

Mass Frontier 8.0 Installation Instructions

Follow these instructions to install and license the Mass Frontier™ software, version 8.0. You must be a system administrator or must have an account with administrator rights on the installation computer to install the application. You may license the application as an administrator or as a standard user.

Contents

- [Installation Options](#)
- [Launching Mass Frontier 8.0 installer](#)
- [Installing as a Local Installation](#)
- [Installing as a Client/Server Installation](#)
- [Technical Support](#)
- [Trademarks](#)

Installation Options

The following options are available for installing Mass Frontier. Choose the option you need and proceed to the installation instructions.

- **Local Installation:** In local installation, Mass Frontier application and the libraries are configured on the same computer. Your computer must have the following services and software:
 - Microsoft SQL Server 2014 Express
 - Mass Frontier Server Manager 3.0
 - Mass Frontier 8.0
- **Client/Server Installation:** In client/server installation, the libraries reside on a dedicated server machine and are shared over the network.
 - **The Server machine must have the following services and software:**
 - Microsoft SQL Server 2014 Express
 - Mass Frontier Server Manager 3.0
 - (Optional) Mass Frontier 8.0
 - **The Client machine must have the following services and software:**
 - Mass Frontier 8.0
 - (Optional) Microsoft SQL Server 2014 Express and Mass Frontier Server Manager 3.0

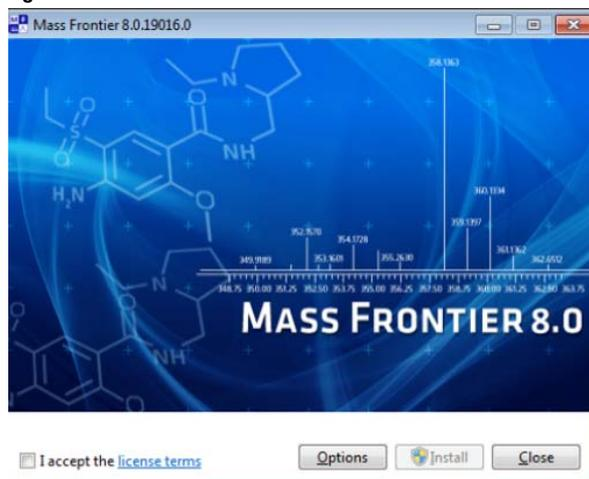
Launching Mass Frontier 8.0 installer

❖ To launch the installer from Mass Frontier 8.0 DVD

1. Insert the Mass Frontier 8.0 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 **MassFrontier80_Setup.exe**.

Installing as a Local Installation

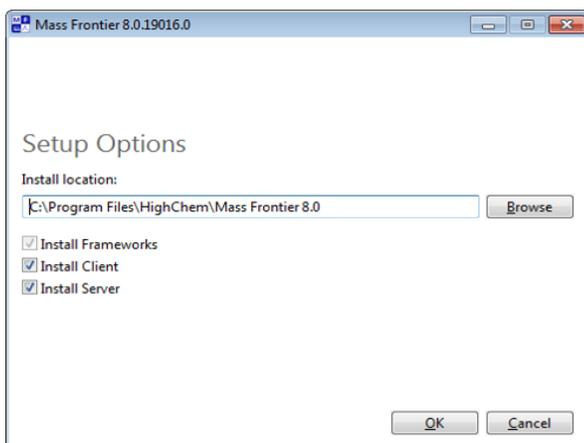
Figure 1. Mass Frontier 8.0 installation



❖ **To install Mass Frontier 8.0 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, Mass Frontier Library Service, and Mass Frontier 8.0 application.

Figure 2. Install Client and Server



4. Click **OK**.
5. Review and accept the license terms.
6. Click **Install**.

Note A User Account Control box displays with the following message:

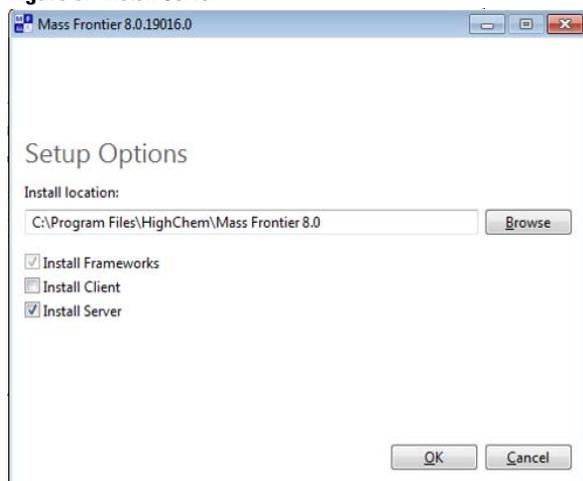
“Do you want to allow the following program to make changes to this computer?” Program Name: Mass Frontier 8.0, Verified Publisher: HighChem s.r.o.
Click **Yes** to start the installation.

Installing as a Client/Server 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 and Mass Frontier Library Service.

Figure 3. Install Server



4. (Optional) To also install the Mass Frontier 8.0 application on the Server computer, select the **Install Client** check box.
5. Click **OK**.
6. Review and accept the license terms.
7. Click **Install**.

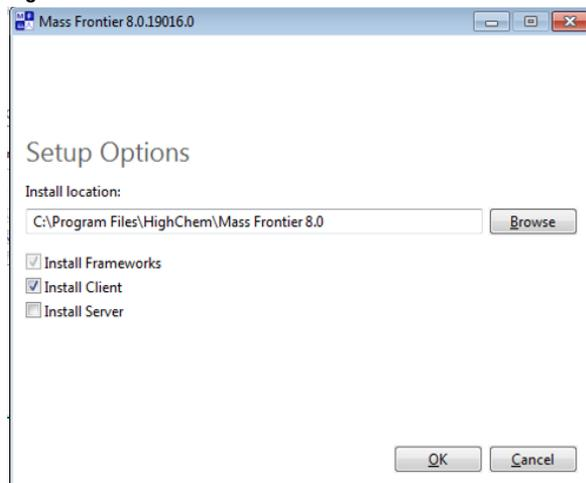
Note A User Account Control box displays with the following message:

“Do you want to allow the following program to make changes to this computer?” Program Name: Mass Frontier 8.0, Verified Publisher: HighChem s.r.o.
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 **Install Client** check box.

Figure 4. Install Client



4. (Optional) To also install the Mass Frontier Server Manager 3.0 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**.

Note A User Account Control box displays with the following message:
 “Do you want to allow the following program to make changes to this computer?” Program Name: Mass Frontier 8.0, Verified Publisher: HighChem s.r.o.
 Click **Yes** to start the installation.

❖ **To activate the Mass Frontier 8.0 license**

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 8.0 application, use the 60-day trial license.

–or–

 - For purchasers of the Mass Frontier 8.0 application, set a permanent license using the activation code provided through your Flexera™ account. You cannot use a license key from an earlier version.
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.

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

Table 1. 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 of HighChem, LLC., Slovakia.

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.