

SIMCA-Q

Embedded solution

Readme SIMCA-Q 17

15 June 2021

Thank you for your interest in SIMCA®-Q 17.

SIMCA-Q 17.0.1 is a minor release that introduces a new licensing system.

New features in SIMCA-Q 17:

- Support for SIMCA 17 models
- CPython updated to version 3.7.9
- Python sample

SIMCA-Q 17 is also compatible with SIMCA 13, SIMCA 14, SIMCA 14.1, SIMCA 15 and SIMCA 16 projects. New vectors from SIMCA 17 have been added, and Unicode (utf-8) is supported.

Supported Versions

SIMCA-Q embedded solution 17 is supported on:

- Windows 8.1 and 10
- Linux CentOS 8

Please contact us for other Linux distributions.

Installation

Administrator permissions are necessary to install SIMCA-Q 17.0.1. For more details on general installation and activation see "SIMCA-Q Installation Instructions.pdf"

Licensing

To run SIMCA-Q 17.0.1, a valid license is required. Due to the change of the licensing system, the license file for SIMCA-Q 17 will not work.

Run the 'RegisterSIMCA-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 Data Analytics Knowledge base](#).

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