

SARTORIUS

Sartoclear Dynamics[®] Lab

Filter hands-free up to 6 × 50 mL cell culture samples in parallel with the Sartolab[®] Multistation!

Facilitate your filtration of small volumes (≤ 50 mL) using the new innovative stand, Sartolab[®] Multistation, specially designed to hold 1 to 6 vacuum filtration units, allowing simultaneous filtration of up to 6 samples.

Discover our Solutions: [sartorius.com/sdl](https://www.sartorius.com/sdl)

Simplifying Progress

