

# SIMCA-Q

**Embedded solution** 

## Readme SIMCA-Q 17

9 February 2021

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

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 and 10
- Linux CentOS 8

Please contact us for other Linux distributions.

#### Installation

Administrator permissions are necessary to install SIMCA-Q 17. For more details on general installation and activation see "SIMCA-Q 17 Installation Instructions.pdf"

#### Licensing

To run SIMCA-Q 17, a valid license is required.

License files for SIMCA-Q 16 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 <u>Sartorius Data Analytics</u> Knowledge base.

For support send an email to  $\underline{umetrics\ support-oem@sartorius.com}$