

"Donations Instead of Gifts": Sartorius supports people in the Congo

Goettingen, December 20, 2019 - For Christmas, Sartorius is supporting an aid project that improves health care for people in the Democratic Republic of Congo. The company's donation of 85,000 euros will go to action medeor e. V., the European medical aid organization, which is supplying 14 health care facilities, among others, in this African country with medical equipment and nutritional supplements. Groups of rebels, refugee influxes and unstable economic development have caused many people in the Congo to depend on humanitarian aid.

"As a life science group, Sartorius contributes to the development of new and better therapies and affordable medicine. Through our donation, people in the Congo can be provided with life-saving medicines, medical personnel can be trained, and access to drinking water can be improved. In this way, we directly support medical care for people in need," says Petra Kirchhoff, Head of Corporate Communications.

As part of the "Donations Instead of Gifts" initiative, Sartorius is supporting an aid organization from the medical and pharmaceutical sectors for the sixth time, foregoing Christmas presents for customers and business partners. Moreover, Sartorius promotes further charitable initiatives at several company sites: For example, the company has been donating to an SOS Children's Village in Bangalore, India, which gives orphaned children access to medicine and education. This year, 46,000 euros were made available for this purpose.

Follow Sartorius on [Twitter](#) @Sartorius_Group and on [LinkedIn](#).

A profile of Sartorius

The Sartorius Group is a leading international partner of life science research and the biopharmaceutical 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 2018, Sartorius earned sales revenue of around 1.57 billion euros. Currently, more than 8,900 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