

Sartorius supports people in Yemen with a Christmas donation

Göttingen, December 12, 2018 – Sartorius is donating 85,000 euros at Christmas time for humanitarian help in Yemen. This year's charitable contribution will be made to the European medical aid organization "action medeor e. V.," which supplies life-saving medications and food supplements to health facilities in Yemen. After around four years of ongoing war and due to a failing economy in Yemen, not all groups of the population have sufficient access to basic health care.

As part of its Christmas Donations Instead of Gifts campaign, Sartorius is supporting a charitable organization in the medical and pharmaceutical sector for the fifth time in a row and, for this special purpose, has again decided to forego giving Christmas gifts to customers and business partners.

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 | www.sartorius.com