

FreeStyle 1.8 SP2 Installation Instructions

This guide provides instructions for installing the FreeStyle™ 1.8 SP2 application to view and analyze raw mass spectrometry data.

Contents

- [System requirements](#)
- [Installing the FreeStyle application](#)
- [Upgrading the FreeStyle application](#)
- [Trademarks](#)

IMPORTANT You must have administrator rights to install the FreeStyle 1.8 SP2 application on your computer.

IMPORTANT The FreeStyle 1.8 SP2 application was tested and supported for US-English Only locale settings. To change the format on your computer, go to Control Panel > Region and specify the language format.

System requirements

These are the minimum hardware and software configurations required for FreeStyle 1.8 SP2 operation.

System	Requirements
Computer	<ul style="list-style-type: none">• 3.6 GHz quad core processor with a minimum of 8 GB RAM (16 GB recommended)• CD/R-ROM or DVD drive• 1 TB hard drive• Video card and monitor capable of 1920 × 1080 resolution
Software	<ul style="list-style-type: none">• Adobe™ Reader™ 10.1 or later• Microsoft™ .NET Framework 4.7.2• Microsoft Office™ 2013 (for exported data)• Microsoft Windows™ 7 SP1 (64-bit), Windows 10 (64-bit)• Thermo Foundation™ 3.1 SP6 or later

Installing the FreeStyle application

Follow one of these procedures:

- [To install or upgrade the FreeStyle application on a computer with Xcalibur installed](#)
- [To install the FreeStyle application on a computer without Xcalibur installed](#)

❖ To install or upgrade the FreeStyle application on a computer with Xcalibur installed

1. Download the FreeStyle 1.8 SP2 software from Flexera (<https://thermo.flexnetoperations.com/control/thmo/login>) to your local processing computer.

If the computer has Xcalibur installed, the FreeStyle application should also be installed.

Before upgrading to FreeStyle 1.8 SP2, verify the Foundation version installed on your system by going to Version Info.

2. Follow the instructions in [Upgrading the FreeStyle application](#).

Upgrading the FreeStyle application

❖ To install the FreeStyle application on a computer without Xcalibur installed

Download the FreeStyle 1.8 SP2 software from Flexera (<https://thermo.flexnetoperations.com/control/thmo/login>) to your local processing computer.

- If the computer has never had the Foundation platform or the FreeStyle application installed before (clean installation), first install Foundation 3.1 SP6 or later and then install the FreeStyle 1.8 SP2 application.
- If the computer has Foundation 3.1 SP6 or later installed, uninstall the previous version of the FreeStyle application from the Control Panel and then install the FreeStyle 1.8 SP2 application.
- If the computer has Foundation 3.1 SP5 or earlier installed, do the following in this order:
 - a. Uninstall the FreeStyle application.
 - b. Uninstall the Foundation platform.
 - c. Install the Foundation 3.1 SP6 or later platform.
 - d. Install the FreeStyle 1.8 SP2 application.

Before upgrading to the FreeStyle 1.8 SP2 application, verify the Foundation platform version installed on your system by going to Version Info.

Follow one of these procedures:

- [To upgrade to the FreeStyle 1.8 SP2 application for a system with Foundation 3.1 SP6 or later](#)
- [To upgrade to the FreeStyle 1.8 SP2 application for a system with Foundation 3.1 SP5 or earlier](#)

❖ To upgrade to the FreeStyle 1.8 SP2 application for a system with Foundation 3.1 SP6 or later

1. Download the FreeStyle 1.8 SP2 software from Flexera (<https://thermo.flexnetoperations.com/control/thmo/login>) to your local processing computer.
2. Open the Windows Control Panel and click **Programs and Features**.
3. Uninstall the previous version of the FreeStyle application.
4. Install the FreeStyle 1.8 SP2 application.

❖ To upgrade to the FreeStyle 1.8 SP2 application for a system with Foundation 3.1 SP5 or earlier

1. Download the FreeStyle 1.8 SP2 software from Flexera (<https://thermo.flexnetoperations.com/control/thmo/login>) to your local processing computer.
2. Open the Windows Control Panel and click **Programs and Features**.
3. Uninstall the FreeStyle application.
4. Uninstall any previous version of the Xcalibur application.
5. Uninstall any previous version of the Foundation platform.
6. Install the Xcalibur 4.2 SP1 application or later, if needed.
7. Install the FreeStyle 1.8 SP2 application.

Trademarks

FreeStyle and Foundation are trademarks; and Xcalibur 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, Windows, and Office are registered trademarks of Microsoft Corporation. Adobe and Reader are registered trademarks of Adobe Systems Inc. in the United States and other countries.

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