

## SIMCA®

Multivariate Data Analysis Solution

### Readme SIMCA® 17.0.2

November 16, 2021

Thank you for your interest in SIMCA® 17.

SIMCA® 17.0.2 is a minor release that addresses some bugs that were found in the previous release of SIMCA® 17.

SIMCA® is a versatile multivariate data analysis software, praised for its accessibility and ease of use and this release has a focus on making analysis of spectroscopy data easier, faster and more intuitive. This involves context-based ribbons, extended library of spectral preprocessing and a calibration wizard that makes creating multivariate calibration models easier. All users of SIMCA® will also benefit from the improvements in performance and plot responsiveness made in this version.

#### Changes in SIMCA® 17.0.2

- Addressed stability issues found in SIMCA® 17

#### New features and changes in SIMCA® 17:

- Integrated spectroscopy project
  - Context based ribbons with spectra focus
  - User defined Spectral ID used in plots
  - New spectra visualization
  - Full range of spectral preprocessing options
  - Preprocessing wizard for single or chained filters
  - Calibration wizard for single Y prediction models
- Batch Data and Database Import
  - Possibility to control Batch alignment
  - BCC from dataset
  - SimAPI:s available at installation
  - Database interface improvements including node hierarchies
- Plot Interactivity
  - Color by Rank – new option for better vector coloring
  - Reverse and log axis control in properties pane
  - Color BCC by Batch condition vectors
- Other Improvements
  - Multiple file import for CSV, TXT and DIF files
  - Import configuration template
  - CPython updated to version 3.7.9
  - NumPy, SciPy, and Pandas installed by default
  - Possible to change name of dataset
  - Dataset merge keeps un-matched content in new dataset
  - New bias vectors (MBE)
  - Possibility to Save as version 16
  - Multiple performance related improvements

Information about this and other functionality in SIMCA® is available at our [website](#).

### Installation

Administrator permissions are necessary to install SIMCA® 17.0.2.



The Omics skin is included in the installation file. For information on how to enable the skin and more details on general installation and activation see “SIMCA 17 Installation instructions.pdf”.

## Licensing and Activation

To run the full version of SIMCA® you need to have a valid license that has been activated. The licensing system was updated in SIMCA® 17.0.1. This means that if you have a license for 17.0.1 it works for 17.0.2 as well. A license from an earlier version does not work for 17.0.2.

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.

## Virtual Machine

To run SIMCA® on virtual machine, your license must specifically allow this. Contact your sales representative for more information.

## Minimum System Requirements

- CPU: Intel Core i5-4460, 4:th generation, (3.20GHz) or AMD FX-6300
- RAM: 4 GB
- Disk space: 700 MB available disk space
- OS: Microsoft Windows 8.1 or 10
- Video card: Nvidia GeForce GTX 460, ATI Radeon HD 4850, or Intel HD Graphics 4400

## Support

The latest information and support questions can be found in [Sartorius Data Analytics Knowledge base](#).

For Technical support, contact [umetrics\\_support@sartorius.com](mailto:umetrics_support@sartorius.com).