

# Compound Discoverer 3.0 Installation Instructions

Thermo Compound Discoverer™ is a licensed application. Thermo Fisher Scientific provides a full version and a demo version of the application. For either version, you must use the license key provided to successfully install the software.

- Full version—Install Compound Discoverer 3.0 from the software media kit that includes a key-shaped USB flash drive with the installation executable and a small ID card with the license key.
- Demo version—Install Compound Discoverer 3.0 software from the DVD. The license key is on the back of the DVD case.

**IMPORTANT** The Compound Discoverer installation requires Microsoft™ Windows™ 7 SP1 64-bit or Windows 10 64-bit operating system. Refer to the release notes for the minimum and recommended data system computer configurations.

The Compound Discoverer application is ONLY supported with US-English locale settings.

The installation process for Compound Discoverer 3.0 requires an Internet connection to validate the software license.

**Note** To install Compound Discoverer 3.0 software, you do not need to uninstall Compound Discoverer version 1.0, 2.0, or 2.1. These software versions can coexist on the same computer.

## Contents

- [Installing Compound Discoverer](#)
- [Trademarks](#)

Follow one of these procedures:

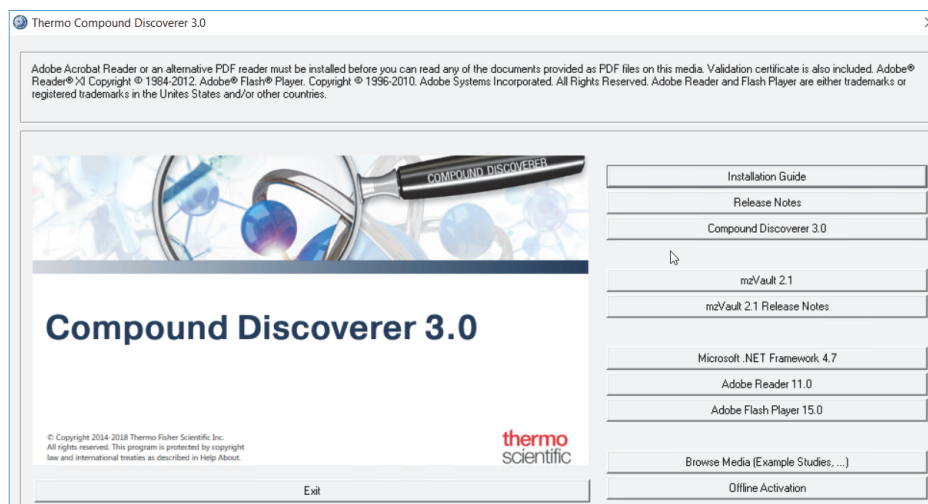
- [To install Compound Discoverer on a computer with Internet access](#)
- [To install Compound Discoverer on a computer without Internet access](#)

### ❖ To install Compound Discoverer on a computer with Internet access

1. Do one of the following:
  - For the full version, insert the Compound Discoverer USB flash drive into a USB port on your computer.
  - For the demo version, insert the Compound Discoverer DVD into the DVD drive on your computer.
2. Open Windows Explorer and select the USB or DVD drive to view its contents.
3. Double-click **XStart\_Compound Discoverer.exe**.  
The installation wizard starts ([Figure 1](#)).

## Installing Compound Discoverer

**Figure 1.** Compound Discoverer installation wizard



4. Click **Compound Discoverer 3.0** and continue the installation process.

The Minimum Product Requirements screen appears with a product requirements summary. If the computer meets the system requirements, the box displays the following:

All the system minimum requirements checks passed.

5. When the License Activation dialog box opens, type the activation code in the Activation Code boxes.
  - For the full version, enter the license key from the small ID card in the Compound Discoverer 3.0 software media kit.
  - For the demo version, enter the license key from the back of the Compound Discoverer 3.0 DVD case.
6. Complete the installation process.

❖ **To install Compound Discoverer on a computer without Internet access**

1. On the computer without Internet access and where you will run the Compound Discoverer application, do the following:
  - a. Follow [step 1](#) through [step 4](#) of “[To install Compound Discoverer on a computer with Internet access](#)” on [page 1](#).

**IMPORTANT** If an Activation Error dialog box opens, temporarily connect the computer to a network switch or hub that is turned on but is not connected to a network. Wait a few minutes before you continue the installation process.

- b. When the License Activation dialog box opens, do the following:
  - i. Type the activation code in the Activation Code boxes.
  - ii. In the area labeled If You Do Not Have an Internet Connection on this Computer, click **Export the License Request**.  
The Export License Request dialog box opens.
  - iii. Select the directory where you want to store the license request file, type a name for the file in the File Name box, and click **Save**.
- c. Eject the flash drive or DVD from the computer and leave the License Activation dialog box open.
- d. Transfer the license request file to a computer with Internet access.

For example, copy the license request file to a read-writable USB flash drive. Insert the USB flash drive into a USB port on the computer that has Internet access and copy the **FileName.licrequest** file to an appropriate directory.

2. On a computer with Internet access, activate the license as follows:
  - a. Verify that you transferred the *FileName.licrequest* file to this computer.

**Note** Ensure that the license request file is not write-protected.

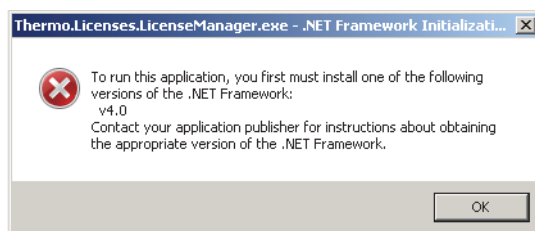
- b. Follow [step 1](#) through [step 3](#) of “[To install Compound Discoverer on a computer with Internet access](#)” on [page 1](#).
- c. Click **Offline Activation**.

The Import License Activation Request dialog box opens with the Thermo License Activation dialog box in the background.

–or–

If the computer does not have the appropriate version of the .NET Framework, a message box appears in front of the installation screen ([Figure 2](#)).

**Figure 2.** .NET Framework initialization message box



**IMPORTANT** The License Activation dialog box requires Microsoft Visual C++™ redistributables and .NET Framework 4.0 or later. If the redistributables are not present on the computer, the program installs them automatically. If the appropriate version of the .NET Framework is not present, click OK to close the message box. Then, click Microsoft .NET Framework 4.7.

- d. Browse to the license request file that you installed on this computer and click **Open**.

The Import License Activation Request dialog box closes, the Thermo License Activation dialog box becomes available, and the Activation Code boxes are populated with the license activation code from the license request file.
  - e. Enter the appropriate information in the Contact Information area.
  - f. Click **Activate**.

The program activates the license and stores it in the license request file.
  - g. With the license activation confirmed, click **OK**.
  - h. Close the installation wizard.
  - i. Transfer the activated license request file back to the computer without Internet access.
3. On the computer without Internet access and where you will run the Compound Discoverer application, do the following:
    - a. Verify that you transferred the activated *FileName.licrequest* file (see [step 2](#)) to this computer.
    - b. In the License Activation dialog box that you left open in [step 1c](#), click **Reimport the Activated License Request to Continue Activation**.

The Import License Activation Request dialog box opens.
    - c. Browse to the activated license request file and click **Open**.

The Activation Success box opens and displays the following message:

Licenses successfully imported.
  4. Complete the installation process.

---

## Trademarks

Compound Discoverer is a trademark of Thermo Fisher Scientific Inc. in the United States.

Microsoft, Windows, and Visual C++ are registered trademarks of Microsoft Corporation in the United States and other countries.