



MODDE®

Design of Experiments
Solution

Simplifying Progress

SARTORIUS

Create the Knowledge You Need More Efficiently

The more efficient your Design of Experiments (DOE) solution, the faster you can get products to market. MODDE® is an innovative DOE solution from Sartorius Stedim Data Analytics. Its straightforward graphical interface and support for data analytics lets you interpret your results with confidence. With over 30 years of experience in data analytics, MODDE® gives you the insight to set up your experiment and get it right from the start.

What does MODDE® offer?

MODDE® is a lot more than just DOE software. It also provides a quality analysis on your decisions and looks at the risks – warning you about critical settings and guiding you towards more robust conclusions. It will:

- Reduce the number of required experiments
- Guide you through the set-up of your experiment
- Provide confidence in your data handling
- Help you make better decisions
- Integrate with your systems
- Meet your quality goals

MODDE® at a Glance

- Automated analysis wizard
- Robust optimum identification
- Interactive setpoint analysis with risk estimate
- Design Space visualization
- Generalized subset designs
- Stability testing design setup

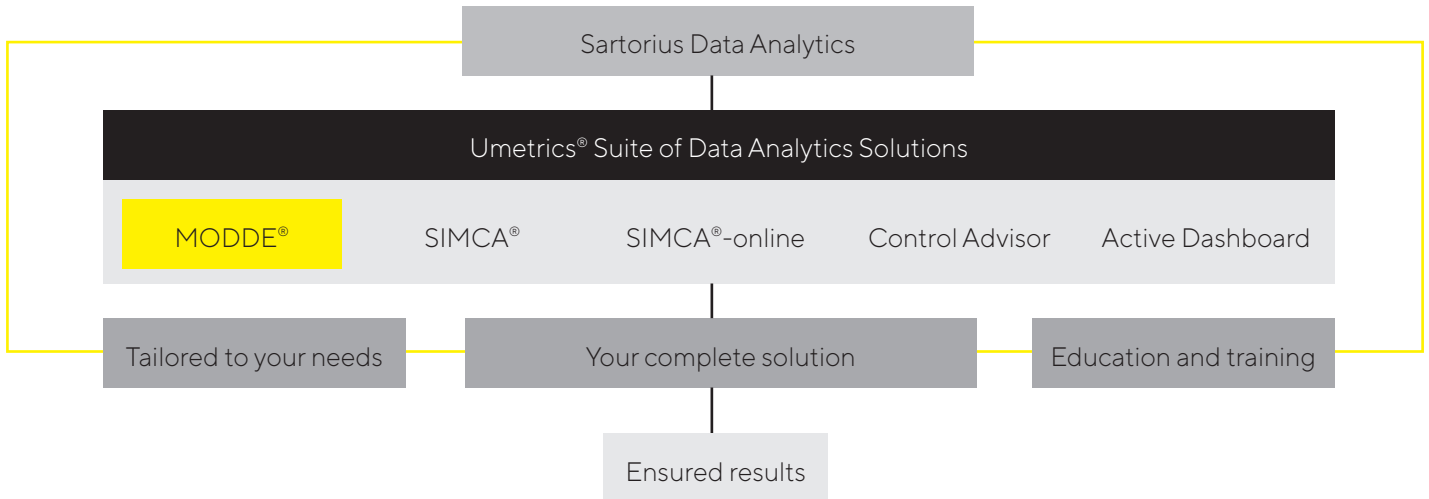


Who is using MODDE®?

MODDE® is helping companies in many different industries to design more effective experiments and create more effective strategies for tackling process problems. For instance:

- In pharma and biopharma, there are a wide variety of aspects to consider when determining the correct tablet formulation. These include how the drug dissolves, its hardness and how it is administered, which can now be addressed with MODDE's advanced formulation DOE toolbox.
- In manufacturing there needs to be a careful balance between productivity and quality. MODDE® helps producers use DOE to maximize production quality through applications like robust optimization – where MODDE® has the best available tools.

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:

SIMCA®

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

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.



“The Umetrics® Suite simplifies the entire process of handling, analyzing and managing our data.”

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