

Sartorius Data Analytics

Technical Note

Manual (Offline) Activation of Umetrics® Suite Products

March 17, 2023

This technical note applies to:

- MODDE® and MODDE®-Q 13.0.2 and later
- SIMCA® 17.0.2 and later
- SIMCA®-Q 17.0.1 and later
- SIMCA®-online 17 and later

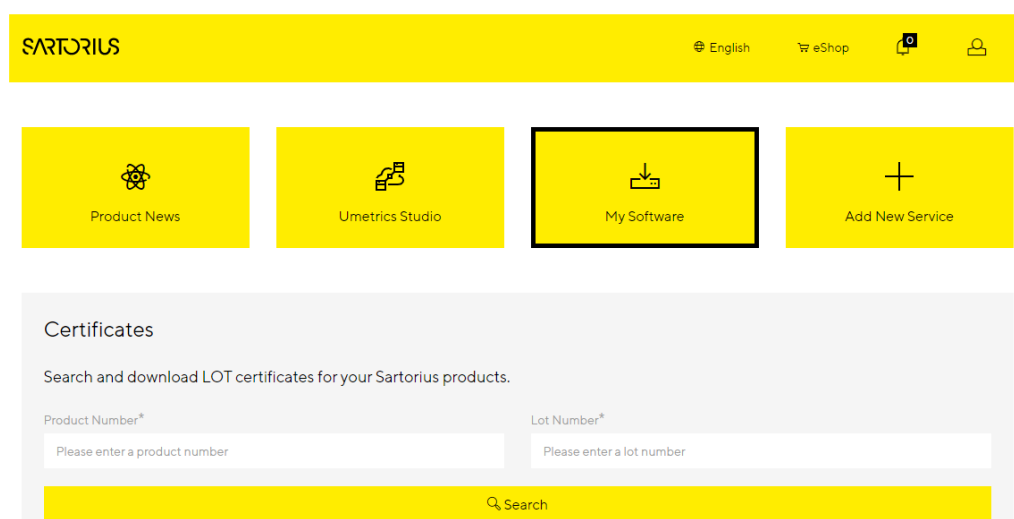
If your computer cannot be activated over the Internet, or if the online activation process isn't working, you can manually download a license file for your product on a different PC where you have Internet access. Here is how:

Download a License from My Sartorius Portal

Your activation ID can either be connected to an individual license or to a bulk license. The difference is how many times the license can be activated.

To activate your license:

1. These instructions assume you have created an **offline activation file** in the software you want to activate. You'll need this file in the steps below.
2. Visit my.sartorius.com and login with the credentials used when downloading the application.
3. Click on **My Software**:

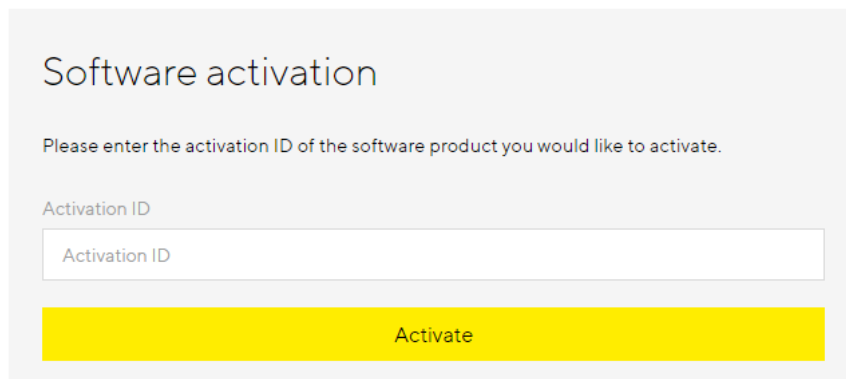


4. Enter Activation ID provided:

Activate Your Software Purchase

You have recently purchased a Sartorius software product and received two separate emails, one containing a link to this site and another one containing the license of your software purchase.

To access the software download, please activate your software.



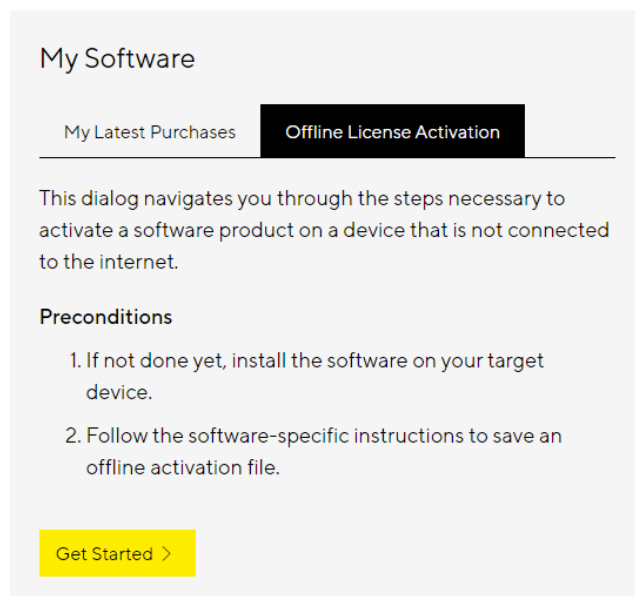
Software activation

Please enter the activation ID of the software product you would like to activate.

Activation ID

Activate

5. In My Sartorius portal, go to **Purchased Software -> Offline License Activation**:



My Software

My Latest Purchases Offline License Activation

This dialog navigates you through the steps necessary to activate a software product on a device that is not connected to the internet.

Preconditions

1. If not done yet, install the software on your target device.
2. Follow the software-specific instructions to save an offline activation file.

Get Started >

6. Choose the offline activation file to upload and click **Request Activated License**:

My Software

My Latest PurchasesOffline License Activation

This dialog navigates you through the steps necessary to activate a software product on a device that is not connected to the internet.

Step 1 of 3:

1. If not done yet, copy the offline activation file to the device you are currently using to display this website.
2. Upload the offline activation file via the below form.

Drop an offline activation file here

or

Select an offline activation file (*.bin) from computer

< Back

Request Activated License >

7. After successful activation, you will have an option to download the activated license:



My Software

My Latest PurchasesOffline License Activation

This dialog navigates you through the steps necessary to activate a software product on a device that is not connected to the internet.

Step 2 of 3:

1. Click "Request Activated License".
2. You will be navigated to the next step automatically once an activated license has been created.

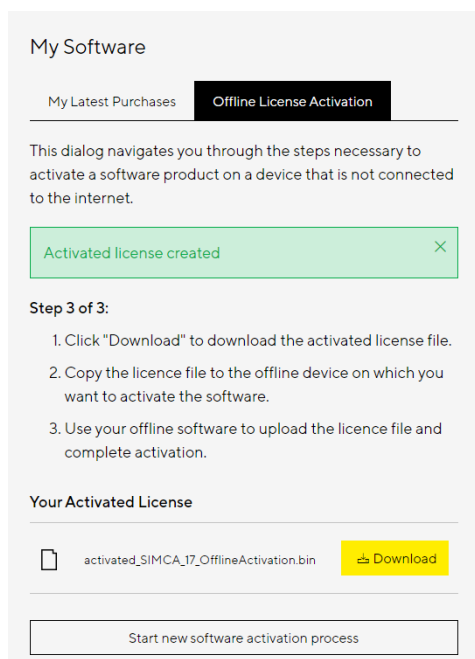
SIMCA_17_OfflineActivation.bin0.5KB

Discard File

< Back

Request Activated License >

8. Click **Download** to download your activated license. Take a note where it was saved. Copy the file to the computer to activate:



9. In the Sartorius Data Analytics product, if the Activation dialog is still open, click **Next** to browse to the downloaded license file and click **OK**. Alternatively, if the dialog isn't open, click **File | Help | Manage license**. Then, in the dialog that opens:
 - a. Click **Activate using activation ID**.
 - b. Select the checkbox **Perform offline activation** and click **Next** to continue the offline activation
 - c. Click **Next** again, browse to the downloaded license file and click **OPEN**.

Contact Sartorius Data Analytics to Receive a License File

If you couldn't download a license as described above, you can contact Sartorius for manual processing. **Please note that this can take up to 10 workdays to process.**

To process your request, we need the following information:

1. The **product name** (MODDE, SIMCA, ...) and **version number**. You can find the version number by clicking **File | Help**.
2. Your **activation ID** and the **host ID**. You find this information in the Activate dialog.
 - a. Your **name** and the **company name**.
 - b. If applicable, the company **department** and **site**.

Send the email to umetrics_delivery@sartorius.com.

After your information has been processed, you will receive an email containing a manually created license file. Go to **File | Help | Manage license** and click **Import license file**.