

Sartorius Stedim Biotech launches new versions of data analytics software solutions

SIMCA® and SIMCA®-online software updates in the Umetrics™ Suite help manufacturing industries make better business decisions and optimize process control

Goettingen, Germany / Malmö, Sweden – December 07, 2017: Sartorius Stedim Biotech (SSB), a leading international supplier for the biopharmaceutical industry, has introduced a new version of its SIMCA® and SIMCA®-online data analytical solutions, which are offered by its subsidiary Sartorius Stedim Data Analytics, formerly known as Umetrics.

Every day, businesses generate a vast array of data derived from a variety of different sources. This data holds the key to better performance. The challenge is to interpret this information in a meaningful way. However, with so many parameters and such vast system complexity to consider, it is hard to find a solution that is both powerful and smart enough. SIMCA®, an established advanced data analytics and visualization program as part of the company's proven Umetrics™ Suite, makes it possible to combine and analyze data from all sources to isolate, understand and act on the hidden gems that hold the secret to better decision-making and greater business success.

SIMCA®'s multivariate data analysis engine enables companies to swiftly detect and analyze deviations from normal operating conditions by modeling an idealized process. Once this model is transferred into SIMCA®-online, it serves as a valuable reference for your current production. The newly enhanced software offers an intuitive graphical interface and the flexibility to handle complex data, such as reworking, splitting and merging, and more. SIMCA® projects can be uploaded directly to an available SIMCA®-online server for real-time visualization of the process from a data point of view.

The real-time monitoring and prediction system, SIMCA®-online, constantly monitors processes to provide a continuous snapshot of the users' operations. It not only helps to identify when set parameters change, but also enables remedial action to be taken before production is affected, ensuring that product quality remains consistent. With this level of control, it is possible to maximize resource efficiency, minimize operational costs and benefit from increased confidence in end-product quality. Among the new features in SIMCA®-online is the self-service analytics capability, which allows anyone to create fundamental process models, regardless of their

background. The new notification system, along with the new web client, gives the user peace of mind about the quality of production anywhere and anytime as this system provides an overview of the production processes on devices such as tablets or mobile phones. SIMCA® and SIMCA®-online have been developed according to GAMP5 and have been extensively tested and validated. These programs are also used by the EMA and FDA for Real-Time Release testing.

See what others don't

As leading experts for analyzing data, Sartorius Stedim Data Analytics helps companies in any industry, which include the pharma, chemical and food sectors, find the growth opportunities they need using the comprehensive Umetrics™ Suite. These solutions enable them to harness the wealth of data within an organization, identifying vital elements to improve the results of research, product development and manufacturing processes. With improved process understanding and more consistent product quality, users will be able to reduce risk, speed up time to market and accelerate business growth.

A complete solution encompasses software, training, support and project management. Also, as part of the Sartorius Group, a global company with approximately 7,500 employees, Sartorius Stedim Data Analytics provides the backup of an international network.

Experience the benefits for your business today

Find out how our solutions can help your business grow, whichever industry you are in. Visit www.umetrics.com to access a trial version of SIMCA® and view the details of our free SIMCA®-online simulation.

Image File:



The real-time monitoring and prediction system, SIMCA®-online, constantly monitors processes to provide a continuous snapshot of the users' operations.

Link for Download: [SIMCA-online photo](#)

A profile of Sartorius Stedim Biotech

Sartorius Stedim Biotech is a leading international supplier of products and services that enable the biopharmaceutical industry to develop and manufacture drugs safely and efficiently. As a total solutions provider, Sartorius Stedim Biotech offers a portfolio covering nearly all steps of biopharmaceutical manufacture. The company focuses on single-use technologies and value-added services to meet the rapidly changing technology requirements of the industry it serves. Headquartered in Aubagne, France, Sartorius Stedim Biotech is quoted on the Eurolist of Euronext Paris. With its own manufacturing and R&D sites in Europe, North America and Asia and an international network of sales companies, Sartorius Stedim Biotech has a global reach. In 2016, the company employed approx. 4,700 people, and earned sales revenue of 1.051,6 million euros.

Contact:

Dominic Grone
Senior Manager Corporate Communications
Sartorius Stedim Biotech
Phone: +49.(0)551.3083324
dominic.grone@sartorius.com