

Göttingen, Germany | September 9, 2024

Next-generation TFF cassette, Vivaflow® SU, delivers intuitive, effortless, and sustainable ultrafiltration to research scientists

- Enhanced ease-of-use for fast and intuitive setup right out of the box
- All-in-one design can be operated with a standard laboratory pump
- Advanced flow path delivers speed without tedious process optimization
- Extended membrane range in single-use format improves molecule recovery and integrity
- Responsibly made with 30.5% less plastic and more recyclable packaging

The life science group Sartorius is launching Vivaflow® SU, setting a new standard for laboratory-dedicated tangential flow filtration (TFF). Designed for enhanced ease of use and flexibility, Vivaflow® SU ensures more efficient and sustainable ultrafiltration and diafiltration processes for feed volumes ranging from 100 to 1,000 mL.

Vivaflow® SU is a unique TFF cassette that reduces cost and complexity, streamlines molecule isolation workflows, and champions sustainability. Each all-in-one cassette is ready to use with a standard laboratory pump, enabling simple set-up, while avoiding the need for expensive systems and process optimization. It allows for faster and more reliable concentration and rebuffering of any molecule, eliminating the need for cleaning, reducing contamination risk, and minimizing waste.

“Vivaflow® SU is a testament to our promise to always simplify progress,” explains John Cashman, Product Manager for Lab Ultrafiltration at Sartorius. “It takes the ease of use that our customers already value about TFF with Vivaflow® and elevates the experience with improved handling and performance, while reducing waste. Our next generation cassette eliminates the need for feed flow rate, TMP, and permeate flux calculations, while also offering sustainability improvements that are truly impressive for a single-use consumable.”

By significantly reducing single-use waste compared to previous TFF cassettes, Vivaflow® SU demonstrates Sartorius' commitment to environmental responsibility. The new housing uses 30.5% less plastic and incorporates several eco-friendly design improvements, including locally sourced raw materials, customized tubing kits for each cassette configuration, a digital user manual, and reduced non-recyclable packaging, all contributing to a lower research footprint.

Vivaflow® SU replaces the long-standing and successful Vivaflow® 50, as the next generation TFF cassette for research scientists who need a fast and reliable solution for concentrating and rebuffering large feed volumes in their molecule isolation workflows.

More information about Vivaflow® SU:

www.sartorius.com/en/products/lab-filtration-purification/ultrafiltration-devices/tangential-crossflow

A profile of Sartorius

Sartorius 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](#).

Contact

Lisa Dabney

Head of Marketing Programs

Lab Products & Services Division

+1 734 474 7117