

Information that Gets Right to the Point

The ProControl@Informer software module from Sartorius evaluates data quickly and flexibly

Sartorius AG
37070 Goettingen, Germany

Phone +49(0)551.308.3684
Fax +49(0)551.308.3572
www.sartorius.com

Göttingen, September 05, 2008 – Complete transparency throughout all areas of a company is more important than ever before in today's industrial processes. Particularly in the areas of processing and quality assurance, fast and clear information are of the essence. This is the only way corrective measures can be taken while the process is still running. For this reason, evaluations and data presentations have to "get right to the point" to be relevant.

Sartorius ProControl for Windows, or SPCfWin for short, has long since proven to be invaluable in quality assurance. The SPCfWin QA system can collect process data on a wide variety of quality attributes in samples with individual values, as well as in condensed form from dynamic checkweighers.

The new Sartorius ProControl@Informer software module, or SPCInformer, is the solution for flexible and fast selection and evaluation of all process data collected by SPCfWin and Sartorius ProControl@Inline (SPCInline). Sartorius ProControl@Informer is both easy to use and easy to adapt to individual user requirements and preferences. Users who are accustomed to spreadsheet and database programs will find the operation of SPCInformer intuitive in many aspects, including in its filter and sorting functions. SPCInformer can be installed on the same computer that runs SPCfWin or SPCInline, or on any other computer on the network that has access to the SPCfWin or SPCInline database.

Complex information made simple

The ease of operation offered by Sartorius ProControl@Informer is crucial for the operator. It is no problem to return to the previous processing stage from any point in the program, making it easy to adjust the selection and sorting of data until the display contains exactly the desired information. This saves both time and printing resources. The SPCInformer program structure is straightforward and logical, so it is easy to create even highly complex data analyses for simple applications in minutes. Sophisticated selection options and the possibilities for generating data presentations as printed reports, PDF files or MS Excel files ensures that the most important information is available at any time, providing the user with reliable data for rapid decision-making.

Getting the right information quickly

The SPCInformer system for SPCfWin operates in logical sequences: The type of evaluation is assigned first; this may consist of sampling, statistics or special evaluations, or even special steps or comments. In the second step, data is selected according to time period, product group, product, machine, batch, characteristics, and other criteria. Sophisticated selection options and the use of wildcards enable a high degree of flexibility. Once a particular set of criteria has been selected, the selection can be saved as a template for later use. The third step is the display of data in the form of a table. You can actively influence grouping and sorting by dragging-and-dropping column headers. In the final step, the data is presented as a report. At this point, too, report layouts can be stored as templates for later use. You can use all of the tools available in the powerful MS Excel spreadsheet program to modify the display of data.

A profile of Sartorius

The Sartorius Group is an internationally leading laboratory and process technology provider covering the segments of biotechnology and mechatronics. In 2007, the technology group earned sales revenue of 622.7 million euros. Founded in 1870, the Goettingen-based company currently employs approximately 4,500 persons. The major areas of activity in its biotechnology segment focus on fermentation, filtration, purification, fluid management and laboratory applications. In the mechatronics segment, the company primarily manufactures equipment and systems featuring weighing, measurement and automation technology for laboratory and industrial applications. Key Sartorius customers are in the pharmaceutical, chemical and food and beverage industries as well as in numerous research and educational institutes in the public sector. Sartorius has its own production facilities in Europe, Asia and America as well as sales subsidiaries and local commercial agencies in more than 110 countries.

Image File/Photo for Press Release:



Illustration: The new Sartorius ProControl@Informer software module.

Media Contact: Dominic Grone | Elke Schild, Group Corporate Communications
 Sartorius Corporate Administration GmbH; 37070 Göttingen, Germany
 Phone: +49(0)551.308.3684; fax: +49(0)551.308.3572
 E-mail: dominic.grone@sartorius.com; elke.schild@sartorius.com
 Web: <http://www.sartorius.com>; Press Service: <http://www.sartorius.com/index.php?id=1036>