



MassHunter BioConfirm Software **Installation Guide**

Notices

Document Identification

D0145962 A.00

May 2026

Copyright

© Agilent Technologies, Inc. 2026

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
www.agilent.com

Software Revision

This Guide is valid for the 14.0 revision or higher of Agilent MassHunter BioConfirm until superseded.

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	In This Guide	4
2	Preparation	5
	Prepare for installation	7
	Upgrading from BioConfirm 11 or 12.1	8
	Uninstall previous versions	10
3	Installation	11
	Install BioConfirm Software	12
	Run Software Verification	23
4	OpenLab Control Panel Initialization	25
	Getting and Installing the License	27
	To get a license from SubscribeNet	29
	To get a license offline	31
	Adding a License	32
	Create Initial Project	35
	Start Application	41



1

In This Guide

Use this guide to install or upgrade the MassHunter BioConfirm software.

This guide applies to the installation of:

- MassHunter BioConfirm 14.0 and later



2 Preparation

Prepare for installation	7
Upgrading from BioConfirm 11 or 12.1	8
Uninstall previous versions	10

MassHunter BioConfirm software has two possible installation configurations which support five possible application topologies:

- **Workstation** (standalone system with no content management)
 - Workstation configuration with CDS Acquisition or MassHunter Acquisition
 - Workstation configuration on a stand-alone DA only PC
- **Networked Workstation** (client computer with content management on a separate server)
 - Client/Server topology on CDS AIC
 - Client/Server topology on CDS Client PC

Prepare the MassHunter Workstation computer the same way, regardless of the configuration.

NOTE

A Networked Workstation or CDS client requires an OpenLab Server/ECM XT with OpenLab Shared Services. These server products provide a central secure file storage for scientific data.

CAUTION

For CDS Workstation as well as CDS Client/Server AICs and Clients, the CDS must be installed before installing MassHunter BioConfirm.

WARNING

Installing MassHunter Acquisition 12.2 and below together with MassHunter BioConfirm 14.0 on the same PC is unsupported and will render the system unusable. Recovery requires a complete PC reimage.

Prepare for installation

To simplify installation of the software, decide on Workstation or Networked Workstation configuration before continuing.

1. Create an account on the PC with local administrative privileges to run the installation.
2. For Networked Workstation, a separate network connection is required to connect to OpenLab Server/ECM XT or ECM 3.x with OpenLab Shared Services Server.
3. Set all required anti-virus exclusions as described in the [GC/MS and LC/MS MassHunter Workstation and Networked Workstation Requirements Guide](#).
4. Ensure that Windows Date & Time is set correctly for the local time zone before installation.

CAUTION

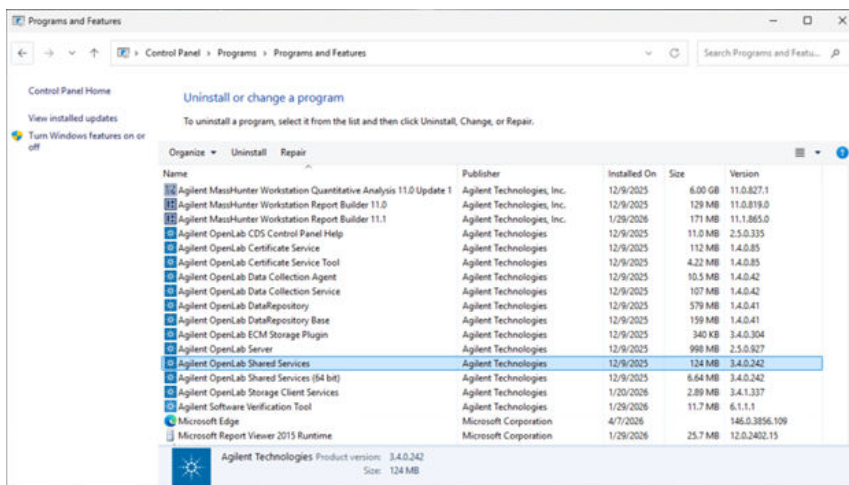
Following installation, it is not possible to modify the Date & time settings to a date and time earlier than the actual installation date and time. Doing so may have adverse effects on OpenLab Shared Services.

5. Check Agilent SubscribeNet (<https://agilent.subscribe.net.com/>) for the latest software updates or revisions for installation packages, software products, and instrument firmware/embedded software (ESW). For additional information, see [Firmware and Embedded Software \(ESW\) Updates](#).

Upgrading from BioConfirm 11 or 12.1

Follow this addendum when upgrading from BioConfirm 11 or BioConfirm 12.1 to BioConfirm 14 to verify the installed OpenLab Shared Services version and to complete required upgrade steps before beginning the BioConfirm 14 installation.

1. On the installation PC, determine the currently installed version of OpenLab Shared Services.
 - a. Open **Windows Control Panel**.
 - b. Select **Programs > Programs and Features**.
 - c. In the list of installed programs, locate OpenLab Shared Services and note the installed version number.



2. If OpenLab Shared Services version 3.4.x.x. is installed, upgrade OpenLab Shared Services to the intermediate 3.7 version before installing BioConfirm 14.0.

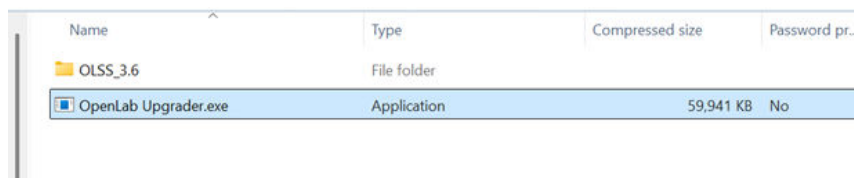
NOTE

If OpenLab Shared Services 3.7 is installed, no Shared Services upgrade is required. Continue with the [BioConfirm 14 installation](#).

- a. From the BioConfirm 14.0 installation media, run **OpenLab Upgrader.exe**.

Preparation

Upgrading from BioConfirm 11 or 12.1



Name	Type	Compressed size	Password pr...
OLSS_3.6	File folder		
OpenLab Upgrader.exe	Application	59,941 KB	No

- b. Complete the upgrade by following the on-screen prompts.
3. After the OpenLab Shared Services upgrade completes, install BioConfirm 14 using the [standard installation procedure](#).

Uninstall previous versions

Before installing software, it is necessary to uninstall previous versions of:

- MassHunter Acquisition for LC/MS TQ
- MassHunter Acquisition for Ultivo LC/MS
- MassHunter Acquisition for LC/MS TOF/Q-TOF
- MassHunter Quantitative Analysis software

Use Programs and Features in Windows Control Panel to remove Agilent MassHunter Workstation BioConfirm Software.

1. Close all MassHunter Workstation programs, including the Data Acquisition, Qualitative Analysis, Quantitative Analysis, and Report Builder programs.
2. Open Windows Control Panel.
3. Click **Programs and Features**.
4. Find and click the desired software to uninstall.
5. Click **Uninstall**.
6. Repeat steps 4 and 5 as needed.
7. Reboot the computer.
The computer is now ready to have the software installed.

WARNING

For Networked Workstations and CDS Client/Server only, ensure the OpenLab Server/ECM XT server is installed and available. If not, reach out to your local Agilent sales and support contact: <https://www.agilent.com/en/support>



3 **Installation**

Install BioConfirm Software	12
Run Software Verification	23

Install BioConfirm Software

CAUTION

If OpenLab CDS is required on the target PC, it must be installed before installing the MassHunter BioConfirm program. OpenLab CDS cannot be installed after BioConfirm; failure to install it first will prevent the use of OpenLab CDS on that system. For more information, reach out to your local Agilent sales and support contact: <https://www.agilent.com/en/support>

NOTE

The screenshots in this guide reflect a clean installation environment with standard instrument installation selections. While minor interface differences may occur depending on system configuration, the installation steps and workflow remain the same. Follow the on-screen installation wizard to complete the process.

NOTE

Before installing the MassHunter BioConfirm software, uninstall any previous versions. See [Uninstall previous versions](#).

CAUTION

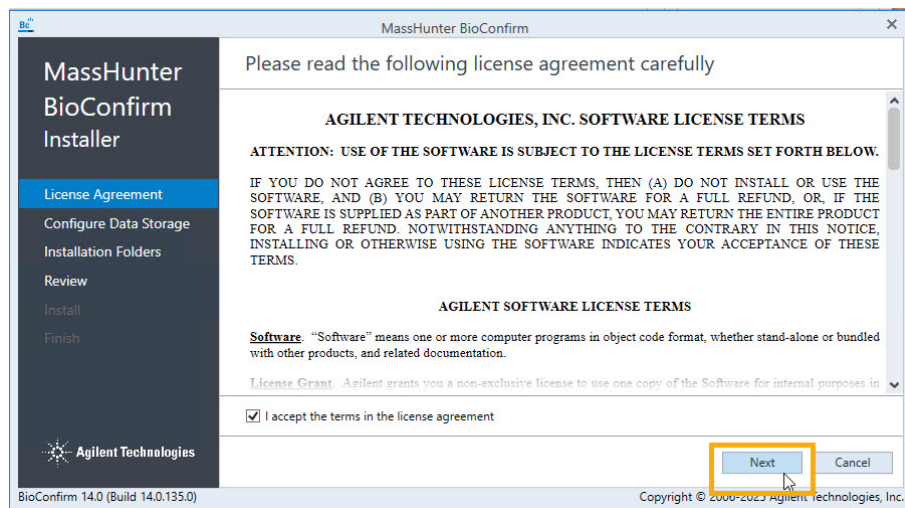
If the installation media was downloaded from SubscribeNet, mount the image file and copy all contents from the virtual drive to a local folder. All installations should be conducted from a local folder or physical media. Any required restarts, as described below, will unmount a virtual drive and result in an error requiring a PC re-image.

1. From the installation media, run the setup program **BioConfirm.exe**.
2. Read and agree to the **License Terms** and select the check box next to **I accept the terms in the license agreement**.

Installation

Install BioConfirm Software

- Click **Next**.



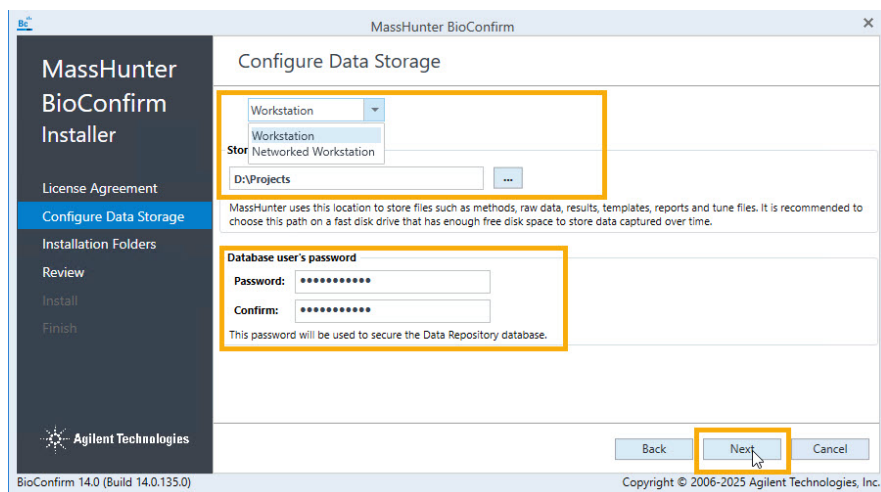
CAUTION

Once the Data Storage Configuration has been set and installed, this setting should only be changed by removing the software and deleting the OLSS databases that have been installed, then reinstalling.

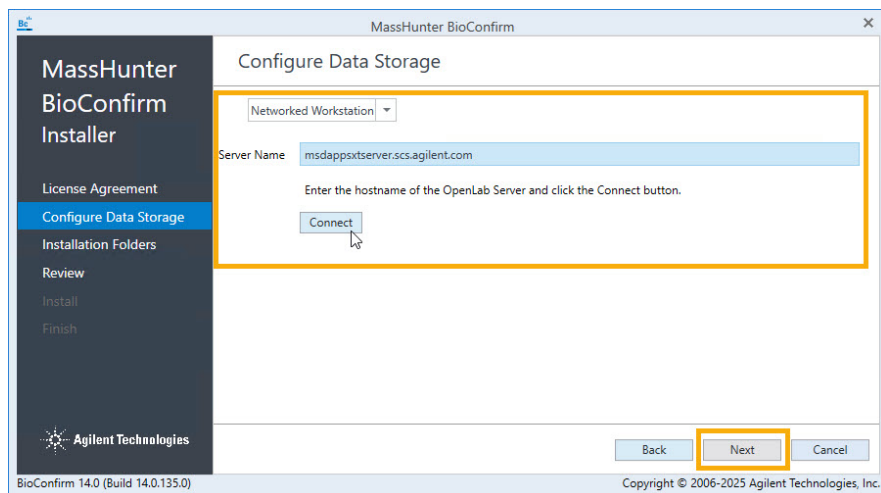
Re-imaging the PC is the recommended way to change the Data Storage Configuration.

- If OpenLab Shared Services has already been installed, skip to Step 5. If OpenLab Shared Services have never been installed on the PC, a prompt for the data storage type displays. Configure the data storage type as appropriate, and then click **Next**.
The choices are:

- **Workstation** - Files are stored on the local file system with no content management. See [Preparation](#) for information on topologies. All MassHunter files (such as data, methods, studies, report templates, and reports) are stored on the local file system. For Storage Path, enter or browse for the folder to store all user-generated data, including methods and reports. Enter a Password for use with the Datastore database that hosts the Compound Database Service, a core component of the Optimizer workflow. Make sure to remember it for future installations.



- **Networked Workstation** - Files are stored on an external OpenLab content management server. See [Preparation](#) for information on topologies. All MassHunter files (such as data, methods, studies, report templates, and reports) are stored on an external OpenLab Server/ECM XT server. For this configuration type, enter the name of the server to connect to, then click **Connect** to verify the connection to this server.

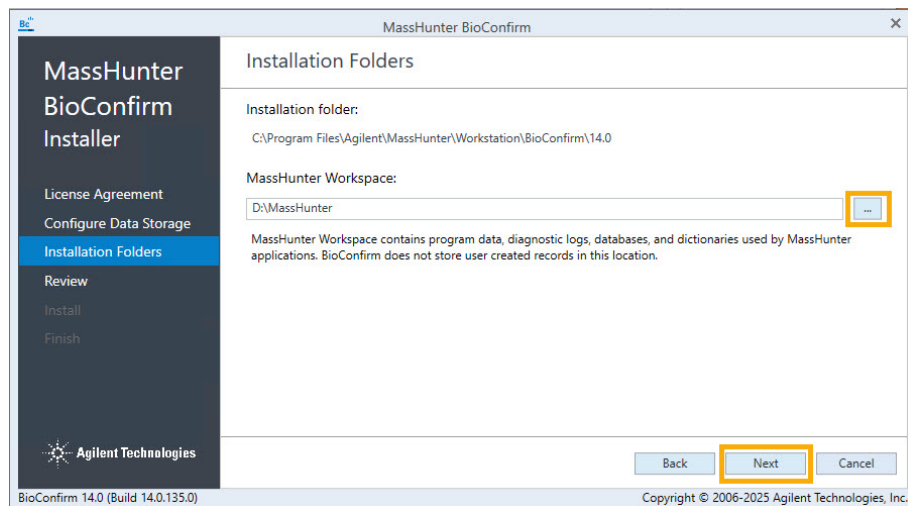


5. In the Installation Folders page, select the folders for the installation files and MassHunter Workspace, and then click **Next**. The workspace is used to store the instrument configuration, log files, the factory default methods, glycan

Installation

Install BioConfirm Software

databases, report templates and the Data Repository to store the Sequences and Chemical Data Dictionary information.



NOTE

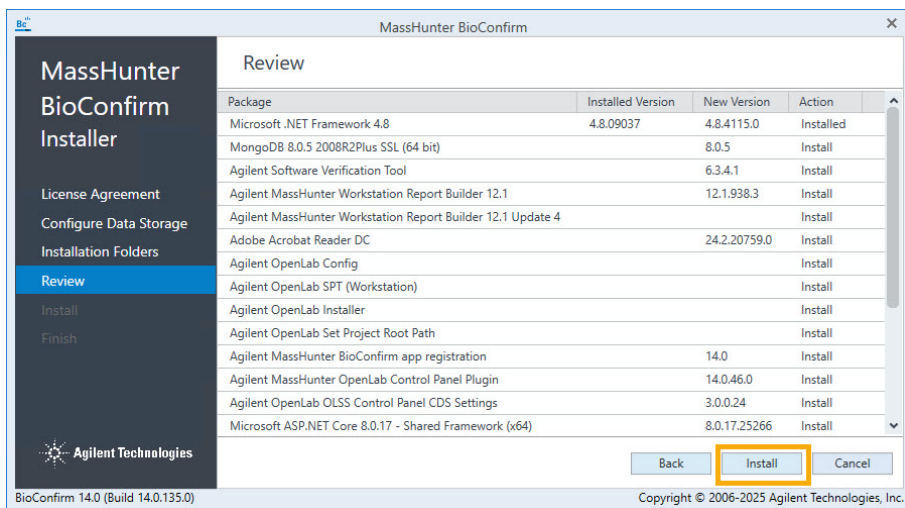
If a MassHunter application has been installed previously, the option to select a different MassHunter Workspace is disabled and cannot be changed.

Installation

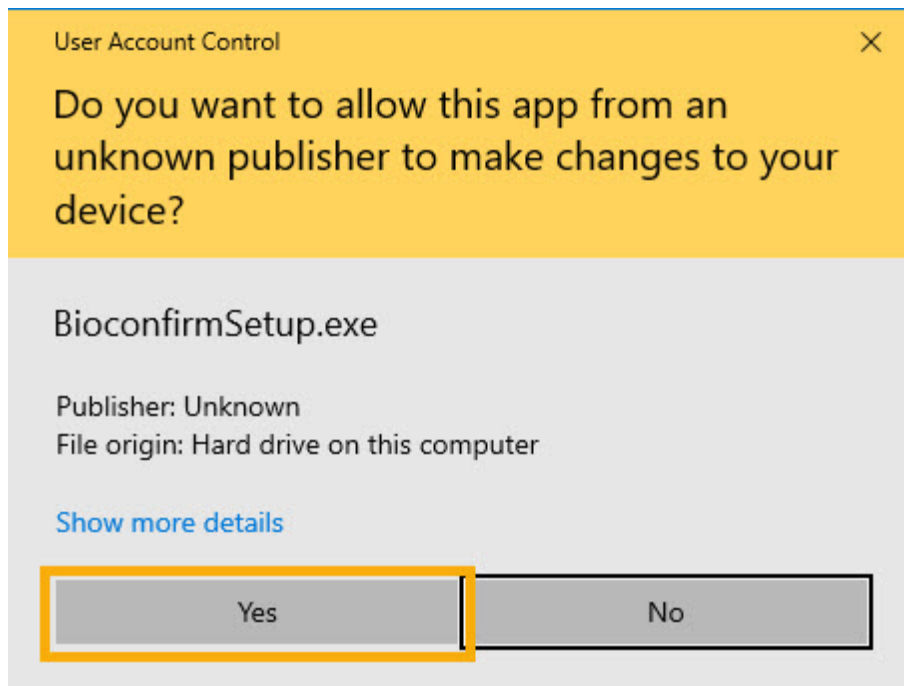
Install BioConfirm Software

- In the **Review** page, review the applications that will be installed, and then click **Install**.

The installation of the required packages starts.



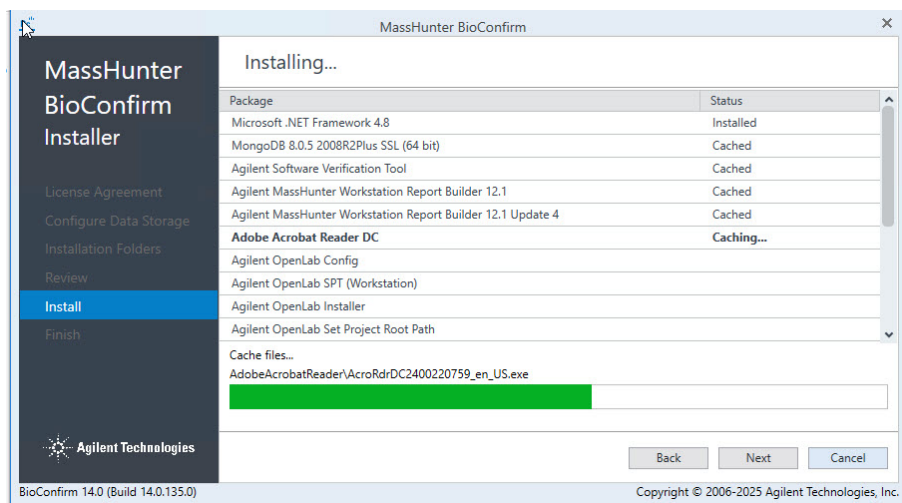
7. If prompted whether to allow the publisher to make changes to the device, click **Yes**.



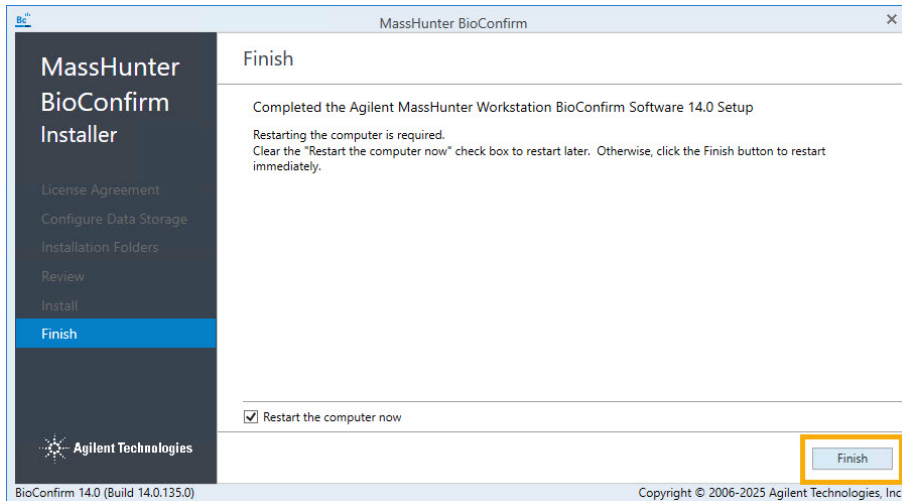
Installation

Install BioConfirm Software

The installation continues, displaying the progress of the installation.



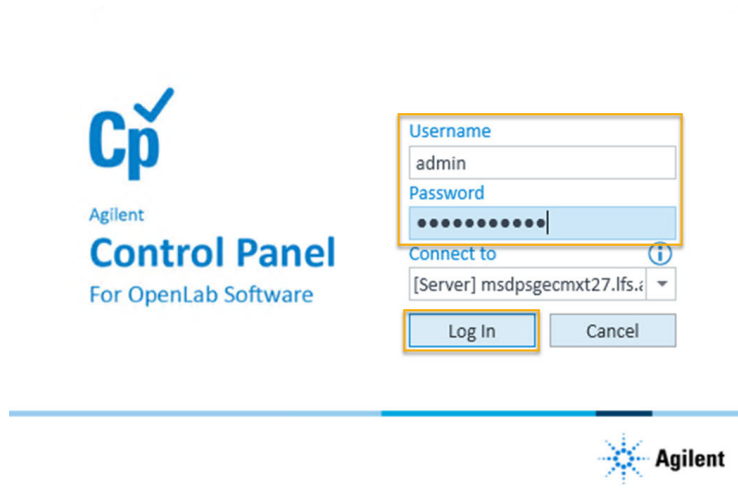
8. Click **Next** to continue the installation as prompted.



Installation

Install BioConfirm Software

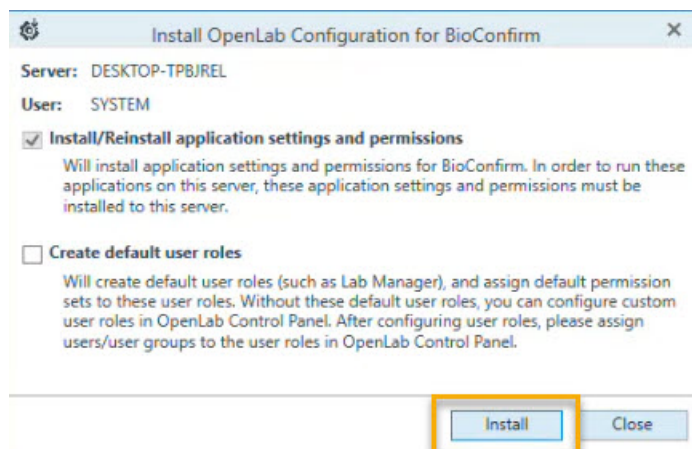
- If prompted, select **Restart the computer now** each time it displays. This restart can happen multiple times. The computer is restarted immediately after clicking **Next** or **Finish**.
- For networked workstations, if prompted during installation, enter the username and password for the user with administrative permissions on the ECM XT server, then click **Log In**.



- Click **Next** or **Finish** to continue the installation for each prompt.
9. Installation of OpenLab Configuration Bio Confirm for continues automatically. When completed, click **Install**.

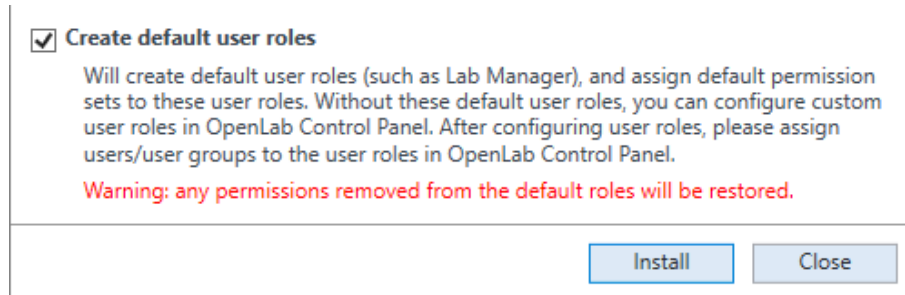
Installation

Install BioConfirm Software

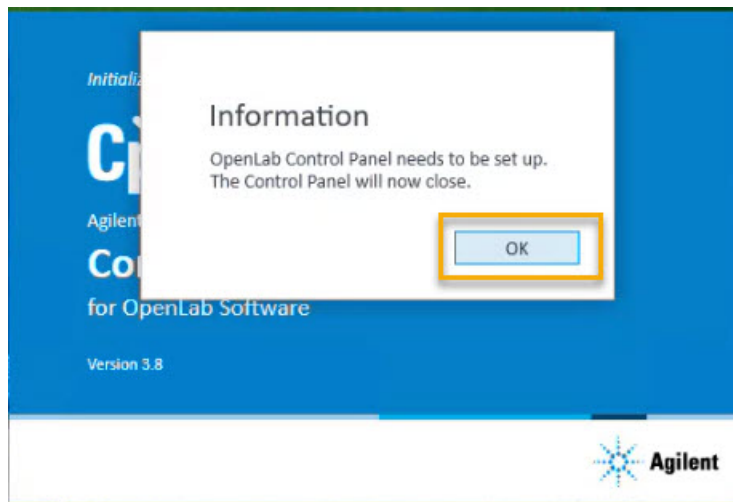


NOTE

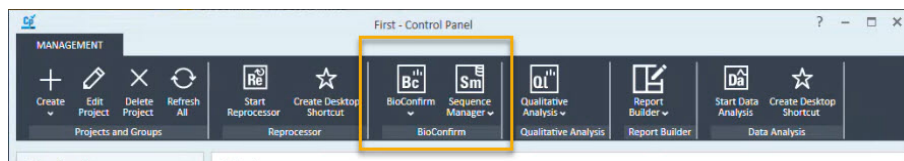
The application registration is only required once per workstation or server (Networked Workstation). Selecting **Create default user roles** will restore any permissions removed from the default roles.



10. Close and reopen OpenLab Control Panel to refresh the toolbar.



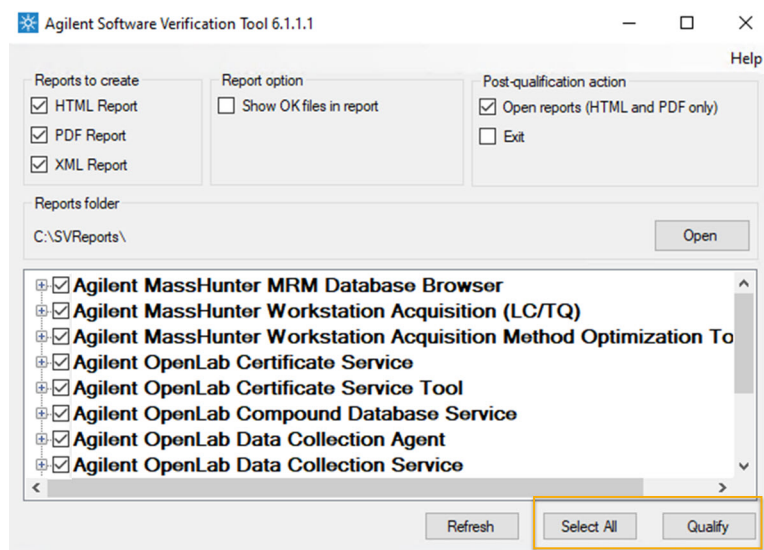
11. The BioConfirm icon appears in the Management tab of Control Panel, in the Project context.



Run Software Verification

The Software Verification Tool (SVT) provides evidence that the system has been installed correctly.

1. From the Windows Start menu, go to **Agilent Technologies > Software Verification Tool**.
2. Select the check boxes for each component that needs to be verified. Click **Select All** to qualify all components.
3. Click **Qualify**.
The system will run the application and generate a Software Verification Report.



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

Installation

Run Software Verification

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

Software Verification Report					
Date:	Wednesday, March 27, 2024	Time:	12:17:17 PM [UTC -07:00:00]	Host Name:	DESKTOP-655PFQM
Windows User Name :	admin	Base Revision Number:	1.7.0	Product Name :	Agilent OpenLab Data Collection Agent
Install Type:	Agilent OpenLab	Additional Packages:	NA		
Base Reference File Name : DCA_IQT_Ref.xml					
Summary :					
Overall Evaluation of Installation Check : PASS					
File Report Summary					



4 **OpenLab Control Panel Initialization**

Getting and Installing the License	27
To get a license from SubscribeNet	29
To get a license offline	31
Adding a License	32
Create Initial Project	35
Start Application	41

This chapter describes the steps needed or required to initialize the OpenLab shared services control panel when running software for the first time. If connecting a Networked Workstation to an existing OpenLab Server/ECM XT server/ECM 3.x server, some steps might already be done.

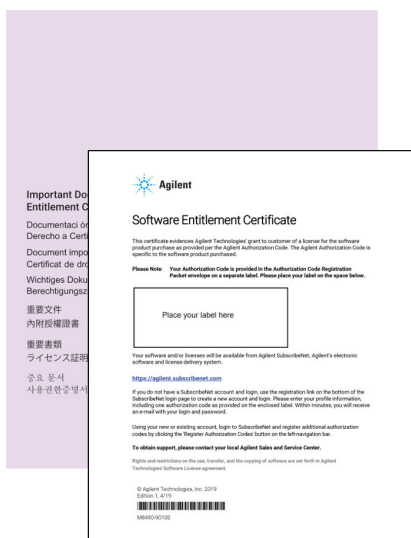
Getting and Installing the License

There are two ways to get a license:

- From SubscribeNet (<https://agilent.subscribe.net/>)
- Offline (from an Agilent representative)

To begin, locate the following information:

- The authorization code from the labels provided in the lavender envelope along with the Software Entitlement Certificate that had additional information on registering an authorization code.



If you did not receive a lavender envelope with the product, contact your vendor or internal support at subscribe.net_support@agilent.com.

OpenLab Control Panel Initialization

Getting and Installing the License

- The host name and MAC address of the license server.
 - For Workstations, this information comes from the PC where MassHunter BioConfirm is installed.
 - For Networked Workstations, this information comes from the server running OpenLab Server/ECM XT.

To retrieve the MAC address, open the Control Panel and browse to the **Administration > Licenses** section. Use the **Copy MAC Address** or **Save MAC Address** function to obtain the MAC address for license generation.

NOTE

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

NOTE

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

NOTE

Licenses can be obtained at a future date. The included startup license enables full use of BioConfirm programs for 60 days.

To get a license from SubscribeNet

An Internet connection is needed to access SubscribeNet. If you do not have one, see ["To get a license offline"](#).

1. If new to SubscribeNet:
 - a. Go to <https://agilent.subscribenet.com/control/agil/AgilRegisterToAccount> to register the product with SubscribeNet.
 - b. On the registration page, enter the authorization code from the label and complete the profile information (required fields are marked with an asterisk *).
The email address that you enter will become your login ID.
 - c. Click **Submit**. The system will generate and display an account name for you. SubscribeNet will send a welcome email with your login ID and password.
2. Go to <https://agilent.subscribenet.com/>.
3. Log in to SubscribeNet using your login ID and password.
After logging in, use the online user manual link for help with any questions.
4. As an existing SubscribeNet user:
 - a. If there is more than one user account, select the SubscribeNet account associated with this authorization code.
 - b. In the navigation pane, select **Register Authorization Code**.
Enter a new authorization code and make available the new license entitlements.
5. In the navigation bar on the left side of the page, select **Generate** or **View licenses**.
6. Follow the prompts to generate a new license.
Enter a name in the prompt for the HOST NAME of the computer. The host name entered must match with the network name of the computer where the Control Panel is running. Do not include any DNS suffix (*domain.com*) references in the entered machine name.

After the system generates the license, view its details, then click **Download License File**. Save the license file to a computer and to a backup location (such as a portable storage device).

Use the login ID and password when revisiting the Agilent SubscribeNet site to regenerate a license file, add new authorization codes, or further configure the license for a system.

To get a license offline

If you are unable to generate a license, contact your nearest Agilent technical support office. Agilent collects the necessary information from you to create a license account on your behalf. For phone support in your region, call the sales and service number for your region.

1. Collect profile information:

Your account name will be your company name and Lab name separated by a comma. Employee information provided here will be used to define the first administrator of your account for future access to the system as required.

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

2. Collect Authorization Code(s).

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

3. Contact your Agilent representative and provide the collected information for Agilent to generate a license through SubscribeNet.

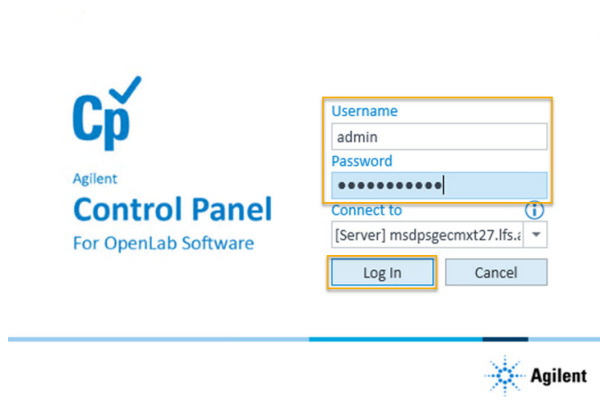
Once the license file is generated, it will either be sent to your shipping address, or your local Agilent-certified service professional will deliver it in person.

Adding a License

1. On the desktop, double-click the Control Panel icon.



For Networked Workstation, the server login dialog box is displayed. Log in as a user with an account on the OpenLab Server/ECM XT server/ECM 3.x server to continue.

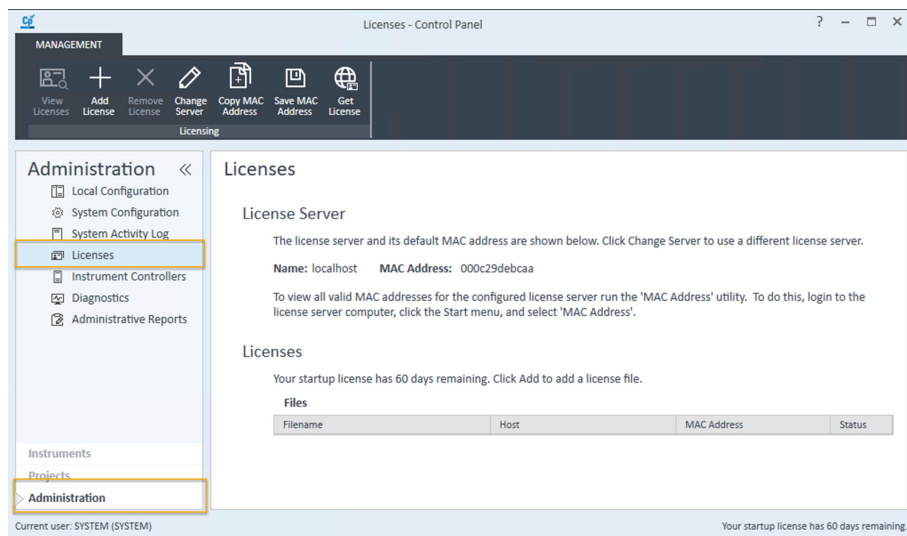


The Control Panel initializes and then opens.

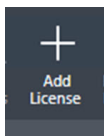
OpenLab Control Panel Initialization

Adding a License

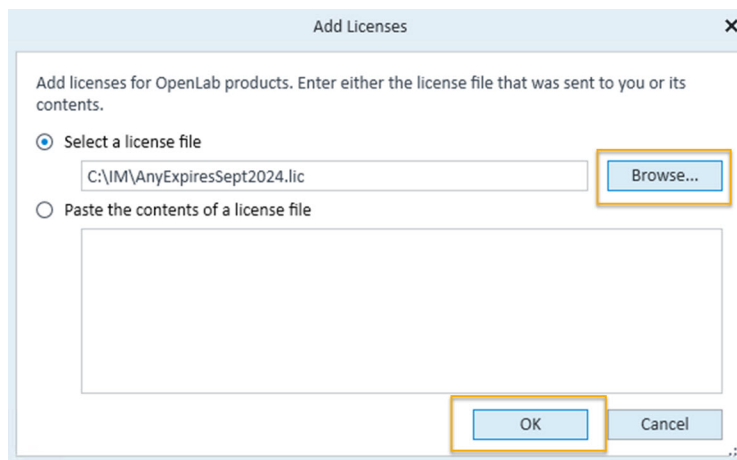
- In the navigation panel, select **Administration > Licenses**.



- Click **Add License**.



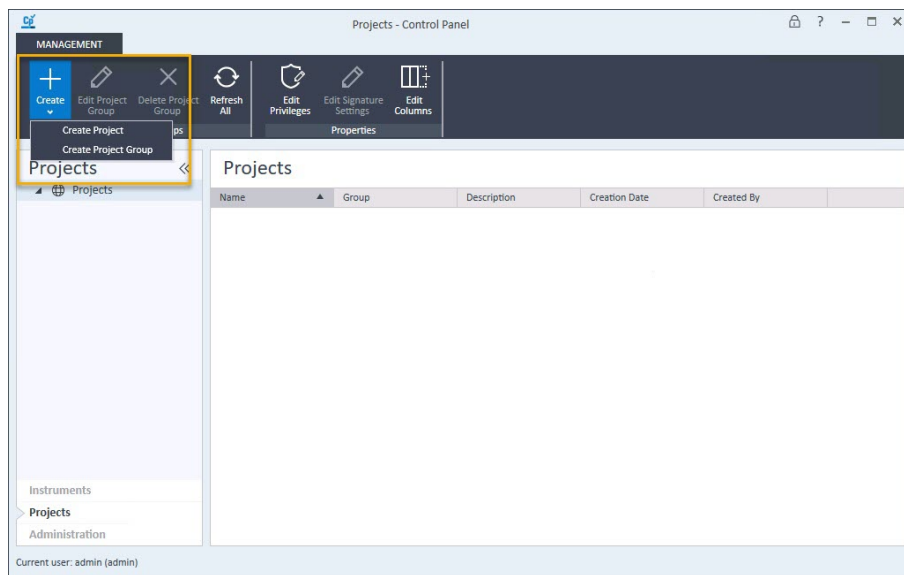
4. Browse to select a license file, then click **OK**.



Create Initial Project

In this example, the Storage path is set to C:\Projects.

1. In the navigation pane, select **Projects**.
2. On the ribbon, click **Create > Create Project**.



3. In the **Properties** tab, specify a **Name**, **Project folder path**, **Description**, and the **Application** (*OpenLab CDS* or *MassHunter Workstation*). Click **OK**.

WARNING

If the Project folder path is C:\ with no set subfolder, do not proceed with the remaining procedures in this section. If the project root is undefined, the installation will not work.

Projects - Control Panel

MANAGEMENT

Create Edit Delete Refresh
Project Project Project All

Properties Edit Privileges Edit Signature Settings

BioConfirm Sequence Manager Quantitative Analysis Report Builder

Projects <<

▲ Projects

Instruments

Projects

Administration

Current user: admin (admin)

Create Project

Properties MassHunter Workstation

Name: Example

Project folder path: C:\Projects\Example Browse

☐ Include project groups in project path

Description:

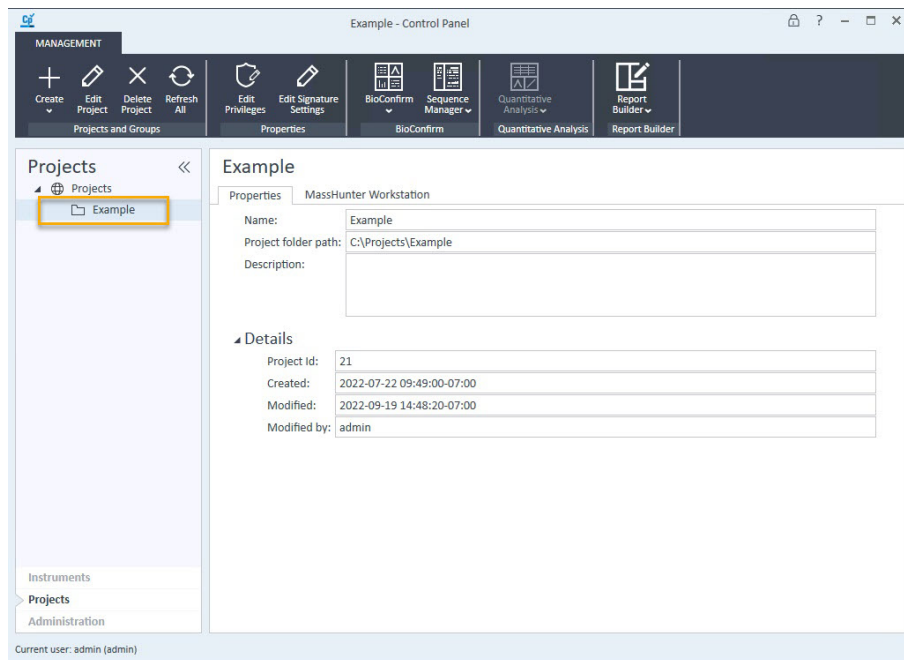
Applications: ☒ MassHunter Workstation

OK Cancel

OpenLab Control Panel Initialization

Create Initial Project

- In the navigation pane, select the **Project** that was just created and verify the values.

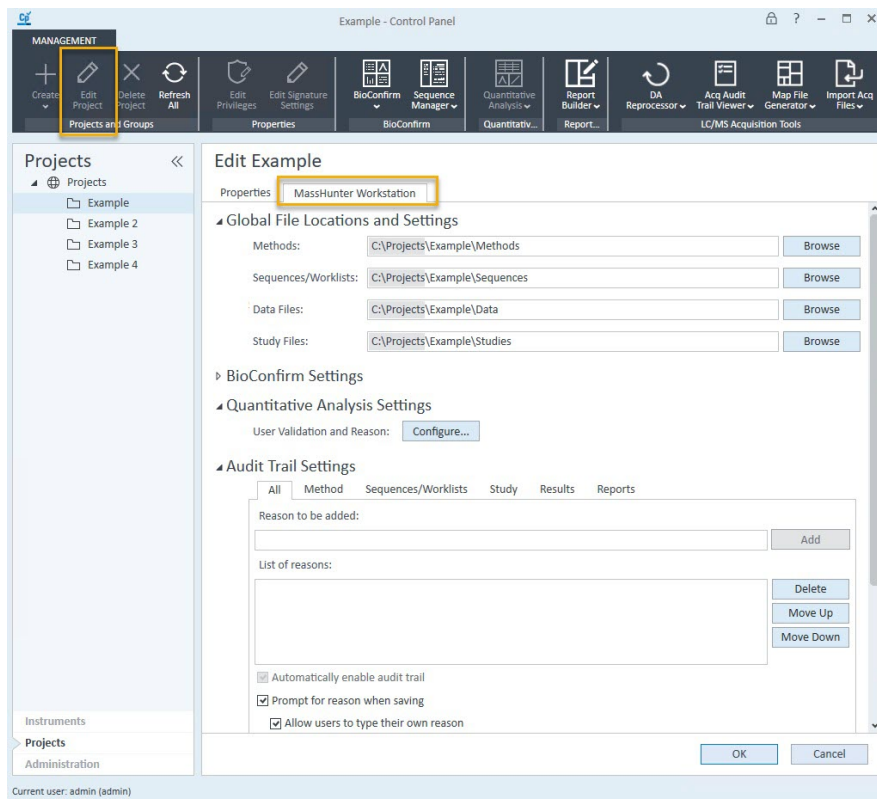


- On the Ribbon click **Edit Project**.
In the **MassHunter Workstation** tab:

OpenLab Control Panel Initialization

Create Initial Project

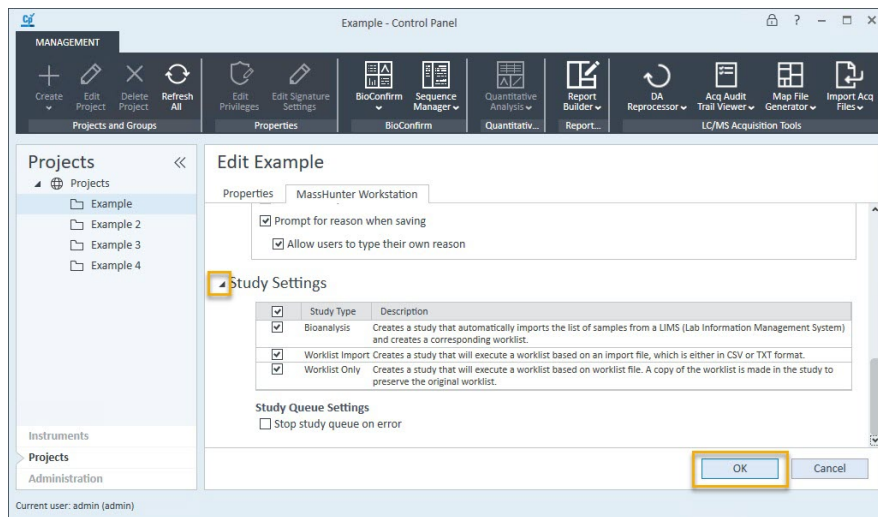
- Confirm the **Global File Locations** and **Settings**.
- In the **All** tab, select **Prompt for reason when saving** to require the user to enter a reason when saving MassHunter objects. The ability to require the user to enter a reason for specific MassHunter objects is available. If **Allow users to type their own reason** is selected, users are allowed to manually enter a reason.
- If needed for this project, under **Audit Trail Settings**, add default reasons for methods, worklists, studies, results, and reports on the appropriate tab. The ability to add general reasons to the **All** tab is available, which will display in all contexts.



OpenLab Control Panel Initialization

Create Initial Project

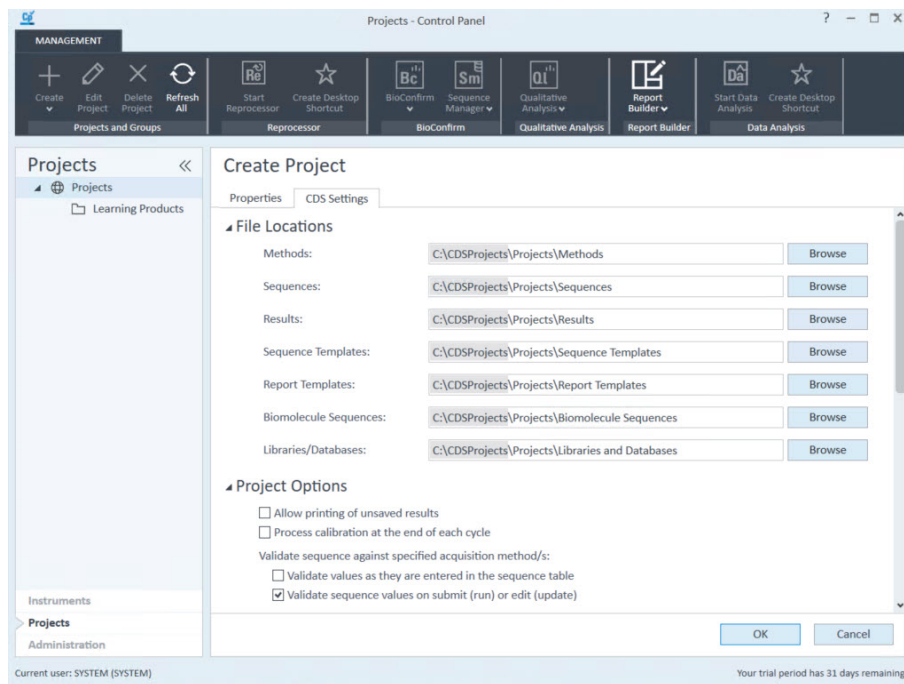
- Expand **Study Settings**. Several different types of studies are enabled by default. Mark or clear the study types that are used with this project.



In the CDS settings tab:

OpenLab Control Panel Initialization

Create Initial Project



- If needed for this project, under Project Options, adjust printing and validation options.
- If needed for this project, under **Audit Trail Settings**, add default reasons for methods, worklists, studies, results, and reports on the appropriate tab. The ability to add general reasons to the **All** tab is available, which displays in all contexts.

NOTE

See *Control Panel Online Help* for more information on creating Projects and field definitions.

6. Click **OK**.

Start Application

To start the BioConfirm program, select the project that you are using. If no projects exist, you need to first create the project.

1. Double-click the OpenLab Control Panel icon on the desktop or click **Start > Agilent Technologies > Control Panel**.
2. Click **Projects** in the left pane.
3. If needed, click **Create > Create Project** in Projects and Groups on the ribbon. All data files, methods, sequences, report templates, and databases are stored in a project. See the online Help for OpenLab Control Panel for more information.
4. Select the project to use in the left pane.
5. On the ribbon in the BioConfirm group, click **BioConfirm > Start BioConfirm** or click **BioConfirm > Create BioConfirm Shortcut** to create a desktop icon to use with the selected project.

www.agilent.com

© Agilent Technologies, Inc. 2026

D0145962 A.00

May 2026

