

Sartorius Stedim Biotech S.A. - 13781 Aubagne France

## **ANNUAL REPORT OF THE LIQUIDITY CONTRACT WITH THE STOCKBROKER COMPANY GILBERT DUPONT**

Aubagne, France | January 2, 2019

Under the liquidity contract concluded between Sartorius Stedim Biotech S.A. and the stockbroker Gilbert Dupont, the following assets appeared on the liquidity account at December 31, 2018:

- Number of shares: 3,416
- Liquidity account cash balance: € 357,282.50

For information, the following assets appeared on the half-year report of the liquidity contract of June 29, 2018:

- Number of shares: 2,794
- Liquidity account cash balance: € 441,850.72

### **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**

Ben Orzelek  
Investor Relations  
+49(0)551.308.1668  
[ben.orzelek@sartorius.com](mailto:ben.orzelek@sartorius.com)

### **Sartorius Stedim Biotech SA**

Z.I. Les Paluds  
Avenue de Jouques, CS 91051  
13781 Aubagne Cedex, France  
Phone: +33.(0)4.42.84.56.00  
Fax: +33.(0)4.42.84.56.19  
[www.sartorius-stedim.com](http://www.sartorius-stedim.com)