



SIMCA[®]-online

Ensuring Manufacturing
Success

Simplifying Progress

SARTORIUS

Predicting Process Problems Boosting Business Growth

Every production process follows a strict set of rules. If those rules are broken, you need to know – and the earlier the better. SIMCA®-online from Sartorius Data Analytics monitors your processes in real-time for a continuous snapshot of your operations. You can identify when set parameters change, fix them before they affect production and keep quality consistent. With this level of control, you can maximize efficiency and minimize costs. You will enjoy the confidence of high quality in your product and a real boost to your business growth.

How Does SIMCA®-online Work?

Instead of monitoring each variable, you can concentrate them into one view that is key to your whole process.

Easy-to-understand graphics make interpretation simple.

- Monitor in real time and swiftly detect deviations. With SIMCA® you can model your ideal process from your collected data. Transferred into SIMCA®-online, the model acts as a valuable reference for your current production.
- Predict with confidence. You can predict final quality from the properties of the raw material and the process parameters as well as forecast the final quality.
- Control at a glance. SIMCA®-online uses an 'ideal process' model to anticipate the effect of changes and recommend immediate adjustments. This will ensure product performance according to specifications and optimize throughput.

SIMCA®-Online at a Glance

- Remote predictive monitoring
- Detect and correct deviations in multi-unit processes with improved access to process data
- Root-cause-analysis
- Predicting final quality attributes
- Soft sensing
- Real-time supervisory control



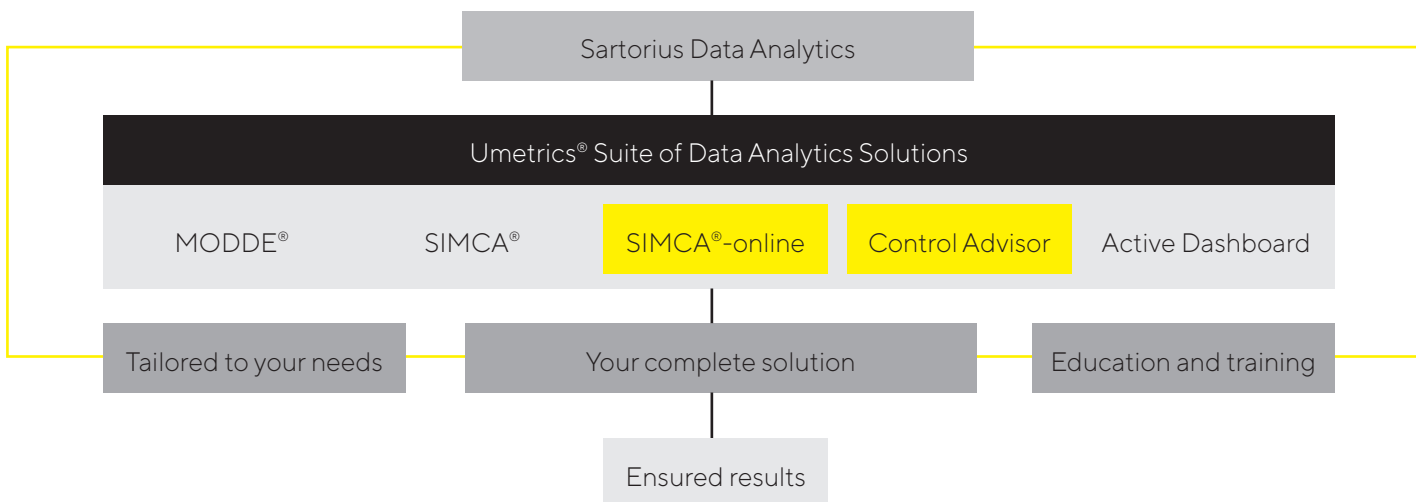
Proven in a Wide Range of Industries

SIMCA®-online has been adopted widely in many different industries, including pharma & chemical, pulp & paper and food & beverage.

Our customers have seen excellent results in cost savings, efficiency and product quality. For instance:

- A multi-national paper company reduced costs, achieved a more consistent product quality and gained a deeper understanding of their data.
- A major food processing company saved over USD 3 million each year in production costs.
- A pharmaceutical company paid for their investment several-fold in recovered batches alone.

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®

- Multivariate Data Analysis Solution to help you see what others don't

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.



“SIMCA®-online improves overall understanding of process and equipment.”



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 demo version of SIMCA®-online.



Sartorius Data Analytics

Phone: +46 40 664 2580

E-mail: umetrics@sartorius.com