



# Active Dashboard

Interactive Performance  
Insight

Simplifying Progress

SARTORIUS

# Better Insight for Better Decisions

Active Dashboard gives you in-depth, at-your-fingertips insight into production performance across all your sites. That means you can achieve significant savings and ensure product quality.

## Compare Performance Across Your Sites

What if you could compare performance across all of your production sites? Active Dashboard let's you do just that. It takes the data from your SIMCA®-online solutions and visualizes it in a series of easy-to-understand interactive charts. Or you can connect Active Dashboard to other data sources like OSIsoft's PI System™ via the Asset Framework infrastructure.

Active Dashboard gives you options for innovative role-based data visualization. For example, easy, self-service visual analytics with cross-filtering make the involvement of data analysis experts less important. You can view information about your final products and the relationship between processes and product performance. You can also view information about the processes themselves and about your raw materials. So you'll be able to see which sites are performing well and which sites are performing less well and then investigate further to find out why – and do something about it. The bottom line for your business is maximized yield, optimized quality, and lower costs across your production. Active Dashboard gives you the improved transparency you need to make the right decisions. If you need further analysis it will also export seamlessly into SIMCA®.

## Active Dashboard at a Glance

- Production transparency
- Interactive charts and maps
- Product quality assurance
- Performance insight
- A summarized real-time view of all your sites and products

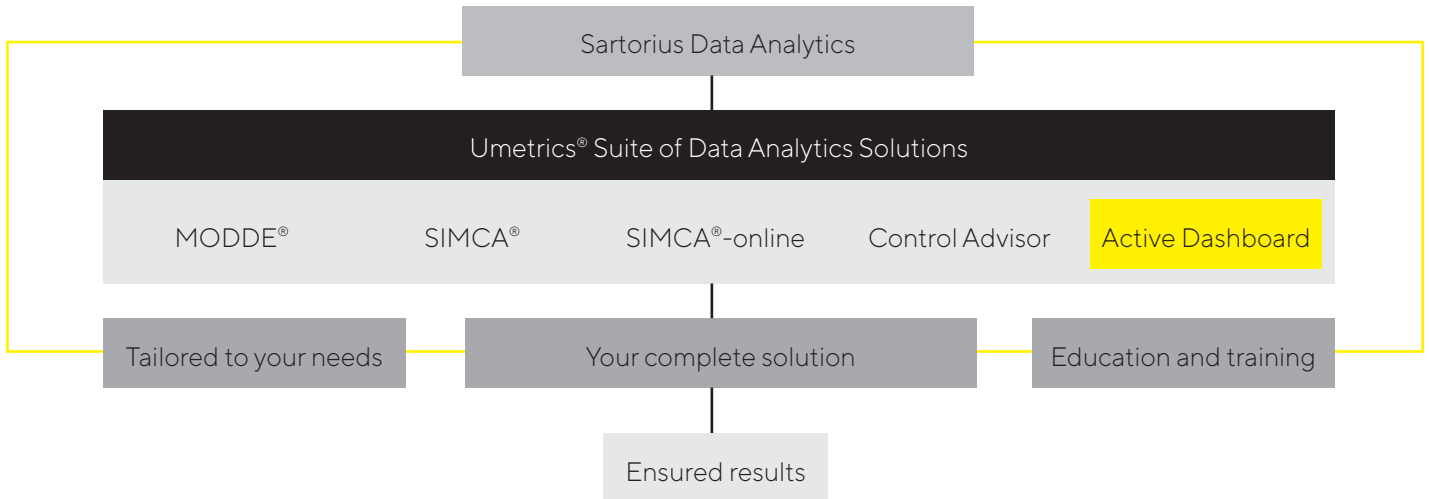


## How does it work?

Active Dashboard allows you to choose the most appropriate from three types of analysis tools.

- View summarized production data across locations, products and time.
- Compare products, processes, and processing performance.
- Use interactive filters to slice and dice data as you prefer.
- Drill down for information about a specific site or sites.
- Take action to improve performance across your sites.

# 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

### SIMCA®-online

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

### Control Advisor

- Predictive capabilities to be able to forecast the output

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.umetrics.com/product/umetrics-active-dashboard](http://www.umetrics.com/product/umetrics-active-dashboard)

# 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](http://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](mailto: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](mailto:umetrics@sartorius.com)

 For further contacts, visit  
[www.sartorius.com/umetrics](http://www.sartorius.com/umetrics)