

Mass Frontier 8.1 Installation Instructions

Follow these instructions to install and license the Mass Frontier™ software, version 8.1. You must be a system administrator or must have an account with administrator rights on the installation computer to install the application. Thermo Fisher Scientific does not limit the licensing permissions, so you may license the application as an administrator or as a standard user.

Contents

- [System Requirements](#)
- [Installation options](#)
- [Launch the installer](#)
- [Install as a local installation](#)
- [Install as a client/server installation](#)
- [Technical support](#)
- [Trademarks](#)

System Requirements

These are the minimum required and recommended hardware and software configurations for Mass Frontier 8.1 application.

Table 1. System requirements

System	Minimum configuration	Recommended configuration
Hardware	<ul style="list-style-type: none">• Intel compatible processor with a minimum speed of 1 GHz.• 8GB RAM• 30GB of disk space• Monitor resolution display of 1024 x 768pixels.• DVD-ROM drive (This is optional and not required if you can download from Flexera).	<ul style="list-style-type: none">• Intel compatible processor with a minimum speed of 2.5 GHz.• 32GB RAM• 100GB of solid-state drive disk space• Dual monitor resolution display of 1920 x 1080pixels. DVD-ROM drive (This is optional and not required if you can download from Flexera).
Software	<ul style="list-style-type: none">• Microsoft Windows™ 10 (English) 64bit.• Microsoft™ .NET Framework 4.7.2• Adobe™ Acrobat Reader DC• Google Chrome version 81.0 or higher.	<ul style="list-style-type: none">• Microsoft Windows™ 10 LTSC 64bit.• Microsoft™ .NET Framework 4.7.2• Adobe™ Acrobat Reader DC• Google Chrome version 81.0 or higher.
Computer	<ul style="list-style-type: none">• To run processing workflows with online mass spectral database searches, the computer must have an unblocked access to the mass spectral data via the internet .• The time and date settings in computer must be synchronized with the internet time. .• Set the Region and Language setting to English (United States).	

Installation options

The following options are available for installing Mass Frontier.

- Local Installation

In a local installation, the Mass Frontier application and the libraries reside on the same computer, and this computer has the following services and software:

- PostgreSQL 15
- Mass Frontier Server Manager 8.1
- Mass Frontier 8.1 application

- Client/Server Installation

In a client/server installation, the libraries reside on a dedicated server machine and are shared over the network.

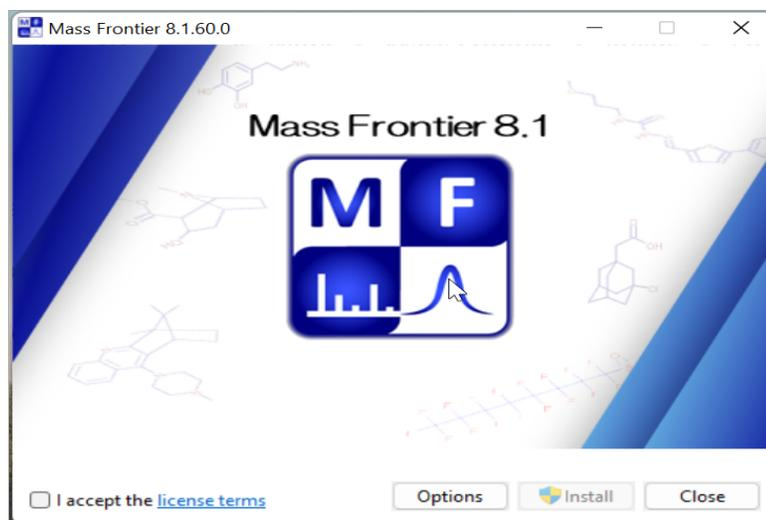
- The server machine has the following services and software:
 - PostgreSQL 15
 - Mass Frontier Server Manager 8.1
 - (Optional) Mass Frontier 8.1 application
- The client machine has the following services and software:
 - Mass Frontier 8.1 application
 - (Optional) PostgreSQL 15 and Mass Frontier Server Manager 8.1

Launch the installer

❖ **To launch the installer from the Mass Frontier 8.1 DVD**

1. Insert the Mass Frontier 8.1 DVD into the DVD/CD-ROM drive.
The AutoRun program automatically launches the installer wizard.
2. If the wizard does not start, navigate back to the DVD in Windows Explorer and double-click **MassFrontier81_Setup.exe**.

Figure 1. Mass Frontier 8.1 installer



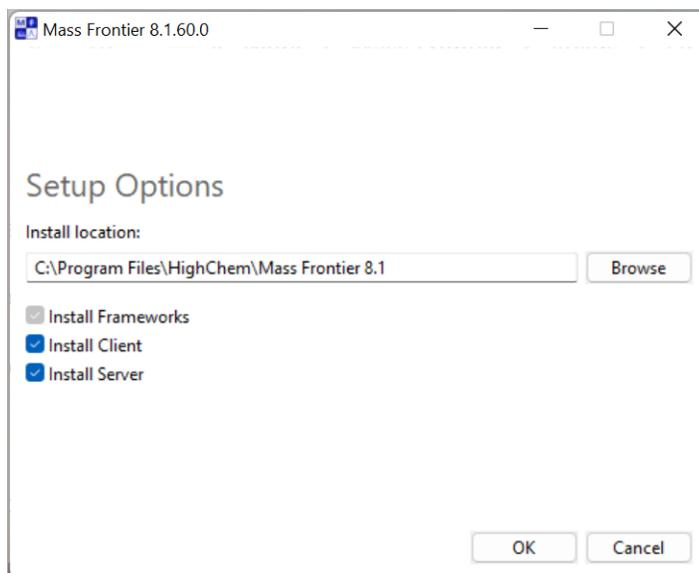
Install as a local installation

❖ **To install Mass Frontier 8.1 on your local computer**

1. Click **Options** after launching the installer.
2. Browse to the location where you want to install the software.
3. Select the **Install Server** and **Install Client** check boxes to install the Server Manager 8.1, Mass Frontier Library Service, and Mass Frontier 8.1 application.

Install as a client/server installation

Figure 2. Setup options for the installation

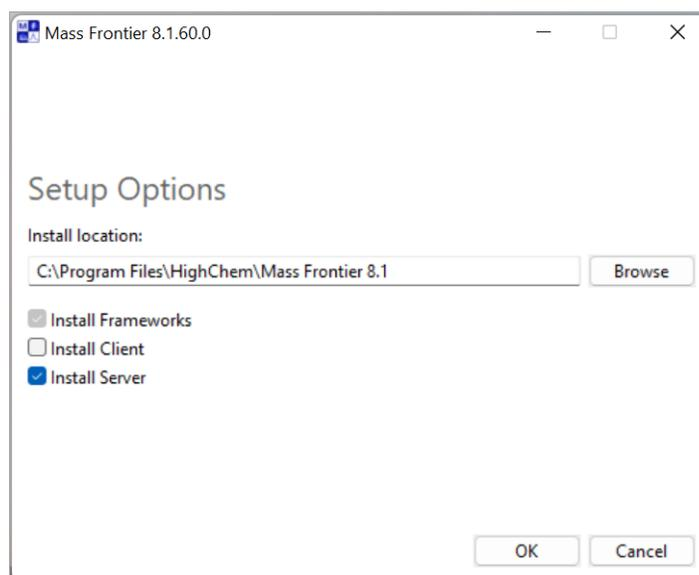


4. Click **OK**.
5. Review and accept the license terms.
6. Click **Install**.
7. At the prompt, click **Yes** to start the installation.

❖ To install components on the server computer

1. Click **Options** after launching the installer.
2. Browse to the location where you want to install the software.
3. Select the **Install Server** check box to install Server Manager 8.1 and Mass Frontier Library Service.

Figure 3. Install Server



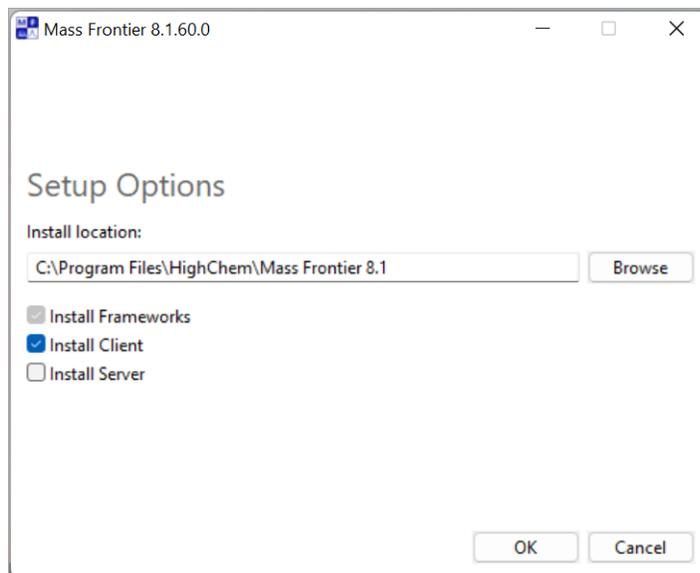
4. (Optional) To also install the Mass Frontier 8.1 application on the server computer, select the **Install Client** check box.
5. Click **OK**.
6. Review and accept the license terms.
7. Click **Install**.

8. At the prompt, click **Yes** to start the installation.

❖ **To install components on the client computer**

1. Click **Options** after launching the installer.
2. Browse to the location where you want to install the software.
3. Select only the **Install Client** check box.

Figure 4. Install Client



4. (Optional) To also install the Mass Frontier Server Manager 8.1 application and library service on the client computer, select the **Install Server** check box.
5. Click **OK**.
6. Review and accept the license terms.
7. Click **Install**.
8. At the prompt, click **Yes** to start the installation.

Mass Frontier 8.1 license activation

To activate the Mass Frontier 8.1 software, do the following:

1. After the installation is completed, start the application from the desktop icon or Start menu.
 2. Choose the appropriate licensing option:
 - For the demo version of the Mass Frontier 8.1 application, use the 60-day trial license that comes preinstalled.
- or–
- For purchasers of the Mass Frontier 8.1 application, set a permanent license using the activation code provided through your Flexera™ account. You cannot use a license key from an earlier version.
- You will receive an e-mail from Thermo MS licensing with subject line “Mass Frontier 8.1 software is now ready for download!” This e-mail contains the information for activation code.
3. Choose **Mass Frontier Start menu > About > License** and click **Activate License** to display the Product Licensing wizard and follow the on-screen instructions.

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

❖ **Locate or request the activation code**

If you have not received the activation code for Mass Frontier 8.1 software, follow this procedure:

1. Check your Junk e-mail folder.

Technical support

- If the e-mail is not in your junk email folder, log in to your account at the following URL: <https://thermo.flexnetoperations.com>
 - In the left navigation pane, under **Software & Services**, click **Order History**.
 - In the list of ordered products, click the order number to see the activation code.
- If you cannot find your account, send an e-mail message to Licensing at ThermoMSLicensing@thermofisher.com with the following information in the body of the email.
 - Software application: Mass Frontier 8.1 software.
 - End user name
 - End user e-mail
 - Sale order or purchase order number

For technical assistance, use the contact information below in your geographic location.

Table 2. Contact information

Country	Contact
North America	us.techsupport.analyze@thermofisher.com
Europe	eu.techsupport.cms@thermofisher.com
Japan	support.jp@thermofisher.com
China and other APAC regions	Local Support
Latin America, Eastern Europe, Middle East, & Africa	Local Dealer

Trademarks

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

Microsoft and Windows are registered trademarks of Microsoft Corporation in the United States and other countries.

All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.