

SIMCA®

Turn Data Into Growth

Simplifying Progress

SVISCISVS

SIMCA[®] Helps You See What Others Don't

Every day your business generates a variety of data from many sources. Leveraging your data can help you enhance performance.

The challenge is finding a way to collect and interpret these high volumes of data. With so many parameters at play, it's hard to find a solution that's powerful and smart enough to turn data into actionable insight.

SIMCA[®] software enables you to combine and analyze data across all sources. You can isolate, understand, and apply the insights needed to optimize your process.

Who is using SIMCA[®]?

Today's typical users include process experts, from the laboratory to manufacturing, as well as QA/QC professionals and data scientists.

Use SIMCA[®] software to perform tasks like:

Exploratory-focused activities

- Confirm or challenge expected relationships
- Interpret groups or clusters
- Identify process improvement opportunities
- Apply root cause analysis to troubleshoot concerns

Deployment-focused activities

- Process analytical technology with model building for soft sensors
- Golden batch modeling
- Multivariate calibration
- Multivariate characterization

Companies across industries have used SIMCA[®] software to grow their business:

- A major bioprocess company improved process yield by 75%, reduced cycle time by 40%, and trebled plant output.
- A global specialty chemical company enabled savings up to \$10M by detecting, analyzing, and correcting process issues.
- In a pilot study, a university decreased expenses and carbon emissions in their heating processes. They're using SIMCA[®] software to conduct a sustainability project with the goal of avoiding around 10,000 metric tons of carbon emissions.



Multivariate Data Analysis

MVDA

ant output.

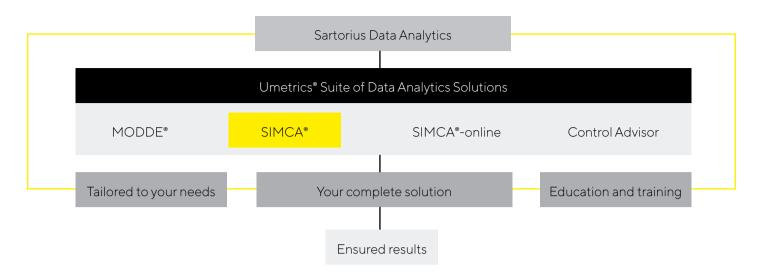
SIMCA® at a Glance

- Spectroscopy features integrated into context-based ribbons
- Guided workflows and interactive graphical interface
- Flexibility to handle complex data in many forms
- Intuitive method to script your workflow
- Manage models and deploy to systems with SIMCA®-Q embedded
- Seamless SIMCA[®]-online integration for model updates





More Than Software



Our complete solution includes everything you need through the whole process and provides results quickly. As our customer, you'll have access to supporting documents, templates, training and consultation to address your specific business challenges. Our courses and webinars help over a thousand people every year develop expertise and confidence in data analytics.

A Complete Suite for Business Growth

The Umetrics[®] Suite is a family of proven data analytics solutions that work seamlessly together. Other software solutions in the Umetrics[®] Suite are:

MODDE[®]

Design of Experiments to get it right from the start

SIMCA[®]-online

• Real-time process monitoring to maintain product quality and reduce risk

Control Advisor

Predictive capabilities to enable forecast of outputs

Web Client

• Process performance insight anywhere at any time

These solutions give you control and confidence in your processes at every stage - from development to manufacturing.

"Using SIMCA" to solve problems has saved us millions of dollars over the years."

Timothy Michaelson, Senior Manufacturing Excellence Consultant at International Paper

Find out more For more information, please visit www.sartorius.com/umetrics



Sartorius Data Analytics

Change a Little - Grow a Lot

We help organizations grow. The Umetrics® Suite of Data Analytics Solutions helps you harness the wealth of data within your organization. Our expertise in data analytics can help you identify vital elements to improve the results of your research, development and manufacturing processes.

With improved process understanding and more consistent product quality, you'll be able to reduce risk, get to market faster, and grow your business.

Our complete solution encompasses software, training, support and project management. And as part of Sartorius, a global company with more than 10,000 employees, we give you the backing of an international presence.

Experience the benefits for your business today

Find out how our solutions can help your business to grow, whatever industry you are in.

Visit **www.sartorius.com/simca** for details or to download a free 30-day trial.



Sartorius Data Analytics

Phone: +46 40 664 2580 E-mail: umetrics@sartorius.com

Germany

Sartorius Stedim Biotech GmbH August-Spindler-Straße 11 37079 Göttingen Phone +49 551 308 0

For further contacts, visit www.sartorius.com/umetrics

Sweden

Sartorius Data Analytics Östra Strandgatan 24 SE-903 33, Umeå Phone: +46 40 664 2580 E-mail: umetrics@sartorius.com