

Sartorius Data Analytics

Technical Note

Manual (offline) Activation of Umetrics Suite's Products

8 June 2020

This technical note applies to MODDE Pro 12.1, MODDE Go 12.1, SIMCA 16, SIMCA-Q 16, and Easy Analytics 1.2.

If your computer cannot be connected to the Internet, or if the online activation process for some reason keeps failing, you can download a license file to your computer. Although manual activation requires the ability to download a license file or to send or receive emails, you can complete this part of the activation process using any computer that can be connected to the Internet, not specifically the one where you installed the product.

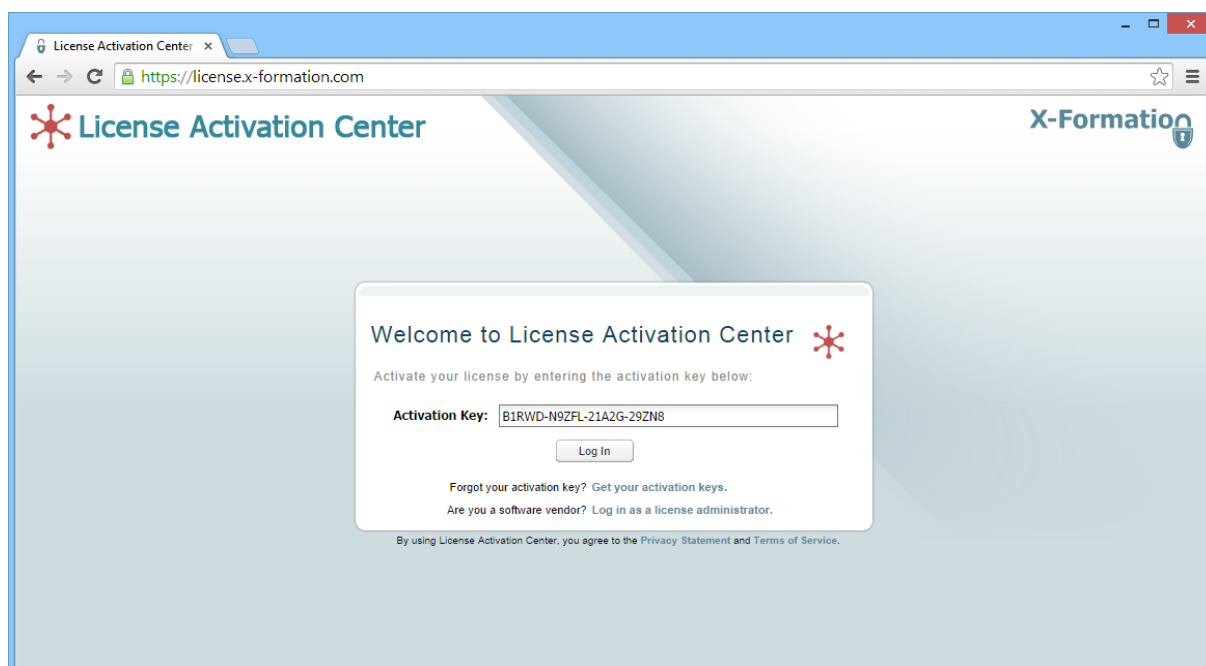
Please follow the instructions below to activate your copy of the program when automatic activation fails:

1. See the "Download a license from License Activation Center" section to manually download a license file.
2. If downloading did not work, see the "Contact Sartorius Data Analytics to receive a license file" section.

If you have installed the trial version and don't have an activation key, please contact Sartorius Data Analytics sales to obtain an activation key. See "Contact Sartorius Data Analytics to receive an activation key".

Download a license from License Activation Center

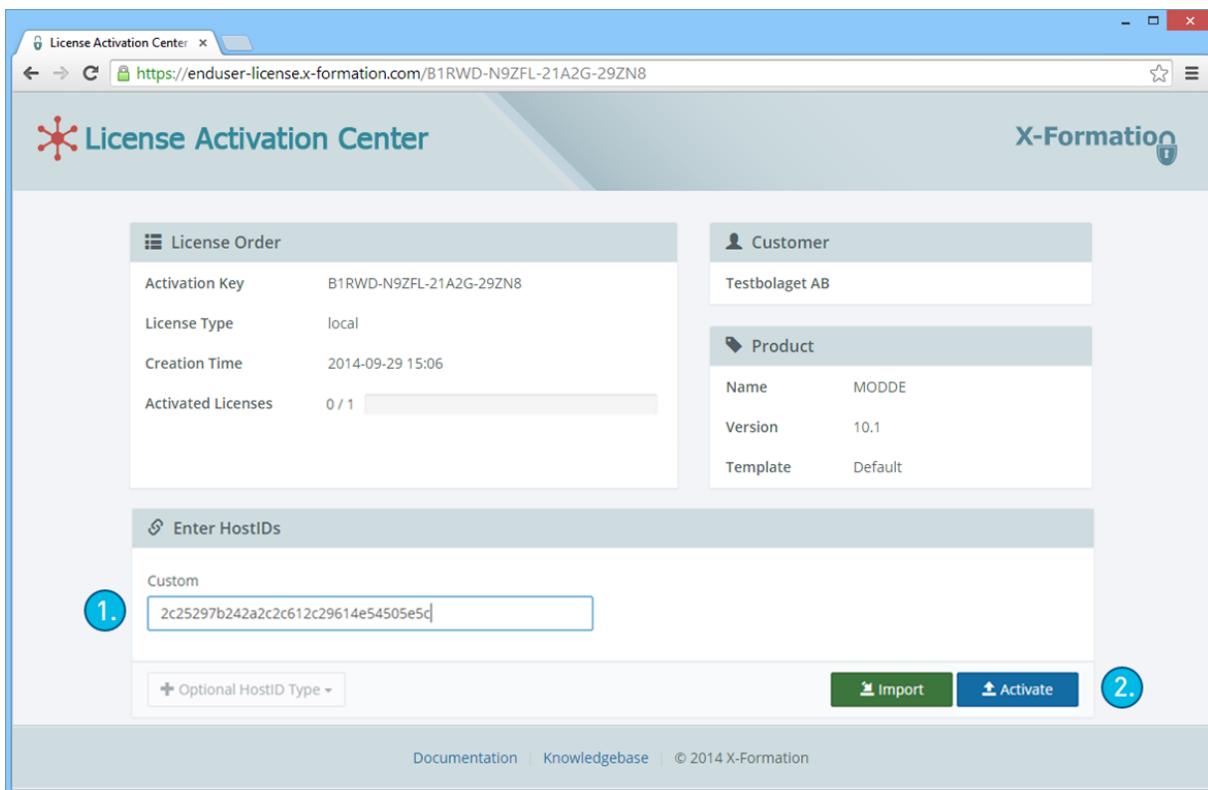
To download a license from License Activation Center (LAC), visit <https://license.x-formation.com> and enter your activation key in the log in screen and then click Log In.



Hint! In the Activation dialog there are links to copy relevant information to clipboard and to open the License Activation Center website with the activation key already filled in.

Activating an individual license

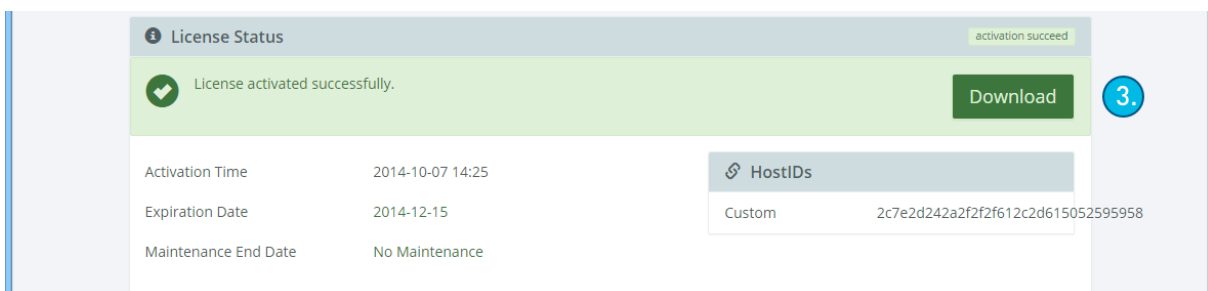
Your activation key can either be connected to an individual license or to a bulk license. The difference is how many times the license can be activated. If you can activate more than one license, see the "Activating a license from a bulk license" section for an initial step you need to complete before proceeding.



To activate your license:

1. Enter your **HostID** in the **Custom** field. Note: Find the HostID in **File | Help** under License information. Click it to copy to Clipboard and paste into the Custom field.
2. Click **Activate** to activate your license.

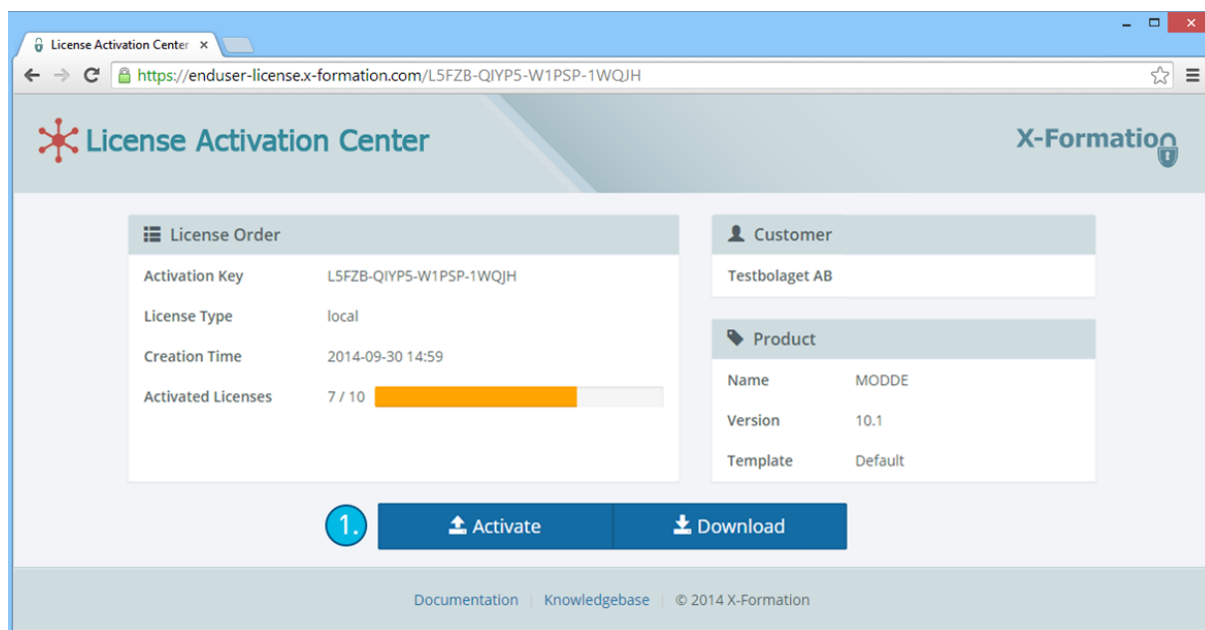
If the activation was successful, the message "License activated successfully" will be displayed.



1. Click **Download** to download your activated license. Take a note where it was saved. If you downloaded the license file on a different computer than the one where the software is installed, copy the file to that computer.
2. 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**. If you have closed the Activation dialog, click **File | Help | Activate**. Then, in the dialog that opens:
 - a. Click **Activate using activation key or license file**.
 - b. Select the **I already have a license file for this computer** check box and click **Next**.
 - c. Browse to the downloaded license file and click **OK**.

Activating a license from a bulk license

The difference between activating an individual license and a bulk license is that you will have to click **Activate** to check out a license before you can enter your **HostID**.



Note that if no licenses have been activated yet, the **Download** button does not appear. The above example shows that 10 activations are allowed and that 7 licenses have been activated, indicating that 3 remain.

1. Click Activate.
2. Proceed to step 1 in the section "Activating an individual license" above. The remaining steps are the same as for activating individual licenses.

Contact Sartorius Data Analytics to receive a license file

Please note that this routine can take up to 10 work days to process.

If, for some reason, downloading a license file through the License Activation Center did not work, please contact Sartorius Data Analytics to receive a license file through email.

To process your request, we'll need the following information:

1. The **product name** (MODDE, SIMCA, SIMCA-online,...) and **version number**. You can find the version number by clicking **File | Help**.
2. Your **activation key** and the **HostID**. 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 the license file.

Contact Sartorius Data Analytics to receive an activation key

If you have installed the program without supplying an activation key, you have 30 days to try it for free. Once you've tried it out and want to buy it, please contact Sartorius Data Analytics sales representatives to purchase the product and receive an activation key. Find Sartorius Data Analytics contact information at: <http://umetrics.com/about-us/contact-us>.

Whether you activate your copy of the product during or after the trial period, you will not need to reinstall. Activate by clicking **File | Help | Activate**, then click **Activate using an activation key...** and enter your activation key.