

## Perfect Popcorn Production and Color-Free Fanta – Night of Knowledge at Sartorius

Göttingen, January 15, 2019 – How special data analysis helps to make the perfect microwave popcorn is one technology Sartorius will be demonstrating on January 26 from 5:00 p.m. to 12:00 a.m. on Sartorius Campus. As part of the 4<sup>th</sup> Göttingen Nacht des Wissens – the Night of Knowledge event – visitors can also use Sartorius pipettes and filters to remove colorants from Fanta Orange sodas, weigh their fingerprints or watch on a live-cell analysis system how tumor cells are destroyed. For the first time, Sartorius will be taking part in the City of Göttingen's Night of Knowledge event series.

Sartorius staff members will hold presentations on the growing importance of the biopharmaceutical industry for society and laboratories of the future. In addition, visitors will receive information about the company's history and on how Sartorius products help in the development and manufacture of new medications. This event is free of charge.

To take one of the tours of the Sartorius filter production facilities that will begin every hour, visitors are required to make reservations at [www.ndw.uni-goettingen.de](http://www.ndw.uni-goettingen.de). Shuttle buses will bring visitors to Sartorius Campus and back to the inner city on a regular schedule for free. Visitors can purchase food and beverages on Campus.

### A Profile of Sartorius

The Sartorius Group is a leading international partner of biopharmaceutical research and the industry. With innovative laboratory instruments and consumables, the Group's Lab Products & Services Division concentrates on serving the needs of laboratories performing research and quality control at pharma and biopharma companies and those of academic research institutes. The Bioprocess Solutions Division with its broad product portfolio focusing on single-use solutions helps customers to manufacture biotech medications and vaccines safely and efficiently. The Group has been annually growing by double digits on average and has been regularly expanding its portfolio by acquisitions of complementary technologies. In fiscal 2017, Sartorius earned sales revenue of more than 1.4 billion euros. Currently, more than 8,000 people work at the Group's approximately 60 manufacturing and sales sites, serving customers around the globe.

---

### Contact

Timo Lindemann | Corporate Communications

+49 (0)551.308.4724 | [timo.lindemann@sartorius.com](mailto:timo.lindemann@sartorius.com) | [www.sartorius.com](http://www.sartorius.com)