



SIMCA®

Turn Data Into Growth

Simplifying Progress

SARTORIUS

SIMCA® Helps You See What Others Don't

Every day your business creates a wide variety of data from many different sources. This data holds the key to better performance.

The challenge is to interpret this information in a meaningful way. But with so many parameters in play, it's hard to find a solution that's both powerful and smart enough. SIMCA® gives you the ability to combine and analyze all these different data sources. It helps you to isolate, understand and act upon the hidden gems that hold the secret to greater business success.

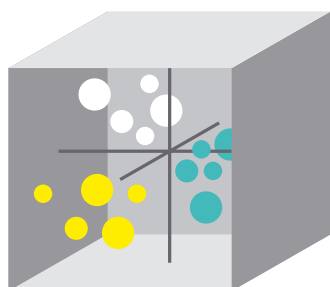
Who is using SIMCA®?

Wherever you create data you can use SIMCA®. That's why companies in many different industries have worked with us to help their business grow.

- A major bio-process company improved process yield by 75%, reduced cycle time by 40% and trebled plant output.
- An international food processing company resolved a logistics issue and saved USD 1 million per year in shipping costs.
- A wastewater treatment company used SIMCA to improve their processes for a cleaner, safer environment.

SIMCA® at a Glance

- Integrated spectroscopy features through context based ribbons
- Interactive graphical interface
- Flexibility to handle complex data in many forms
- An easy way to script your workflow
- Seamless model update integration with SIMCA®-online

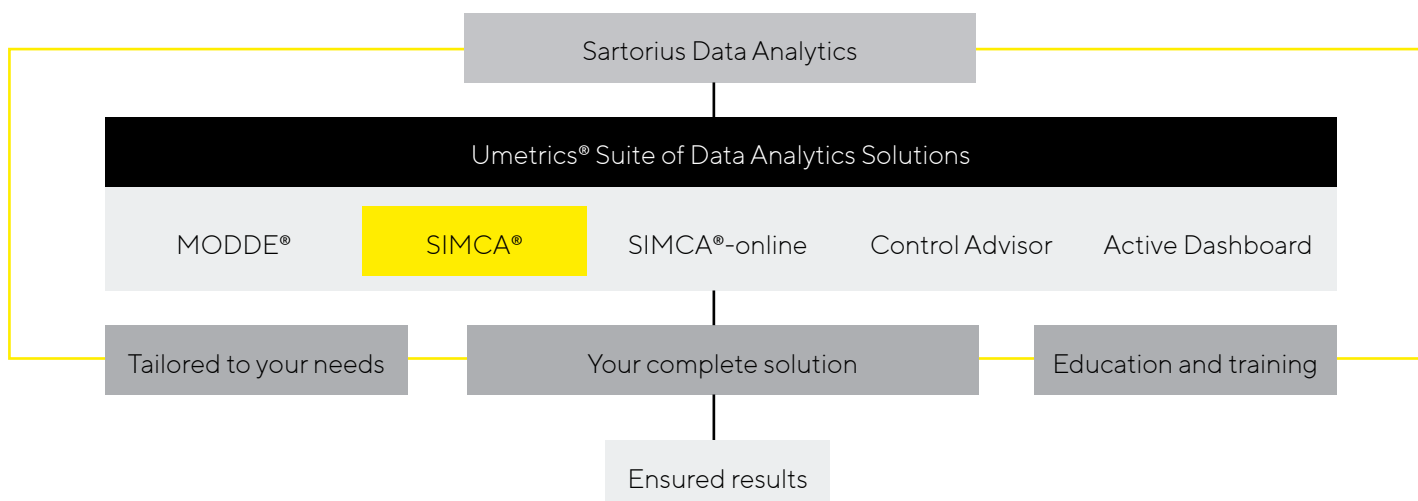


MVDA

Multivariate Data Analysis



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

- Online – Real-time process monitoring to maintain product quality

Control Advisor

- Predictive capabilities to be able to forecast the output

Active Dashboard

- Interactive performance insight

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/umetrics 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

Sweden

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

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

Specifications subject to change without notice.

©2025 Sartorius Stedim Biotech GmbH, August-Spindler-Strasse 11, 37079 Goettingen, Germany

Last updated: 06 | 2025