

Eliminate Centrifugation

Discover the most convenient
way to harvest your
mammalian cell cultures!





Directly harvest your mammalian cell cultures with Sartoclear Dynamics® Lab

The Sartoclear Dynamics® Lab V Kits combine the use of a filter aid for clarification and a vacuum filtration unit with a 0.22 µm polyethersulfone membrane for sterile filtration of up to 1 L mammalian cell cultures. The ready-to-use filter aid is supplied in dust-free and gamma-irradiated pouches. The Sartolab® RF bottle top filters with receiver flasks, sterilized by gamma irradiation are available in volume capacities ranging from 150 mL to 1,000 mL.

How does it work?

- ▶ Attach the Sartolab® RF unit to a vacuum source
- ▶ Pour the filter aid into your cell culture and mix to homogenize
- ▶ Pour the cell culture | filter aid mixture into the Sartolab® RF and filter

Feel free to request a demo at:
Info-LF@Sartorius.com

or visit us and watch our video:

