

## Sartorius Expands Production Site in England

- New facility at the Stonehouse site for the manufacture of laboratory products
- Officially opened by HRH Princess Anne, The Princess Royal

Goettingen, (Germany) | Stonehouse, England, July 18, 2016 – Sartorius, a leading international pharmaceutical and laboratory equipment provider, today opened a new plant at its Stonehouse site (Gloucestershire), doubling its local production capacities. Together with Princess Anne, member of the Royal Family, and nearly 100 guests, Group CEO Dr. Joachim Kreuzburg and Chris Biddell, the managing director of the U.K. subsidiary Sartorius Stedim Lab Ltd., officially opened the new building covering approximately 4,800 square meters (52,000 sq. ft.). Around 110 employees currently work at this facility in production and R&D.

In his opening speech, Joachim Kreuzburg stated: "Sartorius is growing dynamically, particularly with single-use products as are manufactured here in Stonehouse. To promote this positive development, we are currently investing substantially into the expansion of our global infrastructure. Since our acquisition of Vivascience back in 1999, this Stonehouse site continuously contributed to the strong development of the Sartorius Group. Our local team has achieved a very successful track record, so the new, larger plant is the next logical step."

In Stonehouse, Sartorius develops and manufactures a variety of single-use plastic products, such as filtration units and laboratory-scale bioreactors. The approximately 9-million-euro new plant additionally serves as a competence center within the Sartorius Group for designing and manufacturing plastic components. Such products are used in laboratories and production facilities of pharmaceutical and biopharmaceutical companies and in research labs in academia. "The plant incorporates the latest technology and significantly expands our production capacity," said Biddell. "At the same time, we have built it with an eye toward the future so that we will have sufficient space to once more double our production."

Apart from the Stonehouse site, Sartorius also operates other sites within the United Kingdom in Royston, Epsom and Glasgow and employs nearly 500 people in total.

### **A Profile of Sartorius**

The Sartorius Group is a leading international pharmaceutical and laboratory equipment provider with two divisions: Bioprocess Solutions and Lab Products & Services. Bioprocess Solutions with its broad product portfolio focusing on single-use solutions helps customers produce biotech medications and vaccines safely and efficiently. Lab Products & Services, with its premium laboratory instruments, consumables and services, concentrates on serving the needs of laboratories performing research and quality assurance at pharma and biopharma companies and on those of academic research institutes. Founded in 1870, the company earned sales revenue of more than 1.1 billion euros in 2015. More than 6,200 people work at the Group's 50 manufacturing and sales sites, serving customers around the globe.

### Current Image Files

Opening Ceremony in Stonehouse: HRH Princess Anne; Group CEO Dr. Joachim Kreuzburg and Chris Biddell, the managing director of Sartorius Stedim Lab Ltd.



Sartorius\_Stonehouse\_01.jpg



Sartorius\_Stonehouse\_02.jpg

[https://www.sartorius.com/en/detail/sartorius\\_expands\\_production\\_site\\_in\\_england/](https://www.sartorius.com/en/detail/sartorius_expands_production_site_in_england/)

---

### Contact

Petra Kirchhoff, Vice President of Corporate Communications  
+49 (0)551.308.1686 | [petra.kirchhoff@sartorius.com](mailto:petra.kirchhoff@sartorius.com)