



Library Manager for OpenLab CDS

Installation and Configuration

Notices

Document Information

Document No: D0126153 Rev. A.00
Edition: 07/2025

Copyright

© Agilent Technologies, Inc. 2025

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

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

USA

Software Revision

This design specification is valid for revision 1.0 of
Library Manager for OpenLab CDS.

Warranty

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

Technology Licenses

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

Restricted Rights Legend

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

Safety Notices

CAUTION

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

WARNING

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

Contents

- 1 About Library Manager for OpenLab CDS 4**
- 2 Requirements 5**
- 3 Installation 6**
 - Install and Configure the Library Manager Web Application 7
 - Install the Library Manager Add-on for Data Analysis 9
 - Confirm Privileges 10
- 4 Configuration 12**
 - Library Administrator 13
 - Spectra Administrator 14
 - Spectra Contributor 15
 - Library Consumer 16
 - Chemist or Corresponding Default Role 17
- 5 Get Started 18**
 - Run the Library Manager Web Interface 18
 - Create First library 19
 - Configure Access Control 20
 - Use Library Manager 21



1

About Library Manager for OpenLab CDS

Library Manager for OpenLab CDS provides tools to organize and search UV spectra:

- A web interface offers the organization of spectra in libraries. Access control can be configured for each library.
- An Add-on for OpenLab CDS provides tools to search for similar spectra from Data Analysis or during unattended processing.

Spectra from all UV/Vis detectors are supported.



2 Requirements

Supported OpenLab CDS systems:

- Workstation, Workstation Plus, and Client/Server systems.
- OpenLab CDS rev, 2.8 Update 08 or higher

Supported browsers:

- Edge (current version)
- Chrome (current version)

Required .NET framework on the server and on workstations:

- .NET 8

Users and privileges:

- Installation user: Windows user with local administrative privileges
- If user authentication is enabled in the OpenLab CDS system, you need the credentials of the OpenLab CDS administrator.

3 Installation

The installation includes two parts, a web application to host the Library Manager web interface, and an add-on for the Data Analysis client. In a Client/Server system, they are installed on different machines. The following table shows where you need to install them.

Table 1: Installation of Library Manager components

OpenLab CDS system	Web application	Add-on for Data Analysis
Workstation Workstation Plus	X	X
Shared Services server	X	-
Analytical Instrument Controller (AIC)	-	X
Client	-	X

Install and Configure the Library Manager Web Application

- 1 Copy the downloaded zip file to your local hard drive, and unzip its content.
- 2 Run **Agilent.OpenLab.LibraryManager.msi**.
- 3 The installation is performed in the background. No user entries are required.
- 4 From the Start menu, select **Library Manager Configuration Tool**.

The tool checks if the required .NET 8 components are installed. If a component is still missing in your Windows system, you are asked to download and install it. Use the provided link to download the relevant file. After installing it, run the configuration tool again.

The following .NET components are required

- .NET 8.0 Desktop Runtime
- ASP.NET Core 8.0 Runtime

- 5 Click **Test URL** to confirm the connection to your OpenLab system.
The button turns green if the test succeeded.
- 6 If user authentication is enabled in the OpenLab CDS system: Provide the user credentials of the OpenLab CDS administrator created on OpenLab CDS installation.
- 7 Choose the language in which the privileges for the Library Manager shall be shown in the Control Panel.

NOTE

You will not be able to change the language after installation!

In case your server operating system has a different language than used on your clients or instrument controllers: Select the language of your client systems to allow users to manage the new privileges in their language.

- 8 Click **Test Connection** to confirm the settings.
The button turns green if the test succeeded.

Installation

Install and Configure the Library Manager Web Application

- 9 Under **Data Repository**, it is recommended that you keep the default password. It is prefilled with the password that was set during OpenLab CDS installation under **Component credentials**.

If you set a new password, make sure to document it at a secure location.

NOTE

Passwords must adhere to the following general guidelines:

- Minimum 8 characters
- At least one uppercase letter
- At least one lowercase letter
- At least one number
- Only contains allowed characters:
 - Uppercase letters A-Z
 - Lowercase letters a-z
 - Numbers 0-9
 - Special characters: !@#\$\$%^&-+_ =

- 10 Click **Save**.

A success message confirms your configuration.

- 11 Close the configuration tool.

Wait two minutes until the library service is configured in the background, and the required database and other objects are created.

- 12 To confirm successful installation and configuration, select **(All apps) > Agilent Technologies > Library Manager Web UI**. If needed, log in as OpenLab CDS administrator.

The Library Manager opens.

Installation

Install the Library Manager Add-on for Data Analysis

Install the Library Manager Add-on for Data Analysis

- 1 Copy the downloaded zip file to your local hard drive, and unzip its content.
- 2 Run **Agilent OpenLab Data Analysis Library Manager Addon.msi**, and follow the instructions in the setup wizard.
- 3 To confirm successful installation, open Data Analysis. Load data that contains UV spectra. Confirm that the **Spectra Library Search** layout is available.

Confirm Privileges

After installing Library Manager, new OpenLab CDS privileges should be available. If user authentication is enabled in the OpenLab CDS system, follow the steps below to ensure that all privileges listed in table below are available.

- 1 Launch the Control Panel and navigate to **Administration**.
- 2 In the navigation pane, select **Roles**.
- 3 Click **Create Role**.
- 4 Confirm that the Library Manager privileges are available.

Table 2: Library Manager privileges

Privileges	Purpose	Dependencies
Role type Administrative , under Library Manager		
Access all libraries	Access all libraries independent of access control settings. Use for emergency situations.	Login to Library Manager
Archive library	In the web interface, move a library into the archive	Login to Library Manager
Archive library entry	In the web interface, archive a spectrum to exclude it from being searched	Login to Library Manager Create/edit a library
Create new library spectrum entry	In Data Analysis, add a new spectrum to a library	User configured as Editor for the relevant library via web interface
Create/edit a library	In the web interface, create a new library, and edit name and description of existing libraries	Login to Library Manager
Edit library access control	In the web interface, set users who can edit or view a library in Data Analysis	Login to Library Manager Create/edit a library
Edit metadata of an entry	In the web interface, edit the metadata values of a spectrum, compound, description, or structure.	Login to Library Manager Create/edit a library

Installation

Confirm Privileges

Privileges	Purpose	Dependencies
Import a spectrum	In the web interface, import a spectrum from CSV file to a library	Login to Library Manager Create/edit a library
Import/export a library	In the web interface, import or export an entire library	Login to Library Manager
Login to Library Manager	Log in to the Library Manager web interface	
Review activity log	In the web interface, review the Library Manager activity log	Login to Library Manager
View activity log	In the web interface, view the Library Manager activity log	Login to Library Manager
Role type Project , under Processing Method .		
Edit UV library search parameters	In Data Analysis, view and edit the parameters in the Spectral Analysis > UV Library Search section of a method.	Edit content of project
Role type Project , under Data Processing .		
Do manual UV library search	In Data Analysis, search manually for UV spectra in a library.	Edit content of project



4 Configuration

If user authentication is enabled in the OpenLab CDS system, you must configure Library Manager user roles.

Library Manager can be used in different ways. Users can be configured, for example, to only view the content of an existing library, to add spectra to a library, or to create new libraries.

Configuring library users always includes two aspects:

- Users need specific privileges. Privileges are configured in OpenLab Control Panel.
- Users must be added to the relevant libraries (**Access Control** in the Library Manager web interface)

The following sections describe typical library user roles.

Library Administrator

A library administrator is able to create new libraries, archive existing libraries, edit the library access security, and check the library manager activity logs.

Privileges in OpenLab Control Panel

In OpenLab Control Panel, create a role with the following Library Manager privileges. Add the role to the relevant users or user groups.

- Archive library
- Create/edit a library
- Edit library access control
- Import/export a library
- Login to Library Manager
- View activity log
- Review activity log

Access Control in the Web Interface

For each relevant library, library administrators must be given **Editor** access. When creating a new library, the creator automatically has **Editor** access.

Spectra Administrator

A *spectra administrator* is able to modify the content of existing libraries: add spectra from Data Analysis, import spectra from a file via the web interface, archive spectra, and edit spectrum description and compound information.

Privileges in OpenLab Control Panel

In OpenLab Control Panel, create a role with the following Library Manager privileges. Add the role to the relevant users or user groups.

- Archive a library entry
- Create new library spectrum entry
- Edit metadata of an entry
- Import a spectrum
- Login to Library Manager

To enable the user to search spectra within Data Analysis, also add a role that includes the **Project** privilege **Data Processing > Do manual UV library search**. It is included, for example, in the default role **Chemist**.

Access Control in the Web Interface

For each relevant library, spectra administrators must be given **Editor** access. For details, see [Configure Access Control](#) on page 20.

Spectra Contributor

A *spectra contributor* is able to add spectra from Data Analysis, and to search libraries within Data Analysis. He is not able to use the Library Manager web interface.

Privileges in OpenLab Control Panel

In OpenLab Control Panel, create a role with the Library Manager privilege **Create new library spectrum entry**. Add the role to the relevant users or user groups.

In addition, add a role that includes the **Project** privilege **Data Processing > Do manual UV library search**. It is included, for example, in the default role **Chemist**.

Access Control in the Web Interface

For each relevant library, spectra contributors must be given **Editor** access.

For details, see [Configure Access Control](#) on page 20.

Library Consumer

A *library consumer* is able to search libraries within Data Analysis. He is not able to use the Library Manager web interface, and he will not be able to add any spectra.

Privileges in OpenLab Control Panel

No specific Library Manager privileges are required for this user.

Ensure that the library consumer has a role that includes the **Project** privilege **Data Processing > Do manual UV library search**. It is included, for example, in the default role **Chemist**.

Access Control in the Web Interface

For each relevant library, library consumers must be given **Viewer** access.

For details, see [Configure Access Control](#) on page 20.

Chemist or Corresponding Default Role

The **Chemist** role is by default present in OpenLab CDS. This role includes privileges to create and modify the processing method in Data Analysis. Consider updating the **Chemist** role, or roles that you use instead.

Add the project privilege **Processing Method > Edit UV library search parameters**. This privilege allows users to view and edit the parameters in the **Spectral Analysis > UV Library Search** section of a method. Editing these parameters is required to prepare an automated UV library search during unattended processing.



5 Get Started

Run the Library Manager Web Interface

To run the Library Manager web interface, open <https://localhost/openlab/spectralibraries/> (or <https://servername/openlab/spectralibraries/>) in your browser.

Create First library

- 1 Log in to the Library Manager web interface as a library administrator.
- 2 In the toolbar, click **New Library** to create an empty new library. Alternatively, click **Import Library** to import an existing one from ChemStation or EZChrom.
- 3 In the **New Library** dialog, provide a library name. Optionally, add a description.
- 4 Click **Create**.

Configure Access Control

Library users must have access to an individual library to be able to search or edit it. This library access is configured in the Library Manager web interface.

- 1 In the Library Manager web interface, click the information button (i) for the new library.
- 2 Under **Access Control**, click + to add editors or viewers to this library.

Viewer means that users can search this library in either Data Analysis or in the web interface, but cannot modify its content.

Editor means that users can add spectra from Data Analysis or from the web interface, and can modify information on spectra and compounds in the web interface. Editors must have specific privileges (see **Configuration** on page 12).

Use Library Manager

For details on using the Library Manager web interface, please refer to the interactive online help.

For details on using the Library Manager functions in Data Analysis, please refer to OpenLab Help & Learning.



In This Book

This document provides instructions for installation and configuration of Library Manager for OpenLab CDS.

www.agilent.com

© Agilent Technologies Inc. 2025
Edition: 07/2025

Document No: D0126153 Rev. A.00

