

France's ambassador Descôtes visits Sartorius Stedim Biotech

Göttingen, December 5, 2018 – The ambassador of France to Germany, Anne-Marie Descôtes, visited Sartorius Campus in Göttingen to obtain first-hand information about the development of Sartorius Stedim Biotech. She spoke with the Board of Directors of Sartorius Stedim Biotech headed by CEO Joachim Kreuzburg about the company's rapid growth over more than eleven years since it has been under the Sartorius roof.

"The merger of the two market leaders Stedim Biosystems and Sartorius Bioprocess Technology is a French-German success story – and is viewed as such from the perspective of customers, employees and also shareholders. Over the past decade, Sartorius Stedim Biotech has grown by close to 13% per year, and its share price has increased more than ten-fold. In particular, we have also invested in our French site in Aubagne, where we have approximately quadrupled the number of employees to around 1,000 since 2007," said Kreuzburg.

In 2007, Sartorius AG had merged its Biotechnology Division with French-based Stedim Biosystems S.A., becoming the majority shareholder of Sartorius Stedim Biotech S.A. Quoted on the Euronext in Paris, the company is a market leader for single-use aseptic bags for biopharmaceutical applications, among other products. Currently, more than 5,600 people work at Sartorius Stedim Biotech worldwide.

A profile of Sartorius Stedim Biotech

Sartorius Stedim Biotech is a leading international partner of the biopharmaceutical industry. As a total solutions provider, the company helps its customers to manufacture biotech medications safely, rapidly and economically. 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. The Group has been annually growing by double digits on average and has been regularly expanding its portfolio by acquisitions of complementary technologies. In 2017, Sartorius Stedim Biotech employed approx. 5,100 people and earned sales revenue of €1,081.0 million.

Contact

Timo Lindemann | Corporate Communications
+49 (0)551.308.4724 | timo.lindemann@sartorius.com | www.sartorius.com