



SIMCA[®]-Q

Embedded solution

Readme SIMCA[®]-Q 16

15 May 2019

Thank you for your interest in SIMCA-Q 16.

New features in SIMCA-Q 16:

- Support for SIMCA 16 models
- Support for Python preprocessing when used in SIMCA
- Get the confidence interval for loadings

SIMCA-Q 16 is compatible with SIMCA 13, SIMCA 14, SIMCA 14.1, SIMCA 15 and SIMCA 16 projects. New vectors from SIMCA 16 have been added, and Unicode (utf-8) is supported.

For a detailed list of the changes, see SIMCA-Q Changes.txt in the installation directory.

Supported versions

SIMCA[®]-Q embedded solution 16 is supported on:

- Windows 7, 8 and 10
- Red Hat Enterprise Linux 7.

Please contact us for other Linux distributions.

Installation

Administrator permissions are necessary to install SIMCA-Q 16. For more details on general installation and activation see "SIMCA-Q 16 Installation Instructions.pdf"

Licensing

To run SIMCA-Q 16, a valid license is required.

License files for SIMCA-Q 15 and previous versions will not work with this version.

Run the 'Register SIMCA-Q' application found in the installation directory for information of how to request a license file and to import an existing license file.

Support

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

For support send an email to umetrics_support-oem@sartorius-stedim.com

