

## **Sartolab® MultiStation**

### **For a hands-free filtration of up to 6 x 50 mL samples in parallel**

Goettingen, December 1<sup>st</sup> 2019 – After the introduction of the Sartoclear Dynamics® Lab V kits for harvesting mammalian cell cultures in 2018, Sartorius now completes the Sartolab® RF | BT range and thus the Sartoclear Dynamics® Lab product line with a unit for the filtration of ≤ 50 mL samples.

The new Sartolab® RF 50 vacuum filtration units are available in two versions, one with an asymmetric 0.22 µm PES membrane for sterile filtration and another with a 0.45 µm PES membrane for clarification.

The new Sartoclear Dynamics® lab V50 kits for the filtration of ≤ 50 mL cell cultures comprises the Sartolab® RF 50 for the filtration and either 1 or 2 pouches of filter aid for the clarification. Herewith, the Sartoclear Dynamics® lab product family can now be used for samples from 15 mL up to 1 liter.

Many customers working on small volumes of 30 – 50 mL must filter multiple samples at the same time. This creates a critical pain point in handling: Each unit must be attached to a vacuum tube, then stabilized before filtration and then disconnected again before attaching the next unit to the vacuum source.

Sartorius facilitates the filtration of these smaller sample volumes through the introduction of the new Sartolab® MultiStation, a stand that enables hands-free and parallel filtration of up to 6 samples.

The MultiStation, in combination with the 50 mL vacuum filtration units & the DE filter aid, is the perfect tool for customers struggling with processing many small samples of mammalian cell cultures. The Sartoclear Dynamics® lab V50 kits do not only eliminate the centrifugation step but, together with the MultiStation, facilitate the filtration, avoiding extra time typically consumed for connection and stabilization.

This new product provides a rapid, user-friendly solution unique to the market to simplify and expedite your routine filtration work in the lab!

## Photos | Links for Downloading



**Sartolab® MultiStation** – A stand with one vacuum connection, a rotating, multi-directional head and 6 brackets for hands-free parallel filtration of up to 6 x 50 mL samples

Link for downloading Image above: [Sartolab® MultiStation](#)

For more details on the Sartoclear Dynamics® Lab kits and the Sartolab® MultiStation:  
[www.sartorius.com/sdl](http://www.sartorius.com/sdl)

### **A Profile of Sartorius**

The Sartorius Group is a leading international pharmaceutical and laboratory equipment provider with two divisions: Bioprocess Solutions and Lab Products & Services. Bioprocess Solutions with its broad product portfolio focusing on single-use solutions helps customers produce biotech medications and vaccines safely and efficiently. Lab Products & Services, with its premium laboratory instruments, consumables and services, concentrates on serving the needs of laboratories performing research and quality assurance at pharma and biopharma companies and on those of academic research institutes.

---

**Contact:** Catherine Leluan – Product Manager – Sartorius Lab Instrument GmbH & Co. KG,  
Phone: +49(0)1735368536, email: [catherine.leluan@sartorius.com](mailto:catherine.leluan@sartorius.com); [www.sartorius.com](http://www.sartorius.com)