

**Fluid Management Technology:**

**Opta SFT: Sartorius Stedim Biotech presents a new  
disposable sterile connector**

Sartorius Stedim Biotech S.A.  
Z.I. Les Paluds  
Avenue de Jouques  
B.P. 1051  
13781 Aubagne - France

[www.sartorius-stedim.com](http://www.sartorius-stedim.com)

Goettingen/Aubagne, May, 2009 – Sartorius Stedim Biotech has launched the Opta SFT-I sterile connector, a single-use device that allows sterile connection between two separate, presterilized components in biopharmaceutical manufacturing processes.

The Opta SFT-I sterile connector combines well-established barrier system approaches with an enhanced qualification concept and with improved product features. Opta SFT-I connectors consist of a female and a male coupling body that, once assembled, create a sterile fluid path via easy 4-step operation. In addition to offering 1/2" hose barbs, the Opta SFT-I range includes 1/4" and 3/8" hose barbs. All hose barb dimensions have been specifically designed to ensure the stability and the integrity of the coupling between the hose barb and various types of tubing.

The Opta SFT-I connectors are 100% integrity-tested, gamma irradiated or sterilizable by autoclave to assure the highest quality for critical applications, such as transfer of sterile fluid.

The introduction of the Opta SFT-I sterile connector is a further milestone rounding off Sartorius Stedim's single-use manufacturing range, which enables

users in biomanufacturing to efficiently implement platform technology based disposable solutions.

### **A profile of Sartorius Stedim Biotech**

Sartorius Stedim Biotech is a leading provider of cutting-edge equipment and services for the development, quality assurance and production processes of the biopharmaceutical industry. Its integrated solutions covering fermentation, filtration, purification, fluid management and lab technologies are supporting the biopharmaceutical industry around the world to develop and produce drugs safely, timely and economically. For next generation processes, Sartorius Stedim Biotech focuses on single-use technologies and value-added services to meet the rapidly changing technology requirements of the industry it serves. Strongly rooted in the scientific community and closely allied with customers and technology partners, the company is dedicated to its philosophy of "turning science into solutions" on a daily basis.

Headquartered in Aubagne, France, Sartorius Stedim Biotech is listed on the Eurolist of Euronext Paris. With its own manufacturing and R&D sites in Europe, North America and Asia and a global network of sales companies, Sartorius Stedim Biotech enjoys a worldwide presence. Its key manufacturing and R&D site is in Germany. The company employs over 2,300 people, and in 2008 generated sales revenue of 368.0 million euros.

### **Photo | Image File for Press Release:**



**Image | Photo:** OPTA SFT-I connector.

---

#### Contacts:

Dominic Grone | Stephanie Sinagowitz, Group Corporate Communications  
Sartorius Corporate Administration GmbH; 37070 Goettingen, Germany

Phone: +49(0) 0551/ 308-3965 0551/ 308-3572;

E-Mail: [dominic.grone@sartorius.com](mailto:dominic.grone@sartorius.com); [stephanie.sinagowitz@sartorius.com](mailto:stephanie.sinagowitz@sartorius.com)

<http://www.sartorius.com>; Press Service: <http://www.sartorius.de/index.php?id=73>