

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 beta irradiation are available in volume capacities ranging from 50 mL to 1,000 mL.



How Does It Work?

- Attach the Sartolab® RF unit to a vacuum source (on the Multistation or as standalone)
- 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 visit us and request a demo at:

www.sartorius.com/sdl



Or watch our video:




Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906

 For further contacts, visit
www.sartorius.com