

SIMCA® Multivariate Data Analysis Solution

Readme SIMCA® 16

May 7, 2019

Thank you for your interest in SIMCA 16. SIMCA is a versatile multivariate data analysis software, praised for its accessibility and ease of use and this release is aiming at improving this further. In addition to the context based ribbons and side panes we are also adding some useful tools for model interrogation and finally, a brand new analytical technique for multi-block analysis, MOCA.

New features and changes in SIMCA 16:

- Advanced merge of datasets
 - Data Explorer pane which includes
 - Appearance for direct customization of plot visualisation
 - Properties for direct interaction with plots
 - o Quick info
 - o Trimming/Winzorizing
 - o What-If
 - o Score space exploration
 - o Solver
 - o Advisor
- Getting started guided tour for first time users
- New Python possibilities
 - o Updated to CPython 3.7.1
 - Create and use plugins for custom file formats and spectral filters
 - Easy installation of additional Python packages using Pip install
 - MOCA Multiblock Orthogonal Component Analysis
- OPLS has been extended to include all O2PLS
- calculations. O2PLS has therefore been removed
- Workset dialog improvements
 - o Integrated Quick info
 - Wizard mode for improved workflow
- New report generator with report saved inside the project file

Information about this and other functionality in SIMCA is available at our website: http://umetrics.com/product/simca

Installation

Administrator permissions are necessary to install SIMCA 16. Omics and Spectroscopy skins are included in the installation file. For information on how to enable skins and more details on general installation and activation see "SIMCA 16 Installation instructions.pdf"

Licensing and activation

To run the full version of SIMCA 16 you need to have a valid license that has been activated.

For single installations an activation key was delivered in your delivery e-mail.

If you have been instructed to connect to a license server, you may need information from your local IT department about which Host/IP to use.

Trial limitations

SIMCA can be run as a trial for 30 days.

In the trial version copy, paste, print and the report generator are disabled. Scripting is not available in the trial version. All other functionality is available.

Terminal Services and virtual machine

To run SIMCA under Windows Terminal Services or on a virtual machine, your license must specifically allow this. Contact your sales representative for more information.

Minimum system requirements

- Microsoft Windows 7, 8 or 10
- 1024x768 screen resolution color display

Support

The latest information and support questions can be found in <u>Sartorius Stedim Data Analytics</u> Knowledge base.

For Technical support, contact <u>umetrics_support@sartorius-</u> stedim.com.

