

Göttingen, Germany | May 21, 2024

Sartorius Refill Towers earn ACT Ecolabel, leading the way in sustainable pipetting

- Sartorius receives first ACT Ecolabel for its pipette tip Refill Towers
- Refill Towers are precision engineered to reduce plastic and packaging waste
- Innovative manufacturing facility powered by 100% renewable electricity reduces environmental impact through recycling partnerships

The life science group Sartorius has reached a significant milestone in sustainability and transparency by earning its first Accountability, Consistency, and Transparency (ACT) Ecolabel for its Pipette Tip Refill Towers. This internationally recognized ecolabel is offered by My Green Lab®, a non-profit organization dedicated to building a global culture of sustainability in science.

The ACT Ecolabel is a standardized measure of the environmental impact of laboratory products. The evaluation process examines multiple criteria across the product life cycle, resulting in the most comprehensive ecolabel available. Its aim is to provide scientists and procurement specialists with clear, third-party verified information, enabling better decision-making when comparing products from different manufacturers.

"Laboratories are increasingly aware of their environmental footprint and are seeking suppliers that help them reduce that impact. Access to clear information is vital for making informed purchase decisions about essential consumables, such as tips," said Maria Voutilainen, Product Manager of Liquid Handling at Sartorius. "The ACT Ecolabel effectively communicates the sustainability attributes of our Refill Towers, simplifying the process of choosing a more eco-friendly product."

Refill Towers are a convenient and ecological way to refill single trays of Sartorius' purity certified Optifit standard pipette tips. They are specifically designed to reduce plastic and packaging waste, encourage the reuse of tip boxes, and lower transportation emissions due to their space-efficient packaging.

All Sartorius pipette tips and Refill Towers are manufactured at the company's Kajaani facility in Finland, which has implemented substantial energy, water, and waste reduction strategies to reduce its environmental footprint. The facility operates on 100% renewable electricity and participates in local recycling programs to repurpose production waste, demonstrating a deep commitment to sustainability.

More information about the ACT Ecolabel and Refill Towers:

<https://www.sartorius.com/en/products/pipetting/pipetting-and-sustainability>

A profile of Sartorius

Sartorius (XETRA: SRT3 | SRT) is a leading international partner of life sciences research and biopharmaceutical manufacturing. With innovative laboratory instruments and consumables, the Group's Lab Products & Services division focuses on laboratories performing research and quality control at pharmaceutical and biopharmaceutical companies as well as academic research institutes. The Bioprocess Solutions division, with its broad product portfolio focusing on single-use solutions, helps customers manufacture biotech medications, vaccines, and cell and gene therapies safely, rapidly, and economically. The company, based in Göttingen, Germany, has a strong global reach with around 60 production and sales sites worldwide. Sartorius regularly expands its portfolio through the acquisition of complementary technologies. In 2023, the company generated sales revenue of around 3.4 billion euros. Currently, around 14,600 employees are working for customers around the globe.

Visit our [Newsroom](#) or follow us on [LinkedIn](#).

About My Green Lab®

My Green Lab® is a non-profit environmental organization with a mission to build a global culture of sustainability in science. The organization is the world leader in developing internationally recognized sustainability standards for laboratories and laboratory products, bringing sustainability to the community responsible for the world's life-changing medical and technical innovations. Laboratories are one of the most resource-intensive spaces in any industry, but they don't have to be. By introducing both a new perspective and proven best practices within a carefully crafted framework, My Green Lab has inspired tens of thousands of scientists to make a positive change in their labs by reducing the environmental impact of their work.

Contact

Lisa Dabney
Head of Marketing Programs
Lab Products & Services Division
+1 734-474-7117
Lisa.Dabney@Sartorius.com.