

# Agilent OpenLAB Chromatography Data System (CDS)

Guide for Administrators



**Agilent Technologies**

# Notices

© Agilent Technologies, Inc. 2010-2016

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

## Manual Part Number

M8305-90017

## Edition

03/2016

Printed in Germany

Agilent Technologies  
Hewlett-Packard-Strasse 8  
76337 Waldbronn

### Software Revision

This guide is valid for revision A.02.02 SR2 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

If software is for use in the performance of a U.S. Government prime contract or subcontract, Software is delivered and licensed as “Commercial computer software” as defined in DFAR 252.227-7014 (June 1995), or as a “commercial item” as defined in FAR 2.101(a) or as “Restricted computer software” as defined in FAR 52.227-19 (June 1987) or any equivalent agency regulation or contract clause. Use, duplication or disclosure of Software is subject to Agilent Technologies’ standard commercial license terms, and non-DOD Departments and Agencies of the U.S. Government will

receive no greater than Restricted Rights as defined in FAR 52.227-19(c)(1-2) (June 1987). U.S. Government users will receive no greater than Limited Rights as defined in FAR 52.227-14 (June 1987) or DFAR 252.227-7015 (b)(2) (November 1995), as applicable in any technical data.

## Safety Notices

### CAUTION

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

### WARNING

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

## In this Guide ...

This manual describes the concepts of OpenLAB CDS (system architecture, licensing strategy, data integrity) and the administration of OpenLAB CDS with the OpenLAB Control Panel. In addition, it contains specific information on the administration of OpenLAB CDS ChemStation Edition. This Edition also includes information on the OpenLAB Data Analysis Add-on.

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

<b>Term</b>	<b>Description</b>
ChemStation	OpenLAB CDS ChemStation Edition
EZChrom	OpenLAB CDS EZChrom Edition
ECM	OpenLAB Enterprise Content Manager
Data Store	OpenLAB Data Store
AIC	Agilent Instrument Controller

### 1 Architectural Concepts of OpenLAB CDS

This chapter gives you an overview of the general concepts of Agilent OpenLAB CDS architecture. In the following, the terms ChemStation and EZChrom refer to OpenLAB CDS ChemStation Edition and OpenLAB CDS EZChrom Edition.

### 2 OpenLAB Control Panel

Using the OpenLAB Control Panel, you can access OpenLAB Shared Services control features such as security policy, central configuration, or lab status at a glance. These features are described in more detail in this chapter.

### 3 OpenLAB Shared Services Server

This chapter describes the OpenLAB Server Utility Tool.

#### **4 ChemStation-Specific Administration**

This chapter describes various tools that are helpful for diagnosis, support and troubleshooting.

#### **5 EZChrom-Specific Administration**

This chapter describes fail over procedures and tools for AIC administration.

#### **6 Appendix**

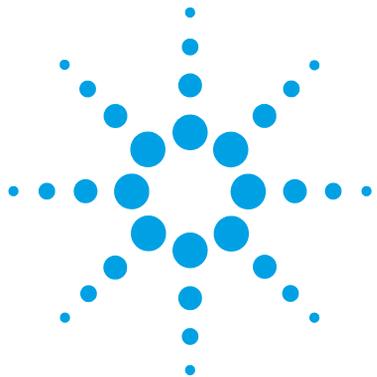
This chapter contains information on privileges used in OpenLAB CDS, and on driver license features for instruments of other vendors than Agilent.

# Contents

<b>1</b>	<b>Architectural Concepts of OpenLAB CDS</b>	<b>7</b>
	Overview	8
	OpenLAB CDS System Architecture	10
	OpenLAB CDS Licensing	23
	Security and Data Integrity	42
<b>2</b>	<b>OpenLAB Control Panel</b>	<b>45</b>
	Project Management	46
	Instrument Management / Lab Status at a Glance	46
	License Management	48
	System Activity Log	50
	Diagnostics	51
	Administrative Reports	52
	Authentication Provider	53
	Security Policy	56
	User Management	58
<b>3</b>	<b>OpenLAB Shared Services Server</b>	<b>63</b>
	OpenLAB Server Utility Tool	64
	Maintenance Procedures	71
<b>4</b>	<b>ChemStation-Specific Administration</b>	<b>75</b>
	ChemStation Administration Tool	76
	Support Reports	89
	OpenLAB CDS Config Checker	90
	Fail Over Procedures for ChemStation Edition	91

## Contents

<b>5 EZChrom-Specific Administration</b>	<b>113</b>
Fail Over Procedures for OpenLAB CDS EZChrom Edition	114
AIC Driver and Install Tool	120
Change the Password of the AFS Administrator	122
<b>6 Appendix</b>	<b>125</b>
Privileges in OpenLAB Control Panel	126



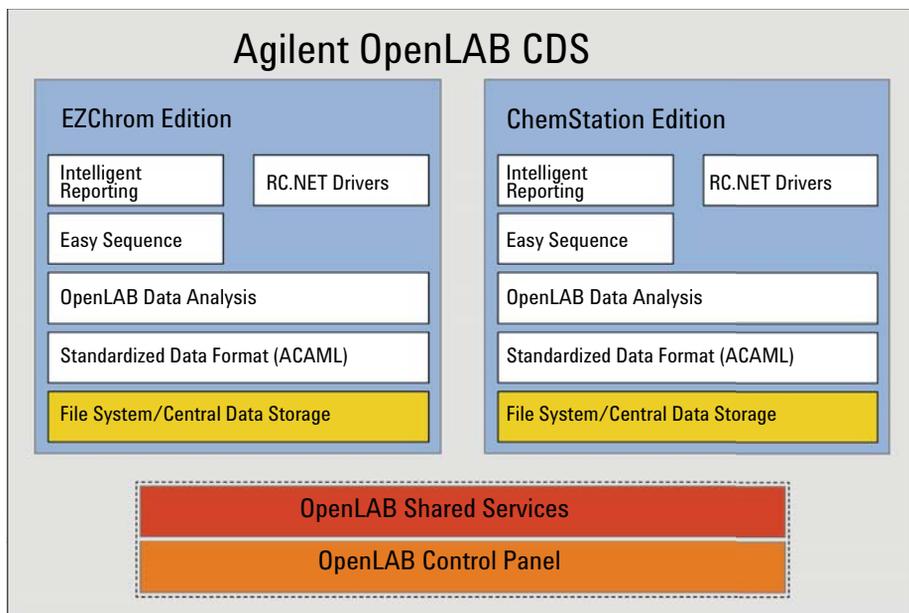
# 1 Architectural Concepts of OpenLAB CDS

Overview	8
OpenLAB CDS System Architecture	10
Workstation	10
Networked Workstation	11
Distributed Systems	14
Networked Workstations in Distributed Systems	20
Data Compatibility Between Different Versions	22
OpenLAB CDS Licensing	23
General Product Structure	23
License Types	24
Licensing Scheme	25
Main Features for Licensing	28
License Features Associated with Agilent OpenLAB CDS Products	31
Licensing Examples	35
VL License Features	38
Flexera License Manager	39
License Upgrade	40
Windows Server Licensing	41
Security and Data Integrity	42
Security Aspects	42
Data Integrity	42

This chapter gives you an overview of the general concepts of Agilent OpenLAB CDS architecture. In the following, the terms ChemStation and EZChrom refer to OpenLAB CDS ChemStation Edition and OpenLAB CDS EZChrom Edition.



## Overview

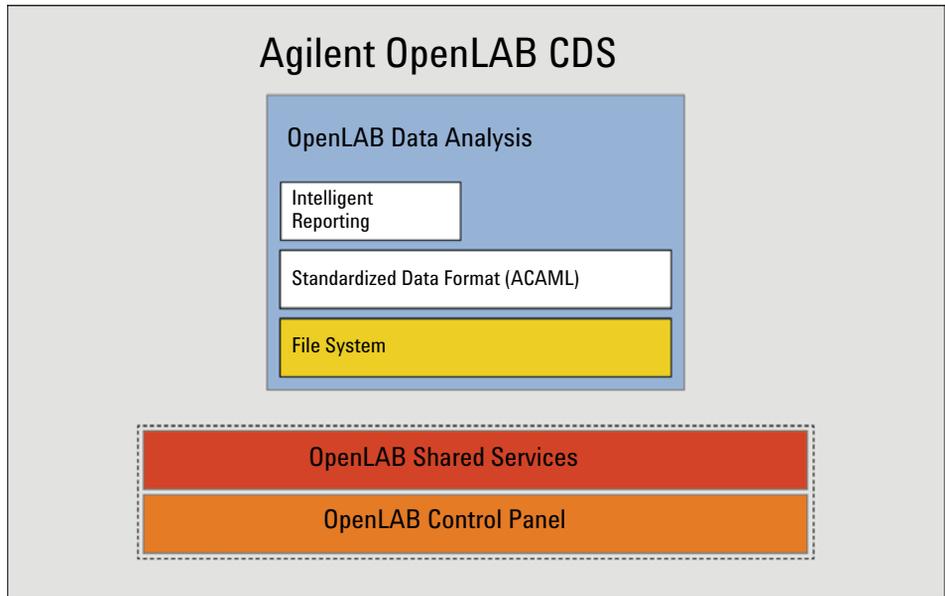


**Figure 1** OpenLAB CDS architecture

OpenLAB CDS contains the following software and interface modules:

- OpenLAB Control Panel  
The OpenLAB Control Panel is the user interface that provides access to OpenLAB CDS components, as well as administrative functions used for managing OpenLAB Shared Services features.
- OpenLAB Shared Services  
These services offer central access, central configuration, lab status at a glance, and remote control of instruments. The central functions can be used by all OpenLAB modules.
- Instrument control, data acquisition and data analysis module (ChemStation/EZChrom)  
This module is available as either ChemStation or EZChrom Edition. Data acquired with previous versions of the respective software can be

processed. ChemStation and EZChrom share several common functions such as Easy Sequence, RC.NET drivers, or Intelligent Reporting. OpenLAB Data Analysis can be installed as an add-on to ChemStation or EZChrom, or as a standalone application for offline data review.



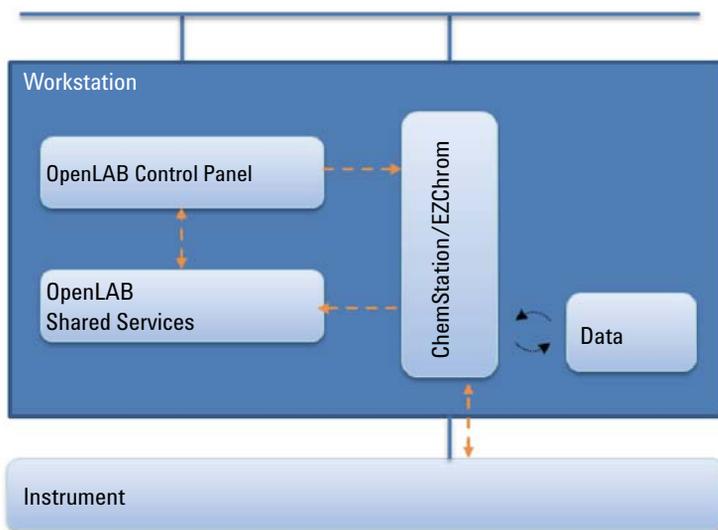
**Figure 2** OpenLAB Data Analysis architecture

## OpenLAB CDS System Architecture

### Workstation

In small laboratories, you can install all components of OpenLAB CDS on one single workstation. As a consequence, OpenLAB Shared Services runs on the same PC as ChemStation or EZChrom.

The following figure shows the configuration of an OpenLAB CDS Workstation. Only one ChemStation/EZChrom instance is shown, but you can configure multiple instances and associated instruments on the workstation.



**Figure 3** OpenLAB CDS Workstation



## OpenLAB Data Analysis with Workstations

With OpenLAB Data Analysis, you can process and analyze the data acquired with ChemStation or EZChrom. OpenLAB Data Analysis is supported on single workstations under specific conditions (see [Table 2](#) on page 11). You can install OpenLAB Data Analysis as an add-on to OpenLAB CDS on a single workstation or as a stand-alone application with OpenLAB Shared Services. For more information on OpenLAB Data Analysis, refer to the *OpenLAB Data Analysis - Getting Started* guide.

**Table 2** Supported Scenarios

Installed component	Storage Type	Supported?
ChemStation Edition with OpenLAB Data Analysis	Data Store ECM	No
EZChrom Edition with OpenLAB Data Analysis	Data Store ECM	No
OpenLAB Data Analysis stand-alone (ChemStation data)	Data Store ECM	Supported using the Agilent OpenLAB Data Analysis Upload/Download tool
OpenLAB Data Analysis stand-alone (EZChrom data)	Data Store ECM	No  Supported using the Agilent OpenLAB Data Analysis Upload/Download tool

## Networked Workstation

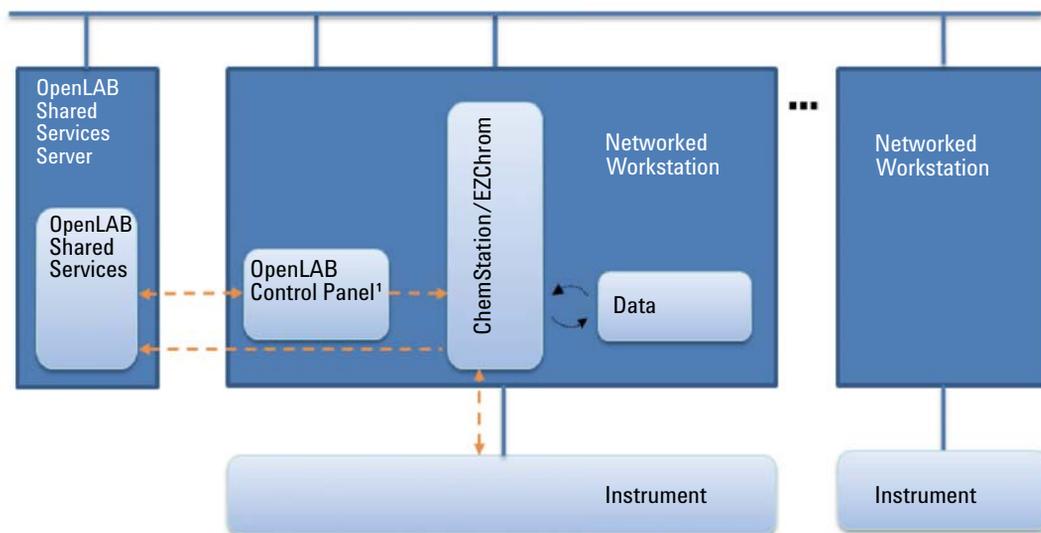
In larger laboratories with many instruments in a network, you can install OpenLAB Shared Services on a separate server that acts as an OpenLAB Shared Services server. On the Networked Workstations, the OpenLAB Control Panel accesses the dedicated OpenLAB Shared Services server. In this scenario, you can access all information provided by OpenLAB Shared Services from any workstation configured as part of the Networked Workstation system. For example, you can see on each workstation which instruments are available and which status (Online, Offline, Error, In Run, Not Ready, etc.) the instruments currently have.

## 1 Architectural Concepts of OpenLAB CDS

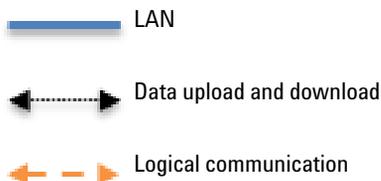
### OpenLAB CDS System Architecture

Since Networked Workstations cannot be remotely controlled, you can launch and configure instruments only from the specific PC on which you configured the instrument.

The following figure shows an OpenLAB CDS Networked Workstation configuration. There can be multiple workstations that are part of the Networked Workstation system. The figure shows only one ChemStation/EZChrom instance, but you can configure multiple instances and associated instruments on the same machine.



**Figure 4** OpenLAB CDS Networked Workstation



You can use the OpenLAB CDS Networked Workstation configuration with or without central data storage. If you are connected to a central repository, the data stored on the workstation is synchronized with the central repository. For more information on ChemStation with central data storage, refer to the *OpenLAB CDS ChemStation Edition with Central Data Storage - User's Guide*.

In case of OpenLAB Data Store systems, OpenLAB Data Store and OpenLAB Shared Services are installed on one server. In case of OpenLAB ECM systems, separate OpenLAB Shared Services servers are required per ECM Account.

With EZChrom Edition, you can also use a network file share (the Enterprise Path) for storing your data. ChemStation Edition does not support networked file shares.

### OpenLAB Data Analysis with Networked Workstations

With OpenLAB Data Analysis, you can process and analyze the data acquired with ChemStation or EZChrom. OpenLAB Data Analysis is supported on Networked Workstations under specific conditions (see [Table 3](#) on page 13). For more information on OpenLAB Data Analysis, refer to the *OpenLAB Data Analysis - Getting Started* guide.

**Table 3** Supported scenarios

Storage type	CDS Edition	OpenLAB Data Analysis supported
Local file system	ChemStation	Yes
	EZChrom	Yes
Network file share	ChemStation	No
	EZChrom (Enterprise Path without Advanced File Security)	Yes
ECM	ChemStation	No
	EZChrom	No
Data Store	ChemStation	No
	EZChrom	No

## Distributed Systems

With OpenLAB CDS installed as a Distributed System, you are able to access and run instruments from any client PC in the system.

As in the Networked Workstation installation, OpenLAB Shared Services provide an overview of all instruments in the system. You can access all information provided by OpenLAB Shared Services from any OpenLAB CDS Client. For example, you can see which instruments are available in which location and which status (Online, Offline, Error, In Run, Not Ready, etc.) the instruments currently have.

In contrast to the Networked Workstation installation, in a Distributed System you can configure, launch and control any instrument from any OpenLAB CDS Client PC.

The Distributed System configuration enables you to work more flexibly with instrument sessions. You can, for example, launch an online instrument, start a sequence, and then disconnect only the CDS Client while the instrument continues to run on the Agilent Instrument Controller machine (AIC). You or another user can later connect to this instrument again from a different OpenLAB CDS Client to finish work on the online instrument or to analyze the data.

### OpenLAB Data Analysis in Distributed Systems

With OpenLAB Data Analysis, you can process and analyze the data acquired with ChemStation or EZChrom. OpenLAB Data Analysis is supported in a Distributed System under specific conditions (see [Table 4](#) on page 15). For more information on OpenLAB Data Analysis, refer to the *OpenLAB Data Analysis - Getting Started* guide.

**Table 4** Supported scenarios

<b>Storage type</b>	<b>CDS Edition</b>	<b>OpenLAB Data Analysis supported</b>
Local file system	ChemStation	Yes
	EZChrom	Yes
Network file share	ChemStation	No
	EZChrom (Enterprise Path without Advanced File Security)	Yes
ECM	ChemStation	No
	EZChrom	No
Data Store	ChemStation	No
	EZChrom	No

## ChemStation Architecture

**NOTE**

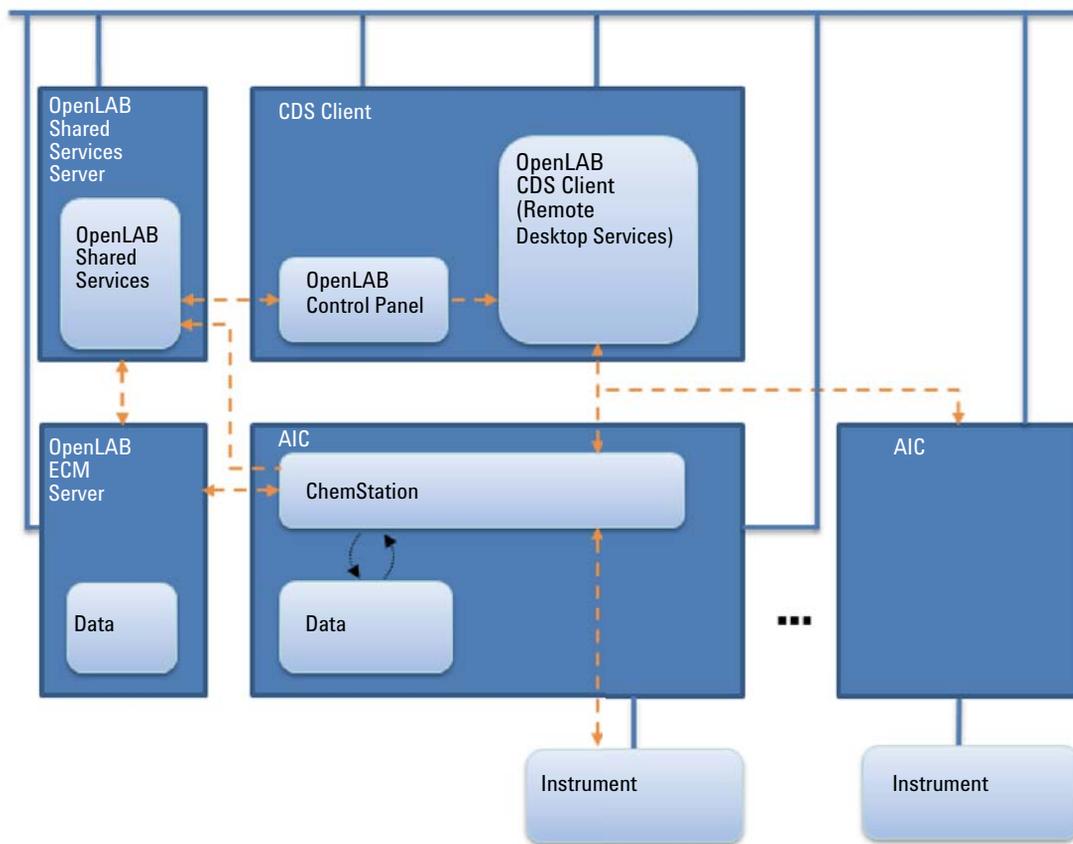
The Distributed System configuration requires either OpenLAB Data Store or OpenLAB ECM.

---

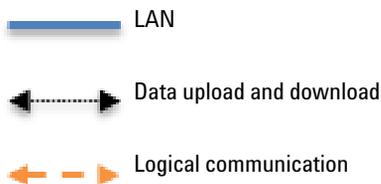
In a ChemStation Distributed System, the instruments are configured on the AIC. The AIC is a Windows server. Each AIC can control up to ten instruments. You can access the ChemStation instance on the AIC machine from any OpenLAB CDS Client. The connection between the CDS Client and the AIC is transparently using Remote Desktop Services. When you disconnect the CDS Client, you disconnect the Remote Desktop Connection. ChemStation continues to run on the AIC. For more information on remote instrument control, session takeover, session disconnect, or force shutdown, refer to the *OpenLAB CDS ChemStation Edition Concepts and Workflows Guide*.

The Distributed System configuration of OpenLAB CDS always includes central data storage. The data stored on each AIC is synchronized with the central repository. For more information on ChemStation with central data storage, refer to the *OpenLAB CDS ChemStation Edition with Central Data Storage User's Guide*.

The following figure shows the system architecture of ChemStation with OpenLAB ECM. In case of OpenLAB Data Store systems, there would be no separate Data Store server, as OpenLAB Data Store and OpenLAB Shared Services are always installed on one server. In case of OpenLAB ECM systems, separate OpenLAB Shared Services servers would be required per ECM Account.



**Figure 5** OpenLAB CDS Distributed System with ChemStation and OpenLAB ECM



## **EZChrom Architecture**

**NOTE**

The Distributed System configuration is supported with Network File Share, OpenLAB Data Store, or OpenLAB ECM.

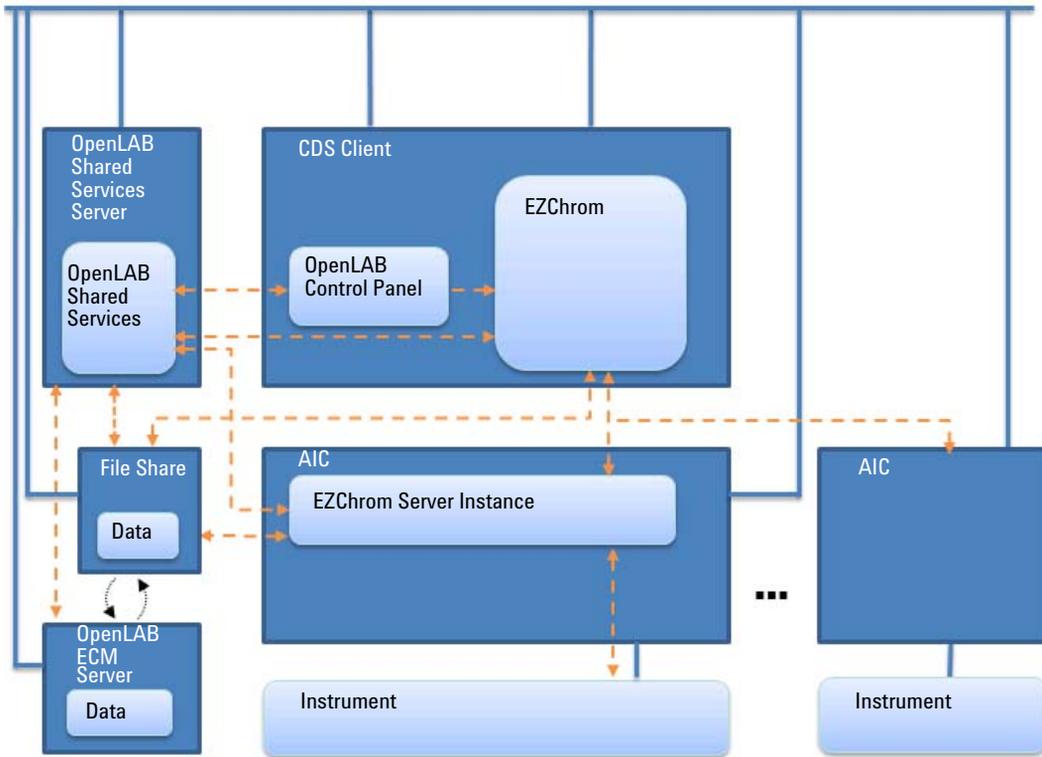
---

In an EZChrom Distributed System, the instruments are configured and controlled from the Client PC through the instrument connection on the AIC. The AIC is a regular PC and can control up to four OpenLAB CDS Clients. You can access the EZChrom server instance on the AIC machine from any OpenLAB CDS Client with EZChrom installed. When you close the EZChrom application on the CDS client, you disconnect from the EZChrom Server instance running on the AIC. The EZChrom server instance continues to run on the AIC until the connection is closed in the OpenLAB Control Panel.

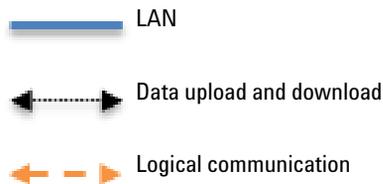
The data is written on a file share. If you use central data storage, the data is synchronized with the central repository. The following figure shows the system architecture of EZChrom with OpenLAB ECM.

In case of OpenLAB Data Store systems, there would be no separate Data Store server, as Data Store is always installed on the OpenLAB Shared Services Server. In case of OpenLAB ECM systems, separate OpenLAB Shared Services servers would be required per ECM Account.

With EZChrom Edition, you can also use a central file-based data storage.



**Figure 6** OpenLAB CDS EZChrom Edition with Distributed System with ECM



## Networked Workstations in Distributed Systems

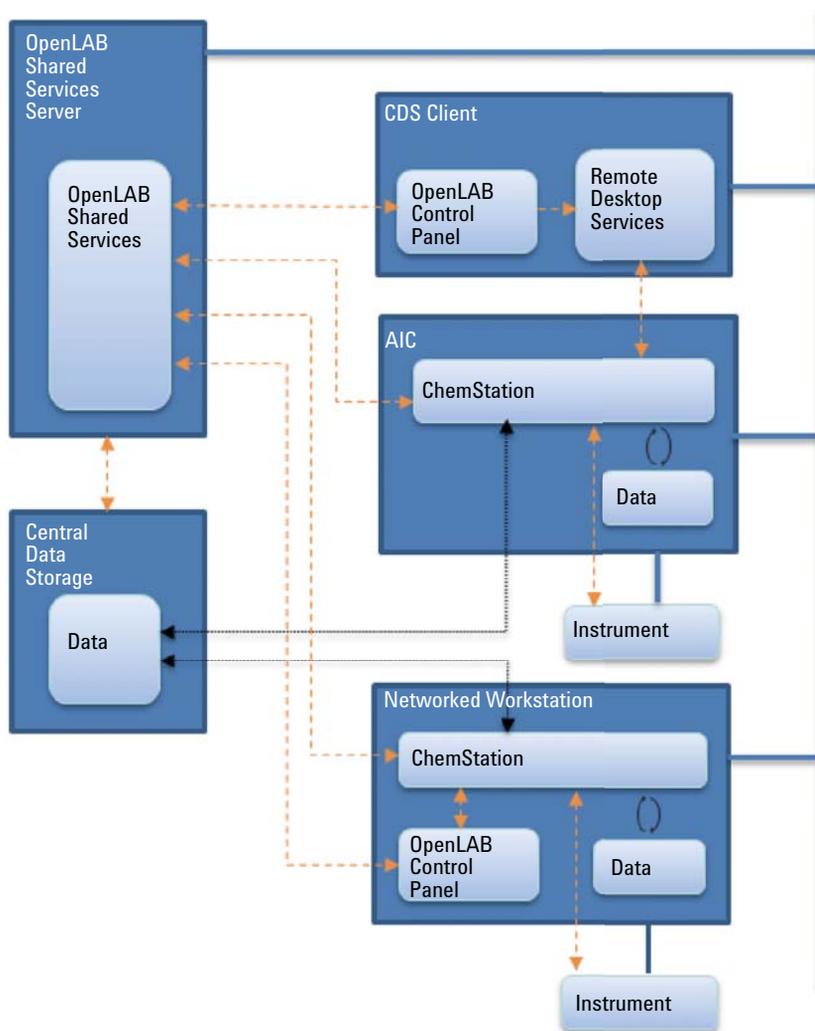
When an instrument cannot be controlled by an AIC or needs to be isolated and controlled by a single computer, you can add one or multiple Networked Workstations to a Distributed environment.

The following figure shows a mixed topology with one Networked Workstation, one CDS Client, one AIC, and central data storage. You can have multiple Networked Workstations, multiple CDS Clients, and multiple AICs in this topology.

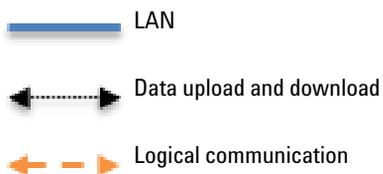
In a mixed environment, you can use the Networked Workstation to access all instruments configured on this workstation, and you can also use it as a CDS Client for all instruments configured on AICs. Instruments configured on a Networked Workstation cannot be controlled from a CDS Client.

On Networked Workstations, OpenLAB CDS Clients, or OpenLAB Shared Services Servers, OpenLAB Data Analysis is only supported with EZChrom Edition; you must use an Enterprise Path without Advanced File Security to store your data.

Systems with both ChemStation Edition and EZChrom Edition being controlled by the same OpenLAB Shared Services Server are not supported.



**Figure 7** OpenLAB CDS mixed topology



## Data Compatibility Between Different Versions

Backward compatibility provided by OpenLAB Shared Services and OpenLAB Data Store was introduced with OpenLAB CDS A.02.01 to allow lab operation (or minimal lab downtime) during an upgrade phase by having the existing and the newer version of the CDS on the same Networked or Distributed environment. The mixed version system is to be used only during the upgrade phase. Procedures must be in place so older clients or workstations are not used to access and reprocess data created by newer clients or workstations.

In the OpenLAB Control Panel, the **Instruments** and **Projects** tabs will be fully functional in the older revisions. The **Administration** tab must not be used with the older revisions, for example: A.02.01, A.01.05 or A.01.04. The OpenLAB Control Panel may be used to start/run instruments that were previously configured and working prior to the upgrade. No other functionality is supported until the system upgrade is completed. Administrative tasks must be performed from an OpenLAB Control Panel running the newer version of the software.

An AIC is configured to run a specific CDS version which is compatible with the instruments drivers connected to the AIC and compatible with the database backend to store the data. The advantage of the CDS solution with AICs is the easy access to the AICs via clients installed somewhere in the lab. Data acquisition, data analysis, and data reprocessing must be used with the compatible version of AIC software and client software.

Typically, newer versions of software are backward compatible with older versions. For this release of OpenLAB CDS, the following specific mixed versions are supported:

- The A.02.02 OpenLAB Shared Services server is compatible with client versions A.02.01, A.01.05, and A.01.04. A hot fix is required for OpenLAB CDS revisions A.01.05 and A.01.04.
- The A.02.02 OpenLAB Data Store server is compatible with the client version A.02.01. A hot fix is required for the OpenLAB CDS revision A.01.05 (or Data Store client A.01.02).

### NOTE

Agilent recommends keeping installations on the same software version. Acquisition and data analysis are supported on the same version of software last used for reprocessing or on a newer version of the software.

## OpenLAB CDS Licensing

This chapter summarizes the main components and features of Agilent OpenLAB CDS, and outlines the licensing strategy for these components and features.

For procedures on how to obtain and install licenses, please refer to the *Software License Installation Guide* (CDS\_LicenseInstallationGuide.pdf on disc 1).

### General Product Structure

OpenLAB CDS has been designed as a combination of software modules for instrument control, data acquisition and data analysis (integration, quantification and reporting), automation and customization.

Technique-specific single instrument products provide control capabilities for a given separation technique, e.g. for a Gas Chromatograph or Liquid Chromatograph. A single instrument configuration may be expanded by additional software modules (Add-ons).

The sections below describe the product modules. For more information on the product licenses, see “[Licensing Scheme](#)” on page 25.

#### Core module

The core module provides data analysis, automation and customization functions (no instrument control) for the following separation techniques:

- Gas chromatography (GC)
- Liquid chromatography (LC)
- Analog data acquisition with external event protocol (A/D)
- ChemStation only: Capillary Electrophoresis (CE)

## Instrument drivers

Technique-specific single instrument drivers provide control capabilities for the separation techniques listed above. By installing multiple instrument drivers, Agilent OpenLAB CDS is capable of controlling more than one analytical system, for example, two LCs, two GCs, or an LC and a GC.

The instrument control capability of OpenLAB CDS may be expanded by purchasing additional modules to allow mixed technique configurations.

## Add-ons

The acquired data is typically two-dimensional ("2D"), that is, by measuring detector response over time. Spectroscopic detectors can deliver three-dimensional ("3D") data, by additionally measuring detector response over a third axis (for example, wavelength or mass range).

For 2D data, OpenLAB Data Analysis provides a single data analysis product for LC and GC data from both ChemStation and EZChrom.

For 3D data, the following modules permit analysis and reporting:

- OpenLAB CDS 3D UV Add-on
- ChemStation only:
  - OpenLAB CDS ChemStation CE 3D MS Add-on
  - OpenLAB CDS ChemStation LC 3D MS Add-on
  - OpenLAB CDS ChemStation LC/MS Deconvolution and Bioanalysis

## License Types

The new licensing strategy introduced with OpenLAB CDS helps you use your licenses more effectively. In contrast to previous ChemStation or EZChrom revisions, licenses for instrument control, drivers and add-ons are *floating licenses*. Any instrument that starts up requests the necessary licenses from license management, and when the instrument is closed, it returns the licenses. Therefore, you only need licenses to cover the maximum number of concurrently running instruments rather than one for each installed instrument. License management is part of the OpenLAB Shared Services.

Licenses are of two types:

- *Counted licenses* are consumed once for each associated software or instrument module.
- *Shared licenses* can be shared per PC or instrument. For example, the Agilent OpenLAB CDS core license is a shared license, which means you need only one license per PC, no matter how many ChemStation or EZChrom instances you run on it.

There is a 60-day Startup License for the entire OpenLAB CDS installation. The expiration period starts with the first launch of an application.

## Licensing Scheme

Figure 8 on page 27 shows the licenses for OpenLAB CDS in the different installation scenarios:

- OpenLAB CDS Workstation
  - One OpenLAB CDS core license
  - Instrument and Add-on licenses as needed; you can run up to four 2D LC or GC instruments on the same workstation.
  - A license for the new OpenLAB Data Analysis is automatically included. The installation is optional.
- Networked Workstation
  - One OpenLAB Shared Services server license
  - One OpenLAB CDS core license for each Networked Workstation; you can connect multiple Networked Workstations to the OpenLAB Shared Services server.
  - Instrument and Add-on licenses as needed; you can run up to four 2D LC or GC instruments on the same Networked Workstation.

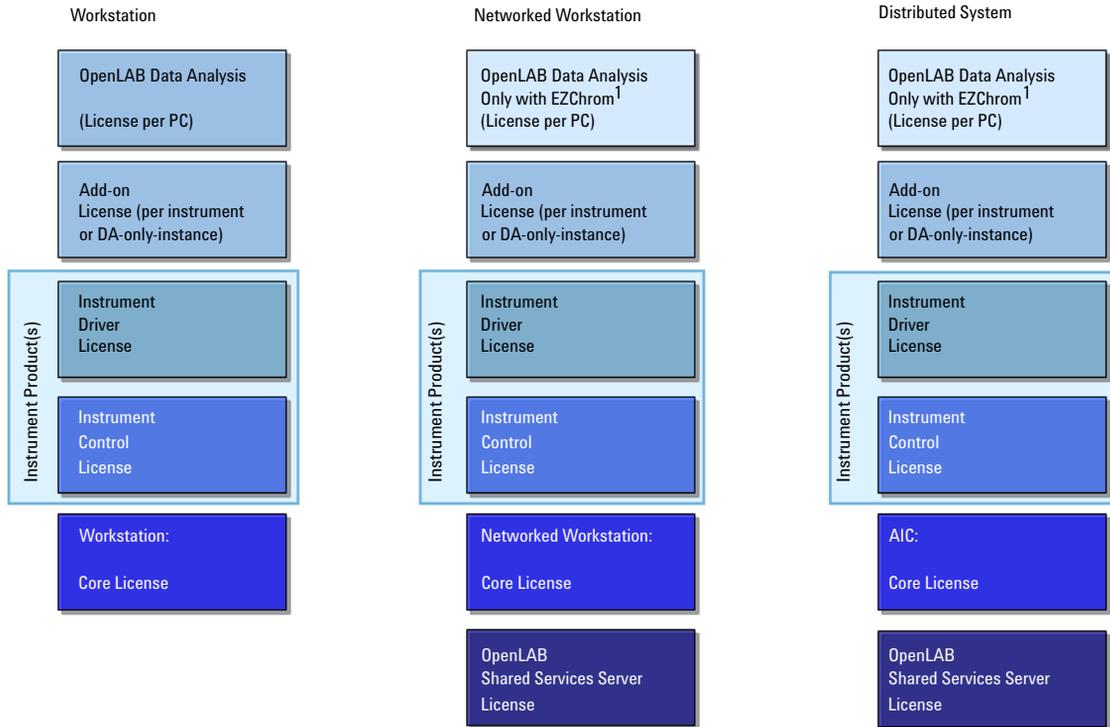
## 1 Architectural Concepts of OpenLAB CDS

### OpenLAB CDS Licensing

- Distributed System
  - One OpenLAB Shared Services server license
  - One OpenLAB CDS core license and AIC-Add-on license for each Agilent Instrument Controller (AIC) machine. You need one AIC license for each AIC connected to the Shared Services server.
  - Instrument and Add-on licenses as needed; you can run up to ten 2D LC or GC instruments on the same AIC.

To control instruments from vendors other than Agilent, an Agilent Instrument Control License is required in addition to the driver license for the other vendor's instrument. A list of driver licenses that are available for instruments from other vendors can be found in the *Supported Instruments and Firmware Guide*.

For Agilent instruments, Agilent Instrument Control licenses and Agilent Driver licenses are always bundled together. They are shown as one product license in the Shared Services License Management. You can see these items as separate lines in the license file itself only.



<sup>1</sup> OpenLAB Data Analysis is supported with an OpenLAB CDS Workstation, or with OpenLAB CDS EZChrom Edition systems using an Enterprise Path without Advanced File Security.

**Figure 8** Licensing schema

## Main Features for Licensing

The following table shows the main features that have to be licensed. If you purchase an Agilent product, the licenses for several features are already included by default. The following tables show the license features associated with Agilent products.

**Table 5** Main license features

License Feature	License Type	Required For	Comment
<b>AgilentOpenLABCDSCHEMStation</b>	Shared per PC	All ChemStation instances	ChemStation full core license Supports up to four instruments with the LC full driver or GC full driver package (including CE, ADC, CE/MS or LC/MS).
<b>AgilentOpenLABCDSEZChrom</b>	Shared per PC	All EZChrom instances	EZChrom full core license Supports up to four instruments with the LC full driver or GC full driver package.
<b>AgilentOpenLABCDSCHEMStationVL</b>	Shared per PC	1220/1260 Infinity LC and 7820 GC	ChemStation VL core license Sufficient to control one 1220/1260 Infinity LC (selected modules only), 7820 GC, 490 micro GC or 790 micro GC.
<b>AgilentOpenLABCDSEZChromVL</b>	Shared per PC	1220/1260 Infinity LC and 7820 GC	EZChrom VL core license Sufficient to control one 1220/1260 Infinity LC (selected modules only), 7820 GC, 490 micro GC or 790 micro GC.
<b>AgilentOpenLABCDSEECOMPACT</b>	Shared per PC	Compact LC, Compact GC	EZChrom compact core license. Sufficient to control up to two compact instruments (1220, 6820/7820, and 490 Micro GC).
<b>AgilentOpenLABCDSAICAddOn</b>	Shared per PC	All AICs with ChemStation or EZChrom instances	This feature turns a core license into an AIC-core license.
<b>AgilentOpenLABCDSCSDAOnly</b>	Shared per PC	ChemStation "classic" Data Analysis	License for the "classic" data analysis in OpenLAB CDS ChemStation Edition (required for "DA Only" workstations).
<b>AgilentOpenLABDataAnalysis</b>	Shared per PC	New OpenLAB Data Analysis	Core license for the new OpenLAB Data Analysis.

**Table 6** Shared Services license features

License Feature	License Type	Required For	Comment
<b>AgilentOpenLABSharedServices</b>	Counted	Only for OpenLAB Shared Services running on a separate server	The OpenLAB Control Panel does not need an extra license. Also OpenLAB Shared Services running on an OpenLAB CDS Workstation do not need an extra license.

**Table 7** Driver license features

License Feature	License Type	Required For	Comment
<b>AgilentInstrumentControl</b>	Counted	Only online instances	The license is consumed whether OpenLAB CDS can connect to the instrument or not. The instrument control license comes as part of the driver product.
<b>AgilentDriversLC AgilentDriversGC AgilentDriversADC</b>	Counted	Only online instances	The license is consumed whether OpenLAB CDS can connect to the instrument or not.
<b>AgilentDriversLCVL AgilentDriversGCVL</b>	Counted	Only online instances	Drivers for OpenLAB CDS VL are not available separately. They are always bundled with an instrument or with the OpenLAB CDS VL core license. See <a href="#">“VL License Features”</a> on page 38. These license features can be combined with a OpenLAB CDS VL core license or a full core license.
<b>AgilentLCCoreVL</b>	Counted	Only online instances	This driver comes with the OpenLAB CDS VL version and allows to control one 1260 Infinity Instrument. See <a href="#">“VL License Features”</a> on page 38.
<b>AgilentDriversCE AgilentDriversMS</b>	Counted	Only ChemStation online instances	The license is consumed whether ChemStation can connect to the instrument or not.

## 1 Architectural Concepts of OpenLAB CDS

### OpenLAB CDS Licensing

**Table 7** Driver license features

License Feature	License Type	Required For	Comment
<b>AgilentDriversLCCompact</b> <b>AgilentDriversGCCCompact</b>	Counted	Only EZChrom online instances	Compact drivers are not available separately. They are bundled with the EZChrom compact core license.
<b>VarianCP_4900</b>	Counted	Only online instances	The license is consumed whether OpenLAB CDS can connect to the instrument or not.
Other drivers	Counted	Only online instances	There are a several driver license features available for instruments of vendors other than Agilent. See Appendix for a complete list.

**Table 8** Add-on license features

License Feature	License Type	Required For	Comment
<b>AddOn3DUV</b>	Shared (per instrument)	Online and offline instances (only if 3D option is selected in the Setup Wizard.)	The license is optional. If the license is not available, spectra analysis is disabled in the software.
<b>AgilentAddOnMSDataAnalysis</b>	Shared (per instrument)	ChemStation online and offline instances.	If MS is configured in ChemStation, the MS Data Analysis Add-on is mandatory.
<b>AgilentAddOnMSDeconvolution</b>	Shared (per instrument)	ChemStation online and offline instances (only if Bioanalysis option is selected in the Setup Wizard).	This license is only mandatory for LC/MS with Deconvolution in ChemStation.
<b>AgilentAddOnSystemSuitability</b>	Shared (per instrument)	EZChrom online and offline instances.	This license is bundled with EZChrom core licenses. If more than the bundled licenses are required, you will need to purchase them separately.

## License Features Associated with Agilent OpenLAB CDS Products

When you purchase an Agilent OpenLAB CDS product, several license features are included by default.

**Table 9** License Feature Associated with Agilent OpenLAB CDS Products

Description	License Feature(s) in OpenLAB CDS
OpenLAB CDS EZChrom Edition Workstation	1 x AgilentOpenLABCDSEZChrom 4 x SystemSuitability 1 x AgilentOpenLABDataAnalysis
OpenLAB CDS EZChrom Edition Workstation Right-to-copy	1 x AgilentOpenLABCDSEZChrom 4 x SystemSuitability 1 x AgilentOpenLABDataAnalysis
OpenLAB CDS EZChrom Edition AIC	1 x AgilentOpenLABCDSEZChrom 1 x AgilentOpenLABCDSAICAddon 8 x SystemSuitability 1 x AgilentOpenLABDataAnalysis
OpenLAB CDS EZChrom Compact	1 x AgilentOpenLABCDSEECCompact 2 x AgilentDriversLCCompact 2 x AgilentDriversGCCompact 2 x VarianCP_4900 2 x AgilentInstrumentControl 2 x SystemSuitability 1 x AgilentOpenLABDataAnalysis
Upgrade EZChrom Elite Compact to OpenLAB CDS EZChrom Compact	1 x AgilentOpenLABCDSEECCompact 2 x AgilentDriversLCCompact 2 x AgilentDriversGCCompact 2 x VarianCP_4900 2 x AgilentInstrumentControl 2 x SystemSuitability 1 x AgilentOpenLABDataAnalysis
OpenLAB CDS EZChrom Edition WorkStation VL	1 x AgilentOpenLABCDSEZChromVL 1 x AgilentOpenLABDataAnalysis 1 x AgilentLCCoreVL
Upgrade OpenLAB CDS EZChrom Compact to OpenLAB CDS EZChrom Edition Workstation	1 x AgilentOpenLABCDSEZChrom 2 x SystemSuitability
OpenLAB CDS EZChrom Edition Agilent GC Headspace	1 x AgilentHeadspaceControl

# 1 Architectural Concepts of OpenLAB CDS

## OpenLAB CDS Licensing

**Table 9** License Feature Associated with Agilent OpenLAB CDS Products

<b>Description</b>	<b>License Feature(s) in OpenLAB CDS</b>
OpenLAB CDS EZChrom Edition VICI Valve Control	1 x VICIValveControl
OpenLAB CDS EZChrom Edition PE Nelson Iface Control	1 x PENelsonInterface
OpenLAB CDS EZChrom Edition Hitachi LaChrom Control	1 x HitachiLC
OpenLAB CDS EZChrom Edition PE LC Series 200	1 x PerkinElmerLC
OpenLAB CDS EZChrom Edition Varian CP38/3900, CP-200X	1 x Varian_3800_3900_200x_GC
OpenLAB CDS Instrument Driver for 490uGC and 790 uGC	1 x VarianCP_4900 1 x AgilentInstrumentControl
OpenLAB CDS EZChrom Edition Waters LC Control	1 x WatersLC
OpenLAB CDS EZChrom Edition Hitachi LC LaChrom PDA	1 x HitachiLCPDA
OpenLAB CDS EZChrom Edition GPC/SEC Subsystem	1 x AddOnGPCSEC
OpenLAB CDS EZChrom Edition System Suitability Lic.	1 x SystemSuitability
OpenLAB CDS ChemStation Edition Workstation	1 x AgilentOpenLABCDSChemStation 1 x AgilentOpenLABDataAnalysis
Right-to-copy OpenLAB CDS ChemStation Edition Workstation	1 x AgilentOpenLABCDSChemStation 1 x AgilentOpenLABDataAnalysis
OpenLAB CDS ChemStation Edition Workstation VL	1 x AgilentOpenLABCDSChemStationVL 1 x AgilentOpenLABDataAnalysis 1 x AgilentLCCoreVL
OpenLAB CDS ChemStation AIC	1 x AgilentOpenLABCDSChemStation 1 x AgilentOpenLABCDSAICAddon
OpenLAB CDS Shared Services Server Software	1 x AgilentOpenLABSharedServices
OpenLAB CDS ChemStation Edition LC Upgrade	1 x AgilentOpenLABCDSChemStation 1 x AgilentInstrumentControl 1 x AgilentDriversLC 1 x AddOn3DUV 1 x AgilentOpenLABDataAnalysis
OpenLAB CDS ChemStation Edition GC Upgrade	1 x AgilentOpenLABCDSChemStation 1 x AgilentInstrumentControl 1 x AgilentDriversGC 1 x AgilentOpenLABDataAnalysis

**Table 9** License Feature Associated with Agilent OpenLAB CDS Products

Description	License Feature(s) in OpenLAB CDS
OpenLAB CDS ChemStation Edition CE Upgrade	1 x AgilentOpenLABCDSChemStation 1 x AgilentInstrumentControl 1 x AgilentDriversCE 1 x AddOn3DUV 1 x AgilentOpenLABDataAnalysis
OpenLAB CDS ChemStation Edition LC-MS Upgrade	1 x AgilentOpenLABCDSChemStation 1 x AgilentInstrumentControl 1 x AgilentDriversLC 1 x AddOn3DUV 1 x AgilentDriversMS 1 x AgilentAddOnMSDataAnalysis 1 x AgilentAddOnMSDeconvolution 1 x AgilentOpenLABDataAnalysis
OpenLAB CDS 3D UV Add-on	1 x AddOn3DUV
OpenLAB CDS ChemStation Edition CE 3D MS Add-on	1 x AgilentDriversMS 1 x AgilentAddOnMSDataAnalysis
OpenLAB CDS ChemStation Edition LC 3D MS SQ Add-on	1 x AgilentDriversMS 1 x AgilentAddOnMSDataAnalysis
OpenLAB CDS ChemStation Edition LC/MS Deconvolution Bioanalysis	1 x AgilentAddOnMSDeconvolution
OpenLAB CDS ChemStation Edition LC Dissolution	1 x AgilentAddOnCSLCDissolution
OpenLAB CDS ChemStation Data Analysis Only	1 x AgilentOpenLABCDSDataOnly
OpenLAB Data Analysis	1 x AgilentOpenLABDataAnalysis
OpenLAB CDS Instrument Control License	1 x AgilentInstrumentControl
OpenLAB CDS Instrument Driver for Agilent GC	1 x AgilentInstrumentControl 1 x AgilentDriversGC
OpenLAB CDS Instrument Driver for Agilent A/D	1 x AgilentInstrumentControl 1 x AgilentDriversADC
OpenLAB CDS Instrument Driver for Agilent GC VL	1 x AgilentInstrumentControl 1 x AgilentDriversGCVL
OpenLAB CDS Instrument Driver for Agilent LC	1 x AgilentInstrumentControl 1 x AgilentDriversLC

## 1 Architectural Concepts of OpenLAB CDS

### OpenLAB CDS Licensing

**Table 9** License Feature Associated with Agilent OpenLAB CDS Products

<b>Description</b>	<b>License Feature(s) in OpenLAB CDS</b>
OpenLAB CDS Instrument Driver for Agilent CE	1 x AgilentInstrumentControl 1 x AgilentDriversCE 1 x AddOn3DUV
OpenLAB CDS Instrument Driver for Agilent LC VL	1 x AgilentInstrumentControl 1 x AgilentDriversLCVL
OpenLAB CDS Waters Acquity LC Driver	1 x AgilentDriversWatersAcquity

## Licensing Examples

### Example 1: Workstation (with ChemStation full core license)

**Table 10** Purchased products and associated license features

Quantity	Product	License Feature
1	OpenLAB CDS ChemStation Workstation	AgilentOpenLABCDSChemStation AgilentOpenLABDataAnalysis
1	OpenLAB CDS Instrument Driver for Agilent LC	AgilentInstrumentControl AgilentDriversLC
1	OpenLAB CDS Instrument Driver for Agilent CE	AgilentInstrumentControl AgilentDriversCE AddOn3DUV

A ChemStation set up with the above set of licenses allows running a 2D LC and a CE instrument. If a 3D LC is started, the 3D-UV license required for the CE is consumed and the CE will not start:

- No license is required for OpenLAB Shared Services with a standalone workstation.
- An LC ChemStation with 3D option enabled is started. The following licenses are consumed: 1x OpenLAB CDS ChemStation; 1x Instrument Control; 1x LC Driver; 1x Add-on 3D UV.
- On the same PC, a CE ChemStation should be started. ChemStation would require additionally 1 Instrument Control, 1 CE Driver and 1 Add-on 3D UV. The ChemStation fails to start, because no Add-on 3D UV license is available.
- OpenLAB Data Analysis can be started multiple times. The license is shared per PC.

## Example 2: Networked Workstation

**Table 11** Purchased products and associated license features

Quantity	Product	License Feature
1	OpenLAB CDS Shared Services server	AgilentOpenLABSharedServices
2	OpenLAB CDS EZChrom Workstation	AgilentOpenLABCDSEZChrom SystemSuitability AgilentOpenLABDataAnalysis
2	OpenLAB CDS Instrument Driver for Agilent LC	AgilentInstrumentControl AgilentDriversLC
1	OpenLAB CDS 3D UV Add-on	AddOn3DUV

- The OpenLAB Shared Services server license is consumed when the OpenLAB Shared Services are started on the Shared Services server.
- On PC1, EZChrom with 3D option is started. The following licenses are consumed: 1x OpenLAB CDS EZChrom; 1x Instrument Control; 1x LC Driver; 1x Add-on 3D UV.
- On PC2, a second EZChrom application is started. The 3D option is not available, as the 3D UV Add-on license has already been consumed on PC1. The following licenses are consumed on PC2: 1x OpenLAB CDS EZChrom; 1x Instrument Control; 1x LC Driver.

### Example 3: Distributed System

**Table 12** Purchased products and associated license features

Quantity	Product	License Feature
1	OpenLAB CDS Shared Services server	AgilentOpenLABSharedServices
1	OpenLAB CDS ChemStation AIC	AgilentOpenLABCDSChemStation AgilentOpenLABCDSAICAddon
6	OpenLAB CDS Instrument Driver for Agilent GC	AgilentInstrumentControl AgilentDriversGC
2	OpenLAB CDS Instrument Driver for Agilent LC	AgilentInstrumentControl AgilentDriversLC
2	Add-on	AddOn3DUV

- The OpenLAB CDS Shared Services server license is consumed when the OpenLAB Shared Services are started on the Shared Services server.
- On an AIC machine, six GC ChemStations are remotely started. The following licenses are consumed: 1x OpenLAB CDS ChemStation; 1x AIC Add-on; 6x Instrument Control; 6x GCDriver.
- On the same AIC machine, two LC ChemStations with 3D option enabled are remotely started. No additional ChemStation or AIC Add-on licenses are required. The following licenses are consumed: 2x Instrument Control; 2x LC Driver; 2x Add-on 3D UV.

## VL License Features

With an OpenLAB CDS VL core license, you can control one LC or GC instrument without purchasing additional instrument driver licenses.

OpenLAB CDS VL licenses are supported only for standalone workstation scenarios. They do not run on Networked Workstations or Distributed Systems. LC-MS configurations are not possible with an OpenLAB CDS VL system, neither is an integration with OpenLAB ECM.

The 3D UV Add-on always requires the full AddOn3DUV (M8360AA) license.

Driver licenses for OpenLAB CDS VL are not available separately. They are either bundled with the respective instruments or with the OpenLAB CDS VL core license. Depending on the instrument type chosen when creating the instrument the following applies:

- *Agilent LC Core*: Driver bundled with the OpenLAB CDS VL core license. You can run one full 1260 Infinity LC Instrument with selected modules (for details, see *Supported Instruments and Firmware Guide*).
- *Agilent 1220 LC System*: Driver bundled with 1220 Infinity LC instruments. You can run one Agilent 1220 Infinity LC system, including individual Agilent LC modules except modular pumps (for details, see *Supported Instruments and Firmware Guide*).
- *Agilent 7820 GC system*: Driver bundled with the 7820 GC instrument

### Example: Pure VL configuration

This example applies to ChemStation as well.

**Table 13** Purchased products and associated license features

Quantity	Product	License Feature
1	OpenLAB CDS EZChrom Workstation VL	AgilentOpenLABCDSEZChromVL
1	Agilent 1220 LC	AgilentInstrumentControl AgilentDriversLCVL
1	Agilent 7820 GC	AgilentInstrumentControl AgilentDriversGCVL
1	OpenLAB CDS 3D UV Add-on	AddOn3DUV

- No license is required for OpenLAB Shared Services with a standalone workstation.
- An Agilent 1220 LC with 3D option enabled is started. The following licenses are consumed: 1x EZChrom VL; 1x Instrument Control; 1x 1220 LC Driver for OpenLAB CDS VL; 1x Add-on 3D UV.
- On the same PC, an Agilent 7820 GC is started. As the EZChrom core license is shared, no additional EZChrom core license is required. The following licenses are consumed: 1x Instrument Control; 1x 7820 GC Driver for OpenLAB CDS VL.

## Flexera License Manager

OpenLAB Shared Services use a 3rd party tool called *FlexNet Producer Suite* from Flexera to manage the licenses. The required components are installed by default together with OpenLAB Shared Services. The license server can be the local PC, a remote OpenLAB Shared Services server, or a server with an already existing Flexera license manager in your environment. If you use an existing Flexera license manager, you can provide the hostname or IP address of the license server in the OpenLAB Control Panel.

License Management in OpenLAB Shared Services requires an additional Windows service to be running. This Windows service is called *Agilent*

*OpenLAB License Server.* This service must be running on the server where you manage your licenses. Each time you start an instrument, the instrument requests licenses from the License Server service; therefore, you can only start an instrument if this service is running.

## License Upgrade

If you are upgrading from an older version, 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 which are upgraded to the new core software version, without a new workstation license, will not work until the new workstation licenses are added to the OpenLAB Control Panel.

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

- 1** Log into the Agilent Electronic Software and License Delivery (<https://agilent.subscribenet.com/>).
- 2** Navigate to **Manage Licenses by Host**, and search for the machine hosting your workstation. Alternatively, select the machine from the drop-down list.
- 3** Click **Upgrade All** to upgrade all licenses on this host to the most recent version. Confirm the upgrade on the next screen.

This upgrades the license file to the most current version. SubscribeNet will send you an email with a new license file.

- 4** Put the new license file on your system.

If you have multiple standalone Workstations, repeat this step for each individual workstation.

If you use a Networked System, you only need to provide the license file once, as licenses are managed on the central OpenLAB Shared Services server.

## Windows Server Licensing

OpenLAB CDS ChemStation Edition Distributed Systems are taking advantage of Microsoft Remote Desktop Services. In addition to the Client Access Licenses, this technology requires additional licenses called Remote Desktop Services Client Access License (RDS CAL).

### Client Access License (CAL)

A Client Access License permits client computers to connect to Microsoft server hosting shared resources.

### Remote Desktop Services CAL (RDS CAL)

The OpenLAB CDS ChemStation Instrument Controller (AIC) software uses the Remote Desktop Services (RDS) role in Windows Server. Using this role requires a Windows Server Remote Desktop Services CAL (RDS CAL), in addition to the Windows Server Client Access License (CAL), to access any application or graphical user interface remotely hosted by Windows Server.

You need at least one Remote Desktop Services Licensing Server deployed and activated in your environment. During a grace period of 120 days no license server is required. At the end of the grace period, remote connections will be refused.

Both the Client Access License (CAL) and Remote Desktop Services CAL (RDS CAL) can apply to a device or to a user.

For more information about CAL and RDS CAL requirements, see:

- Client Access Licenses and Management Licenses  
(<http://www.microsoft.com/licensing/about-licensing/client-access-license.aspx>)
- Understanding Remote Desktop Licensing  
(<http://technet.microsoft.com/en-us/library/cc772298.aspx>)
- Licensing Windows Server 2012 R2 Remote Desktop Services  
([http://download.microsoft.com/download/3/D/4/3D42BDC2-6725-4B29-B75A-A5B04179958B/WindowsServerRDS\\_VLBrief.pdf](http://download.microsoft.com/download/3/D/4/3D42BDC2-6725-4B29-B75A-A5B04179958B/WindowsServerRDS_VLBrief.pdf))

## Security and Data Integrity

This chapter explains the built-in security and how it complies with the FDA 21 CFR Part 11. It also explains the system security features provided by OpenLAB Shared Services.

### Security Aspects

In OpenLAB CDS, security aspects are mainly controlled by OpenLAB Shared Services. In addition, some aspects that are only relevant for ChemStation are controlled by the ChemStation Administration Tool.

The OpenLAB Shared Services functionality related to security includes the following (see [“OpenLAB Control Panel”](#) on page 45 for details):

- System Activity Log
- Selection of authentication provider
- User Management
- Security Policy

### Data Integrity

The result data is stored either locally or in a central data repository, depending on the installed OpenLAB CDS configuration. If you store the data in the local file system, you must manually back up the data. With ChemStation, you can reach full compliance with 21 CFR Part 11 together with OpenLAB ECM or OpenLAB Data Store only. These systems store data in compliance with 21 CFR Part 11. They provide secure data storage with access control and an audit trail. Data files are versioned to ensure data integrity and traceability. In addition, OpenLAB ECM and OpenLAB Data Store provide electronic signatures allowing users to sign off on data. The systems can be configured to automatically back-up and archive data on a regular basis.

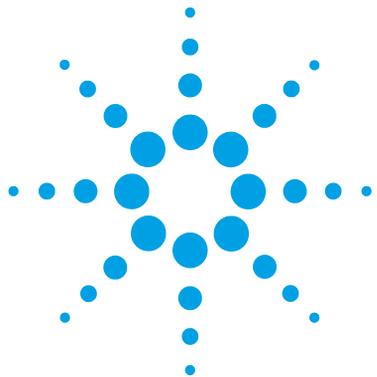
For more information on ChemStation with central data storage, refer to the *OpenLAB CDS ChemStation Edition with Central Data Storage User's Guide*.

With EZChrom, you can also use the central data storage system as described above. In addition, with EZChrom networked systems you can also use a network file share for storing your data. In this case, you must configure Advanced File Security (AFS) to be compliant with 21 CFR Part 11. AFS provides enhanced security on the enterprise path in order to prevent any unauthorized access to project data outside of the data system. This configuration sets the appropriate Windows sharing and security settings to allow only a defined group to access the enterprise data from Windows Explorer. This may only be configured if your system uses Windows Domain as the Shared Services authentication provider.

For more information on configuring AFS, refer to the *Networked and Distributed System Installation and Configuration* guide.

# 1 Architectural Concepts of OpenLAB CDS

## Security and Data Integrity



## 2 OpenLAB Control Panel

Project Management	46
Instrument Management / Lab Status at a Glance	46
License Management	48
System Activity Log	50
Diagnostics	51
Administrative Reports	52
Authentication Provider	53
Setting the Authentication Provider and Storage Location	54
Security Policy	56
User Management	58
Users	58
Groups	60
Roles and Privileges	60
Specific Roles for Individual Instruments or Projects	62

Using the OpenLAB Control Panel, you can access OpenLAB Shared Services control features such as security policy, central configuration, or lab status at a glance. These features are described in more detail in this chapter.



## Project Management

A project is the set of directories that store related methods, data, sequences, and templates. Individual access rights can be applied to each project.

Project Groups can be used to arrange projects to match your organization.

The Projects pane is automatically available when the Control Panel is used with EZChrom Edition. The Projects pane is not automatically available when it is used with ChemStation Edition.

## Instrument Management / Lab Status at a Glance

The **Instruments** view in OpenLAB 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 (including color-coded state, current run, and run status)

The status is stored and constantly updated in the OpenLAB Shared Services database in which the instrument is stored. It can be monitored by every OpenLAB Control Panel connected the corresponding server.

- Instrument Name
- Instrument location
- Instrument type
- Last change of configuration

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

You can create a tree of different locations in the OpenLAB 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.

The type of instrument is preconfigured. When you install OpenLAB CDS, you can choose between ChemStation Edition or EZChrom Edition. The selected edition determines the type of instrument you can use in the Instrument Management.

**NOTE**

Mixed scenarios with ChemStation and EZChrom instruments are not supported with OpenLAB CDS.

---

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 on the local PC or AIC, but you access the configuration tool from the OpenLAB Control Panel.

- Launch the instrument online or offline.

**Workstation and Networked Workstation:** As the instrument configuration is stored on the local PC, you only launch instruments that are configured on this PC.

**Distributed System:** As the instrument configuration is stored on an AIC, you can launch all 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 62).

## License Management

This service includes the administration of all licenses that are required for your instrument modules and Add-ons. When you start an instrument, OpenLAB CDS automatically checks whether the required licenses are available in the license pool, and reserves the licenses needed to operate the instrument. When you stop the instrument, the freed licenses can be used by other instruments.

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, refer to the *Software License Installation Guide*.

License Management in OpenLAB 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 get a license.

For more information on adding license files and viewing the license properties, refer to the OpenLAB Control Panel online help.

The following properties are shown for installed licenses:

- **Feature:** This indicates the type of license used, for example, AgilentOpenLABCDSChemStation, AgilentInstrumentControl, or AgilentDriversLC.
- **Version:** If a license is versioned, you can see the version number, for example 1.1 for Agilent OpenLAB CDS C.01.01. 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 24).
- **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.

## System Activity Log

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

The following types of events are recorded:

- System
- Instrument Management
- Instrument
- Project Management (only applicable to EZChrom)
- 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. ChemStation records these events in its own environment but also sends the events to the System Activity Log. 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.

### NOTE

By default, activity logging is disabled. To enable it in OpenLAB Control Panel, you must have the **Edit activity log properties** privilege. Once enabled, activity logging cannot be disabled again.

## Diagnostics

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

- Ping the OpenLAB Shared Services server.
- Create a report, either for the local system or for the OpenLAB Shared Services 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 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 (EZChrom Only)**

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.

### **Users and Groups Report**

This report provides an overview of all users and groups access rights to instruments and projects on the system. Note that users and groups that have not been granted access to instruments or projects are not included in this report.

## Authentication Provider

Authentication providers are used to prove the identity of users that log in to the system. OpenLAB Shared Services support the following Authentication providers:

- **None**

In this mode, no login screen is shown when you access the OpenLAB Control Panel. The user is automatically logged in to the application with security disabled. All log entries record the user as "Anonymous". With the authentication provider **None**, the Security Policy and User Management nodes are unavailable in OpenLAB Control Panel.

**NOTE**

With the authentication provider **None**, any activity logs will display a generic **System** operator with no additional identification. This is not recommended for compliant setups.

- **Internal**

In this mode, the user's credentials are stored in the OpenLAB Shared Services database. You are asked to create an administrator account for OpenLAB Shared Services 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 Local** or **Windows Domain**

You import existing Windows users into OpenLAB Shared Services. The authentication is done either by a local Windows user management, Windows Active Directory domain or NT 4.0 Domain within the Enterprise. OpenLAB Shared Services only use the identity and password of the mapped users; roles and privileges for OpenLAB CDS are still configured with OpenLAB Shared Services.

- **ECM**

In this mode, an OpenLAB ECM system is responsible for authentication. When you start the OpenLAB Control Panel, the application will prompt for ECM credentials to validate a user. You must choose an existing ECM user as an administrator for OpenLAB Shared Services. The Search function helps you to find specific ECM users. OpenLAB Shared Services only use the identity and password of the mapped users; roles and privileges for OpenLAB CDS are still configured with OpenLAB Shared Services.

## Setting the Authentication Provider and Storage Location

- 1 Open OpenLAB Control Panel from the OpenLAB Control Panel shortcut on the desktop, or go to **Start > All Programs > Agilent Technologies > OpenLAB > OpenLAB Control Panel**.
- 2 From the navigation pane, select **Administration > System Configuration**.
- 3 In the **System Configuration** toolbar, select **Edit System Settings**.

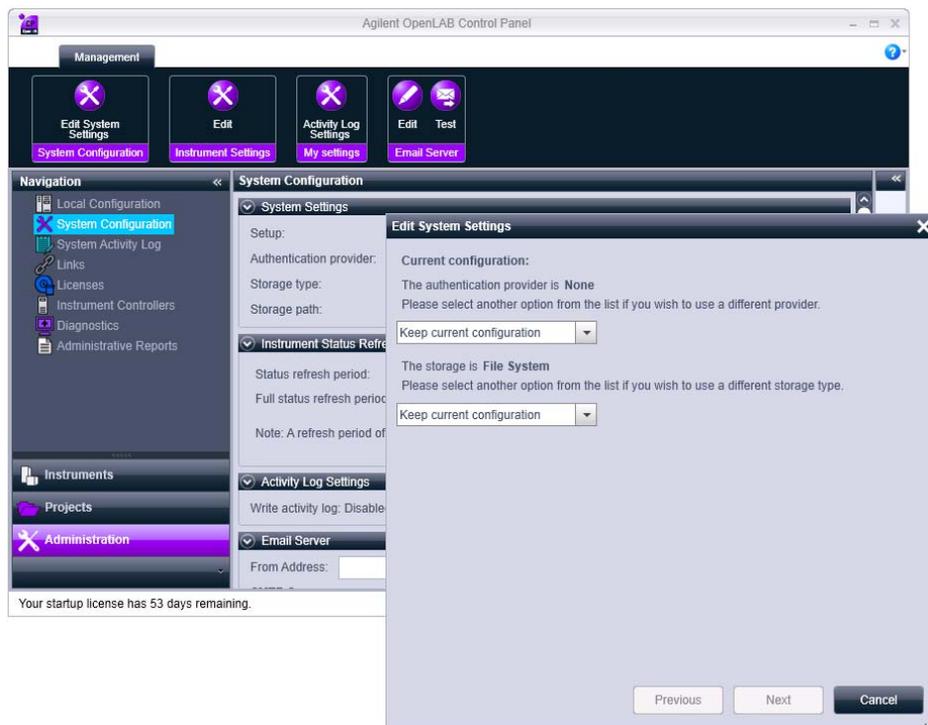


Figure 9 Edit System Settings dialog in the OpenLAB Control Panel

- 4 In the **Edit System Settings** window, select your authentication provider from the drop-down list.

If you installed a Data Store server, the **Internal** authentication provider is automatically configured and a user called 'admin' with password 'openlab' is created during the installation of the Data Store server. If required, you can change the authentication provider to **Windows domain**.

- 5 Select your storage provider from the drop-down list.  
The storage type **Data Store** is available only if Data Store has been deployed.
- 6 Select **Next**.
- 7 Select a user to administrate the system.
  - a If you selected **Internal** as an authentication provider:
    - Select **Create Account**.
    - In the **Create Administrator Account** dialog box, enter a **Name** and a **Password**.
  - b If you selected **Windows Local** as an authentication provider:
    - Select **Select Account**.
    - Enter a search string.
    - In the **Search Users** dialog box, select **Search** to view a list of users.
    - Select a user.
  - c If you selected **Windows Domain** as an authentication provider:
    - Select the check box to activate the input fields.
    - Enter a **Domain**, **User**, and **Password**.
    - Select **Select Account**.
    - Enter a search string.
    - In the **Search Users** dialog box, select **Search** to view a list of users.
    - Select a user.
  - d If you selected **ECM** as an authentication provider:
    - Provide the **ECM Server URL** and enter the ECM user credentials.
    - Select **Select Account**.
    - Enter a search string.
    - In the **Search Users** dialog box, select **Search** to view a list of users.
    - Select a user.
- 8 Select **OK**.
- 9 Select **Next**.
- 10 Review your settings and select **Apply**.

## Security Policy

The Security Policy is only available if you select an authentication provider other than **None**.

With the authentication provider **Internal**, you can set all of the parameters described below in the OpenLAB Control Panel. With an external authentication provider (e.g. Windows Domain), you can only set the inactivity time in the OpenLAB Control Panel; all other parameters are defined by the external system.

For more information on 21 CFR Part 11 requirements for ChemStation, refer to the *OpenLAB CDS ChemStation Edition with Central Data Storage User's Guide*.

**Table 14** Security Policy settings

Setting	Description	21 CFR Part 11 Requirements
<b>Minimum password length</b>	If users change their passwords, they must choose a password with at least the given number of characters. The default setting is 5. Only available for authentication provider <b>Internal</b> .	You should require a minimum password length of at least 5 characters.
<b>Password expiration period (days)</b>	The default value is 30 days. When the user tries to log in after this period of time, the system will ask him to change the password. The expiration period starts with the last password change or with the creation of a user with a new default password. Only available for authentication provider <b>Internal</b> .	You should use an expiration period of 180 days or less.
<b>Maximum unsuccessful login attempts before locking account</b>	If a user tries to log in with invalid user credentials a number of times, the user is locked out of the system for a certain period of time ( <b>Account lock time</b> , see below). Login is impossible, even with valid user credentials. You can define the number of allowed login attempts. The default setting is 3. Only available for authentication provider <b>Internal</b> .	You should limit the number of allowed login attempts to three.

**Table 14** Security Policy settings

Setting	Description	21 CFR Part 11 Requirements
<b>Account lock time (minutes)</b>	Once a user has exceeded the maximum number of allowed unsuccessful login attempts, this is the amount of time that must pass before he can try again. The default setting is <i>5 min</i> . Only available for authentication provider <b>Internal</b> .	
<b>Inactivity time before locking the application</b>	If the OpenLAB Control Panel is inactive for this amount of time, the user interface will be locked. This setting is also used to set the time-based session lock in ChemStation. The default setting is <i>10 min</i> . Set the value to zero to never lock.	
<b>Single Sign-On</b>	With Single Sign-On enabled, the user will not see the OpenLAB Control Panel login screen. Only available for authentication provider <b>Windows Domain</b> .	

## User Management

OpenLAB Shared Services allow you to assign specific roles to users or user groups. If you manage your users within an external system (for example, OpenLAB ECM or a Windows domain), you can map those existing users into OpenLAB Shared Services.

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.

The roles are equipped with numerous specific privileges which define what the users are allowed to view or do in OpenLAB Control Panel and in OpenLAB CDS.

## Users

The following information is required if you create a new internal user:

**Table 15** User Credentials

Value	Description	Mandatory
<b>Name</b>	Username to log in to the system.	Yes
<b>NOTE</b> The following characters are not allowed as part of a username: < > : " / \   % * ? ' °		
<b>Description</b>	Additional information about the user (e.g. department, function etc.)	No
<b>Password</b>	Password for the user; minimum password length is defined in the Security Policy.	Yes
<b>Email address</b>	Email address of the user.	No
<b>Full name</b>	The full (long) name of the user.	No

**Table 15** User Credentials

Value	Description	Mandatory
<b>Contact Information</b>	General contact information (e.g. telephone number, pager etc.)	No
<b>Password never expires</b>	The password will never expire, even if a password expiration period is set in the security policy.	No
<b>Account is disabled</b>	Select the check box to disable an account. Users with disabled accounts cannot log in any more. Accounts may be automatically disabled after too many failed login attempts. If an account is disabled, a corresponding message is displayed instead of the check box. After a given time (see <b>Account lock time</b> in the <b>Security Policy</b> settings), the account is automatically enabled again.	No
<b>User cannot change password</b>	Flag that indicates whether the user can change his own password. The flag is false by default (that is, users CAN change their passwords).	No
<b>User must change password at next login</b>	If set to true, the user has to change his password at the next login. The flag is automatically set to false after the user has changed the password successfully. The flag is true by default for new users.	No
<b>Group Membership</b>	Assign the user to the relevant groups.	
<b>Role Membership</b>	Assign roles directly to the user.	

If you use an external authentication provider (Windows domain), you cannot create new users, but must import users that exist in the authentication systems. A search function helps you find specific users in the authentication system. In the Control Panel, you can manage the roles for those external users, but not the actual user credentials such as user name and password. If you want to remove an external user, you unmap the user in the Control Panel. The user continues to exist in the external authentication system.

## Groups

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 OpenLAB Control Panel. If you need additional user assignments that are relevant only for OpenLAB CDS, create them in the OpenLAB Control Panel. Otherwise it is sufficient to only import the groups and assign the required roles to the groups.

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 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 *ChemStation Operator* will always need another role such as *Instrument User* with the privilege to run an instrument.

You can create a tree of different locations in the OpenLAB 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 62). For example, a user can have the role **Instrument Administrator** for one instrument, and **Instrument User** for another instrument.

With EZChrom or with Data Store, you can also create a tree of different projects or project groups in the OpenLAB Control Panel, and assign different Project roles for different projects (see also “[Specific Roles for Individual Instruments or Projects](#)” on page 62). For example, a user can have the role **Project Administrator** in one project, so that he can manage the settings in the OpenLAB 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.

**Table 16** 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, Manage security, Manage printers</b> etc.
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 OpenLAB Control Panel.
Project privileges	Privileges for accessing or modifying different levels of data. <ul style="list-style-type: none"><li>• With EZChrom, you can assign these privileges on project level.</li><li>• With ChemStation, these privileges are globally assigned to all ChemStation users</li></ul>

For a detailed list of privileges, see the Appendix.

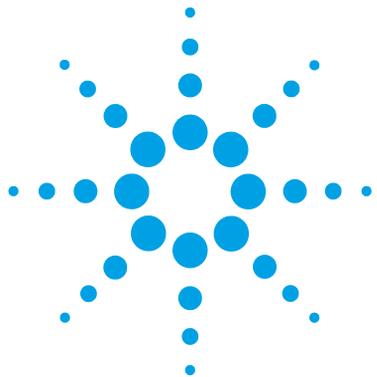
## Specific Roles for Individual Instruments or Projects

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 node, you can deselect the **Inherit privileges from parent** check box in the **Edit Privileges** dialog for the required node. 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.



## 3 OpenLAB Shared Services Server

OpenLAB Server Utility Tool	64
Exporting and Purging Log Entries	65
Using Backup and Restore Functions	66
Configuring Access to Windows Domain	69
Managing Server Settings	70
Maintenance Procedures	71
Update Database Statistics	71
Procedures for PostgreSQL Database	71
Procedures for SQL Server	72
Monitor Resource Usage on OpenLAB Shared Services Server	73
Purging the Activity Logs	73
Domain or Server Name Change	73
Additional Best Practices	74

This chapter describes the OpenLAB Server Utility Tool.



## OpenLAB Server Utility Tool

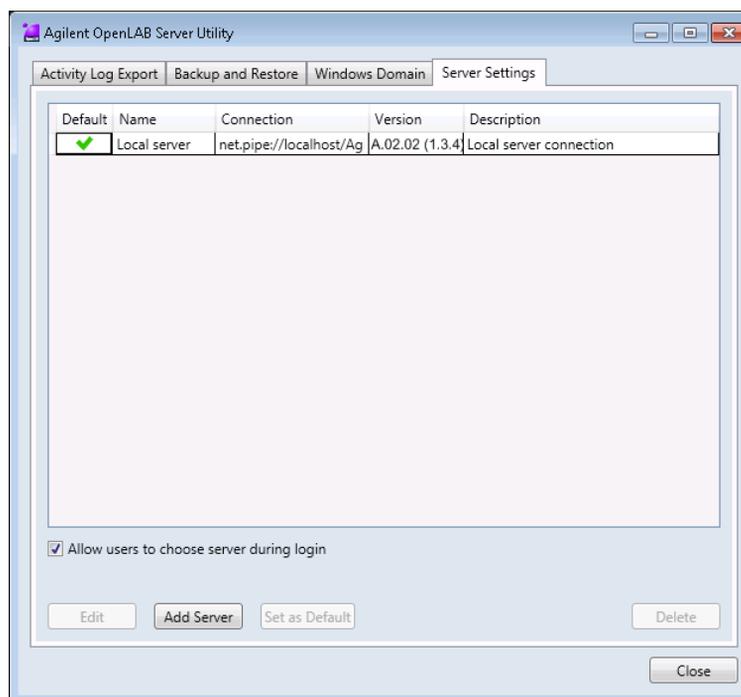
The Agilent OpenLAB Server Utility Tool helps managing the system. It is automatically installed with your OpenLAB software.

The settings made in this utility affect all users of workstations, AICs, or servers.

### NOTE

The OpenLAB Server Utility Tool can only be started by administrators.

- 1 To start the application, go to **Start > All Programs > Agilent Technologies > OpenLAB > OpenLAB Server Utility**.

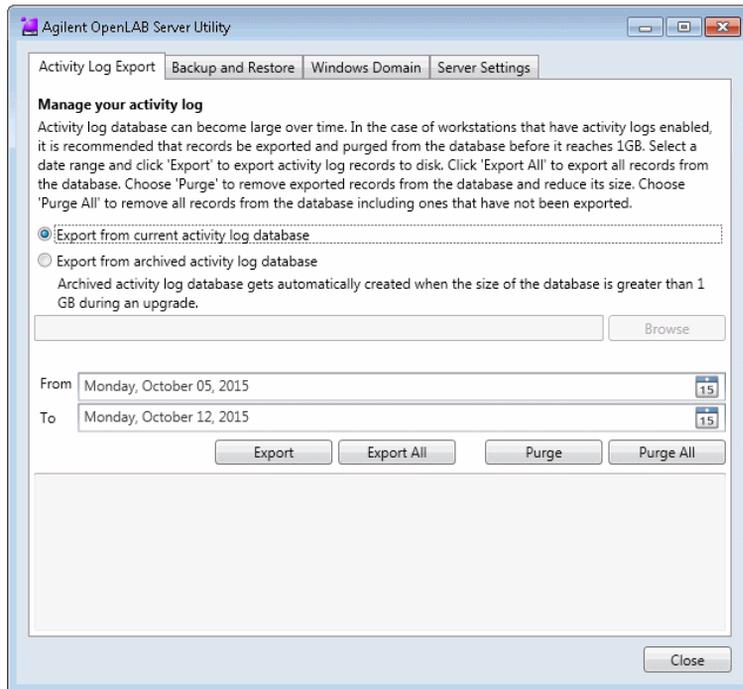


**Figure 10** OpenLAB Server Utility (for example with **Server Settings** tab selected)

## Exporting and Purging Log Entries

Activity log databases can become large over time and affect performance of activity log related operations. Use the **Export** and **Purge** functions on the **Activity Log Export** tab to export the activity log entries to an XML file and purge them from the Activity Log database.

This export can only access logs that are stored on the computer where you are using the **Server Utility** program.



**Figure 11** Activity Log Export tab (here: using an SQL CE database, therefore including the **Export from archived log database** option)

### **Export current activity log for a Networked or Distributed System**

- 1 Access the **Server Utility** program that is installed on the server.
- 2 Specify a date range, and click **Export**.

### **Export current activity log for a workstation**

- 1 Access the **Server Utility** program that is installed on the server.
- 2 Specify a date range, and click **Export**.

### **Export an archived activity log for a workstation**

You can create an archived activity log only during an upgrade from a system using *SQL CE* (primarily associated with a workstation solution) with a database larger than 1 GB. The **Export from archived activity log database** option is not available for other database types.

- 1 Access the **Server Utility** program that is installed on the workstation.
- 2 Select **Export from archived activity log database**.
- 3 Browse for and select the archived database.
- 4 Click **Export**.

During the export or purge, the OpenLAB Control Panel is disconnected from the server. Agilent recommends that you notify all users before beginning an export.

## **Using Backup and Restore Functions**

In order to simplify backup and restore tasks for the OpenLAB Shared Services database, the **Backup and Restore** tab of the OpenLAB Server Utility provides a simple interface for performing these tasks.

These functions are supported by the following database types:

- SQL Server Express or Compact
- PostgreSQL

For details on the maintenance of OpenLAB Data Store and supported database types, refer to the Data Store documentation on the OpenLAB Data Store DVD.

## Backup

Backup operations behave slightly different depending on what version of SQL Server you are using.

Backup operations are:

- supported with PostgreSQL databases
- supported with SQL Server databases (Workstation and Distributed System installations)
- supported with SQL Server databases installed by the OpenLAB CDS Master Installer
- not supported when creating a new database in an existing SQL Server.

### To perform a backup:

- 1 Specify the backup directory and retention time.

When a new backup is performed, the currently set retention time is used to delete any files older than specified.

- 2 Click **Backup**.

The backup is placed in the specified backup directory. Backups older than the retention time are deleted.

#### NOTE

The tool automatically generates filenames for the backup files. Never change these filenames, as the tool relies on a specific naming convention.

---

#### NOTE

When using SQL Server Compact (installed on workstations), you are only allowed to perform *full* database backups. This is a direct file copy of the existing database files used for the workstation.

---

- 3 When using SQL Server Express (installed with a Distributed System), you have the option to perform both full and incremental (transaction log) backups. Select the type of backup.

#### NOTE

For an SQL Server, backups are grouped into related sets, where each set contains a single full backup and all the incremental backups based on it.

---

For an SQL Server, a new backup file is created for each month regardless of the method selected.

Retention time applies on a per transaction basis so that no month's backup file is deleted unless every transaction in that backup is younger than the set time. This includes the full and incremental log transactions.

## Restore

Restore operations behave slightly different depending on what version of SQL Server you are using.

- 1 Specify the backup directory and click **Restore**.

### NOTE

The database is taken offline while this process executes a restore.

---

The restore function operates on backup sets, which include a full backup and all related incremental backups.

When restoring SQL Server Express, Standard or Enterprise database backups for client or server systems again, you will restore to the selected backup in the list.

- 2 Verify that all connections to the system are shut down before performing a restore.
- 3 If you have selected the most recent backup, and if additional transactions have been executed against that database, the tool will also ask if you would like to save those changes prior to restoring the database. Click **Yes** to effectively perform a transaction log backup prior to the restore.

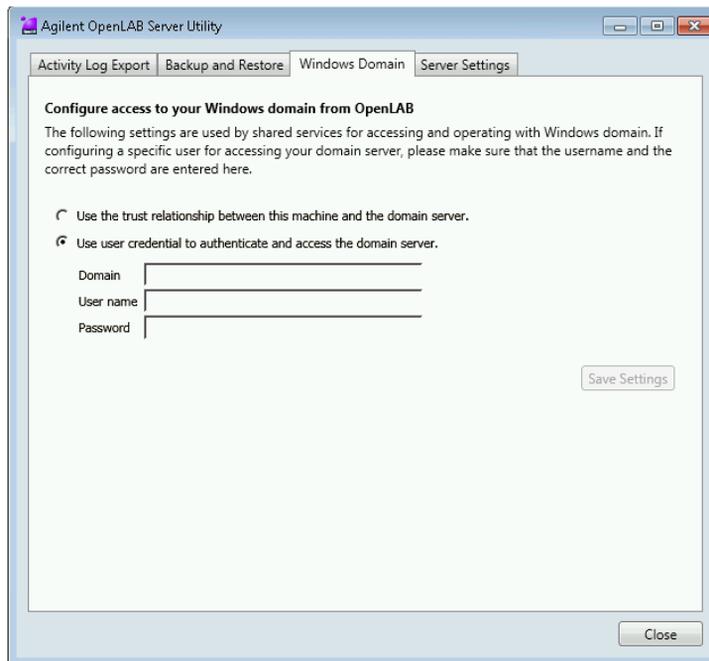
This message will not display for restore operations with an SQL Server Compact (installed on workstations).

## Configuring Access to Windows Domain

If you use Windows domain authentication to identify your OpenLAB users, OpenLAB must be given access to the server where these credentials are stored.

Normally, if a machine has been joined to the domain, the trust relationship between this machine and the domain server is sufficient to grant this access. If a problem occurs while selecting an account: Use the **Windows Domain** tab in the Server Utility program to specify or change the credentials that OpenLAB will use to access your Windows domain server. The user specified here must have the privilege to obtain user and group information from the domain.

This feature can only access credentials that are stored on the computer where you opened the Server Utility program.



**Figure 12** Windows Domain tab with specific user credentials

- 1 Choose the **Windows Domain** tab.
- 2 Enter the user credentials and click **Save Settings**.

## Managing Server Settings

In a client/server configuration, use **Server Settings** to manage server connections for your local system. The list of servers here determines which servers users may choose to connect to when they log into OpenLAB. Administrators can limit users from switching to a non-default server from this tab.

This feature manages server connections for the computer where you are using the **Server Utility** program.

The server connections for each client in a client/server system are managed through each client, therefore to change the server connections for a client, access the **Server Utility** program installed on that client.

In a workstation configuration, there is typically one server connection so this feature is not used.

To move your workstation from a domain to a workgroup, or from one domain to another domain, the SQL Server must be configured to a local account (not a domain account). See SQL Server help for more information.

**1** Select the **Server Settings** tab.

A table shows all server entries and the default server connection.

For workstation installations, by default there is one entry for the local OpenLAB Shared Services server. For distributed or networked workstation installations, there is a second entry for the central OpenLAB Shared Services server (default).

**2** You can add more servers by clicking **Add Server**.

You can switch the default server by selecting a server and clicking **Set as default**.

**3** By default, the **Allow users to choose server during login** check box is selected. To forbid users to connect to non-default servers, clear this check box.

Users will need to contact their administrator in order to connect to another server.

All connections provided here will be listed under **Local Configuration** in the OpenLAB Control Panel.

## Maintenance Procedures

### Update Database Statistics

To maintain optimal database performance, periodically update the OpenLAB Shared Services server database statistics. These statistics are used by the database engine to determine the most optimal way to execute queries.

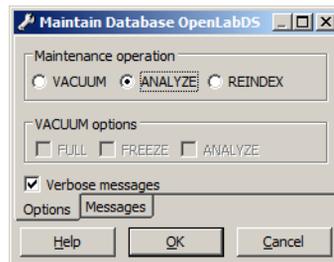
You need to update statistics for the OLSharedServices database. If a custom database name was chosen during installation, use the correct name from your installation notes.

### Procedures for PostgreSQL Database

For PostgreSQL database, these procedures must be performed on a regular basis. The frequency depends on the usage of the system. As a guideline, you should at least do this every time a full backup is taken.

#### Updating Statistics Using the Maintenance Wizard

- 1 Start PostgreSQL pgAdmin, connect as the database administrator, and select the database you want to update the statistics. The default database administrator user name is 'postgres' and the default password is an empty string (no characters).
- 2 Right-click the database and select **Maintenance....** The following form is displayed.



**Figure 13** Maintain Database

- 3 Choose **ANALYZE**, and click **OK** to analyze the database.

### Additional Maintenance for PostgreSQL Database

PostgreSQL supports some additional maintenance commands that can be beneficial to helping keep your database system running smoothly. These include VACUUM and REINDEX. See the PostgreSQL documentation for additional details about these commands.

<http://www.postgresql.org/docs/9.0/static/maintenance.html>

#### CAUTION

PostgreSQL Service packs or Hotfixes

→ Only apply Agilent provided service packs or Hotfixes to your OpenLAB PostgreSQL server.

---

## Procedures for SQL Server

For MS SQL Server database the procedure to update statistics can be easily automated using the SQL Server Management Studio.

### Updating Statistics Using Maintenance Plan Wizard

- 1 Start SQL Server Management Studio and connect as the database administrator.
- 2 Expand the server.
- 3 Expand the Management folder.
- 4 Right-click **Maintenance Plans** and select **Maintenance Plan Wizard**. Use the wizard to create a plan customized to meet your maintenance requirements.
  - a Select a **Weekly Schedule** to be executed at a time when there may be minimal activity (for example, Sunday, 12:00 noon).
  - b Select Update Statistics as the maintenance task.
  - c Choose the Shared Services database (OLSharedServices) as the database against which the task will be executed.

## Monitor Resource Usage on OpenLAB Shared Services Server

Administrators of the system must regularly monitor disk space usage on all disks where data is stored. When the disks get close to 80% full, consider increasing disk space.

CPU, memory and network usage must be monitored to check for performance bottlenecks on the server.

### Recommended Best Practices for Monitoring Resource Usage

- 1 Monitor the disk usage of OpenLAB Shared Services server at least weekly.
- 2 Optionally, implement automated disk space monitoring tools that send e-mail alerts when disk usage exceeds the thresholds. Examples of such tools are: Monit, Munin, Cacti, and Nagios.
- 3 Monitor system resource usage such as memory, CPU, and network throughput. Windows Performance Monitor can be used for this purpose.

## Purging the Activity Logs

Activity Log database can become large over time and affect performance of activity log related operations. Use the OpenLAB Server Utility tool to archive the Activity Log databases (see [“Exporting and Purging Log Entries”](#) on page 65). This utility is automatically installed with your OpenLAB software.

## Domain or Server Name Change

Contact Agilent before changing the domain membership or renaming the OpenLAB Shared Services server.

## Additional Best Practices

- 1 Apply 3rd party updates and patches on the OpenLAB Shared Services server.

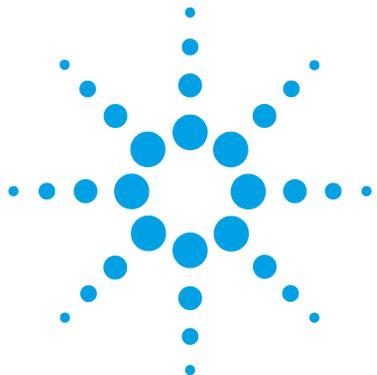
On the Agilent Customer Care Portal, Agilent regularly posts information on 3rd party updates and patches that have been validated for use with the OpenLAB software suite. These include OS security patches and updates, database updates, and application updates.

The Customer Care Portal is available at:

<http://www.ccportal.chem.agilent.com/PortalHome>

- 2 Apply Agilent software updates.

Apply software updates for Agilent OpenLAB Shared Services on your Data Store server. When you receive notification of an update, please take note and read the information to determine if the update is applicable, and its urgency.



## 4 ChemStation-Specific Administration

ChemStation Administration Tool	76
About the ChemStation Administration Tool	76
Enabling users to start the ChemStation Administration Tool	77
Machine-dependent encryption	79
ChemStation Session Locks	79
Automatic Data Transfer Settings	81
Transfer Management Settings	82
Audit Trail Settings	83
Configuration Profiles	84
Folder Protection with Secure File IO	85
Support Reports	89
OpenLAB CDS Config Checker	90
Fail Over Procedures for ChemStation Edition	91
Introduction	91
Scenarios	94
Preparations for Networked Workstations	97
Fail Over Instructions for Networked Workstations	104
Preparations for Distributed Systems	107
Fail Over Instructions for Distributed Systems	111

This chapter describes various tools that are helpful for diagnosis, support and troubleshooting.



## ChemStation Administration Tool

### About the ChemStation Administration Tool

The ChemStation Administration Tool offers a number of functions related to the ChemStation configuration. One of these functions is to configure the data handling and audit trail settings per ChemStation instance (that is, per instrument and per online or offline instance). Alternatively, this function allows using the same settings for all instances running on a workstation, networked workstation, or Agilent Instrument Controller (AIC). Access to the ChemStation Administration Tool is therefore strictly limited:

- The ChemStation Administration Tool can only be opened directly on the ChemStation PC. In Distributed System installations, you must open the tool on the relevant AIC.
- The ChemStation Administration Tool can only be started by users who are a member of the local user group **CSAdministrators** (see [“Enabling users to start the ChemStation Administration Tool”](#) on page 77).

#### To start the ChemStation Administration Tool:

- 1 From the Start menu in the Task Bar, select **Start > All Programs > Agilent Technologies > OpenLAB CDS ChemStation Edition > ChemStation Administration Tool**.

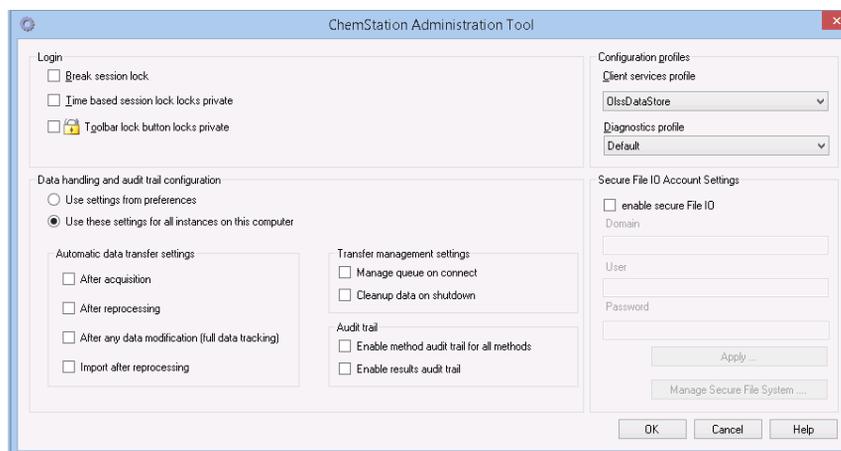


Figure 14 ChemStation Administration Tool

## Enabling users to start the ChemStation Administration Tool

During the installation of OpenLAB CDS ChemStation Edition, the local user group **CSAdministrators** is automatically created. Only members of this group are allowed to run the ChemStation Administration Tool. The user who installs ChemStation is automatically added to the **CSAdministrators** group. Additionally, the Windows group **Administrators** and the user who installs ChemStation are granted Full Control privileges on the Administration Tool program executable (Agilent.ChemStation.ECM.ECMAdmin.exe) itself, thus they are able to run the tool.

To add a Windows user to the **CSAdministrators** group:

- 1 Go to **Start > Control Panel > Administrative Tools**<sup>1</sup>.

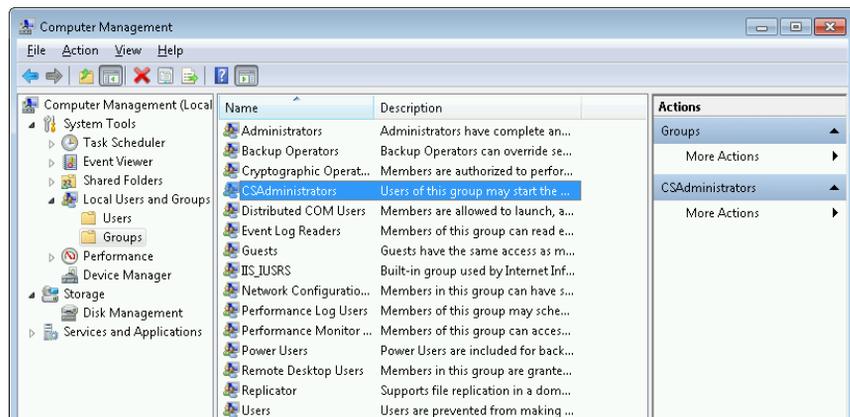


Figure 15 Computer Management window

<sup>1</sup> View the items by icon to see a list of all items

## 2 Select Computer Management

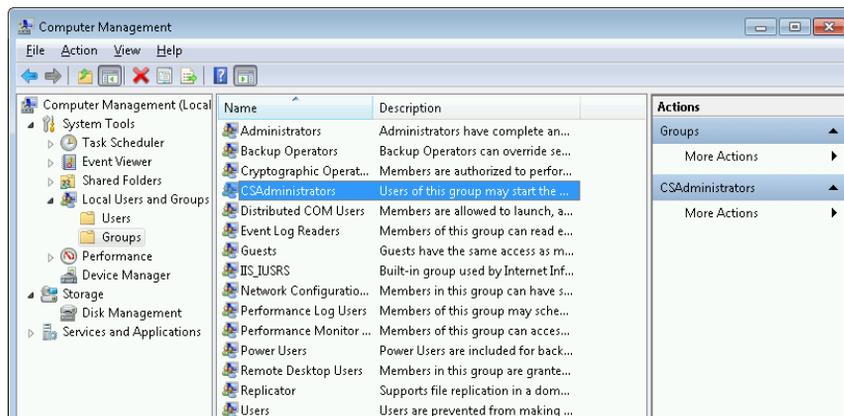
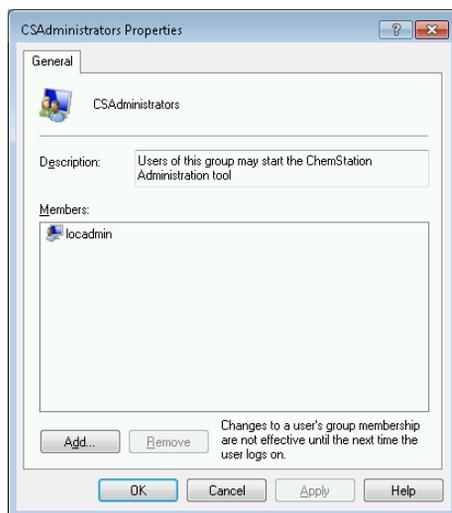


Figure 16 Computer Management window

- 3 Under **Groups**, right-click the group **CSAdministrator** and select **Add to Group...** from the context menu.

The **Properties** dialog shows the users who are currently members of the group.



- 4 Use the **Add** button to add the required users.

After confirming with **OK**, the **Properties** dialog also contains the newly added users.

## Machine-dependent encryption

When you call the ChemStation Administration Tool, it encrypts parts of the ChemStation application configuration file on the system. This encryption is machine-dependent.

### NOTE

If you want to create an image of a ChemStation installation that can be cloned and distributed to other PCs, you must not call the ChemStation Administration Tool prior to creating this image. ChemStation will not work on images with an encrypted ChemStation application configuration file.

---

## ChemStation Session Locks

### Concept of Session Locks

If you leave the ChemStation computer for a certain period of time, you can lock ChemStation so that no other user can access the application. This is a safety feature to ensure that there is no unauthorized access to ChemStation. When you activate the session lock, you or another user must first provide a valid login before continuing to work with ChemStation.

In ChemStation, there are the following options to activate the session lock:

- *Privately* (**User > Lock Session > privately**): Only the user who activated the session lock, or a user with the **ChemStation: Break session lock** privilege, can log in. You can set this privilege in the OpenLAB Control Panel. It is part of the project privileges (see “Roles and Privileges” on page 60).
- *Non-privately* (**User > Lock Session > non privately**): Any valid user can log in. This is useful, for example, if there is a shift change, and the personnel ending their shift secure ChemStation until the personnel of the new shift start work.

## 4 ChemStation-Specific Administration

### ChemStation Administration Tool

- *Toolbar lock button:* The toolbar lock button can be configured to lock the ChemStation session privately or non privately.
- *Time-based:* Depending on the configuration in OpenLAB Control Panel, ChemStation is automatically locked after a given period of time without any user interaction (see **Inactivity Timeout** under “[Security Policy](#)” on page 56).

The time-based session lock can be configured to lock the ChemStation session either privately or non-privately (see [Figure 17](#) on page 80).

### Session Lock Settings in the Administration Tool



**Figure 17** ChemStation Administration Tool

In the ChemStation Administration Tool, you can set the following options for creating and breaking session locks:

- **Break session lock:** If you select this check box, a locked ChemStation can be accessed by any user just by clicking **Cancel** in the **Login** dialog. If a ChemStation is currently locked and the authentication provider is unavailable, selecting this check box is the only way to regain access to the current ChemStation session.

#### NOTE

Be aware that, as a consequence, the user who regained access to ChemStation now has unlimited access to all ChemStation functions.

- **Time based session lock locks private:** If ChemStation has been locked by a session time out, only the current user or a user with the required privileges can unlock this session.
- **Toolbar lock button locks private:** If ChemStation has been locked using the Lock button on the ChemStation toolbar, only the current user or a user with the required privileges can unlock this session.

## Automatic Data Transfer Settings

You can apply the following options either individually in each ChemStation instance, or globally across all instances of ChemStation on a workstation or AIC.

### After Acquisition

If you select this check box the data is automatically uploaded to the central repository after an acquisition. The raw data files are written to the local ChemStation file system while the sequence is still running. When the complete sequence is finished, the raw data files are packed in an SSIZip file which is then uploaded to the central repository.

### After Any Data Modification

If you select this check box, the result set is automatically uploaded to the central repository after you changed the data analysis parameters for a sample. The data is uploaded even if you do not reprocess the sequence.

### After Reprocessing

If you select this check box, the result set is automatically uploaded to the central repository each time after you reprocessed the sequence.

### Import after Reprocessing

If you select this check box, a sequence that is only stored locally is automatically uploaded to the central repository after reprocessing. This setting is useful, for example, if you reprocess data from an older version of ChemStation.

## Transfer Management Settings

You can apply the following options either individually in each ChemStation instance, or globally across all instances of ChemStation on a workstation or AIC.

### Manage queue on connect

If the connection to the central data storage is interrupted for any reason, a running data upload might be interrupted. In this case, the remaining data is written to an internal queue on the workstation or AIC.

If you select the **Manage queue on connect** check box, ChemStation tries to upload the remaining data once the connection to the central data storage is established (that is, each time a user logs on to the central data storage system from ChemStation).

If you select this check box, ChemStation checks the local file system at each shutdown. It deletes all local data and sequence files that have been stored in the central repository. Methods and sequence templates remain on the local file system.

#### NOTE

Consider selecting the **Cleanup data on shutdown** check box when the system shall be compliant to 21 CFR Part 11. Using this function prevents unauthorized access to the local data files.

#### CAUTION

Inadequate check box selection

Loss of data

→ If you select the **Cleanup data on shutdown** check box, you should also select the Automatic data transfer settings **After acquisition**, **After reprocessing**, and **After any data modification**. Otherwise, if the users forget to upload the data to the central data storage before closing ChemStation, data might be lost.

## Audit Trail Settings

### Audit Trail

In the **Audit Trail** group, specify the conditions for the automatic audit trails for methods and results.

<b>Enable Method Audit Trail for all Methods</b>	Mark this check box to switch on the Method Audit Trail for all methods.
<b>Enable Results Audit Trail</b>	Mark this check box to switch on the Results Audit Trail for all results, and the Sequence Audit Trail for all sequences created with <b>Unique folder Creation ON</b> .

These settings override the settings in the **Audit Trail** tab of the **Preferences** dialog box in the ChemStation.

## Configuration Profiles

### Client Services Profile

By providing specific client services profiles, you activate specific functions and behaviors in ChemStation. The following profiles are relevant if you use OpenLAB Shared Services:

- **OIss**

ChemStation communicates with OpenLAB Shared Services; it sends status information to OpenLAB Shared Services, and uses the settings defined in OpenLAB Shared Services (for example, user authentication, roles and privileges, configuration settings, audit trail settings). Use this profile for ChemStation instances that are connected to OpenLAB Shared Services, but have no connection to a central data storage system.

- **OIssEcm**

In addition to communication with OpenLAB Shared Services, ChemStation allows transferring data files to and from OpenLAB ECM. Use this profile for ChemStation instances that are connected to both OpenLAB Shared Services and OpenLAB ECM.

In a fail over scenario, where the connection between the workstation and OpenLAB ECM is interrupted, you can set the authentication provider in OpenLAB Shared Service to **None**. This allows you to log on to ChemStation in the absence of OpenLAB ECM.

In this case, using the **OIssEcm** profile allows ChemStation to send the data files to the queue. Once the authentication provider in OpenLAB Shared Services is set back to ECM, the upload can be resumed.

- **OIssDataStore**

In addition to communication with OpenLAB Shared Services, ChemStation allows transferring data files to and from OpenLAB Data Store. Use this profile for ChemStation instances that are connected to both OpenLAB Shared Services and OpenLAB Data Store.

## Diagnostics Profile

For each ChemStation instrument, the system creates two log files with information for diagnostic purposes:

- Review.svclog for the offline instrument
- Acquisition.svclog for the online instrument

With the **Diagnostics profile** setting, you can define the level of logging information that is written to these files. You can select a level of logging between **None** and **Maximum**; the normal amount is provided by the **Default** selection. With the **ECM Troubleshoot** selection, more of the ECM specific information is added to the log files. With the **Troubleshoot** selection, the system not only adds the information to the log files but also creates small memory dump files in case of critical errors.

### NOTE

Only change the **Diagnostics profile** setting if instructed to do so by an Agilent Support representative.

## Folder Protection with Secure File IO

ChemStation metadata such as methods or sequences are stored in various local folders. By default, the content of these folders is not protected against modification or deletion from local file browsers or from file dialogs within ChemStation.

To ensure data integrity, enable the secure file IO function for ChemStation. All relevant paths will then be protected. The relevant paths are set by default, but can also be configured. The content of protected paths can no longer be modified by Windows file operations because members of the Windows user group **Interactive**<sup>1</sup> are denied access. The protected folders can only be used inside ChemStation. If audit trails are enabled, all modifications inside ChemStation will be documented by the corresponding audit trail entries.

<sup>1</sup> In Windows, all users who log on interactively are automatically assigned to the *Interactive* group. The membership persists for the time the users remain logged in.

#### Prepare Secure File IO in Windows:

- 1 Ensure that all relevant ChemStation folders are on an NTFS file system.
- 2 Do not use ChemStation folders on removable storage devices (for example, external disks or USB sticks).
- 3 Create a standard user account (local or domain) with the same or fewer privileges than a normal ChemStation user. Do not grant administration or power user privileges.

This account will be used by the ChemStation for internal file operations. As the user is not supposed to log on interactively, do not use the option **Change password at next logon**.

#### NOTE

Consider using a password that never expires. If this is not possible due to your security policy, plan a downtime for ChemStation while updating the password.

As long as the password for this user account is expired, do not use the ChemStation. ChemStation will not work properly until you have updated the password in the ChemStation Administration Tool.

---

#### Enable Secure File IO in ChemStation:

- 1 Click **Start > All Programs > Agilent Technologies > OpenLAB CDS ChemStation Edition > ChemStation Administration Tool** to open the ChemStation Administration Tool.
- 2 In the ChemStation Administration Tool, select the **enable secure file IO** check box.
- 3 Provide the credentials of the special user account that you have prepared.
- 4 Click **Apply**.
- 5 A dialog lists all directories that are going to be protected. Click **OK** to confirm.

A protection information file is stored inside the folder to document the date, time and user who has enabled the protection. The file can be used to demonstrate the uninterrupted protection status of a folder.

- 6 For administrative tasks, you can temporarily disable the protection of specific folders:
  - a Shut down all ChemStation sessions.
  - b In the ChemStation Administration Tool, click **Manage Secure File System ...**
  - c In the list of directories, find the relevant path and click **Unprotect**.  
The protection information file is deleted to document the interruption of the protection state.
  - d After finishing your tasks, click **Protect All** to restore protection for all relevant folders.  
New protection information files are created.
  - e Start ChemStation again.  
ChemStation will only start if all folders are protected.

#### **Check privileges for Chemstation users**

Users can load data, methods, or sequences from paths that are set up in the ChemStation Preferences. New paths can be added to the Preferences. Adding paths to directories that are not empty requires specific privileges which are granted by default.

To remove the corresponding privileges, edit the user's role in OpenLAB Control Panel and clear the following check boxes:

- **ChemStation: Data > Load data using not configured path**
- **ChemStation: Method > Load/Save method using not configured path**
- **ChemStation: Sequence > Load/Save sequence template using not configured path**

### **Important Notes for Working with Secure File IO**

When using Secure File IO to protect your ChemStation folders:

- Do not share protected folders. Sharing a folder will break the protection.
- Do not log in interactively with the user account provided in the ChemStation Administration Tool.

If this user logs in interactively, you will not be able to save any changed data in ChemStation until the user logs out again.

- If the password for this user account has expired, do not use the ChemStation until you updated the password in the ChemStation Administration Tool. ChemStation will not work properly as long as the password is outdated.
- When uninstalling ChemStation, it is recommended that you first disable Secure File IO and start the uninstallation afterwards.

## Support Reports

ChemStation allows you to create a support report. This report contains status information on the ChemStation and its environment, for example, the serial number and firmware revision of configured instruments.

### To create a support report:

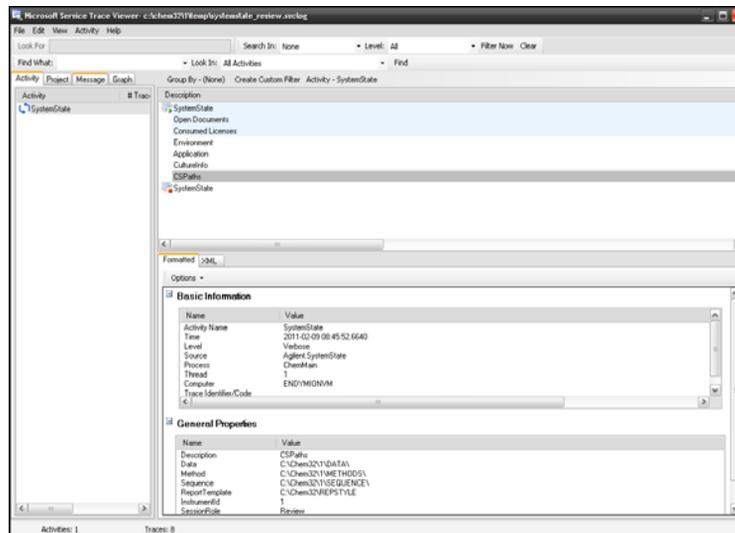
- 1 Click **Help > Generate Agilent Support Information** in ChemStation.

The file SystemState\_Review.svclog is created in a temporary ChemStation directory, typically chem32\X\temp with x as the instrument number.

**NOTE**

Fine granular device information can only be provided for RC.NET and is not supported by classic drivers.

- 2 You can view this file with the Microsoft Service Trace Viewer.



**Figure 18** Support Report, opened with Microsoft Service Trace Viewer

## OpenLAB CDS Config Checker

### NOTE

This tool is only available for OpenLAB CDS ChemStation Edition on Windows 7.

---

The OpenLAB CDS Configuration Checker helps to troubleshoot the Operating System configuration and to prevent computer problems. It checks and repairs all *mandatory* settings. It does not check optional settings or settings that improve the performance.

The OpenLAB CDS Configuration Checker comes as .diagcab file, which is a file format used with the Microsoft Windows Troubleshooting Platform (WTP) program. The Microsoft Windows Troubleshooting Platform (WTP) is a platform to locate and fix hardware and software settings in Windows. It is used specifically for diagnosing and repairing computer settings.

In general, .diagcab files are useful for deploying troubleshooting packs because they are self-contained and require no installation. They can be deployed onto web sites, network shares, or copied to USB keys. The .diagcab file name extension is a registered file name extension that can be executed by WTP.

To start the OpenLAB CDS Configuration Checker, call the file Agilent.Wtp.ChemStation.WindowsConfiguration.diagcab. This file is located on the OpenLAB CDS disc 6 under Tools\OpenLAB CDS ChemStation Edition\Diagnostics\WindowsConfiguration\Cab\

### NOTE

While using the Configuration Checker:

- Ensure that this computer is not turned off by another user.
  - Ensure that the menu bar is enabled.
-

# Fail Over Procedures for ChemStation Edition

## Introduction

OpenLAB CDS is supported in a number of different configurations. All these configurations may require network communication beyond instrument connectivity. The networking infrastructure in different laboratories can have different levels of stability, reliability, and robustness.

OpenLAB CDS provides a number of features to make it resilient against network or server outages. In case of a network or server outage, all running or queued sequences continue to run, and the data acquired is stored in the OpenLAB ECM or Data Store queue. The data can be uploaded to the central repository once the systems have been restored to normal operation. However, there will also be situations where users will want to start a ChemStation session during a network or server outage.

This chapter describes fail over and fail back concepts and procedures that ensure continued operation of OpenLAB CDS ChemStation Edition in case of network or instrument failures or server outages.

For the OpenLAB CDS Distributed System, this concept uses a workstation-based approach to remediate server unavailability. It requires dedicated failover workstation licenses that are commercially available.

Networking is involved in multiple areas:

- Instrument communication: Transfer of commands, methods and data
- Authentication: Startup of the OpenLAB Shared Services Control Panel and startup of ChemStation (from OpenLAB Control Panel)
- Status reporting: ChemStation to OpenLAB Shared Services server
- Data upload: ChemStation to central data storage
- Licensing: Licenses are retrieved from the License server (OpenLAB Shared Services server by default)

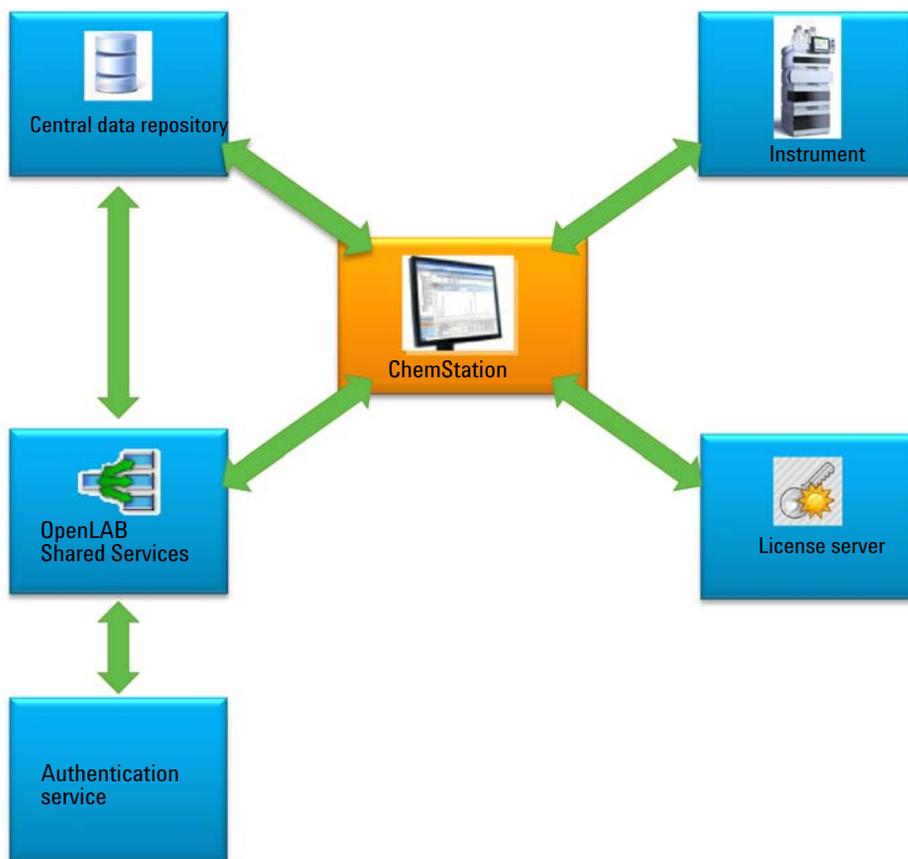
## 4 ChemStation-Specific Administration

### Fail Over Procedures for ChemStation Edition

Communication can be impacted if either the network or one of the following backend services is not available:

- OpenLAB Shared Services server
- OpenLAB ECM or Data Store server
- Windows Domain Controller, DNS server etc.
- License server

Depending on the system configuration, ChemStation has a number of connections to other systems,



**Figure 19** Connections to other systems depending on the system configuration

**Table 17** Purpose of the various communication channels and backend systems

<b>System</b>	<b>Purpose</b>	<b>Startup</b>	<b>Runtime</b>	<b>Comment</b>
OpenLAB Shared Services server	<p>Always required.</p> <ul style="list-style-type: none"> <li>• Transfers authentication information.</li> <li>• Provides privileges for current user (authorization).</li> <li>• Provides information about license server location.</li> <li>• Receives and distributes instrument status.</li> <li>• Receives activity log.</li> </ul>	Authentication/ authorization/ license server	Instrument status and activity log	<p>Runtime communication is not mandatory - status information is discarded in case the OpenLAB Shared Services server is not available. The activity log is buffered.</p> <p>Authentication may need a backend authentication service (e.g. Windows Domain Controller).</p>
License server	<p>Always required.</p> <p>Provides licenses for OpenLAB CDS features.</p>	Provides licenses	N/A	<p>At runtime no further license checks are done. Licenses are mandatory.</p> <p>License server is installed on the OpenLAB Shared Services server host machine.</p>
OpenLAB ECM or Data Store server	Optional data storage backend.	ECM: Authentication (through OpenLAB Shared Services)	Data upload and download	<p>In an OpenLAB ECM-based system the OpenLAB ECM server has two roles - authentication and backend storage.</p> <p>The authentication for Data Store is handled in OpenLAB Shared Services. For ChemStation, the runtime dependency is not mandatory since ChemStation places uploads into a queue if the central repository is not available.</p>

## Scenarios

There are two main scenarios in case of a failure of a backend system or a network outage. In the first scenario the application (OpenLAB Control Panel or ChemStation) is already running and continuous operation must be ensured. In the second scenario the application has not been started and has not yet established connections to one or more of the backend systems.

### Scenario 1: ChemStation is Running

ChemStation acquires all of the required resources at startup:

- Licenses
- Privileges
- Connection token for central data storage (if applicable)

Once the ChemStation is started, the OpenLAB Shared Services server and/or central data storage server may become unavailable:

- If the OpenLAB Shared Services server is not available, then instrument status is no longer reported and activity logs are buffered for deferred upload. Data acquisition and upload to the central repository are still possible.
- If the central data storage is not available, then data upload to the central repository is no longer possible. Instead data is put into the queue and can be uploaded once the central data storage is available again.
- If the authentication provider is unavailable, re-authentication is not possible, i.e. it is not possible to switch to a different user or to log in in case of a session lock. To allow users to break the session lock, select the **Break Session Lock** check box in the ChemStation Administration Tool “[Session Lock Settings in the Administration Tool](#)” on page 80.

**NOTE**

Data is automatically queued if the central data storage is unavailable. Once it is available again, queued data needs to be manually uploaded to the central repository (see the section titled *Troubleshooting - ECM or Data Store Server is not Available After Login* in the *Agilent OpenLAB CDS ChemStation Edition with Central Data Storage User's Guide*).

---

**NOTE**

In the Distributed System, the existing ChemStation sessions on the AIC continue to run and acquire data. However, it is not possible to connect to these sessions from the remote client and to see their status in OpenLAB Control Panel.

---

## Scenario 2: ChemStation Startup

For ChemStation startup two main services are required:

- Licenses: Depending on the instrument configuration some licenses are mandatory in order to start up ChemStation. A core license is always required. This requires a connection to the license server.
- Authentication: When an authentication mode other than **None** is specified, the user must authenticate with his username and password. This requires a connection to the authentication backend (OpenLAB Shared Services server, OpenLAB ECM server, or Domain Controller).

### Licensing

OpenLAB CDS ChemStation Edition needs to connect to a license server that holds valid licenses. If the license server is unavailable or ChemStation cannot connect to it, ChemStation cannot start up. If the central license server is not available it is possible to fall back onto a local licensing service. In this case failover licenses are required and must be configured on the fail over workstation. For details on populating the local licensing service with failover licenses, see [“Preparing Local Failover Licenses \(In Case the Central License Server Becomes Unavailable\)”](#) on page 101.

For a system with central data storage, please ensure that you maintain a local copy on your workstation of all methods and sequences that are required if the central repository is unavailable. This may require periodic updates of the local methods and sequences from master methods and sequences kept in the central repository (for details, see [“Preparing Local Methods and Sequences”](#) on page 104).

## 4 ChemStation-Specific Administration

### Fail Over Procedures for ChemStation Edition

**Authentication Service** If any of the required authentication backends (OpenLAB Shared Services server, OpenLAB ECM server, or Domain Controller) is not available, users can connect to a local instance of OpenLAB Shared Services instead. This allows users to startup ChemStation. Instruments need to be set up in the local OpenLAB Shared Services database manually (for details, see [“Preparing Local Instrument Configuration”](#) on page 100).

Since OpenLAB CDS ChemStation Edition workstations run a local copy of OpenLAB Shared Services and provide their own license server, traditional failover licenses would not apply. Instead, if the workstation failed, there could be a replacement workstation as part of a business continuity plan.

If OpenLAB CDS ChemStation Edition is integrated with and using OpenLAB ECM authentication, if the connection to OpenLAB ECM is lost, then the workstation can no longer be started.

## Preparations for Networked Workstations

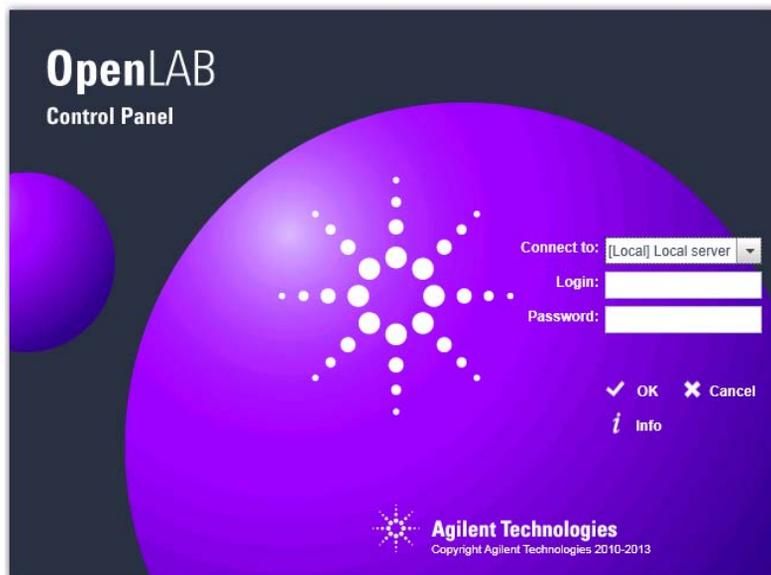
### Connecting to the Local Server

OpenLAB CDS has an option to connect to another instance of OpenLAB Shared Services. This gives you the ability to switch to local OpenLAB Shared Services when the central OpenLAB Shared Services server in a Networked Workstation or Distributed System becomes unavailable.

- 1 To connect to the local OpenLAB Shared Services instances, select **[Local] Local server** under **Connect to** when starting the OpenLAB Control Panel.

#### NOTE

If you are not allowed to switch servers, select the corresponding check box in the OpenLAB Server Utility (see “[Managing Server Settings](#)” on page 70).

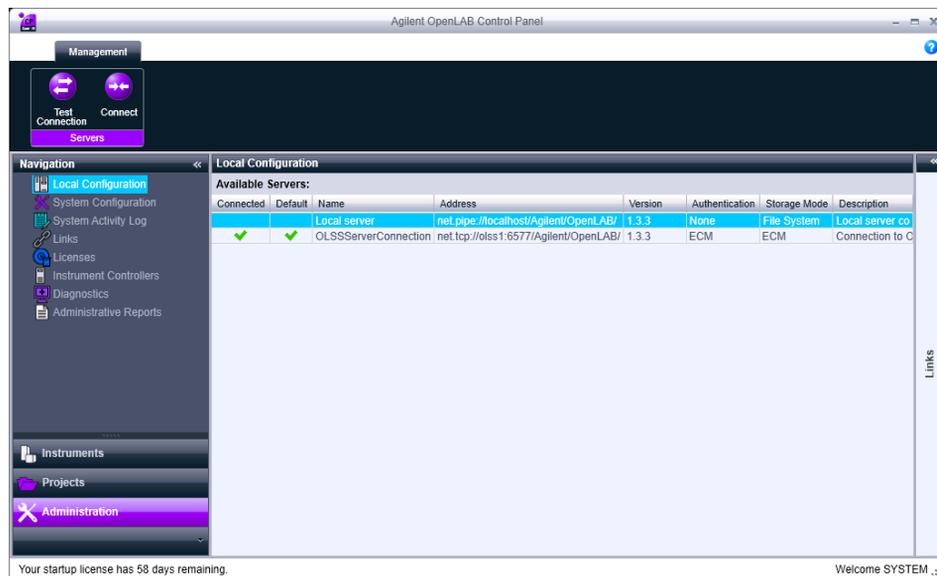


## 4 ChemStation-Specific Administration

### Fail Over Procedures for ChemStation Edition

OR

Alternatively, you can connect to the local OpenLAB Shared Services instance under **Administration > Local Configuration**. Select the local server, and click **Connect**.



You can adjust the list of available servers from the Agilent OpenLAB Server Utility program (see [“Managing Server Settings”](#) on page 70).

## Preparing Local Authentication Mode

If you are connecting to a local OpenLAB Shared Services instance, the default authentication mode is **None**. This allows users to work with the ChemStation without authentication. All users have all privileges. If user authentication is desired, use the authentication mode **Internal**. The other authentication modes (Windows Domain and OpenLAB ECM) are not adequate for a fail over because they introduce dependencies to external systems.

The authentication mode **None** is the default and nothing needs to be done in preparation for the authentication mode **None**. If you wish to use the authentication mode **Internal** for a fail over configuration, you will need to switch to the local OpenLAB Shared Services instance, configure it to use the authentication mode **Internal**, and setup users and assign roles as required.

### Prerequisites

On the central OpenLAB Shared Services server, the **Allow users to choose server during login** check box must be selected in the OpenLAB Server Utility. See [“OpenLAB Server Utility Tool”](#) on page 64.

- 1 Connect to the local instance of OpenLAB Shared Services (see [“Connecting to the Local Server”](#) on page 97).
- 2 Set up the authentication mode **Internal**.
- 3 Set up users, and set up and assign roles as desired.
- 4 When finished, connect to the central OpenLAB Shared Services server again.

## Preparing Local Instrument Configuration

In order to use the instruments that are configured on the central OpenLAB Shared Services server and connected to a given workstation, they need to be made available in the local OpenLAB Shared Services instance on that workstation.

With ChemStation, you can use a batch file to make the instruments available on the workstation PC:

- Open a command prompt on the workstation and navigate to the OpenLAB CDS ChemStation core directory in the ChemStation installation folder (e.g., c:\Chem32\Core).
- Run the OlssFailover.bat script file.

This script copies instruments configured on the central OpenLAB Shared Services to the local OpenLAB Shared Services database.

If the local OpenLAB Shared Services instance has been configured to use an authentication other than **None**, provide the username and password of an administrative user as parameters in the syntax `OlssFailover.bat [username] [password]`.

If you have made instruments available on the local OpenLAB Shared Services and then configured additional instruments, or removed instruments on the central OpenLAB Shared Services server, repeat the above.

You can omit this preparation step and execute the batch file once the fail over has occurred. This has the advantage that all instrument configurations are current. However, we recommend executing the batch file ahead of time because a user in the laboratory may not be able to execute a batch file and will not have the required user credentials available.

Preferably, put a procedure in place to ensure that the batch file is executed every time instruments are added or removed.

### NOTE

All instruments that are part of a business continuity plan using the fail over licenses, must first be configured on the central OpenLAB Shared Services.

## Preparing Local Failover Licenses (In Case the Central License Server Becomes Unavailable)

If the central license server becomes unavailable, the OpenLAB CDS application on a workstation PC can fall back to a local license service.

To prepare for this scenario, you can install failover licenses on the local license service. Failover licenses are additional workstation licenses that are installed on a workstation PC.

**1** Obtain a failover license product.

For example, obtain product number M8205AA *Option 002: OpenLAB CDS ChemStation Failover Workstation License*). This failover license product includes a workstation core license, drivers, and add-ons for a workstation with up to four instruments.

**2** Add the failover license product to the license pool in SubscribeNet.

**NOTE**

You can also use the licenses already available in your license pool. However, this is going to remove them from the pool of concurrent licenses.

**3** In SubscribeNet, register the authorization codes, generate the failover licenses for the workstations, and configure the instruments for fail over mode.

For example, a failover license for a workstation with one GC- and one 3D LC instrument includes:

- one OpenLAB CDS core license
- one OpenLAB CDS Instrument Driver for Agilent GC license
- one OpenLAB CDS Instrument Driver for Agilent LC license
- one OpenLAB CDS 3D UV Add-on license

Configure a license file with the required licenses, and enter the MAC Address of the workstation PC that the license file is to be put on. Store it locally on the workstation PC where you want to install it.

**4** Connect to the local instance of OpenLAB Shared Services (see [“Preparing Local Authentication Mode”](#) on page 99).

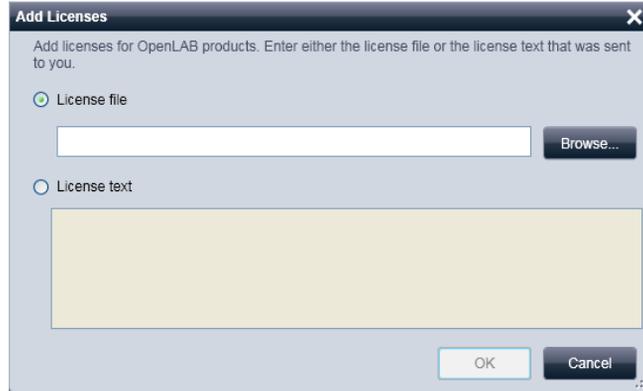
**5** In OpenLAB Control Panel, go to the **Administration** page and click the **Licenses** node.

**6** In the ribbon, click **Add**.

## 4 ChemStation-Specific Administration

### Fail Over Procedures for ChemStation Edition

- 7 In the **Add Licenses** dialog browse to the location of the license file and add it.



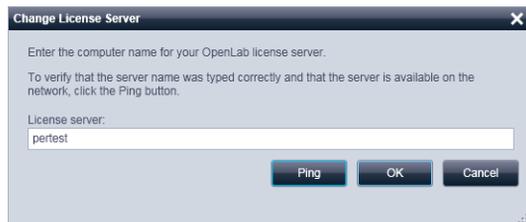
**Figure 20** Entering the failover license for the local license service in the **Add Licenses** dialog

## Preparing Declaration of License Server (In Case the Central License Server is Still Available)

If the central data storage and/or Domain Controller are unavailable, and the central license server (on the OpenLAB Shared Services server host machine) is still available, the central license server can still be used. However, it is necessary to use the local OpenLAB Shared Services instance with the authentication mode **None** or **Internal**. Example: OpenLAB ECM server is unavailable, license server is available. Using local authentication is necessary since authentication cannot be against OpenLAB ECM, but licenses can still be retrieved from the central license server.

The central license server needs to be declared in the local OpenLAB Shared Services instance.

- 1 Connect to the local instance of OpenLAB Shared Services (see [“Preparing Local Authentication Mode”](#) on page 99).
- 2 Go to **Administration > Licenses**.
- 3 Click **Change Server**. In the **Change License Server** dialog, enter the name of the central license server to be used.



**Figure 21** Change License Server dialog

- 4 Once the central license server has been declared in the local OpenLAB Shared Services instance, reconnect to the central OpenLAB Shared Services server.

### **Preparing Local Methods and Sequences**

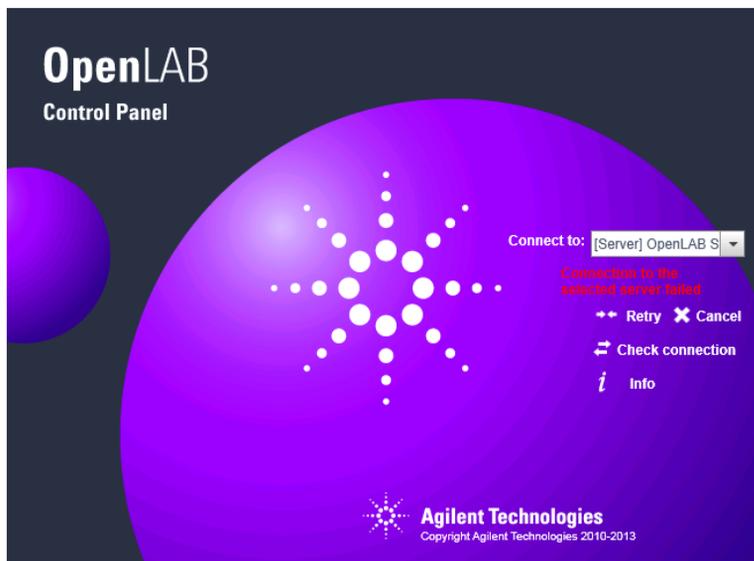
For a system with central data storage, please ensure that you maintain a local copy on your workstation of all methods, sequence templates and report templates that are required if the central repository is unavailable. As mentioned before, this may require periodic updates of the local files from the master data kept in the central repository (see [“Preparing Local Instrument Configuration”](#) on page 100).

## **Fail Over Instructions for Networked Workstations**

### **Connecting to the Local Server in Case of a Fail Over**

If the OpenLAB Control Panel cannot communicate with the OpenLAB Shared Services server or with its authentication backend, the start-up screen will show a corresponding message (see [Figure 22](#) on page 105). You can connect to the local server instead.

If OpenLAB Control Panel cannot communicate with its authentication backend (OpenLAB ECM server and/or Domain Controller), the OpenLAB Control Panel application will not start. If this happens, disconnect the ChemStation workstation PC from the network while starting the OpenLAB Control Panel. When starting a ChemStation workstation PC disconnected from the network, OpenLAB Control Panel shows a dialog box that allows users to connect to the local server.



**Figure 22** OpenLAB Control Panel starting up without connection to the OpenLAB Shared Services server

- 1 Select the **[Server]** entry, and click **Info** button. This will give you some diagnostics information. Click **Check connection** to ping the server.
- 2 If it has become clear that the network or server outage is going to be longer, login to the **[Local]** machine. If authentication is setup on the local OpenLAB Shared Services instance you will be asked for your credentials.
- 3 Since the instruments and licenses have been made available in the preparation steps, instruments can immediately be used. All instrument-related data, methods and configurations are available at their original locations. In addition, the instrument configuration is also available at its original location.

**WARNING**

**Inconsistencies between the local OpenLAB Shared Services configuration and the central OpenLAB Shared Services configuration**

- Do not create new instruments or delete existing ones while connected to the local OpenLAB Shared Services instance.

## Restoring the Connection

Once the connection to the central OpenLAB Shared Services server or authentication backend is restored, the OpenLAB Control Panel can be reconnected to the central OpenLAB Shared Services server.

### Preparations

Shut down ChemStation instances running on the local OpenLAB Shared Services instance before reconnecting to the central OpenLAB Shared Services server.

- 1 When starting the OpenLAB Control Panel, the default connection configured in the OpenLAB Server utility will be used. If this is not the central OpenLAB Shared Services server, select the **[Server]** entry under **Connect to**, and log in with the required user credentials.

OR

Alternatively, you can connect to the central OpenLAB Shared Services instance in OpenLAB Control Panel under **Administration > Local Configuration**.

- 2 OpenLAB Control Panel will connect to the remote OpenLAB Shared Services server.

### NOTE

Instruments are already available in the remote OpenLAB Shared Services server and do not need to be reconfigured.

---

## Preparations for Distributed Systems

If any of the required authentication backends (OpenLAB Shared Services server, OpenLAB ECM server or Domain Controller) is not available, or in case of a network outage, it is not possible to work directly on the CDS client or the AIC.

There are different approaches for the OpenLAB CDS Distributed System. In general, the system offers a workstation-based approach to remediate server unavailability. It requires dedicated failover workstation licenses that are commercially available. Different scenarios that have to be considered include:

- No authentication provider
- No central OpenLAB Shared Services Server
- No network

For instruments that are required to run continuously (24/7), Agilent suggests to set up a Networked Workstation on a separate PC using failover licenses, and to use this workstation in case of a fail over. This *fail over workstation* allows controlling instruments in case of a failure of any of the authentication back-ends as well as a network outage.

## Recommendations for the Fail Over Workstation

The fail over workstation will be used in case ChemStation remote desktop sessions on the AIC are no longer accessible.

### Prepare a Fail Over Workstation

**1** Determine required workstations and licenses.

As part of a business continuity plan, assess how many and which workstations and instruments need to be available as part of the fail over plan. Based on the assessment determine the required number of failover licenses.

An option to reduce hardware costs is to use fail over workstations as client PCs in a Distributed System.

#### NOTE

The OpenLAB CDS hardware and software requirements for Networked Workstations apply. It is recommended to use a PC with two LAN cards:

- One to connect directly to the instrument.
- One to connect to the laboratory network: This allows easy retrieval of data once systems have been restored to normal operation.

---

**2** Determine IP address needs.

The fail over workstation may need a different IP address to connect to the instrument to support fail over mode.

Ensure common network settings for the LAN cards of the fail over workstation and the AIC. Thus you can reuse the instrument's IP address.

If the instrument is set up on the LAN (that is, not directly connected to the AIC), the IP-address of the instrument needs to be adapted to the subnet of the fail over workstation's network card.

After the fail over has occurred, and systems have been restored, the instrument's configuration needs to be restored before it is reconnected to the AIC.

**3** Determine additional equipment needs

You may require additional hardware such as switches, network cabling, etc.

**4** Install an OpenLAB CDS ChemStation Edition Networked Workstation on each fail over workstation. Add central data storage and configure the same database server as for the AIC. Register in the central OpenLAB Shared Services database.

- 5 Execute IQ/OQ procedure if required.
- 6 While connected to the central OpenLAB Shared Services server, configure all instruments for which you require continued operation in case of an anomalous event.

Use a different instrument name, e.g. "<instrument name>\_failover".

Start a session for each instrument (online or offline) and configure the upload path in the preferences.
- 7 Lock all failover instruments using the **Lock** function in OpenLAB Control Panel.

During normal operation these instruments remain locked.
- 8 Close OpenLAB Control Panel, and start it again. This time, connect to the *local* Shared Services database.
- 9 Recreate the instruments on the local server using the script `OlssFailover.bat` developed for networked workstations (see [“Preparing Local Instrument Configuration”](#) on page 100).
- 10 If required, prepare the authentication mode on the local server.

On the fail over workstation, the local OpenLAB Shared Services must use the authentication mode **None** or **Internal**.

In case of internal authentication, you must set up fail over users and roles (see [“Preparing Local Authentication Mode”](#) on page 99).
- 11 Install failover licenses (see [“Preparing Local Failover Licenses \(In Case the Central License Server Becomes Unavailable\)”](#) on page 101).
  - If the license server remains available (that is, only the authentication provider becomes unavailable):

The central license server needs to be declared in the local OpenLAB Shared Services (see *Change the license server* in the OpenLAB Control Panel Help).
  - If the license server becomes unavailable, too:

Failover licenses need to be available in the license pool. Generate the corresponding license file in `SubscribeNet`, and add it to the fail over workstation using the local OpenLAB Shared Services instance.

This licenses file is then bound to the MAC address of the fail over workstation.

## 4 ChemStation-Specific Administration

### Fail Over Procedures for ChemStation Edition

**12** Make all required methods, sequence and report templates available locally.

Consider using acquisition only methods for the anomalous event, and to postpone processing until the system has been fully recovered.

**13** When the anomalous event is over:

- If you want to use the workstation as a normal client or Networked Workstation, switch back to the central OpenLAB Shared Services database.
- If the PC is a dedicated fail over workstation, turn the workstation off.

In your Standard Operating Procedures, define that the usage or changed purpose of the machine needs to be documented.

#### NOTE

Fail over workstations need regular maintenance. They require the same level of operating system updates and OpenLAB CDS updates as the Distributed System.

---

## Fail Over Instructions for Distributed Systems

### Using the Fail Over Workstation

During an anomalous event, the OpenLAB CDS ChemStation Edition instances on the AIC continue to run even if the OpenLAB ECM server, OpenLAB Shared Services server or license server are unavailable.

To make an instrument available during an anomalous event, follow the instructions below.

- 1 Check whether the sequences running on the AIC are completed and the instrument is idle. To confirm this, open the sequence folder and review the log files.
- 2 Reconfigure the instrument. This step depends on the topology used.
  - The AIC has two network cards; the instrument is directly connected to one of them.  
Disconnect the instruments from the AIC (by removing the network cable from the instrument) and connect the instrument directly to the fail over workstation.
  - The instrument is connected to the laboratory network.  
You may need to change the instrument's IP address.
- 3 On the AIC, close the ChemStation sessions using the Windows Task Manager.
- 4 Power on the fail over workstation, or use the dedicated Networked Workstation, and connect to the local server.  
  
Since the instruments and licenses have been made available in the preparation steps, instruments can immediately be used on the fail over workstation.
- 5 Start acquiring data.  
  
The instrument name now includes the term *failover*. The data is securely buffered in the queue.

### Restoring the Connection

Once connection to the OpenLAB Shared Services server or central data storage is reestablished, proceed as follows:

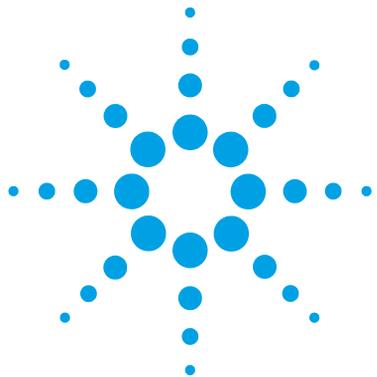
#### NOTE

Once systems have been restored to normal operation, instruments may need to be reconfigured to their original IP-addresses before connecting them to the AIC again.

- 1 On the fail over workstation, connect to the central OpenLAB Shared Services server.
- 2 Start the online or offline ChemStation instrument sessions and resume the queue.  

Acquired data will be uploaded to the central repository.  
The data acquired during the anomalous event will have the user configured in the failover mode as the Acquisition User.

  - If authentication mode **None** was configured in the local Shared Services database, the user **System** will be the Acquisition User.
  - If a fail over user was set up for the failover scenario, this fail over user will be the Acquisition User.
- 3 Close the instrument sessions.
- 4 Power off the fail over workstation.
- 5 Reconnect the instruments to the AIC by unplugging the network cable from the fail over workstation and connecting the LAN cabling with the AIC. If applicable, reassign IP addresses to the instruments.
- 6 On the CDS Client, launch OpenLAB Control Panel and connect to the central OpenLAB Shared Services server. Check the status of the individual instruments in the **Instruments** panel to ensure they are connected and available.



## 5 EZChrom-Specific Administration

Fail Over Procedures for OpenLAB CDS EZChrom Edition	114
Connecting to the Local Server	115
Switching Back to Remote Server	118
AIC Driver and Install Tool	120
Use the AIC Driver and Install Tool	120
Change the Password of the Instrument Service Account	121
Change the Password of the AFS Administrator	122

This chapter describes fail over procedures and tools for AIC administration.



## Fail Over Procedures for OpenLAB CDS EZChrom Edition

OpenLAB CDS EZChrom Edition has an option to connect to another instance of OpenLAB Shared Services. This gives you the ability to switch to local OpenLAB Shared Services on an AIC machine when the central OpenLAB Shared Services server in a Networked Workstation or Distributed System environment becomes unavailable. This will put you in a Workstation mode. You will then need to create projects, instruments, and users (if authentication is setup) on the local OpenLAB Shared Services, so you can connect to instruments and acquire data. When the central OpenLAB Shared Services server becomes available you can then reconnect your system to the central OpenLAB Shared Services server and move over your data.

If an AIC is acquiring data when the OpenLAB Shared Services server and Enterprise Path become unavailable, it will continue to acquire data until the sequence is finished. Once the sequence is finished the AIC will keep trying to put it on the Enterprise Path until it becomes available. When in fail over mode, follow a specific naming convention which distinguishes fail over file names from the file names on the distributed system. For instance, use this convention to prevent issues from copying files from the fail over workstation to the distributed system.

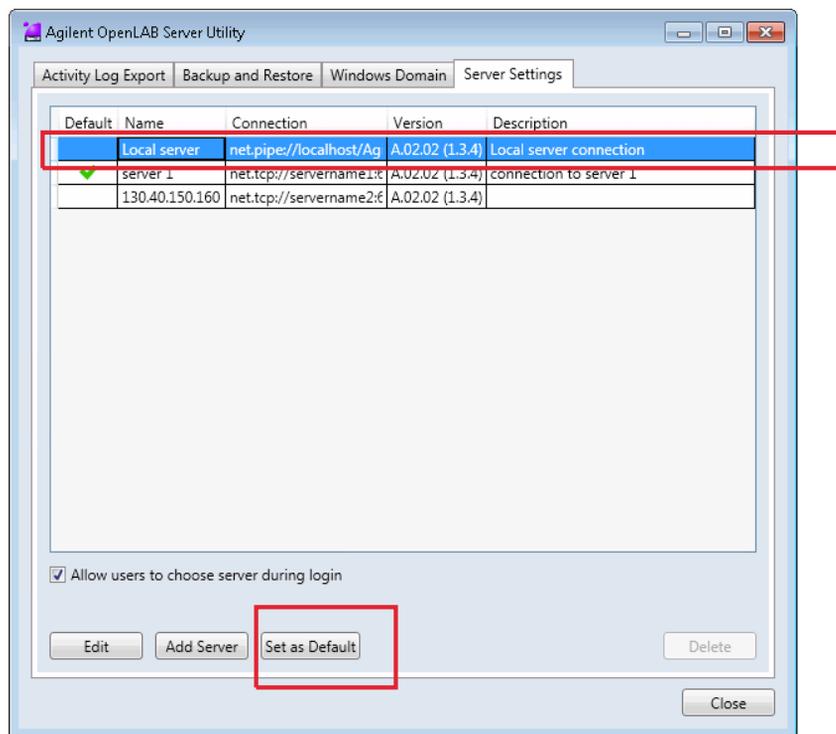
Do not apply electronic signatures while being connected to the local server. The signatures may become invalid after switching back to the central OpenLAB Shared Services server.

## Connecting to the Local Server

**Prerequisites** You have installed an EZChrom AIC.

### Convert an AIC to a local server

- 1 Log in as an administrator to the operating system of the machine that you want to convert.
- 2 If OpenLAB Control Panel is open, close it.
- 3 Launch the **OpenLAB Server Utility** on the AIC, and select the **Server Settings** tab.
- 4 Select the local server, and click **Set as Default**.



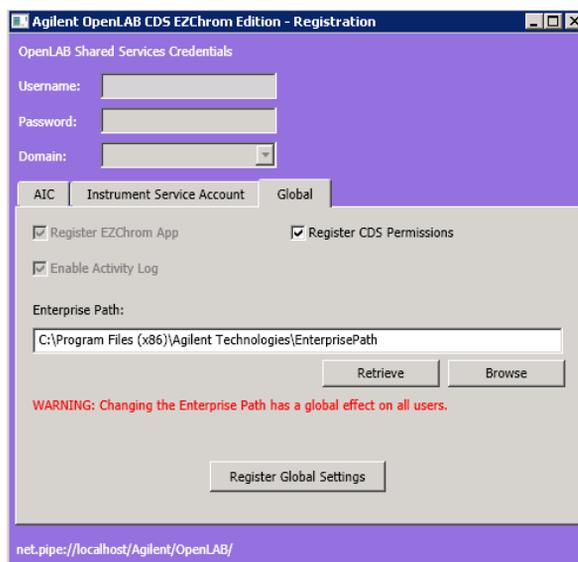
## 5 EZChrom-Specific Administration

### Fail Over Procedures for OpenLAB CDS EZChrom Edition

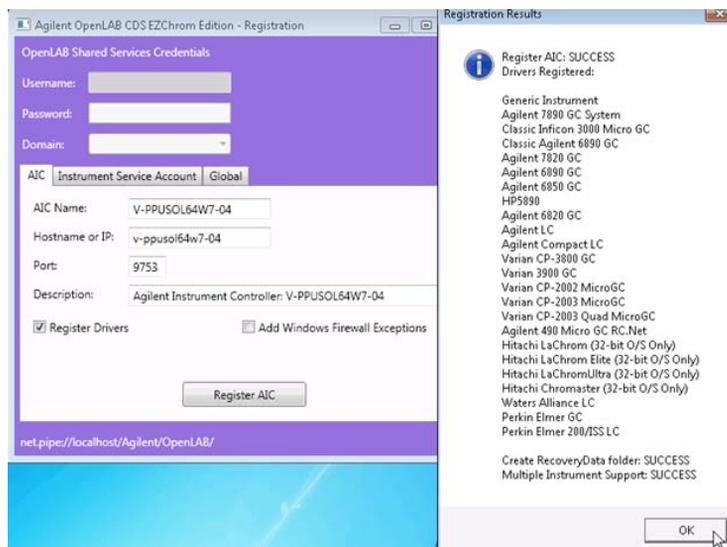
- 5 Close the OpenLAB Server Utility.
- 6 To confirm that the OpenLAB Control Panel connects to the local server, launch the OpenLAB Control Panel.  
If it is connected to the local server, you will only see the **Administration** tab.  
Under **Administration > Local Configuration**, verify that you are connected to the local server.
- 7 Close the OpenLAB Control Panel.
- 8 Launch the **AIC Driver and Install Tool**, and select the **Global** tab.
- 9 Under **Enterprise Path**, enter a local path, and click **Register Global Settings**.

#### NOTE

Proceed with the next step only after you have registered the global settings.



10 Select the **AIC** tab, and click **Register AIC**.



11 Reboot the machine.

12 Launch the OpenLAB Control Panel.

The **Projects** and **Instruments** tabs appear.

### Add failover license and configure instruments

- 1 In the OpenLAB Control Panel, add a failover license.
- 2 Set the security as needed. The recommendation for fail over workstations is to use internal authentication.
- 3 Create new projects, instruments, and users (if authentication is set) on the local OpenLAB Shared Services server.
- 4 Launch the instrument and create a new method. Save the newly created method.
- 5 Create a new sequence. Save the newly created sequence file.

## Switching Back to Remote Server

- 1 Log on to the fail over workstation as a local administrator.
- 2 Launch the **OpenLAB Server Utility**, and select the **Server Settings** tab.
- 3 Select the Distributed System server, and click **Set as Default**.
- 4 Launch the **AIC Driver and Install Tool**.
- 5 On the **AIC** tab, re-register the AIC.
- 6 Reboot the machine.
- 7 Once the remote server is up and running, you need to copy data from the Enterprise folder on the local server to the remote server.
  - a *For a File System* - Manually copy all files from the local path to the storage path on the remote server.
  - b *For ECM and Data Store* - Use the procedure described below.

### To copy data to the remote ECM or Data Store server

- 1 Manually upload files from the Method, Sequence, Data, and Sample Prep folders from the fail over workstation to the Method, Sequence, Data, and Sample Prep folders on the ECM or Data Store machine
- 2 The following changes to the method file on the ECM or Data Store machine may be needed depending on how your methods were created.
  - **Method > Advanced > Files tab** – Make sure the baseline file path is changed to the ECM or Data Store path
  - **Method > Advanced > Advanced Report** – Make sure the report template name field is changed to the ECM or Data Store path.
- 3 The following changes to the sequence file on the ECM or Data Store machine may be needed:
  - a Open your sequence and browse to **Sequence > Properties**. Verify the paths are changed to point to the ECM or Data Store path.
  - b In the Sequence Table:
    - Browse to the run type column. If the run type column has options selected which point to report template paths, these paths will need to be changed to the Data Store or ECM Path
    - Look at the Sample Prep column. If this field is used, the path for the Sample Prep file will need to be changed to the Data Store or ECM path. This does not need to be changed if you are using the default path.

- Look at filename column. Change this to the Data Store or ECM path if necessary. This does not need to be changed if you are using the default path.
  - Look at the Method column. Check the method path to see if it needs to be changed to the Data Store or ECM path. This does not need to be changed if you are using the default path.
- 4** Create a result set using the information from the files above:
- a** Open a sequence that was modified according to the procedures listed in the previous section.
  - b** Reprocess the sequence. After reprocessing a result set will be created in Data Store or ECM. When the result set is created in Data Store, a folder in the result folder will be created containing individual files. In ECM the result set will get zipped up in the Result folder. The result set package options selected in your project will determine which files are placed into the result set.

## AIC Driver and Install Tool

The AIC Driver and Install Tool offers a number of functions related to the EZChrom configuration.

### Use the AIC Driver and Install Tool

- 1 From the Start menu, select **Start > All Programs > Agilent Technologies > OpenLAB CDS EZChrom Edition > AIC Driver and Install Tool**.

The screenshot shows a registration window titled "Agilent OpenLAB CDS EZChrom Edition - Registration". The window has a purple header and contains the following elements:

- OpenLAB Shared Services Credentials:** Fields for Username, Password, and Domain.
- Tabs:** AIC, Instrument Service Account, and Global.
- AIC Tab Fields:** AIC Name (AIC1), Hostname or IP (AIC1), Port (9753), and Description (Agilent Instrument Controller: AIC1).
- Checkboxes:**  Register Drivers and  Add Windows Firewall Exceptions.
- Button:** Register AIC.
- Footer:** net.pipe://localhost/Agilent/OpenLAB/

- 2 In the top area, enter the user name, password, and domain of the OpenLAB Control Panel Administrator. This is required to change the settings on any of the tabs.
  - On the **AIC** tab, you can register a server as an AIC, for example, in case of a fail over. For details, see [“Fail Over Procedures for OpenLAB CDS EZChrom Edition”](#) on page 114.
  - On the **Instrument Service Account** tab, you can retrieve or store the user who runs the instrument service. The user is recommended to

be a domain user with a password that never expires. The user must be part of the local administrator group on the AIC. If you want to store your acquisition data on a network share, this user must have access to the shared folder.

After adding or changing the instrument service account, reboot all AICs.

**NOTE**

An instrument service account can only support up to 64 AICs due to a Microsoft limitation.

- On the **Global** tab, you can retrieve the current Enterprise Path or register a specific Enterprise Path. This is relevant, for example, in the case of a fail over. For details, see [“Fail Over Procedures for OpenLAB CDS EZChrom Edition”](#) on page 114.

## Change the Password of the Instrument Service Account

It is recommended to use a password that never expires. However, if you need to change the password, follow the procedure below.

- 1 Stop all acquisitions and close any OpenLAB Control Panels.
- 2 Power off all Agilent Instrument Controllers (AICs) or Networked Workstations.
- 3 Change the password for the corresponding account in the customer domain.
- 4 From the Start menu, select **Start > All Programs > Agilent Technologies > OpenLAB CDS EZChrom Edition > AIC Driver and Install Tool**.
- 5 In the top area, enter the user name, password, and domain of the OpenLAB Control Panel Administrator. This is required to change the settings on any of the tabs.
- 6 Select the **Instrument Service Account** tab.
- 7 Enter the username, the new password, and the domain, then click **Store**.
- 8 Power on AICs or Networked Workstations.

## Change the Password of the AFS Administrator

Advanced File Security (AFS) is an optional configuration for OpenLAB CDS EZChrom Edition networked systems. It provides enhanced security on the Enterprise Path in order to prevent any unauthorized access to project data outside of the data system. This configuration sets the appropriate Windows sharing and security settings to allow only a defined group to access the enterprise data from Windows Explorer. This may only be configured if your system is configured to use Windows Domain as the Shared Services authentication provider.

It is recommended to use a password that never expires. However, if you need to change the password, follow the procedure below.

- 1** Stop all acquisitions and close any OpenLAB Control Panels.
- 2** Change the password for the corresponding account in the customer domain.
- 3** On any OpenLAB CDS EZChrom Edition client, browse to the directory where the software was installed.  
(by default: C:\Program Files\Agilent Technologies\EZChrom)

#### 4 Launch EnterpriseConfig.exe.

Enterprise Setup Login

WARNING: Before proceeding, make sure there are no CDS components running on the Enterprise machine, instrument controllers or any client PCs.

OpenLAB Control Panel Login

Server: net.pipe://localhost/Agilent/OpenLAB/

User name:

Password:

Domain:

Windows User Information

User name:

Password:

Logon from: Windows domain controller

Domain:

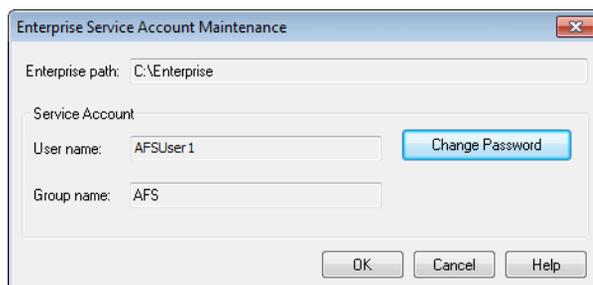
OK Cancel Help

- 5 In the **OpenLAB Control Panel Login** section enter the user name, password, and domain of the OpenLAB Control Panel Administrator.
- 6 In the **Windows User Information** section enter the user name and password of a user with edit permission to the enterprise path.
  - a If this is a domain user account, select **Logon from Windows Domain** and enter the domain name (recommended).
  - b If this is a local PC account, select **Windows Local PC**. This may only be an account local to the machine where the EnterpriseConfig.exe program is being run.
- 7 Click **OK**.

## 5 EZChrom-Specific Administration

### Change the Password of the AFS Administrator

8 In the **Enterprise Service Account Maintenance** dialog, click **Change Password**.



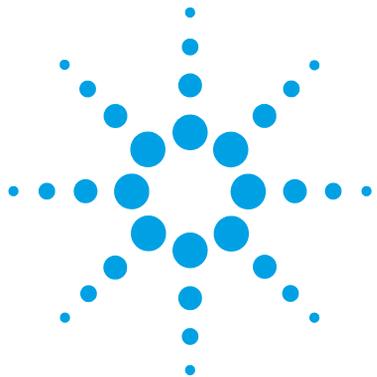
9 Confirm the warning.



10 Enter the *new* password.



11 Click **OK**.



## 6 Appendix

Privileges in OpenLAB Control Panel [126](#)

Project Privileges [127](#)

Instrument Privileges [136](#)

Administrative Privileges [136](#)

This chapter contains information on privileges used in OpenLAB CDS, and on driver license features for instruments of other vendors than Agilent.



## Privileges in OpenLAB Control Panel

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

- Everything
- System Administrator
- Instrument Administrator
- Project Administrator
- Instrument User

Additional default roles are available depending on the installed components. With an OpenLAB CDS ChemStation Edition installation, the following additional default roles are available:

- ChemStation Administrator
- ChemStation Lab Manager
- ChemStation Analyst
- ChemStation Operator

With an OpenLAB CDS EZChrom Edition installation, one additional default role **EZChrom Analyst** is available.

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

### NOTE

Roles are only visible if user authentication is required.

## Project Privileges

**Table 18** Project Management

Name	Description	Relevant for ...
View project or project group	User can see a project and project details but cannot edit. With Data Store: User can view the content of the Data Store repository. <i>Note:</i> This privilege is required for all users.	ChemStation EZChrom OpenLAB Data Analysis Data Store
Manage project or project group	User can create or edit project properties and can move the project but cannot access settings.	EZChrom OpenLAB Data Analysis Data Store
Edit content of project	User can create new versions of documents in Data Store.	Data Store
Manage project or project group access	User can view and edit the project access settings.	EZChrom OpenLAB Data Analysis

**Table 19** E-Signature

Name	Description
E-Signature Sign Data Files	User can sign data files (with ChemStation, this privilege is only relevant if you use OpenLAB Data Store).
E-Signature Revoke	User can revoke the e-signature (EZChrom only).

**Table 20** Intelligent Reporter templates

Privilege	Description
Edit Intelligent Reporter report template	Only relevant for OpenLAB ECM Intelligent Reporter: This privilege is required to edit report templates in the Reporter Client

**Table 21** Agilent OpenLAB Data Analysis

<b>Privilege</b>	<b>Description</b>
Edit sample information	User can edit information in the <b>Injection List</b> window.
Re-process data	User can reprocess injections.
Create new method	User can create a new processing method or save a method under a new name.
Save processing method	User can save an existing processing method.
View/edit integration parameters	User can view and edit the parameters in the <b>Integration Events</b> section of a method.
Edit identification parameters	User can edit the parameters in the <b>Identification</b> section of a method.
Edit calibration parameters	User can edit the parameters in the <b>Calibration</b> section of a method.
Edit report templates	User can create or edit report templates in the <b>Reporting</b> view.
Lock/unlock report items	User can lock or unlock items in report templates.
Do manual compound identification	User can manually assign a compound to a peak.
Do manual integration	User can activate manual integration in the <b>Chromatograms</b> window.

### ChemStation-Specific Privileges

**Table 22** ChemStation: Control

<b>Privilege</b>	<b>Description</b>
Run Acquisition	Start acquisition (single sample or sequence).

**Table 23** ChemStation: Data

<b>Privilege</b>	<b>Description</b>
Delete Data	User can delete data files in ChemStation Explorer.
Load data using not configured path	If Secure File IO is enabled, this privilege is required for ... <ul style="list-style-type: none"> <li>• loading data from unprotected paths</li> <li>• adding paths that are not empty to the ChemStation (<b>Preferences</b> dialog)</li> </ul>
Manual Integration	User can perform manual integration.
Save data to storage	Interactive saving of data to the central data storage system.

**Table 24** ChemStation: Instrument

<b>Privilege</b>	<b>Description</b>
Modify instrument configuration	User can modify the instrument configuration parameters.

**Table 25** ChemStation: Logbook

<b>Privilege</b>	<b>Description</b>
Clear Logbook	Clear the current logbook.
Save Logbook	Save the current logbook.

**Table 26** ChemStation: Method

<b>Privilege</b>	<b>Description</b>
Edit calibration table	Create and modify the calibration table; change calibration settings.
Delete method	Delete a method in ChemStation Explorer.
Edit integration events	Modify integration events and perform Auto Integration.
Edit ion labels	Edit options for ion labels (LC/MS only).
Edit system suitability	Edit noise ranges and performance limits.
Enable audit trail	Enable the audit trail for a specific method.
Load/Save method using not configured path	If Secure File IO is enabled, this privilege is required for ... <ul style="list-style-type: none"> <li>loading a method from an unprotected path</li> <li>saving a method to an unprotected path</li> <li>adding paths that are not empty to the ChemStation (<b>Preferences</b> dialog)</li> </ul>
Modify instrument method	Modify instrument method parameters.
Modify method properties	Modify Run Time Checklist and Method Information.
Perform method recalibration	Perform interactive recalibration.
Save method changes	Save method changes (includes Update Sequence/Master Method in Data Analysis view).

**Table 27** ChemStation: Report

<b>Privilege</b>	<b>Description</b>
Preview/print report	User can preview/print a report.
Modify report	User can modify report calculation/print style and can edit Instrument Curves dialog.
Lock/unlock report template items	Only relevant with Intelligent Reporting: User can lock or unlock report items and composite groups in a report template.

**Table 28** ChemStation: Security

<b>Privilege</b>	<b>Description</b>
Break session lock	Unlock a ChemStation session locked by other users.
Command line	Turn on / off the command line
Manage transfer queue	Access to the Transfer Queue and the Queue Management.
Modify storage transfer preferences	Modify transfer settings for data upload to the central data storage system.
Take over ChemStation Remote Session	Only relevant for ChemStation in a Distributed System: User can take over a running remote session.

**Table 29** ChemStation: Sequence

<b>Name</b>	<b>Description</b>
Change priority in run queue	User can submit samples or sequences to the front of the queue, and change the order of items in the run queue.
Delete entries from run queue	User can remove queued samples or sequences from the run queue.
Delete sequence	User can delete sequences in ChemStation Explorer.
Edit sequence summary	User can modify sequence summary report and extended statistics settings.
Load/Save sequence template using not configured path	<p>If Secure File IO is enabled, this privilege is required for ...</p> <ul style="list-style-type: none"> <li>• loading sequence templates from unprotected paths</li> <li>• saving sequence templates to unprotected paths</li> <li>• adding paths that are not empty to the ChemStation (<b>P</b>references dialog)</li> </ul>

## 6 Appendix

### Privileges in OpenLAB Control Panel

**Table 29** ChemStation: Sequence

Name	Description
Reprocess	User can reprocess a sequence.
Save sequence template	User can save sequence templates locally (on workstation or AIC).

**Table 30** ChemStation: View Access

Privilege	Description
Access Data Analysis view	User has access to Data Analysis view.
Access Diagnostic view	User has access to Diagnostic view.
Access Method & Run Control view	User has access to Method and Run Control view.
Access retention time lock	User has access to the Retention Time Lock menu (GC only).
Access retention time search	User has access to Retention Time Search menu (GC only).
Access Review view	User has access to Review view
Access Tune view	User has access to Tune view (LC-MSD ChemStation only).
Access Verification view	User has access to Verification (OQ/PV) view.
Access Report Layout view	User has access to Report Layout View, ability to create/edit/save report templates.
Enable Batch view	Enables all operations in Batch view.

## EZChrom-Specific Privileges

**Table 31** EZChrom: Method

<b>Privilege</b>	<b>Description</b>
Save Method	Save a method file. Create a new method (requires one or more method privileges listed below.)
Edit Method Properties	View or change the method properties (Description, Options, Calibration, Audit Trail)
Open Method	Open a method file.
Instrument Setup	View or change the method instrument setup parameters.
Integration Events	View or change method integration events.
Peaks Groups	View or change the Peaks and Groups table
Advanced Parameters	View or change advanced method parameters (Export, Custom Parameters, Column/Performance, Files, Advanced Reports)
Custom Reports	Open, edit, save custom method reports
System Suitability	View or change system suitability parameters
Review Calibration	Allows access to the review calibration window
Calibrate	Run a calibration sample to update the method calibration. In order to create or modify the calibration parameters in a method, the user must also have the "Save Method" privilege assigned
Edit Running Method	Allows to edit a running method
Update Master Method	Allows update of the original master method when reviewing results in a container

**Table 32** EZChrom: Data

<b>Privilege</b>	<b>Description</b>
Save Data	Allows to save a data file
Open Data	Allows to open a data file
Edit Properties	View or change the data properties description
Manual Integration Fixes	Allows to add or change manual integration fixes

**Table 33** EZChrom: Sequence

<b>Privilege</b>	<b>Description</b>
Open Sequence	Allows to open a sequence
Sequence Save	Allows to save a sequence file
Process Sequence	Allows to process a sequence
Sequence Properties	Allows to view and edit sequence properties
Edit Any Running Sequence	Allows to edit any running sequence
Edit Own Running Sequence	Allows to edit own running sequence

**Table 34** EZChrom: Sample Prep

<b>Privilege</b>	<b>Description</b>
Open Sample Prep	Allows to open an autosampler sample prep file
Save Sample Prep	Allows to save an autosampler sample prep file
Edit Sample Prep Properties	Allows to view or edit autosampler sample prep properties
Edit Running Sample Prep	Allows to edit a sample prep currently used (running)
Update Master Sample Prep	Allows update of the original master sample prep

**Table 35** EZChrom: Control

<b>Privilege</b>	<b>Description</b>
Preview Run	Allows to preview a run
Single Run	Allows to perform single run acquisition
Sequence Run	Allows to perform sequence acquisition
Print Setup	Allows to set up a printer for the current instrument
Manual Control (idle only)	Allows to access manual control functions while the instrument is idle
Manual Control	Allows access to manual control functions whether or not instrument is idle
Monitor All Data	Allows to monitor real time plot from any running instrument
Abort Any Run	Allows to abort any running sequence or single run

**Table 36** EZChrom: Standard Reports

<b>Privilege</b>	<b>Description</b>
Add Standard Report Template	Allows to add a standard report template

**Table 37** EZChrom: Advanced/Intelligent Report

<b>Privilege</b>	<b>Description</b>
Open Advanced/Intelligent Report	Allows to open an advanced/intelligent report template
Save Advanced/Intelligent Report	Allows to save an advanced/intelligent report template
Lock/unlock report template items	Allows to lock and unlock report template items (tables, chromatograms, groups of items,...) to control who is allowed to modify those

## Instrument Privileges

**Table 38** 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 39** System Administration

Name	Description
Manage printers	Can add/remove printers and print server.
Edit activity log properties	Can change the Activity log Settings in OpenLAB 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).
Manage Security	Can change security settings. Can edit (add, change etc) users, groups and roles. <i>Note:</i> A user with this privilege can grant himself access to all settings in OpenLAB Shared Services. Be careful who you grant the Manage Security privilege.

**Table 39** System Administration

Name	Description
Manage instrument controllers	Can edit AIC configuration, manage the AICs in the Agilent OpenLAB Server Utility tool.
Unlock any locked UI	Can login into any locked portal or instrument session (will be a re-login), even if privately locked.
Manage Data Store admin settings	User can access the <b>Admin</b> sub folder of the <b>Applications</b> root note in Data Store.

**Table 40** Data Store

Name	Description
Archive content	User can archive the content of the Data Store repository.

# Index

## A

add-ons 24  
 administration tool 76, 77  
 administrative privileges 61  
 after acquisition 81  
 after any data modification 81  
 after reprocessing 81  
 architecture 10  
 archiving 42  
 audit trail 42, 83  
 authentication provider 53

## B

back-up 42  
 break session lock 80

## C

ChemStation administration tool 77  
 ChemStation  
   administration tool 76  
   product structure 23  
 cleanup data 82  
 client services profile 84  
 compliance 85  
 Configuration Checker 90  
 configuration profiles 84  
 Control Panel 8  
 core module 23  
 counted licenses 25  
 CSAdministrators 77

## D

Data Analysis 8, 11, 13

data integrity 85  
 Data Store 42  
 diagcab file 90  
 diagnostics profile 85  
 diagnostics 51  
 Distributed System 14

## E

ECM authentication 53  
 ECM 42  
 electronic signature 42  
 encryption 79  
 Enterprise Path 121  
 EZChrom 8

## F

fail over workstation 108  
 fail over  
   ChemStation 91  
   EZChrom 114  
 Flexera 39  
 floating licenses 24  
 folder protection 85

## G

group 77  
 groups 60

## I

image 79  
 import after reprocessing 81  
 instrument management 46  
 instrument

drivers 24  
 privileges 61  
 service account 120  
 internal authentication 53

## L

lab status at a glance 46  
 license file  
   add 48  
   create 48  
   get 48  
   monitor 48  
   remove 48  
 license server 48  
 license  
   failover 101  
   features 28  
   management 48  
   scheme 25  
   server 39  
   types 24  
 local configuration 98  
 locations tree 46  
 lockout  
   break session lock 80  
   inactivity time 57  
   lock button 80  
   lock time 57  
   non-private 79  
   private 79  
   time-based 80, 80  
 login  
   maximum unsuccessful attempts 56

**M**

- manage queue 82
- mapping
  - groups 60
  - users 59
- mixed scenario 47

**N**

- Networked Workstation 12
- non privately locked 79

**O**

- OpenLAB Control Panel 8
- OpenLAB Data Analysis 8, 11, 13
- OpenLAB Shared Services 8

**P**

- password
  - expiration date 56
  - maximum unsuccessful login attempts 56
  - minimum length 56
- privately locked 79
- privileges
  - administrative 61
  - for individual nodes 62
  - instrument 61
  - project 61
  - roles and p. 60
- product structure 23
- projects 46

**R**

- role
  - Everything 60
  - type 60

**S**

- scalability 10
- Secure File IO 85
- security policy 56
- security 42
- session lock 79
- shared licenses 25
- Shared Services server 11
- Shared Services 8
- startup license 25
- storage 42
- SubscribeNet 48
- system activity log 50

**T**

- time based lock 80

**U**

- user
  - credentials 58
  - management 58

**V**

- VL 38

**W**

- Windows
  - authentication 53
  - service 39
- Workstation 10
- WTP 90

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

## In This Book

This manual contains information on the concepts of OpenLAB CDS, and the administration of OpenLAB CDS with the OpenLAB Control Panel. This Edition also includes information on the OpenLAB Data Analysis Add-on.

General concepts:

- System Architecture
- Licensing Strategy
- Security and Data Integrity

OpenLAB Control Panel

- Instrument Management
- License Management
- Log Files and Diagnostics
- Authentication Provider
- Security Policy
- User Management

© Agilent Technologies 2010-2016

Printed in Germany  
03/2016



M8305-90017