

Göttingen, Germany | December 3, 2025

Donations instead of gifts: Sartorius supports laboratory diagnostics in East Africa

- Sartorius supports the “Jambo Labs” project run by action medeor with a donation of 60,000 euros
- The project improves access to vital laboratory services in Tanzania and Zanzibar
- Modern laboratory equipment, training, and quality management systems lay the foundation for reliable diagnoses

Sartorius is continuing its tradition of “donations instead of gifts” and is supporting the “Jambo Labs” project run by the German medical aid organization action medeor with a donation of 60,000 euros this Christmas. The aim of the project is to improve access to vital laboratory services in six hospitals in Tanzania and Zanzibar through equipment, training, and quality management. This will give around 430,000 patients direct access to high-quality diagnostic services each year. In addition, action medeor plans to train laboratory technicians and medical staff.

Many clinics in Tanzania and Zanzibar lack modern laboratory technology and qualified medical staff. This results in delayed or incorrect diagnoses – with sometimes life-threatening consequences for diseases such as malaria, tuberculosis, or diabetes. The “Jambo Labs” project aims to create structures that will improve medical care in the long term through a combination of modern equipment, improved maintenance, training, and quality management. Already, 50 percent more laboratory equipment is in use and 240 additional specialists have been trained for the laboratory teams.

"action medeor and Sartorius as a life science company share the common goal of contributing to better health for more people. We have been supporting the work of the aid organization for around ten years. This has enabled us to promote various projects around the world," said Michael Grosse, CEO of Sartorius. "We also want to thank our customers, partners, and suppliers who support the idea of 'donations instead of gifts'."

About action medeor

For more than 60 years, action medeor has been working as the world's emergency pharmacy to provide sustainable and better health care for people, especially in poorer regions. The non-profit organization is headquartered in Tönisvorst in the Lower Rhine region. From here, medical aid is provided to health stations in around 70 countries in Africa, Latin America, and Asia every year. In emergencies and disasters, action medeor is able to send medicines, medical supplies, and instruments to the disaster area within a

short time. In addition, action medeor carries out health projects in the areas of prevention, diagnostics, and therapy in cooperation with local partners. As part of its pharmaceutical consulting services, action medeor trains health personnel, sets up research laboratories, and supports the establishment of local production facilities for medicines.

Sartorius at a glance

Sartorius is a leading international partner to the biopharmaceutical research and manufacturing industries. The Lab Products & Services Division focuses on innovative laboratory instruments and consumables for research and quality assurance laboratories in pharmaceutical and biopharmaceutical companies as well as academic research institutions. The Bioprocess Solutions Division supports customers with a broad product portfolio focused on single-use solutions for the safer, faster, and more sustainable production of biotech drugs, vaccines, and cell and gene therapies. With around 60 production and sales locations worldwide, the Göttingen-based company has a strong global presence. Sartorius regularly supplements its portfolio with acquisitions of complementary technologies. In 2024, the company generated sales revenue of around 3.4 billion euros. More than 13,500 employees serve customers around the globe.

Visit our [newsroom](#) and follow us on [LinkedIn](#).

Contact

Leona Malorny
Head of External Communications
+49 551 308 4067
leona.malorny@sartorius.com