **Documentation**.
- 4 Click **Install OpenLab Help and Learning Only**.
- 5 Select your language, and click **Next**.
- 6 Accept the terms on the **License agreement** page, and click **Next**.
- 7 On the welcome screen, click **Next**.
- 8 Review the installation directory. If desired, click **Change...** to specify a different directory.
- 9 Click **Install**.
- 10 When the installation is complete, click **Finish**.
- 11 If you plan to use Internet Explorer to view the content, make the following settings to ensure that OpenLab Help and Learning is opened without showing a confirmation prompt. No configuration is required for Google Chrome or Edge.
  - a In Internet Explorer, click **Tools > Internet Options**.
  - b Select the **Advanced** tab.
  - c Under **Security**, select **Allow active content to run in files on My Computer**.
  - d Confirm your settings.
  - e Reboot the computer to make the settings effective.

Without these settings, you will need to click **Allow blocked content** when opening the help.

You can uninstall or repair OpenLab Help and Learning from the same link in the installer (see “[Uninstall OpenLab Help and Learning Only](#)” on page 112).



## Improve Performance on Offline Machines

Computers running OpenLab CDS may exhibit slow performance when they are not connected to the Internet.

Use the following system settings on all workstations, clients, AICs, and servers to remedy this problem.

- 1 Open Internet Explorer and select **Tools > Internet Options**. In the **Advanced** tab, clear the following check boxes:
  - **Security > Check for publisher's certificate revocation**
  - **Security > Check for server certificate revocation**
- 2 Change or add the following registry keys:
  - [HKEY\_LOCAL\_MACHINE\SOFTWARE\Policies\Microsoft\SystemCertificates\AuthRoot]  
"DisableRootAutoUpdate"=dword:00000001
  - [HKEY\_LOCAL\_MACHINE\SOFTWARE\Wow6432Node\Policies\Microsoft\SystemCertificates\AuthRoot]  
"DisableRootAutoUpdate"=dword:00000001
- 3 Document that you turned off the Root Certificates, as this can prevent users from installing other applications.

# 5

## Licensing

About OpenLab CDS Licensing 51

License Types 51

License File 51

Get a License 52

Obtain a License with SubscribeNet 52

Other Ways to Obtain a License 55

Install Your License 57

This chapter provides basic information on OpenLab licensing and describes how you generate a license file with SubscribeNet.

## About OpenLab CDS Licensing

### License Types

The license file is a collection of Product, Instruments and Add-on licenses (or activation keys), and is installed to your OpenLab CDS System.

The licenses or activation keys in the license file can either be Shared or Counted:

- Shared licenses — system computers and other components can have shared, or add-on, licenses — because they share a core license.
- Counted licenses — these licenses are part of the OpenLab CDS floating licensing strategy. They are not permanently assigned to any one component. Instead they are automatically assigned to components, such as AICs and instruments, while the components are starting up. The licenses are automatically returned when the component is closed. The license management program controls license issuance and retrieval.

In this case, the only requirement is that a component is licensed while running. You only need enough licenses for all components running concurrently, rather than for each installed component.

A startup license for the system allows you to run OpenLab CDS for 60 days after the installation. In order to run the data system software after the 60-day period, you must install your license file.

### License File

A license file will contain your software license. This file is installed to the license server, that is, the workstation computer, or the server to which your product was installed in a client/server system. The license file is bound to this server address, and cannot be moved to another server.

Information in the license file defines the number of instruments and other options that may be used concurrently with your system.

The most efficient way to manage and maintain your licensing is through the Internet.

If you have not received a lavender envelope for your product, contact your vendor or internal support.

## Get a License

### Obtain a License with SubscribeNet

Use the following procedure to generate and download your license. In case you do not have internet access, skip to the section [“Other Ways to Obtain a License”](#) on page 55.

If you are a new user who has not registered with SubscribeNet, continue with the section *New Users*.

If you have registered with SubscribeNet, skip to the section *Users registered with SubscribeNet*.

#### Prerequisites

To generate, download, and install a final license for your product, you will need:

- The authorization code label provided in the lavender envelope containing your Software Entitlement Certificate.

If you have not received a lavender envelope for your product, contact your vendor or internal support.

- The URL for SubscribeNet from the Software Entitlement Certificate.
- The host name of the computer where the Control Panel is running.
- The MAC address.

To retrieve your MAC address from a computer where OpenLab CDS is already installed, open the Control Panel and browse to the **Administration > Licenses** section. Use the **Copy MAC Address** or **Save MAC Address** function to obtain the MAC address for license generation.

During this process you will have to enter the MAC address of your license server. For workstations, this is the local computer. For client/server systems, this is the server.

#### NOTE

If any changes are made to the computer name or domain reference after the license is installed, remove the license. A new license will need to be created in SubscribeNet, downloaded, and installed.

#### NOTE

If the network adapter that provides the MAC address used during license creation is removed from the machine, your license will no longer be valid. A new license will need to be generated with a currently available MAC on the license server.

## New Users

- 1 Go to <https://agilent.subscribe.net/control/agil/AgilRegisterToAccount> to register the product with SubscribeNet.
- 2 On the registration page, enter the authorization code from the label and complete the profile information (required fields are marked with an asterisk \*).  
The email address you enter will become your login ID.
- 3 Click **Submit**. The system will generate and display an account name for you. SubscribeNet will send a welcome email with your login ID and password.
- 4 Log in to SubscribeNet using your login ID and password.  
Once you log in, you can use the online user manual link for help with any questions you have.
- 5 Select **Generate or View licenses** from the left navigation bar.
- 6 Follow the prompts to generate your new license.  
You will be prompted for the HOST NAME of the computer. The host name you enter must match with the network name of the computer where the Control Panel is running. Do not include any DNS suffix (*domain.com*) references in the entered machine name.
- 7 When the system generates the license, view its details, then click **Download License File**. Save the license file to your computer and to a backup location (such as a portable storage device).  
Use your login ID and password when you revisit the Agilent SubscribeNet site to regenerate a license file, add new authorization codes, or further configure the license for your system.

**Users registered with SubscribeNet**

- 1 If you already have a SubscribeNet account, use <https://agilent.subscribenet.com/>.  
Lost your SubscribeNet password? Use <https://agilent.subscribenet.com/control/agil/password> to have it emailed to you.
- 2 Select the SubscribeNet account associated with this authorization code, if you have more than one account.
- 3 From the SubscribeNet navigation pane, select **Register Authorization Code**.  
This will allow you to enter your new authorization code and make available the new license entitlements.
- 4 Follow steps 5 through 7 in the previous procedure, *New Users*, to *generate or view* your new licenses.

## Other Ways to Obtain a License

If you are unable to generate a license, contact your nearest Agilent technical support office. A representative will tell you how to submit an OpenLab CDS License Generation Form in your location.

### Offline Licensing

If an internet connection is not available in your laboratory:

You or your local on-site service engineer will collect the necessary information from you to allow Agilent to create a license account on your behalf. For phone support in your region, call the sales and service number for your region. See the Appendix for contact information.

#### *Required Customer Information for Agilent License Support:*

The following information must be provided to Agilent in order to enable us to create a licensing account on your behalf.

##### **1** Collect Account Information:

Your account name will be your company name and Lab name separated by a comma. Employee information provided here will be used to define the first administrator of your account for future access to the system as required. Please prepare the following pieces of information prior to contacting your local Agilent sales and service center in order to expedite service:

- Company Name
- Lab/Department Name
- First Name
- Last Name
- E-mail address
- Job Title
- Phone #
- Address, City, State/Province, Postal Code, Country

##### **2** Collect Authorization Code(s):

The authorization code is an alpha-numeric code provided on a label which is enclosed in a lavender envelope. If you have received more than one code you must provide all codes to ensure that all ordered licenses are granted to your account.


**3** Receiving your license:

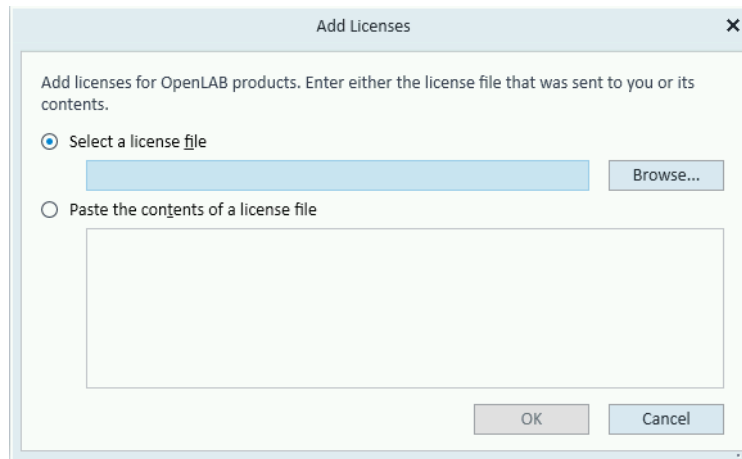
Once the above information is provided Agilent will then work on your behalf to generate a license file through SubscribeNet. The license file will either be sent to your shipping address (on a CD), or your local FSE will deliver it in person (usually on USB media). Once your license is received follow the below section on "Install your License" to finish installing your license on your CDS system(s).



## Install Your License

The license must be added to your system using the Control Panel.

- 1 Start the Control Panel from any machine connected to the system you want to install the license for.
- 2 Navigate to **Administration > Licenses**.
- 3 In the ribbon, click **Add License** .



- 4 Choose to install the license by:
  - Using the license file option to browse to and open the license file (.lic) saved from the license generation process in SubscribeNet.
  - Selecting the License Text option and copying the license text from a text file received into the provided field.
- 5 Click **OK**.

The **Administration** interface in the Control Panel will now display the status of installed licenses.

### NOTE

A full restart is required in order for any license to have an immediate effect.

## 6

# Configure OpenLab CDS

Configure Authentication	59
Configure Security Policy	60
Configure users and roles	61
Create or import users	61
Groups	62
Roles and Privileges	63
Add users or groups to a role	64
Specific Roles for Individual Instruments or Projects	65
Configure Initial Project	66
Configure Initial Instrument	66
Other settings in the Control Panel	67

This chapter describes the initial configuration steps after installing the software. All configuration tasks are performed in the Control Panel. For more details, refer to the Control Panel section in OpenLab Help & Learning.

## Configure Authentication

OpenLab CDS supports the following authentication providers:


- **Internal**

In this mode, the user's credentials are stored in the OpenLab CDS system. You are asked to create an administrator account for OpenLab CDS before setting up other users. This is the only mode in which you can create new users within the system; in all other modes you can only map to users that exist in a different system.

- **Windows Domain**

You import existing Windows users into OpenLab CDS system. The authentication is done either by Windows Active Directory domain or NT 4.0 Domain within the Enterprise. OpenLab CDS only uses the identity and password of the mapped users; roles and privileges for OpenLab CDS are still configured in the Control Panel.

Prior to Client or AIC installation, authentication has been configured on your OpenLab Server. If it is internal authentication, and you are fine with it, nothing needs to be done.

- 1 Launch the Control Panel. Log in with the internal administrator's credentials.
- 2 Navigate to **Administration**.
- 3 In the navigation pane, select **System Configuration**.
- 4 In the ribbon, click **Edit System Settings** .
- 5 Select the authentication provider **Windows Domain** from the drop-down list, then click **Next**.

### NOTE

Do not change the storage type.

- 6 Select the check box to use a domain user, and provide user credentials with the rights to obtain user and group information. Then click **Select Account** to open the **Search Users** dialog and select an administrator account.
- 7 Confirm your settings. When complete, the Control Panel will restart.

## Configure Security Policy

If you need to comply with specific standards (for example, 21 CFR Part 11), adjust the security policy as required.

With the authentication provider **Internal**, you can set all parameters in the Control Panel.

With an external authentication provider (Windows Domain), you can set the inactivity time or enable single sign-on; all other parameters are defined by the external system.

- 1 Launch the Control Panel and navigate to **Administration**.
- 2 In the navigation pane, select **Security Policy**.
- 3 In the ribbon, click **Edit Security Policy**.
- 4 Set the parameters as required, and confirm your changes.  
Restart the Control Panel to apply your changes.

### NOTE

To meet 21 CFR Part 11 requirements, set the **Password expiration period** to 180 days or less. Do not change the other default values, they comply with 21 CFR Part 11.

## Configure users and roles

With internal authentication, you create the required users in the Control Panel. With Windows domain as an external authentication system, you import the Windows domain users.

To define what users are allowed to view or do, OpenLab CDS offers predefined roles and allows you to define your own specific roles.

Roles are equipped with numerous specific privileges. When you set up a customized role, it is therefore recommended starting from a lower role (that is, with fewer privileges) and adding specific required privileges, rather than removing privileges from a higher role.

Each user can be member of multiple groups. You must assign one or more specific roles to each group. You can also assign roles to single users; however, for the sake of clarity, it is strongly recommended to assign roles only on the group level. Every member of a group automatically has all roles of this group.

- 1 Launch the Control Panel and navigate to **Administration**.
- 2 In the navigation pane, select **Users, Groups**, or **Roles**.
- 3 Create new items, or edit the existing ones.

## Create or import users

### Add users (Internal Authentication only)

- 1 From the navigation pane, click **Administration > Users**.
- 2 In the ribbon, click **Create User**.
- 3 In the **Create User** dialog, provide the relevant parameters:
  - Enter the name and password for the new user.
  - By default, the new user will need to change the password at next logon. If this is not required, clear the **User must change password at next logon** check box.
  - In the **Role Membership** tab, assign the user to an appropriate role. You can use the default roles, or prepare your own roles in the Control Panel under **Administration > Roles**.
- 4 Click **OK**.

### Import users (Windows Domain Authentication only)

To add users to your system, you must have privileges to obtain user and group information from the domain.

- 1 From the navigation pane, click **Administration > Users**.
- 2 In the ribbon, click **Import User**.
- 3 In the **Search Users** dialog box, enter search string for the Windows domain username.
- 4 From the **Search Results** list, select the user you want to import, and click **Add**. The user is added to the **Selected Users** list.
- 5 Repeat steps 2 to 4 until you have added all the user names that you want to import to the **Selected Users** list, then click **OK**.

## Groups

Each user can be member of multiple groups. You must assign a specific role to each group. You can also assign roles to single users; however, for the sake of clarity, it is strongly recommended to assign roles only on the group level.

If you use an external authentication provider, you can either import the names of groups that exist in the external system, or create new internal groups. There is no limit on the number of groups that can be mapped or created.

Assign users to groups either in the external system or in the Control Panel. If you need additional user assignments that are relevant only for OpenLab CDS, create them in the Control Panel.

If you delete or unmap a group, the users who were members in this group remain unchanged.

## Roles and Privileges

Roles are used to assign privileges to a user or a user group globally or for a specific instrument, project, project group or location. The system contains a list of predefined roles which are installed as part of the system installation (for example, **Instrument Administrator**, **Instrument User**, or **Everything**). Each role has certain privileges assigned.

Privileges are grouped according to the three main role types (Project role, Instrument role, and Administrative role). When you assign privileges to a role, you first select the required role type and then select the privileges related to this role type. Each role can only have privileges of one specific role type; the only exception is the predefined role **Everything**, which has all privileges of all role types. Users or groups may require multiple roles to perform system functions. For example, a user with the role *Chemist* may need another role such as *Instrument User* with the privilege to run an instrument.

You can create a tree of different locations in the Control Panel, and add instruments to the relevant locations. For each instrument or instrument group, you can assign different instrument roles (see also “[Specific Roles for Individual Instruments or Projects](#)” on page 65). For example, a user can have the role **Instrument Administrator** for one instrument, and **Instrument User** for another instrument.

You can also create a tree of different projects or project groups in the Control Panel, and assign different Project roles for different projects (see also “[Specific Roles for Individual Instruments or Projects](#)” on page 65). For example, a user can have the role **Project Administrator** in one project, so that he can manage the settings in the Control Panel. In a second project, he may have a role that allows him to edit the content of a project, but not to change the project settings.

For more information on privileges, see the Appendix.

**Table 4** Description of role types

Role Type	Description
Administrative privileges	These privileges are globally assigned to a user or group and cannot be changed on the instrument/location level. They are the typical administration privileges such as <b>Backup and restore</b> , <b>Manage security</b> , <b>Manage printers</b> , etc.

Table 4 Description of role types

Role Type	Description
Instrument privileges	These privileges can be assigned globally or on the instrument/location level. Privileges for instruments are, for example, <b>View instrument or location</b> and <b>Run instrument</b> . Users need the <b>View instrument or location</b> privilege on the global level to see the locations and instruments tree in the Control Panel.
Project privileges	Privileges for accessing or modifying different levels of data. You can assign these privileges globally or on project or project group level.

## Add users or groups to a role

- 1 From the navigation pane, click **Administration > Roles**.
- 2 In the **Roles** window, select the role you want to assign to users or groups.
- 3 In the ribbon, click **Edit Role**.
- 4 In the **Edit Role** dialog box, click the **Members** tab.
- 5 Click **Add user or group**.
- 6 In the **Search Users and Groups** dialog box, enter the name of a user or group, and click **Search** to view a list of all users and groups that meet the search criteria.
- 7 Under **Search Results**, select a user or group, and click **Add**.
- 8 Click **OK**.



## Specific Roles for Individual Instruments or Projects

By default, the roles of users or groups are globally set for all locations, instruments, project groups, or projects. The role settings are inherited from the root node **Instruments** or **Projects**, respectively. In order to assign a different role to a user or group for one specific project:

- 1 Select the relevant project.
- 2 In the ribbon, click **Edit Privileges**.
- 3 In the **Edit Privileges** dialog, clear the **Inherit privileges from parent** check box.

Afterwards, you can assign a different role that will be valid only for the specific node.

You can assign **Instrument** roles to individual locations or instruments.

If you use projects, you can assign **Project** roles to individual project groups or projects.

**Administrative** roles are always set globally.

## Configure Initial Project

- 1 Launch the Control Panel and navigate to **Projects**.
- 2 Click **Create** in the ribbon to create a new project.  
On the **CDS Settings** tab:
  - Enter the locations for Methods, Sequences, Results, Sequence Templates and Report Templates.
  - Consider the required audit trail settings for this project.

For more details, refer to the Control Panel section in OpenLab Help & Learning.

## Configure Initial Instrument

- 1 Launch the Control Panel and navigate to **Instruments**.
- 2 Click **Create** in the ribbon to create a new instrument.
- 3 Select the new instrument, and click **Configure Instrument** in the ribbon.
- 4 It is recommended that you use Auto Configuration (if available) to configure your instruments: Click **Auto Configuration**, and provide the instrument's IP address or hostname.

For more details, refer to the Control Panel section in OpenLab Help & Learning.

## Other settings in the Control Panel

Consider also other settings in the Control Panel, such as:

- Changing the instrument status reporting frequency
- Changing audit trail settings for a project
- Editing signature levels for a project

For more details, refer to the Control Panel section in OpenLab Help & Learning.



## 7

# Set Up NIST Library Search

Install Software and Library Components 69

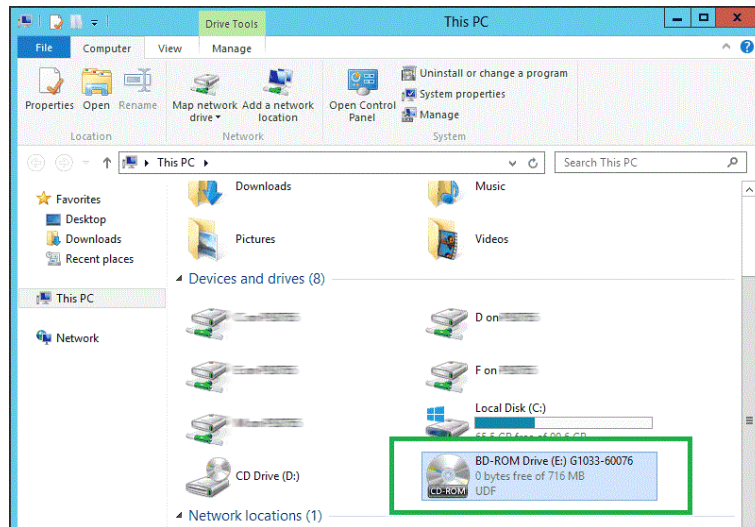
Configure MS Library in OpenLab CDS 79

This chapter describes the installation of NIST software components, and the configuration steps in OpenLab CDS that are required for using the NIST MS library search.

# Install Software and Library Components

Follow these settings to ensure that all the same NIST libraries are used in every session, even if there are two or more Citrix or Terminal Servers.

- 1 Mount the ISO image using a tool such as Daemon.

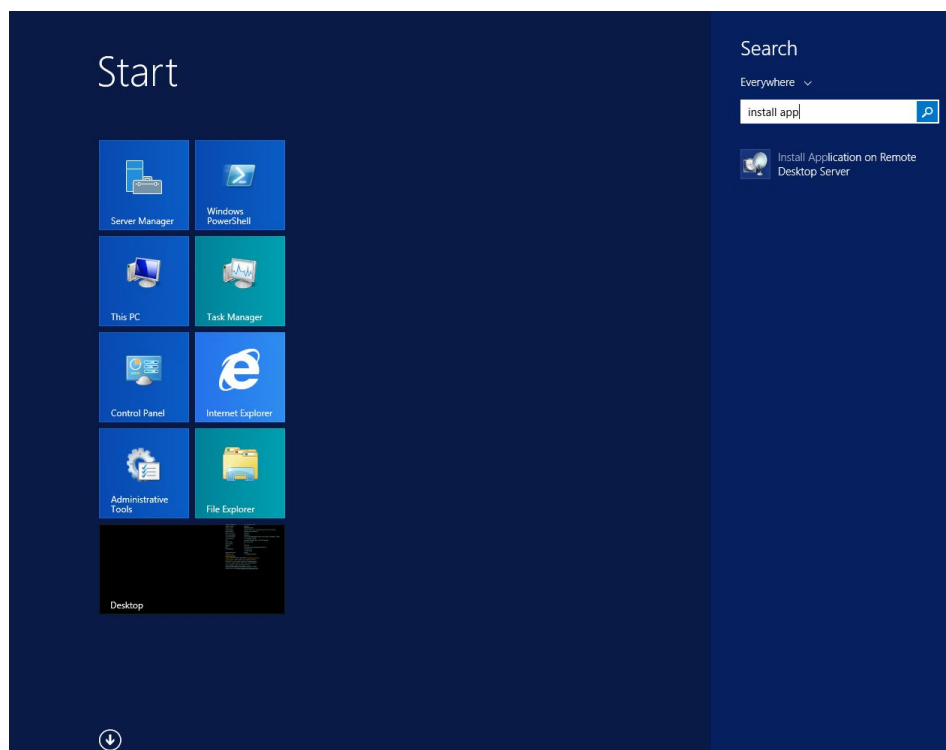


**Figure 2** Example: ISO image mounted as drive E:

## Set Up NIST Library Search

### Install Software and Library Components

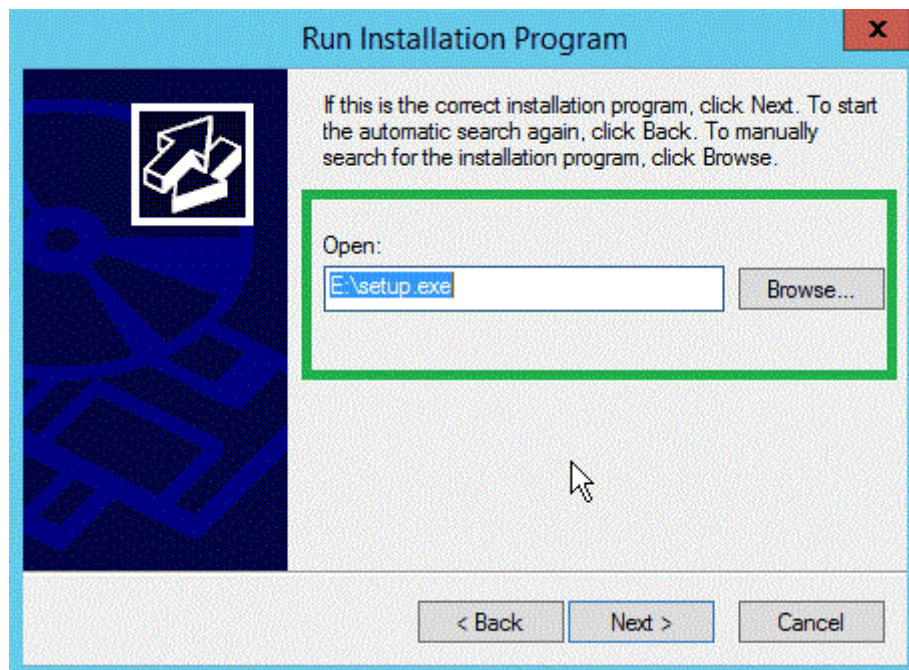
- 2 Start the installation via **Install Application on Remote Desktop...**



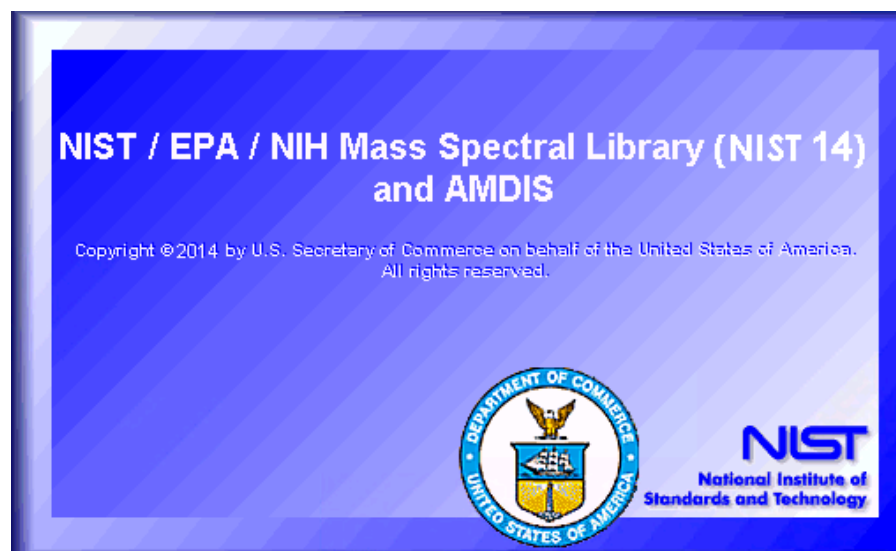
## Set Up NIST Library Search

### Install Software and Library Components

- 3 Run the installation program from the mounted ISO image.



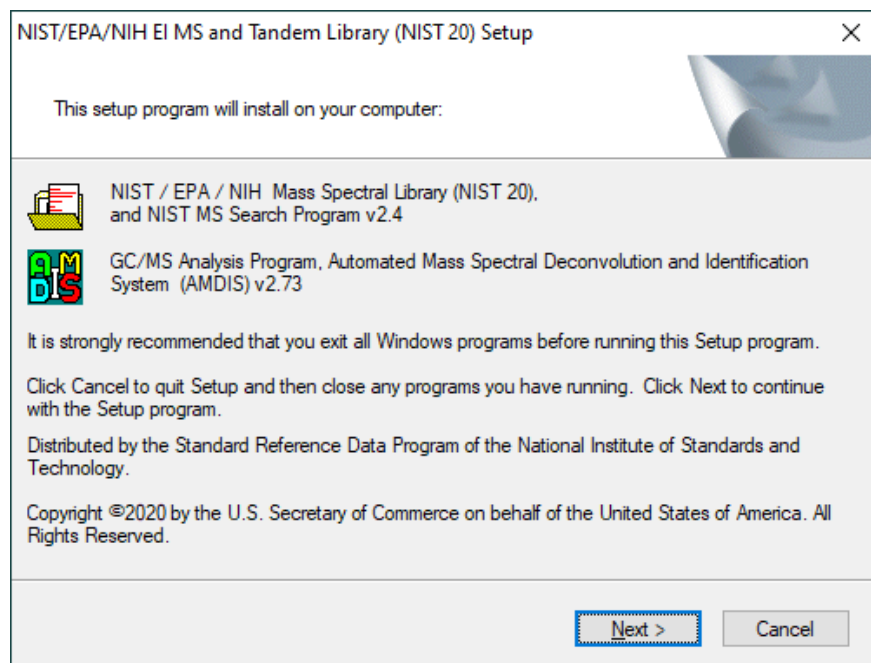
The NIST installation wizard opens.



## Set Up NIST Library Search

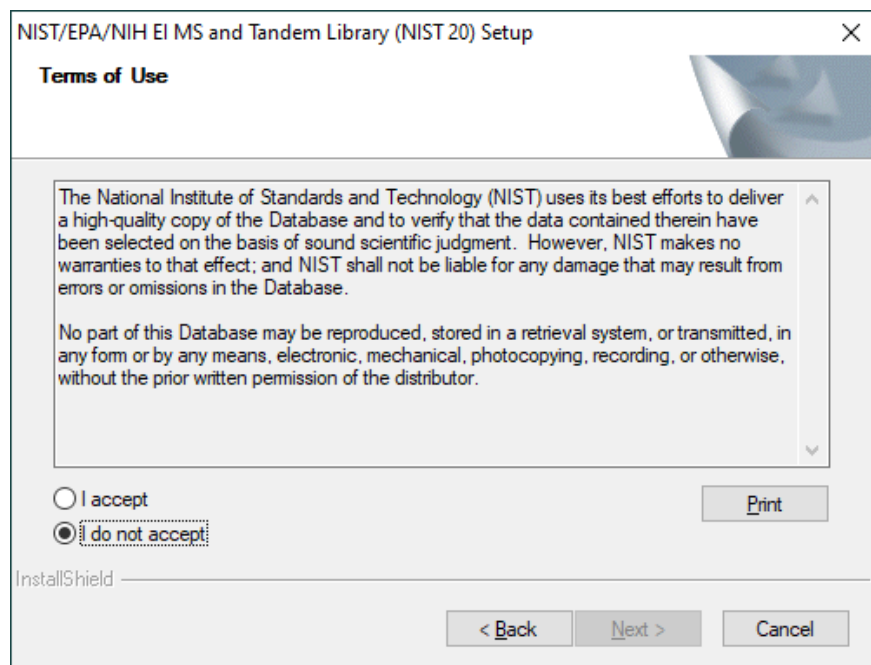
### Install Software and Library Components

#### 4 Click **Next**.

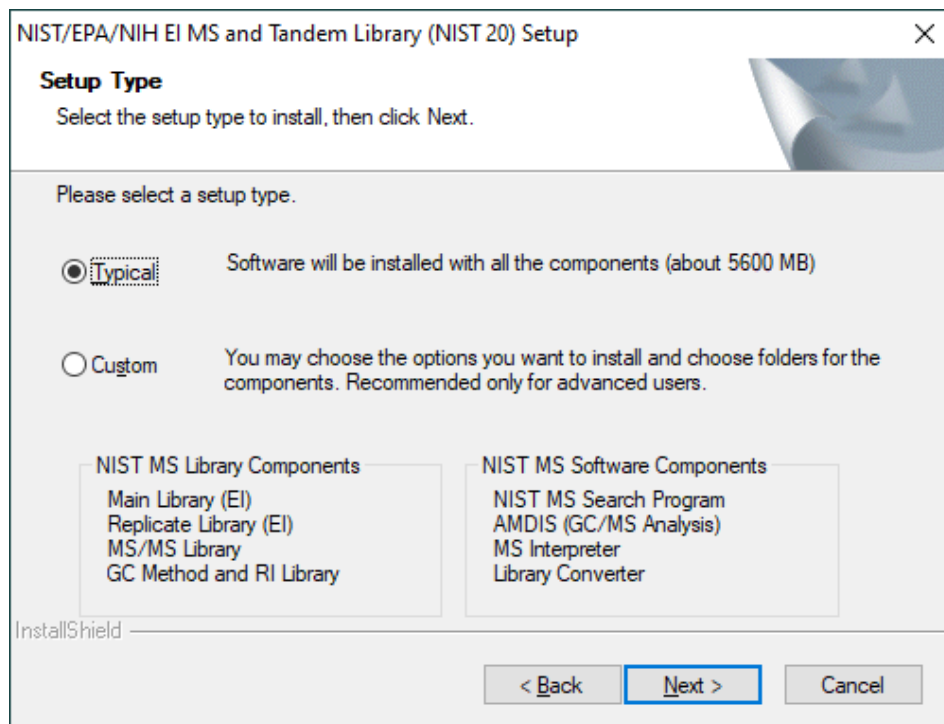




- 5 Accept the license.



- 6 Choose the setup type **Custom**.

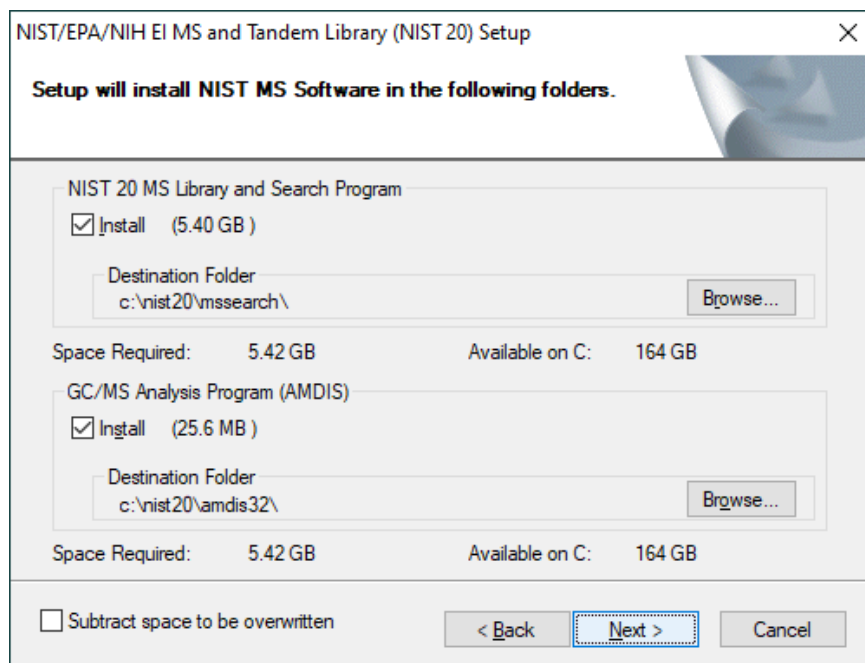


This will install the software with all the components of the NIST software and NIST libraries.

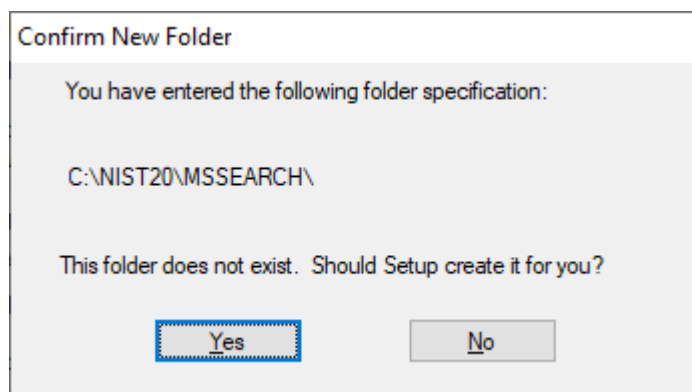
## Set Up NIST Library Search

### Install Software and Library Components

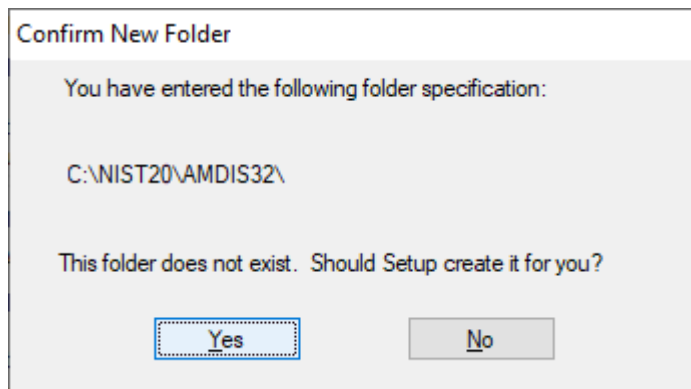
- 7 Specify the folder to install the NIST MS Software (by default it is your local C drive). Make sure the check box has been selected. Click **Next** to continue.



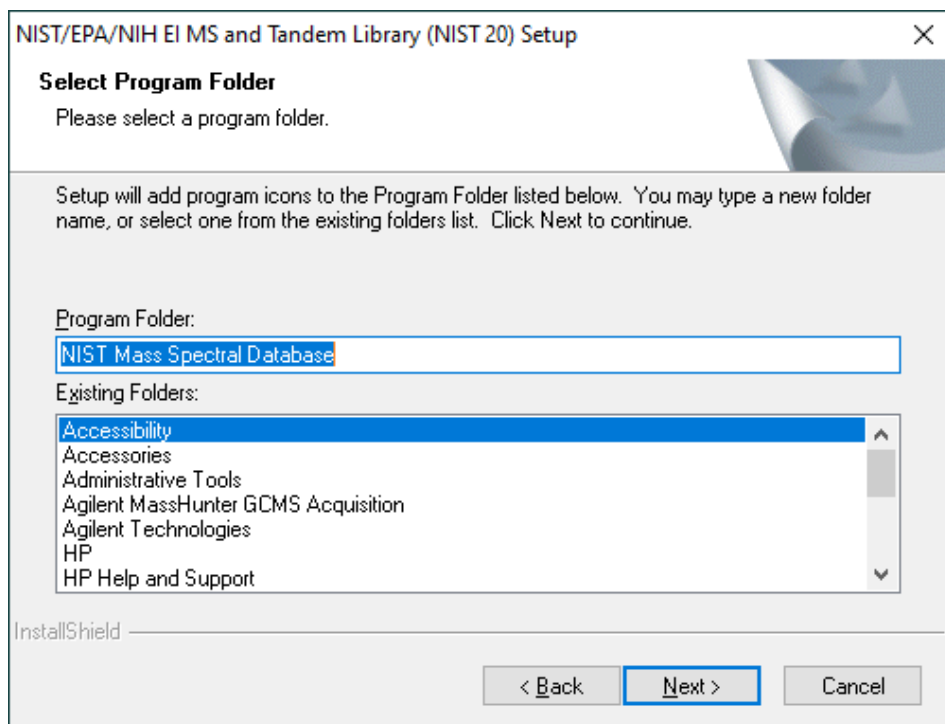
- 8 Click **Yes** to confirm the creation of the new folder for the NIST MS Search Program.



- 9 Click **Yes** to confirm the creation of the new folder for the GC/MS Analysis Program (AMDIS).



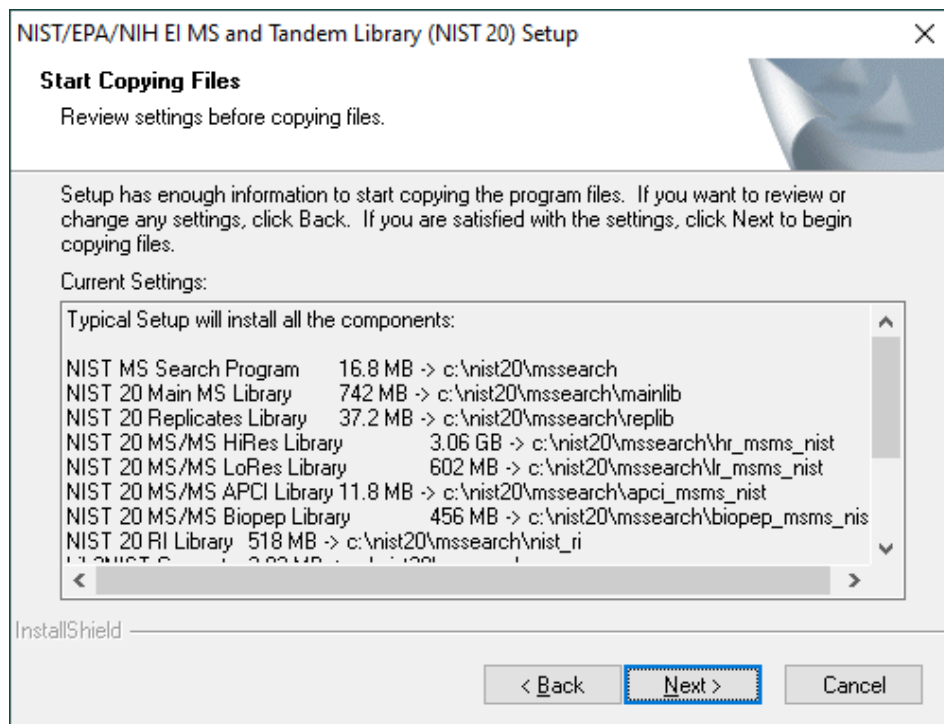
- 10 Select the program folder where the NIST shortcuts will be placed, then click **Next**.



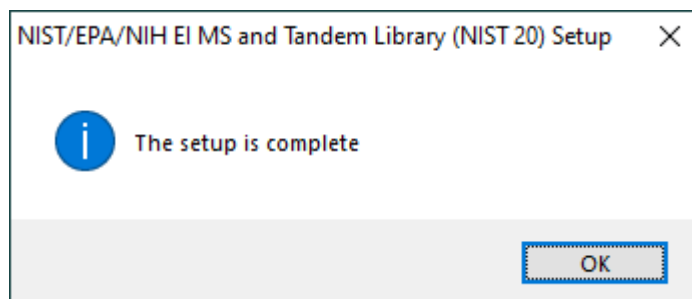
## Set Up NIST Library Search

### Install Software and Library Components

11 Click **Next** to confirm the installation summary.



12 After installation, click **OK** to close the wizard.



13 After installation is done, verify the installation locations.

- a Go to **C:\NIST 20\MSSEARCH\mainlib** and find two files, ALIAS.MSD and LIBSIGN.MSD, and open the ALIAS.MSD file.

You should see the server location (for example, \\nameofserver.site.company.com\subfolder\NIST 20\MSSEARCH\mainlib).

## Set Up NIST Library Search

### Install Software and Library Components

- b** Go to the server location. In the NIST 20 directory.  
There are two folders, AMDIS32 and MSSEARCH. The MSSEARCH folder contains all the NIST library components that are available in Data Analysis.
- c** Check whether the C:\Windows\win.ini file contains the following information. If it does not: Copy the win.ini file to another place for editing. Manually insert the text below to the bottom of win.ini file, save the file and replace C:\Windows\win.ini with the updated win.ini file.

```
[NISTMS]  
Amdis32Path=C:\NIST20\AMDIS32\  
AmdisMSPath=C:\NIST20\AMDIS32\  
Path32=C:\NIST20\MSSEARCH\  
WorkDir32=C:\NIST20\MSSEARCH\
```

#### 14 For Citrix/Terminal Server:

- a** Publish the application from Citrix/Terminal Server setup.
- b** From a thin client, access the Citrix/Terminal Server published link.

On all PCs where NIST is to be installed, select the **Custom** setup type and use the same location for NIST software components (i.e. local C drive) and for NIST MS library components (i.e. shared network folder).

## Configure MS Library in OpenLab CDS

In order to use the common shared NIST MS libraries, you must reference these libraries in a processing method.

- 1 From any client, log in to OpenLab CDS and start Data Analysis.
- 2 Create a new **MS** processing method.
- 3 Under **MS Library Search > Properties**, enter the common shared library paths, and click the + icon to add the libraries. For details, search for *Use the MS library search* in Openlab Help and Learning.
- 4 Save the processing method.
- 5 Load MS data and link it to the prepared processing method.

As long as you use this method, an MS library search will return the same search results, regardless of the client on which it is used.

Alternatively, use any other processing method that references the same NIST MS libraries.

For details on performing an MS Library search, refer to OpenLab Help and Learning.

## 8

# About the OpenLab CDS Software

Software Architecture	81
Components Used in a Client/Server System	83
OpenLab Platform Services and Components	84
21 CFR Part 11 Compliance	85
Data Security	85
Data Integrity	85
Audit traceability	86
Customization	87
Customization via custom calculations	87
Customization via report templates	87
Customize application to start external programs	88
Export raw data and results	88

This chapter contains an overview of the basic software features.



## Software Architecture

OpenLab CDS is a data system solution for analytical workflows that controls a wide variety of instruments including the industry-leading GC and GC/MS-SQ instruments, along with best-in-class LC and LC/MS-SQ. By combining chromatography and single-quad mass spectrometry into a single scalable solution with centralized system administration, you can streamline your laboratory workflows and maximize productivity. A tailored and simplified user interface with a new state-of-the-art user experience, along with e-learning tools, to help you to get up to speed and productive as fast as possible.

The Agilent OpenLab CDS software is provided on read-only USB media that contain all required installables and documents. This includes:

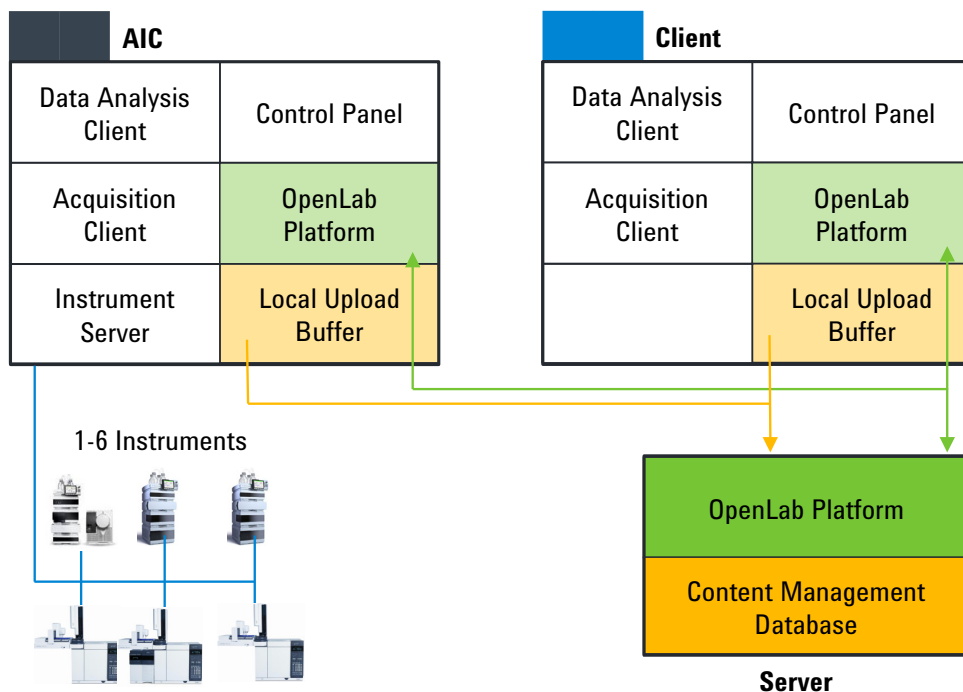
- Acquisition
- Data Analysis and Reporting
- Shared Services
- Content Management
- Custom Calculation Editor
- Help and Learning Platform
- User documentation
- Instrument driver software for Agilent LC, GC, LC/MS, GC/MS, or A/D
- Instrument driver software for non-Agilent instrument drivers
- Instrument driver software for virtual instruments (Data Player)
- Agilent Parts Finder
- Third party tools

With *Client/Server* systems, Content Management and Shared Services are located on a dedicated server. The system supports up to six instrument connections per Analytical Instrument Controller (AIC).



Figure 3 Client/Server System, 1-server system

## Components Used in a Client/Server System



**Figure 4** Components on a Client/Server system

Acquisition client	Acquire raw data with a configured instrument.
Data Analysis Client	Analyze acquired data.
Instrument Server	Control connections to all configured instruments.
Control Panel	User interface to access Shared Services control features.
OpenLab Platform	Provides services and tools that run in the background.
Content Management	Secure storage for your analytical data.

## OpenLab Platform Services and Components

The following services and components are installed as part of the OpenLab Platform. It depends on the type of installation (client, AIC or server) which services are installed.

- **Audit Trail Service**  
Service that encapsulates the retrieval and display of Audit Trail content. Currently not used by OpenLab CDS.
- **Certificate Service and Tool**  
Serves as a solution to secure network traffic in case commercial certificates are not available.
- **Data Collection Service and Agent**  
Service to collect system and topology related information in distributed systems.
- **Data Repository**  
Storage infrastructure to store diagnostic and topology related information.
- **Shared Services**  
Set of services and tools that offer system configuration, security, application configuration, licensing, logging, and notification capabilities.
- **Data Analysis Service**  
Manages Data Analysis instances for improved performance.
- **Distributed Transaction Coordinator Service**  
Handles transaction coordination between microservices. Ensures that all connected actions succeed or fail as one unit.
- **ECM XT Backup services**  
For systems with Content Management or ECM XT: Services to monitor the backup on a schedule, send notifications, and customize the notification process.
- **Electronic Signature Service**  
Provides document signing as a service. To achieve this, ES is based on a centralized storage of electronic signatures.
- **Proxy Configuration Service**  
Provides access to OpenLab components to software applications through an Apache reverse proxy.
- **Test Services**  
Former QualA. Communicates with the Data Repository.

## 21 CFR Part 11 Compliance

To fulfill the FDA rules and guidelines for compliant electronic records and computerized systems, it is important to understand the basic aspects of secure data handling.

- *Data security*: physical protection of data by limiting access to the system and preventing unauthorized access.
- *Data integrity*: protecting raw data and metadata and preventing these from unauthorized modification, and linking raw data and results to reproduce the original results at any time, for example, in an audit situation, and document each new result copy.
- *Audit traceability*: documenting who did what to the results and when, and tracing the user adding new reanalyzed versions to the original raw data.

### Data Security

The Shared Services functionality related to security includes the following (see “[Control Panel](#)” on page 90 for details):

- System Activity Log
- Selection of authentication provider
- Management of users, groups, roles, and privileges
- Security Policy

### Data Integrity

OpenLab CDS stores data in a manner that ensures compliance with 21 CFR Part 11. It provides secure data storage with access control and an audit trail. Data files are versioned to ensure data integrity and traceability. In addition, OpenLab CDS provides electronic signatures allowing users to sign off on data.

## Audit traceability

There are different types of audit trails:

- The *sequence audit trail* is a record of changes made to the sequence when acquiring the data.
- The *method audit trail* provides a detailed list of modifications to a sample preparation method, an acquisition method, or a processing method.
- The *injection audit trail* is the record of a single injection that lists all modifications during the run and in Data Analysis.
- The *result set audit trail* is a superset of injection audit trails for all injections that are contained in a sequence/result set.

The specific behavior of an audit trail depends on the project settings in the Control Panel.

## Customization

OpenLab CDS can be customized to support various workflows and applications. Customizing capabilities are available via different approaches.

For more information on how to use the different tools, refer to OpenLab Help & Learning.

### Customization via custom calculations

Data Analysis can be enriched by calculating additional values. The calculations are done with the Custom Calculation Editor and referenced by or embedded in a processing method.

These calculations can be quite complex. The calculation results are directly visible in Data Analysis, no report generation is needed.

Custom calculations are processed on result set level. They are only computed if all injections of the result set are processed together.

### Customization via report templates

In a report template you can call calculation results from a method-specific custom calculation, or define additional, template-specific calculation expressions. The template-specific values are only visible in the report preview or the final report.

Reports are generated on either injection level, result set level, or across multiple result sets. Reports can be used for automated result evaluation on all the mentioned levels.

Example report templates for typical petrochemical or pharmaceutical applications are provided with the application and can be imported in Data Analysis (see *Import default templates* in OpenLab Help & Learning).

## Customize application to start external programs

The customization capabilities allow to add ribbon groups and icons in the **Data Selection** and **Data Processing** views of OpenLab Data Analysis.

It is possible to start an external program via an icon and to hand over the project data path and the path of the current injection as parameters to the program.

The customization is based on a file CustomToolsConfiguration.xml at C:\ProgramData\AgilentTechnologies\OpenLab DataAnalysis\ that needs to be created by the user. An example CustomToolsConfiguration.xml file is included on the media at **Setup > Tools > Support > UCL > Customization** folder.

For more information, refer to OpenLab Help & Learning.

## Export raw data and results

Data Analysis offers the possibility to execute post processing plug-ins as part of the processing method. These post processing plug-ins allow to export raw data or results as part of the processing routine of single runs as well as sequence runs, also in unattended mode. Scripts are available for the export in the following formats:

- ChemStation Export (.D/.ch format)
- AIA Export (OpenLab CDS 2.x raw data and peak results as netCDF format (revision 3.4) according to AIA Chromatography Data Standard Specification V1.0)
- ASR Export (OpenLab CDS 2.x raw data as an ASR (Analytical Studio Reviewer) file format)
- OpenLab CDS 2.x raw data export (native .dx files for ACE)
- CSV export (OpenLab CDS 2.x raw data as comma separated values in .csv file format)

Other plug-ins such as the Allotrope Data Format (ADF) plug-in are available via SubscribeNet.



## 9

# System Setup and Maintenance

Control Panel	90
Instrument Management / Lab Status at a Glance	90
License Management	91
System Activity Log	92
Diagnostics	93
Administrative Reports	93
Authentication Provider	94
Security Policy	94
User Management	94
Other Maintenance Procedures	95

This chapter contains information on the Control Panel and Shared Services Maintenance. In addition, it contains information on maintenance procedures.

## Control Panel

Using the Control Panel, you can access Shared Services control features such as security policy, central configuration, or lab status at a glance.

### Instrument Management / Lab Status at a Glance

The **Instruments** view in the Control Panel offers an overview of all instruments in the network or on the workstation. You can see the following information for all instruments, summarized on one page:

- Status of the instrument with related color code
- Instrument name
- Instrument location
- Instrument type
- Last change of configuration

Depending on the configuration, this information may be accessed from a single workstation PC or from multiple clients in a network.

You can create a tree of different locations in the Control Panel, and add instruments to these locations. Using locations, you can organize your instruments for example by department, by laboratory, or by lab bench. For each instrument, you can provide basic information such as the name, description, and instrument type.

Depending on your privileges in OpenLab CDS, you can perform several operations on the instruments:

- View instrument information (instrument status, instrument details, activity log)
- View the locations and instruments tree
- Edit the instrument information
- Configure the instrument

The instrument configuration is stored in the Shared Services database. You access the configuration tool from the Control Panel.

- Launch the instrument

On a Workstation, you can only launch instruments that are configured on this PC.

With a Client/Server system, you can launch instruments remotely from any OpenLab CDS client in the network.

Your privileges can differ for the different locations and instruments (see [“Specific Roles for Individual Instruments or Projects”](#) on page 65).

## License Management

This service includes the administration of all licenses that are required for your system.

Before adding a license file, you must first purchase the license and generate the license file using SubscribeNet. For more information on generating new license files, see [“Obtain a License with SubscribeNet”](#) on page 52.

License Management in the Control Panel provides the following functions:

- You can add license files to the license server.
- You can navigate to the license monitor and view the properties of all licenses installed on a given license server.
- You can remove license files from the license server. This may be useful if an invalid license file has been added.
- You can view or change the license server.
- You can view, copy, or save the MAC Address of the license server.
- You can navigate to the Agilent Electronic Software and License Delivery web page to obtain a license.

The following properties are shown for installed licenses:

- **Feature:** This indicates the type of license used.
- **Version:** If a license is versioned, you can see the version number. For licenses that are not versioned, the version is always shown as 1.0.
- **In Use (Available):** This indicates the number of licenses that are currently in use and, in brackets, the total number of licenses. With the OpenLab CDS licensing strategy, a license is only in use as long as a software instance is running (see [“License Types”](#) on page 51).
- **Expiration:** If the license is only valid for a certain period of time, the expiration date is displayed.

In the **Alerts** pane, you are informed if the number of available licenses has gone down to zero for a specific feature, or if you have started a software instance which requires a license that is unavailable.

For more information on adding license files and viewing the license properties, refer to the Control Panel section in OpenLab Help & Learning.

## System Activity Log

The System Activity Log allows you to centrally access all system activities. It contains information on the various events associated with Shared Services or with specific instruments. You can filter the list in order to view only events of a specific type, in a specific time range, created by a specific user, or containing a specific description.

The following types of events are recorded:

- System
- Instrument Management
- Instrument
- Project Management
- Instrument Controller
- User
- Group
- Security
- Printer
- License

The messages can come from other components, such as the user management, or from an instrument module. Instrument messages include error messages, system messages, or event messages. The System Activity Log records these events irrespective of whether you have been alerted to them or not. To get more information on an event, expand the line of interest in the activity logbook viewer.

## Diagnostics

The **Diagnostics** view allows you to access several reports and tools for diagnostic purposes:

- Ping the server.
- Create a report, either for the local system or for the server, with information on the operation system, processors, disk drives, processes, network, and connections.
- Centrally access and download all the log files, trace files, etc. that are created by the registered modules.

## Administrative Reports

In the **Administrative Reports** view, you can additionally create and export various XML or PDF reports related to the system configuration:

### Instrument Controllers Report

Detailed information of all Instrument Controllers. When this report is generated on a Workstation, the information presented relates to the local system. When this report is generated on a client-server system, all Instrument Controllers are included.

### Instruments Report

Provides detailed information about configuration and access privileges for all instruments on the system. On client-server systems, this report includes all instruments on all Instrument Controllers.

### Projects Report

Provides detailed information about configuration and access privileges for all projects on the system.

### Roles and Privileges Report

Describes all roles defined on the system, including details of all privileges included in each role.

### System Report

This report provides a consolidated view of the system, which includes all information about instrument controllers, instruments, projects, roles, users, and groups.

### User's and Group's Role Assignment Report

This report provides an overview of all users and groups with their assigned roles.

## Authentication Provider

The authentication provider is described under *Configure OpenLab CDS*. For details, see ["Configure Authentication"](#) on page 59.

## Security Policy

The security policy is described under *Configure OpenLab CDS*. For details, see ["Configure Security Policy"](#) on page 60.

## User Management

The user management is described under *Configure OpenLab CDS*. For details, see ["Configure users and roles"](#) on page 61.

## Other Maintenance Procedures

### Shared Services and Secure Storage

The Shared Services and secure storage databases are running on the OpenLab Server.

Please refer to the *Agilent OpenLab Server and OpenLab ECM XT Administration Guide* (ECM\_XT\_AdministrationGuide.pdf) for information on:

- FTP server protocol settings
- Database statistics
- Resource monitoring
- Disaster recovery planning

### Backup and Restore

It is mandatory that every OpenLab Server/ECM XT server is backed up regularly. Periodic full backups and differential backups between the full backups are created by OpenLab Server/ECM XT server administrators. These backups are the only way to restore an OpenLab Server/ECM XT server if a hardware or software failure occurs.

The backup only reduces the amount of data loss if a catastrophic system failure occurs. Performing backups guarantees that any data that was committed at the time of the backup can be restored. Data that was queued for upload and not yet committed or was added or updated in the system after the backup was performed will not be recoverable by restoring a backup.

It is also mandatory that the restore procedures are tested to ensure that the backups are performed properly, and can be used for a restore. To do an effective restore, a disaster recovery plan must be created. For details, refer to the *Agilent OpenLab Server and OpenLab ECM XT Administration Guide* guide (ECM\_XT\_AdministrationGuide.pdf).

### Other Products or Databases

For backup procedures of other products, such as the GC Column Database, please refer to the documentation of the respective product.

License Upgrade	97
Get Upgraded License File	97
Add Upgraded License File to the System	98
Upgrade Process	99
Planning the Upgrade	99
Temporary Configuration During Upgrade	99
Step 1: Upgrade the OpenLab Server	101
Step 2: Upgrade an Analytical Instrument Controller	101
Step 3: Upgrade an OpenLab CDS Client	103
Step 4: Upgrade Drivers	105
Register a Client or Instrument Controller on the Server	106
Sample Scheduler for OpenLab	109

This chapter describes the upgrade of the OpenLab CDS software in a Client/Server environment.



## License Upgrade

### Get Upgraded License File

You will need to upgrade your licenses in SubscribeNet prior to upgrading to the next version of OpenLab CDS. We strongly recommend upgrading your licenses *before* upgrading the core software. Standalone workstations or AICs which are upgraded to the new core software version, without a new license, will not work until the new licenses are added to the OpenLab Control Panel.

If you are under SMA subscription, proceed as follows to upgrade your licenses:

- 1 During the following process, you will be prompted in SubscribeNet for the host name or MAC address of the server where OpenLab CDS is already installed.  
  
To retrieve this hostname and MAC address, open the Control Panel and browse to the **Administration > Licenses** section. Note down the host name and use the **Copy MAC Address** or **Save MAC Address** function to obtain the MAC address.
- 2 Log into the Agilent Electronic Software and License Delivery (<https://agilent.subscribenet.com/>).
- 3 Navigate to **Manage Licenses by Host**. In the **Host ID** field, enter the previously noted MAC address, and click **Search**.



<b>Software</b>	<b>Electronic Software and License Delivery</b>
Product List	<b>Manage Licenses by Host</b>
Product Search	Manage Licenses by Host provides a convenient way to upgrade or return all licenses for a license server. For nodelocked license Hosts, licenses can only be returned as upgrades do not apply.
Entitlements	Please select a host by entering the Host ID or Host Name in its search field, or select a host from the pull-down list. You can use * as a wildcard value.
<b>Licenses</b>	
Register Authorization Code	Host ID <input type="text"/> What is my Host ID?
Generate or View Licenses	Host Name <input type="text"/>
View Licenses by Host	<input type="button" value="Search"/>
Manage Licenses by Host	
<b>Administration</b>	
Account Members	
Change Password	

If the relevant host name does not appear, you may be managing your licenses in multiple SubscribeNet accounts. You will need to log into those accounts to upgrade those licenses.

- 4 If your license(s) are eligible for an upgrade, you will see the **Upgrade All** button. Otherwise you will need to contact your Agilent Sales Representative to renew your Software Maintenance Agreement (see [“Sales and Support Assistance”](#) on page 123). To proceed with generating your upgrade license, click the button.
- 5 On the **Upgrade All Licenses for License Host** page, review the data, and confirm by clicking **Upgrade All**.  
This upgrades the license file to the most current version. SubscribeNet will send you an email with a new license file.
- 6 Put the new license file on your system (see [“Add Upgraded License File to the System”](#) on page 98).

## Add Upgraded License File to the System

If you have purchased new options, such as additional instrument controls or client license and regenerated your license in SubscribeNet, the upgraded license file must be re-applied to the system.

- 1 Start the Control Panel from any machine connected to the system you want to install the license for.
- 2 Navigate to **Administration > Licenses**.
- 3 Select a license, then click **Remove License** .
- 4 In the ribbon, click **Add License** .
- 5 Browse to and open the license file saved from the license generation process in SubscribeNet.

### NOTE

A full restart is required in order for any license to have an immediate effect.

## Upgrade Process

### Planning the Upgrade

**NOTE**

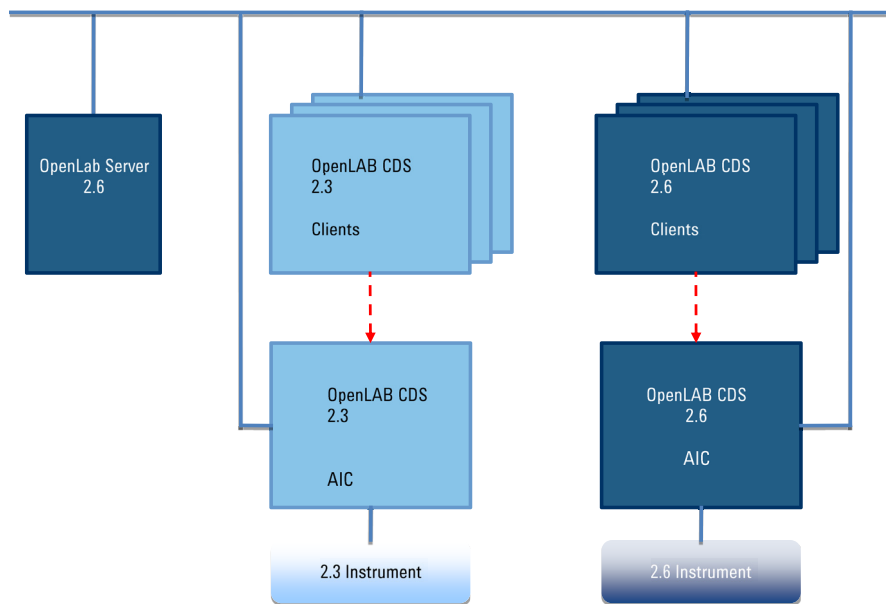
Ensure that all active acquisitions are halted prior to the upgrade.

The components must be upgraded in the following order:

- 1 OpenLab Server
- 2 Analytical Instrument Controller (AIC)
- 3 OpenLab CDS Clients
- 4 If required, upgrade instrument drivers

### Temporary Configuration During Upgrade

During an upgrade project there might be different versions of OpenLab CDS in your environment. An environment with an OpenLab Server 2.6 supports using clients and AICs with OpenLab CDS versions 2.3 or higher. Analytical Instrument Controller and OpenLab CDS Clients must always use the same version of OpenLab CDS.



**Figure 5** Mixed configuration during upgrade

Use the 2.6 clients to access instruments that are controlled by a 2.6 AIC, and use older clients to access instruments controlled by a corresponding older AIC. The data storage is provided by OpenLab Server 2.6. Both environments share the same storage. Use an instrument structure in the Control Panel that helps you differentiating the instruments.

#### NOTE

Always launch and close instruments from the same system. Do not launch an instrument from version 2.6 and close it from version 2.3/2.4/2.5, or vice versa.

#### NOTE

The analysis and reprocessing of data is supported only on versions same as or higher than the one used for the acquisition or last reprocessing.

## Step 1: Upgrade the OpenLab Server

For more information on upgrading an OpenLab Server, refer to the *OpenLab Server and OpenLab ECM XT* installation guide (ECM\_XT\_InstallationGuide.pdf).

## Step 2: Upgrade an Analytical Instrument Controller

### Prerequisites

You have upgraded the OpenLab Server to version 2.6.

### CAUTION

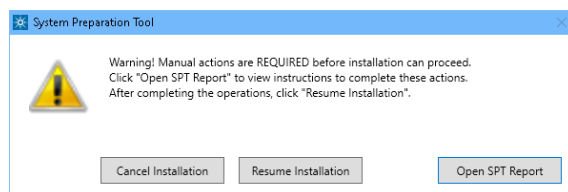
**Uninstallation of ADFExport for OpenLab 1.2 after an upgrade may cause the Data Analysis application to fail to start**

Rev. 1.2 of ADFExport for OpenLab is not compatible to OpenLab CDS 2.6 or higher. Uninstallation of ADFExport for OpenLab 1.2 after an upgrade (e.g. during its own upgrade) may remove vital files from OpenLab CDS. This may cause the Data Analysis application to fail to start. Automated processing and data analysis during acquisition may fail. A re-installation of OpenLab CDS might become necessary.

- ✓ If Rev. 1.2 of ADFExport for OpenLab is installed with your OpenLab CDS version, uninstall it *before* upgrading to OpenLab CDS to Rev. 2.6, or upgrade to the next revision of ADFExport for OpenLab.

- 1 Run the setup.exe file from the installation media as a user with administrative rights.
- 2 Select **OpenLab CDS**.
- 3 In the OpenLab CDS Installer, select the **Installation** screen.
- 4 Click **Install/Upgrade**.
- 5 In the **Storage Configuration** screen, click **Next**.
- 6 In the **Server Information** screen, click **Test Connection** to check the availability of the server. If the server is available, provide the OpenLab CDS credentials and click **Next**.

- 7 In the **System Preparation** screen, the upgrade wizard shows the list of recommended settings for the system. You may unselect items that you do not want to apply on the system. Other mandatory settings will be applied automatically during the upgrade. For an overview of both mandatory and recommended settings, see [Table 2](#) on page 17.  
Click **Next** to proceed to the **Review** page. The settings will be applied as part of the installation.
- 8 In the **Review** screen, click **Upgrade** to start the upgrade.
- 9 If there are still manual actions required before installation, a warning is shown.



Options to proceed

- *Recommended:* Click **Open SPT Report** to view instructions to complete these actions. After completing the operations, click **Resume Installation**.
- **Cancel Installation:** Installation is aborted. Make necessary updates and restart installation.
- **Resume Installation:** The dialog is closed, and installation continues with the installation of the CDS components even if a setting was not applied, or the Operating System does not meet requirements.

- 10 On the **Finish** page, click **Run Software Verification**.
- 11 On the **Finish** page, keep the **Reboot** check box selected, and click **Finish**.

## Step 3: Upgrade an OpenLab CDS Client

### Prerequisites

You have upgraded the Analytical Instrument Controller to version 2.6.

### CAUTION

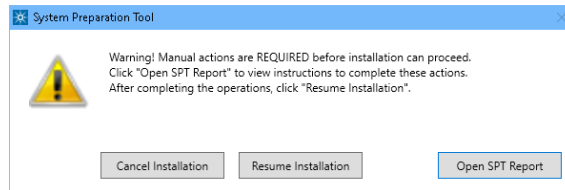
**Uninstallation of ADExport for OpenLab 1.2 after an upgrade may cause the Data Analysis application to fail to start**

Rev. 1.2 of ADExport for OpenLab is not compatible to OpenLab CDS 2.6 or higher. Uninstallation of ADExport for OpenLab 1.2 after an upgrade (e.g. during its own upgrade) may remove vital files from OpenLab CDS. This may cause the Data Analysis application to fail to start. Automated processing and data analysis during acquisition may fail. A re-installation of OpenLab CDS might become necessary.

- ✓ If Rev. 1.2 of ADExport for OpenLab is installed with your OpenLab CDS version, uninstall it *before* upgrading to OpenLab CDS to Rev. 2.6, or upgrade to the next revision of ADExport for OpenLab.

- 1 Run the setup.exe file from the installation media as a user with administrative rights.
- 2 Select **OpenLab CDS**.
- 3 In the OpenLab CDS Installer, select the **Installation** screen.
- 4 Click **Install/Upgrade**.
- 5 In the **Storage Configuration** screen, click **Next**.
- 6 In the **Server Information** screen, click **Test Connection** to check the availability of the server. If the server is available, provide the OpenLab CDS credentials and click **Next**.
- 7 In the **System Preparation** screen, the upgrade wizard shows the list of recommended settings for the system. You may unselect items that you do not want to apply on the system. Other mandatory settings will be applied automatically during the upgrade. For an overview of both mandatory and recommended settings, see [Table 2](#) on page 17.  
Click **Next** to proceed to the **Review** page. The settings will be applied as part of the installation.

- 8 In the **Review** screen, click **Upgrade** to start the upgrade.
- 9 If there are still manual actions required before installation, a warning is shown.



Options to proceed

- *Recommended:* Click **Open SPT Report** to view instructions to complete these actions. After completing the operations, click **Resume Installation**.
- **Cancel Installation:** Installation is aborted. Make necessary updates and restart installation.
- **Resume Installation:** The dialog is closed, and installation continues with the installation of the CDS components even if a setting was not applied, or the Operating System does not meet requirements.

- 10 On the **Finish** page, click **Run Software Verification**.
- 11 On the **Finish** page, keep the **Reboot** check box selected, and click **Finish**.



## Step 4: Upgrade Drivers

If you use a GC/MS instrument: Upgrade the GC/MS firmware to the latest revision. The upgrade is required for the GC/MS software to work properly with OpenLab CDS 2.6.

It is recommended to reconfigure the instrument in the Control Panel.

For more information on upgrading drivers, see ["Install or Upgrade Instrument Driver Software"](#) on page 38.

## Register a Client or Instrument Controller on the Server

This procedure describes how to add an existing OpenLab CDS PC to a distributed environment as a client or Instrument Controller.

Carry out the procedure also after manually installing or upgrading instrument driver software to make sure that all instrument types will be available in the Control Panel.

Ensure that the OpenLab CDS version of your PC is compatible with the OpenLab Server version. See [“Temporary Configuration During Upgrade”](#) on page 99.

- 1 In Windows, select **Start > All Programs > Agilent Technologies > OpenLab Configuration**.
- 2 In the **OpenLab Configuration** tool, enter the hostname (without \\) of the OpenLab server. Click **Connect** to enable the input boxes for the user credentials.

### NOTE

If you register an *Instrument Controller*: Be sure to provide the correct server name. Random switching of Instrument Controllers from one server to another is not supported. Servers can only be switched by following a defined process in the context of a fallback solution in case of network failures.

The image shows a software configuration window titled "OpenLAB Configuration". It is divided into three main sections on the left: "Step 1 - Server", "Step 2 - Authentication", and "Step 3 - Additional Options".

- Step 1 - Server:** Contains a text input field for "Hostname" and a "Connect" button.
- Step 2 - Authentication:** Contains three input fields: "Username", "Password", and "Domain".
- Step 3 - Additional Options:** Contains a checked checkbox labeled "Register as Instrument Controller".

On the right side of the window, there is a blue informational panel with the following sections:

- Server:** Text: "Enter the hostname of the OpenLAB Server that you would like to connect to and click the Connect button."
- Text: "Changes will be applied to the server specified."
- Authentication:** Text: "Enter your login credentials for the server specified."
- Additional Options:** Text: "Choose any additional options and click the Register button."

At the bottom right of the window are two buttons: "Cancel" and "Register".

- 3 Under **Authentication**, enter OpenLab user credentials for the specified server.
- 4 For Instrument Controllers, select the **Register as Instrument Controller** check box. For clients, clear this check box.
- 5 Click **Register**.  
A Registration Results dialog will be shown.
- 6 Reboot the client or AIC.

### Register root certificate

When using self-signed certificates issued by Certificate Service, registering a new server for a client or Instrument Controller requires additional configuration steps. These steps have to be executed once so that the root certificate of the new server is available on the respective machine. Otherwise, connecting to an instrument might fail. Follow these steps to download the root certificate and register it in the local certificate store of a client of Instrument Controller.

- 1 Open the Windows Command Prompt as an administrator.
- 2 Navigate to the Certificate Service Tool folder: C:\Program Files (x86)\Agilent Technologies\OpenLab Certificate Service Tool\Bin.
- 3 Run the following command (replace <servername> by the name of your server):

```
Agilent.OpenLab.CertService.CertTool.exe  
RegisterRootCert -h <servername>
```

## Sample Scheduler for OpenLab

As of OpenLab CDS version 2.6, Sample Scheduler for OpenLab is installed by default with an OpenLab CDS Client/Server installation. The following components are available in the Client/Server topology after upgrading to 2.6:

- Server: Sample Scheduler Database as part of the Data Repository, LIMS Agent, Web Server
- Instrument Controller: Sample Scheduler Agent
- Client: Sample Scheduler Client, Sample Scheduler Configuration

Sample Schedule services are disabled after installation. To try out Sample Scheduler for OpenLab select **Agilent Technologies > Agilent Sample Scheduler** from the Start menu. Click **Access free dashboard**, and follow the instructions.

If you decide to regularly use Sample Scheduler or want to evaluate more capabilities please activate it first. For details on activation see information page that is launched.

# 11

## Uninstall OpenLab CDS

Uninstall a Client or Analytical Instrument Controller (AIC) 111

Uninstall the OpenLab Server 111

Uninstall OpenLab Help and Learning Only 112

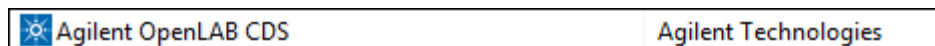
This chapter describes the uninstallation of the OpenLab CDS software.

To uninstall OpenLab CDS:

- 1 Uninstall OpenLab CDS from clients and Analytical Instrument Controllers.
- 2 Uninstall the OpenLab Server.

## Uninstall a Client or Analytical Instrument Controller (AIC)

- 1 Log in as an administrator.
- 2 In the Windows Settings, open **Apps > Apps and features**.
- 3 Uninstall OpenLab CDS add-ons that you installed separately (for example, **ADFExport for OpenLab**).
- 4 Select **Agilent OpenLab CDS**, and confirm uninstallation.



The OpenLab CDS Uninstallation Wizard opens<sup>1</sup>. In the wizard, click **Uninstall**, and follow the wizard.

- 5 In the Windows Settings under **Apps > Apps and features**, select **Agilent Software Verification Tool** and confirm uninstallation.
- 6 Drivers installed by OpenLab CDS are uninstalled automatically. If you installed other drivers, they are still listed under **Apps and features**. Uninstall them manually.
- 7 Reboot.

## Uninstall the OpenLab Server

For details on the uninstallation, please refer to the *Agilent OpenLab Server and OpenLab ECM XT Installation Guide* (ECM\_XT\_InstallationGuide.pdf).

<sup>1</sup> The Uninstallation wizard can also be started from the Installation Wizard used to install the software.

## Uninstall OpenLab Help and Learning Only

If you installed OpenLab Help and Learning only, follow this procedure to uninstall it.

- 1 From the installation media, right-click the setup.exe file, and run it as administrator.
- 2 On the start screen, select **OpenLab CDS**, and click **OK**.
- 3 On the **Documentation** page, select **Install OpenLab Help and Learning Only**.  
The **Agilent OpenLab CDS Help and Learning** wizard opens.
- 4 Click **Remove**.  
The wizard removes OpenLab Help and Learning from your system.
- 5 After completion of the uninstallation, click **Finish** to close the wizard.



Privileges in the Control Panel	114
Project Privileges	114
Instrument Privileges	121
Administrative Privileges	121
Sales and Support Assistance	123

## Privileges in the Control Panel

The privileges described in the following can be associated with different roles in the Control Panel. The following roles are available:

- Everything
- System Administrator
- Instrument Administrator
- Project Administrator
- Project Content Deletion
- Instrument User
- Technician
- Chemist
- Archivist
- Content Management Approver
- Content Management Contributor
- Content Management Reader
- Content Management PDF Template Manager
- Content Management Administrator

In the Control Panel under **Administration > Roles**, you can view or change the associated privileges, or create your own roles.

## Project Privileges

**Table 5      Acquisition Method**

Name	Description
Create and modify acquisition method	Create, edit and save an acquisition method file (*.amx)
Use generic acquisition methods	Use acquisition methods that have the status <b>Generic</b> .

**Table 6      Audit Trail**

Name	Description
Change method audit trail settings	Edit and save method audit trail settings (project properties in the Control Panel).
Review audit trail	Confirm that you reviewed a changed audit trail.
Add manual audit trail entry	Add a manual entry to document your own actions in the audit trail.

**Table 7      Control**

Name	Description
Abort any running sample	Abort any running sequence or single run.
Manual control (in run)	Access manual control functions while the instrument is running.
Manual control (only when instrument idle)	Access manual control functions while the instrument is idle.
MS autotune and manual tuning	Access all MS tune and maintenance functionality, including manual tune, autotune, and check tune.
MS autotune	Perform MS autotune and check tune.
Delete any pending run	Delete pending runs in the run queue submitted by any user. No privilege is required for a user to delete own pending runs
Reorder pending runs	Reorder pending run queue items in the run queue. Moving items around in the run queue.
Run priority sample	Submit <i>Priority</i> single samples.

**Table 8      Custom Tools**

Name	Description
Access Custom Tools section	Start external programs that were added to the application via the customization tool

**Table 9      Data**

Name	Description
Export data	Export data into an OpenLab archive (*.olax).
Import data	Import data from OpenLab archives (*.olax) into the OpenLab system.
Save reports to disk	Save or export a report to a location on a disk or network share.
Edit sample information	Edit information in the <b>Injection List</b> window.

**Table 10 Data Processing**

Name	Description
Create new result set	Combine single samples or sequences from different sources in a new, self-assembled result set.
Disable/enable calibration points	Disable or enable individual points of a calibration curve.
Do manual compound identification	Manually assign a compound to a peak.
Do manual chromatogram extraction	Manually extract MS (TIC-SIM/TIC-SCAN) chromatograms from your data.
Do manual integration	Activate manual integration in the <b>Chromatograms</b> window.
Do manual MS library search	Manually search for matches in an MS library.
Do manual spectrum extraction	Manually extract UV or MS spectra from your data.
Integration Optimizer: Display peak results	Show the retention time, area, and area % for each integrated peak in the Integration Optimizer.
Launch Custom Calculation Editor	Start the Custom Calculation Editor from Data Analysis.
Update master processing method	Save changes from a result set method to the corresponding master processing method in the Methods folder.
Use Integration Optimizer	Edit parameters for the integration optimizer, step through the wizard, and use functions offered by the wizard.
Reprocess data	Reprocess injections or result sets.
Print results reports	Create reports for your methods or results.

**Table 11 E-Signature**

Name	Description
E-Signature Sign Data Files	User can sign data files

Table 11 E-Signature

Name	Description
Revoke E-Signature	User can revoke the e-signature.
Change method status	User can change the status (Generic, Approved, Obsoleted) of sample prep methods, acquisition methods and processing methods.

Table 12 File and Folder Operations

Name	Description
Delete report templates	Delete report templates (*.rdl) in the <b>Data Selection</b> view of Data Analysis (requires also the <b>Project Management &gt; Delete content of project</b> privilege).
Delete sequence templates	Delete sequence templates (*.stx) files in the <b>Data Selection</b> view of Data Analysis (requires also the <b>Project Management &gt; Delete content of project</b> privilege).
Delete methods	Delete processing methods (*.pmx) or acquisition methods (*.amx) in the <b>Data Selection</b> view of Data Analysis (requires also the <b>Project Management &gt; Delete content of project</b> privilege).

Table 13 Lock

Name	Description
Lock Results	Lock a result set to protect it from being changed.
Unlock Results	Unlock a locked result set.

Table 14 Processing Method

Name	Description
Create processing method	Create a new processing method (*.pmx), or save a method under a new name.
Save master method	Save changes to a processing method in the Methods folder.

Table 14 Processing Method

Name	Description
Save result set method	Save changes to a processing method in the result set folder.
Edit integration parameters	View and edit the parameters in the <b>Integration Events</b> section of a method.
Edit identification parameters	View and edit the parameters in the <b>Compounds &gt; Identification</b> section of a method.
Edit chromatogram extraction parameters	View and edit the parameters in the <b>Extraction &gt; Chromatogram</b> section of a method.
Edit spectrum extraction parameters	View and edit the parameters in the <b>Extraction &gt; Spectrum</b> section of a method.
Edit MS library search parameters	View and edit the parameters in the <b>MS Library Search &gt; Properties</b> section of a method.
Edit calibration parameters	View and edit the parameters in the <b>Compounds &gt; Calibration</b> section of method.
Edit spectra parameters	View and edit the parameters in the <b>Compounds &gt; Spectra</b> section of method.
Edit system suitability parameters	View and edit the parameters in the <b>Compounds &gt; System Suitability</b> section of method.
Edit custom calculation parameters	View and edit the parameters in the <b>Tools &gt; Custom Calculation</b> section of a method.
Edit signal parameters	View and edit the parameters in the <b>General &gt; Signals</b> section of a method.
Edit sample purity parameters	View and edit the parameters in the <b>MS Sample Purity</b> section of a method.
Edit reporting parameters	View and edit the parameters in the <b>Reports &gt; Injection Report</b> section of a method.
Edit general parameters	View and edit the parameters in the <b>General &gt; Properties</b> section of a method.
Load older master method	With Content Management, load an older version of a master method.

Table 14 Processing Method

Name	Description
Edit Post Processing Plugins parameters	View and edit the parameters in the <b>Post Processing Plugins</b> section of a method.
Use generic processing methods	Use processing methods that have the status <b>Generic</b> .

Table 15 Project Management

Name	Description
Manage project or project group	User can create or edit project properties and can move the project but cannot view or edit the project access settings.
Manage project or project group access	User can view and edit the project access settings.
View project or project group	User can see a project and project details but cannot edit. <i>Note:</i> This privilege is required for all users.
Access content using web client	User can view the data via the Content Management web interface.
Edit content of project	User can create new versions of documents (e.g. data, methods, or templates).
Delete content of project	<p>If using OpenLab CDS with Content Management:</p> <ul style="list-style-type: none"> <li>• Required to move and delete project files or folders (e.g. data, methods, or templates) via the Content Management web interface.</li> <li>• Required to delete templates and methods in the <b>Data Selection</b> view of Data Analysis.</li> </ul> <p><i>Note:</i> To move or delete files and folders in the Content Management database that are <i>outside</i> of a project, this privilege is not sufficient; the administrative privilege <b>Manage security</b> must be granted!</p>

**Table 16 Report Template**

Privilege	Description
Unlock/lock report template items	Lock and unlock report template items (tables, chromatograms, groups of items, ...) to control who is allowed to modify those.
Validate report template	Confirm usage of report templates that have been modified outside OpenLab CDS.
Create report template	Create and edit report templates in the Reporting view.

**Table 17 Sample Prep**

Name	Description
Create and modify sample prep	View, edit, and save an autosampler sample prep file
Use generic sample prep methods	Use sample prep methods that have the status <b>Generic</b> .

**Table 18 Sequence Template**

Name	Description
Create and modify sequence template	Create, edit and save sequence creation templates (*.stx).

**Table 19 Sequence**

Name	Description
Edit any users running sequence	Edit any user's running sequence (status <b>Acquiring</b> in the Run Queue).
Create and modify sequence	Create, edit and save sequences (*.sqx)
Edit users own running sequences	Edit your own running sequences (status <b>Acquiring</b> in the Run Queue).
Edit method override parameters	Override parameters in a predefined acquisition method.



Table 20 Snapshot

Name	Description
Review snapshot results	From Acquisition, open a currently running sample in Data Analysis.

## Instrument Privileges

Table 21 Instrument Management

Name	Description
View instrument or location	User can view and access a location in the tree, but not edit access security, can view properties.
Manage Instrument or location	User can create and move locations and edit properties (name, description, etc).
Manage instrument or location access	User can view and edit the location access settings.
Run instrument	User can start an instrument session.
Service instrument	User can lock or unlock an instrument (to service it).

## Administrative Privileges

Table 22 System Administration

Name	Description
Manage printers	Can add/remove printers and print server.
Edit activity log properties	Can change the Activity log Settings in the Control Panel (that is, can turn logging on for the System Activity Log).
Create administrative reports	Can create any of the system admin reports.
Manage system components	Can install/remove components (applications).

**Table 22    System Administration**

Name	Description
Manage security	<p>Can change security settings and assign security roles.</p> <p>Can edit (add, change etc) users, groups and roles.</p> <p>In the Content Management web interface: Can move and delete files and folders outside of a project.</p> <p><i>Note:</i> A user with this privilege can grant himself access to all settings in Shared Services. Be careful who you grant the Manage Security privilege.</p>
Manage instrument controllers	Can edit Instrument Controllers in the Control Panel.
Unlock any locked UI	Can log in to another user's locked session.

**Table 23    Content Management**

Name	Description
Archive content	User can archive the content of the Content Management data repository.
Manage Templates	View, create, update and delete PDF templates.

## Sales and Support Assistance

Please check the following web site for your local sales and support contact:

<https://www.agilent.com/en/support>

Agilent Community

<https://community.agilent.com/> To get answers to your questions, join over 10,000 users in the Agilent Community. Review curated support materials organized by platform technology. Ask questions to industry colleagues and collaborators. Get notifications on new videos, documents, tools, and webinars relevant to your work.

## In This Book

This document provides instructions for installation, configuration, administration, and maintenance of OpenLab CDS Clients and Analytical Instrument Controllers. It includes information on the license generation with SubscribeNet and operating system configuration.

The manual describes the following:

- Install an OpenLab CDS Client or AIC
- Generating and Downloading Your Software License
- Configure OpenLab CDS
- Optional Procedures
- Customization
- About the OpenLab CDS software
- System Setup and Maintenance

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

© Agilent Technologies Inc. 2015-2021  
Edition: 04/2021

Document No: D0007072 Rev. A

