

Thermo BioPharma Finder 4.0 QF1 Installation Instructions

Pub. No. XCALI-98250 Rev. B

Follow the instructions in this document to install and license the Thermo BioPharma Finder™ software, version 4.0. You must be a system administrator on the installation computer to install the software. You may license the application as either an administrator or as a standard user.

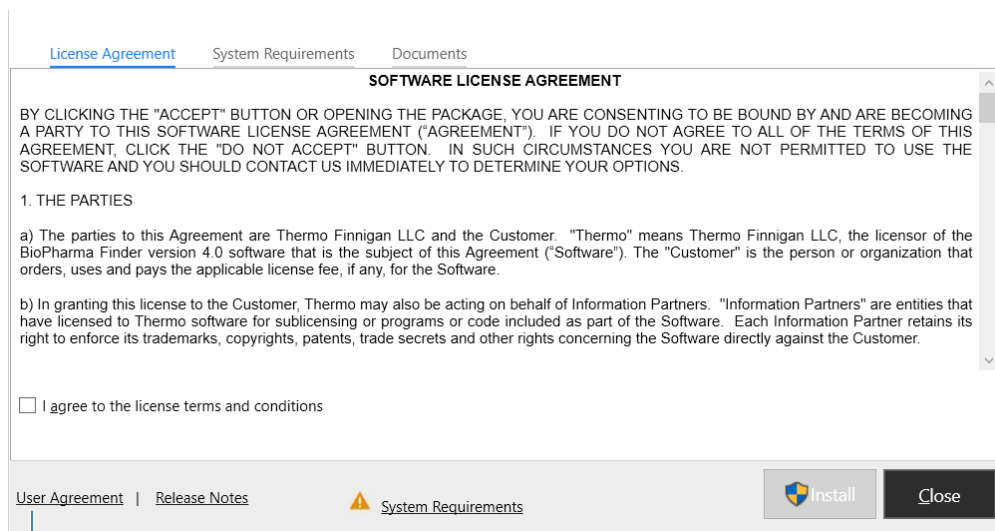
IMPORTANT! If you are upgrading from a previous version and currently have submitted jobs in the run queue, complete these experiments before you upgrade to BioPharma Finder 4.0.

Installing the BioPharma Finder software

The following procedure explains how to install the BioPharma Finder 4.0 software.

To install the BioPharma Finder 4.0 software

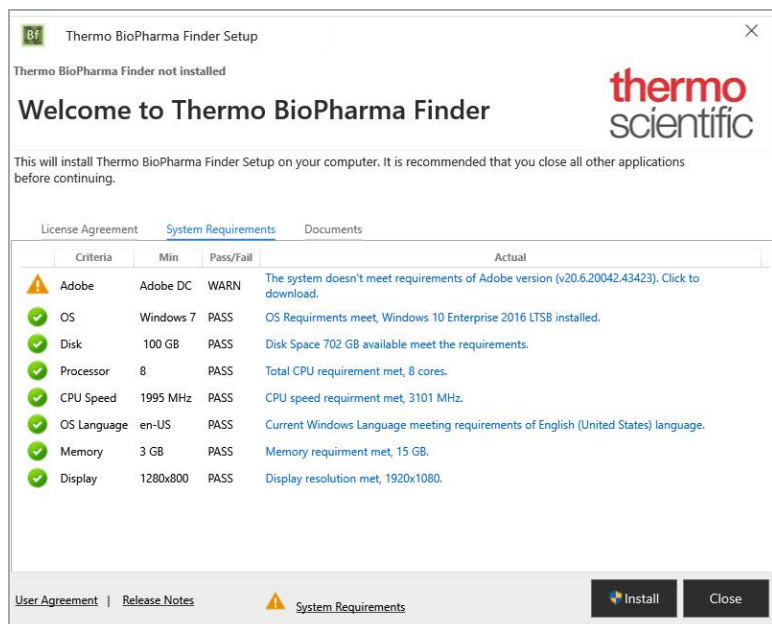
1. Close all the open Thermo Scientific applications.
2. Do one of the following:
 - Insert the BioPharma Finder 4.0 DVD into the DVD/CD-ROM drive. The Autorun program automatically launches the installer wizard. If the wizard does not start, navigate back to the DVD in Windows Explorer and double-click **BioPharma Finder 4.0QF1 - Installer.zip**.
 - Download and unzip the **BioPharma Finder 4.0QF1 - Installer.zip** from the Flexera Software download site: <https://thermo.flexnetoperations.com/control/thmo/index>. Then, double-click the **BioPharmaFinderSetup.exe** to launch the installer wizard.
3. Select the **User Agreement** link to view the terms and conditions in the end-user license agreement.



Opens the Software License Agreement

4. Select the **I agree to the license terms and conditions** check box.
5. In the installer wizard, click the **Documents** tab and click the links to view the Installation Instructions, Release Notes, User Guide, example data, and other documents.

6. In the installer wizard, click the **System Requirements** tab to check the system requirements.



The wizard checks for the minimum system requirements. If any of the specifications do not match, the system requirements list shows an error or warning and displays a summary report of any missing items. You can click the Adobe link to download and install the required software.

Note: If the Adobe Acrobat Reader DC application is not already installed on your system, click the link to install the Acrobat Reader software.

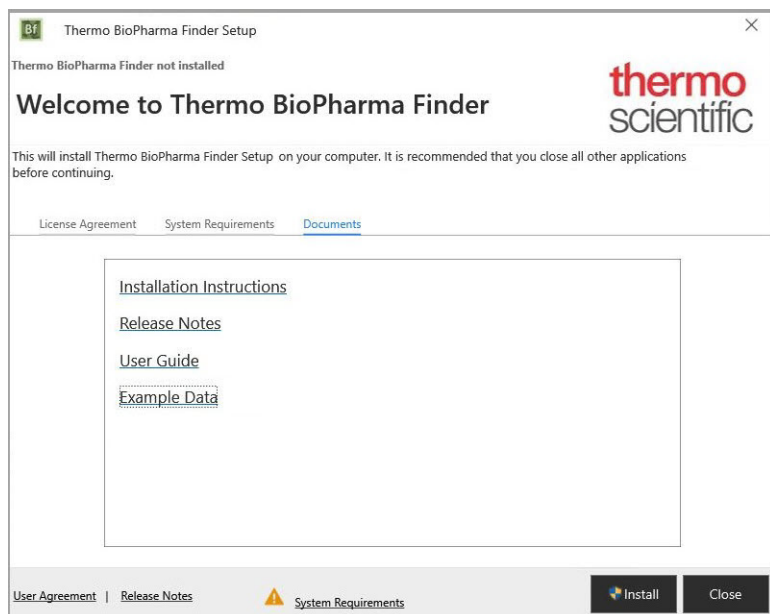
7. If you have previous versions of the BioPharma Finder application on your system, click **Yes** at the prompt to uninstall all legacy applications and convert your legacy data. If the installer determines that there is not enough space for the automatic backup, a message prompts you to delete files on your system to free storage space. Delete files as necessary, and then click **Retry** to retry the automatic backup.

During the installation, the BioPharma Finder installer automatically backs up your current database and other files to the `c:\ProgramData\ThermoScientific\` folder for BioPharma Finder-related files. If this is an upgrade from the previous version, the installer automatically creates a new BioPharma Finder **build version_backup_date-and-time** folder under the **ThermoScientific** folder (for example: 3.1.56.0_backup_2018_12_20_1418), and then moves the BioPharma and Databases folders down one level below the new backup folder.

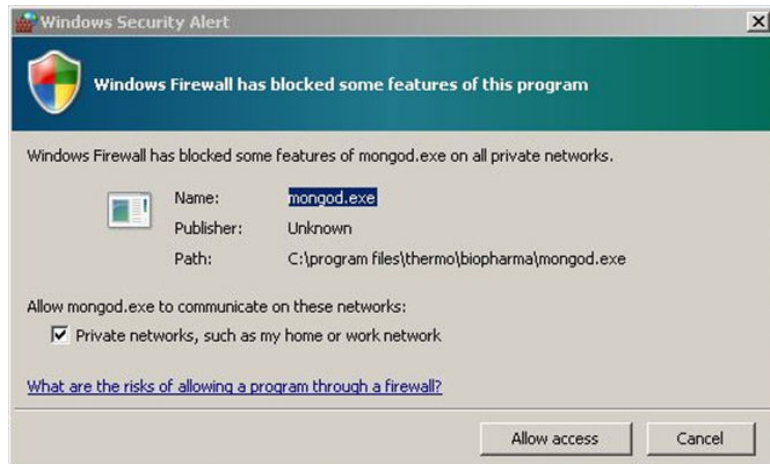
Note: If you have the Protein Deconvolution application, upgrade the software using the BioPharma Finder 3.2 or earlier version. You cannot upgrade the Protein Deconvolution application using the BioPharma Finder 4.0 version.

For more information about the data conversion process, refer to the topic “Converting Data from Legacy Applications” in the *BioPharma Finder User Guide*.

8. After you agree to the terms and conditions in the license agreement, click **Install** or **Upgrade** to install the BioPharma Finder 4.0 software.



If the following Windows Security Alert dialog box appears, click **Allow Access**.



If all the system requirements criteria are satisfied, the installer welcome page appears.

9. Click **Launch** after the installation completes to launch the application or close the installer wizard without launching the application and activate the license later.

Activating the license

To activate the BioPharma Finder 4.0 license

1. After installation is complete, start the application from the desktop icon or Start menu.
2. If you have a trial license and want to install the permanent license, choose **Help ► About BioPharma Finder**. Then, click **Activate** to display the Product Licensing wizard.

Follow the on-screen instructions in the Product Licensing wizard to install the permanent license.

For more information about both trial and permanent licenses, click **Help** on the License Activation page of the wizard.

3. Choose the appropriate licensing option:
 - For the demo version of the BioPharma Finder application, use the 60-day trial license.
 - For purchasers of the BioPharma Finder application, set a permanent license using the activation code provided through your Flexera account. You cannot use a license key from an earlier version.

Checking the language, keyboard, and location settings

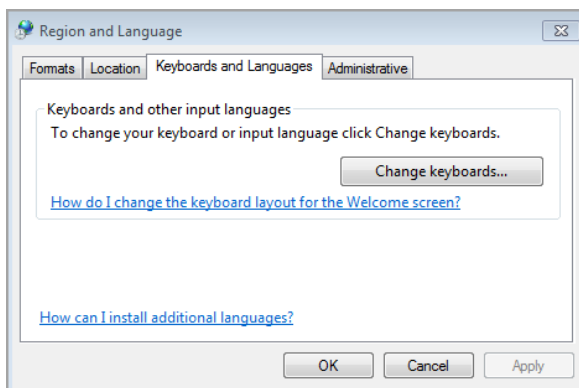
Before installing the software, check your language, keyboard, and location settings on your system and change them as necessary.

To check and change the settings

1. Choose **Start ▶ Control Panel** to open the Control Panel window.
2. Under the Clock, Language, and Region area, click **Change keyboards or other input methods**.



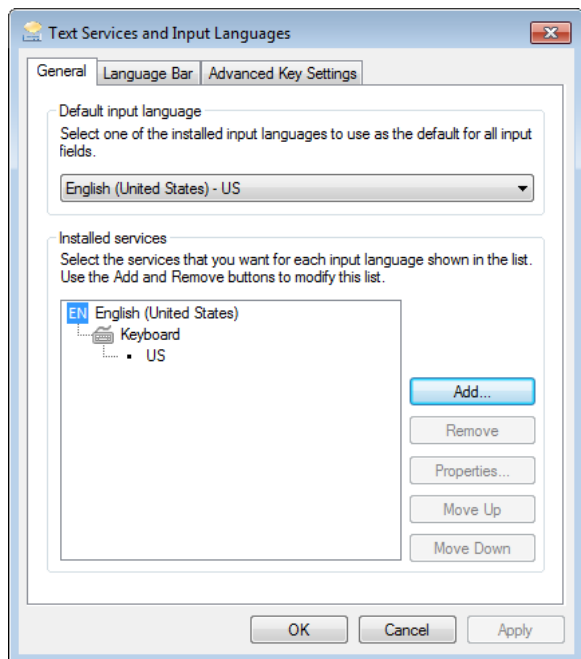
3. In the Region and Language dialog box, click the **Keyboards and Languages** tab, and then click **Change Keyboards**.



4. In the Text Services and Input Languages dialog box, click the **General** tab.

Make sure that the Default Input Language area shows *English (United States)- US*. Otherwise, select this setting from the list.

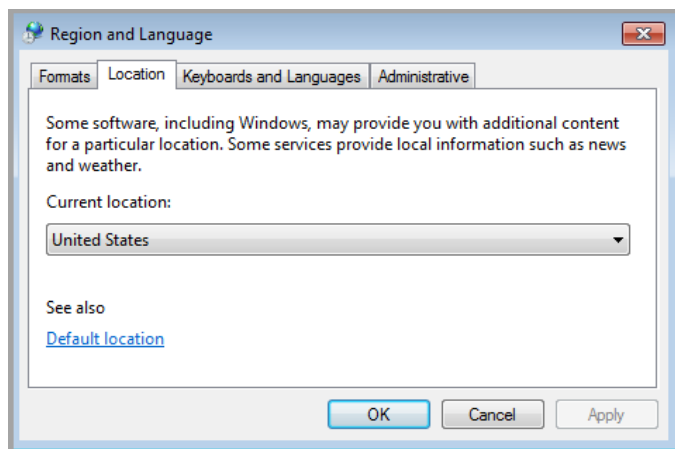
Make sure that the Installed Services area shows *English (United States) > Keyboard > US*. Otherwise, click **Add** and select this setting from the list, and then click **OK**.



5. Click **OK** to close the Text Services and Input Languages dialog box.

6. In the Region and Language dialog box, click the **Location** tab.

Make sure that the Current Location shows *United States*. Otherwise, select this setting from the list.



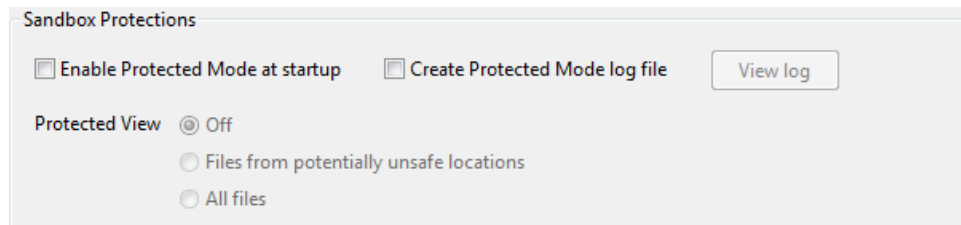
7. Click **OK** to close the Region and Language dialog box.

8. Close the Control Panel window.

Clearing the security setting to view the installation report

(For systems running Microsoft Windows 10) To view the report at the end of the installation, you must clear a security setting in the Adobe Acrobat Reader application before you start the BioPharma Finder software installation. Do one of the following:

- (For systems with Reader) Open the Reader application and choose **Edit ▶ Preferences ▶ Security (Enhanced)**. Clear the Enable Protected Mode at Startup check box, and then click OK.



- (For systems that do not have Reader) Follow “Installing the BioPharma Finder software” on page 1. Then, exit the BioPharma Finder installer wizard, follow the instructions for systems with Reader, and then return to the wizard to complete the BioPharma Finder installation.

Trademarks

BioPharma Finder is a trademark and Thermo Scientific is a registered trademark of Thermo Fisher Scientific Inc. in the United States.

The following are registered trademarks in the United States and other countries: Microsoft and Windows are registered trademarks of Microsoft Corporation. Adobe, Acrobat, and Reader are registered trademarks of Adobe Systems Incorporated.

Flexera Software is a registered trademark of Flexera Software, Inc. in the United States.

©2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.