

SARTORIUS

Free Samples

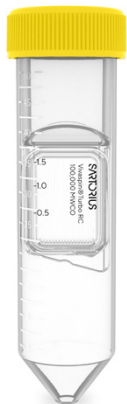
Improve Your Protein Recoveries

Choose the best membrane for your sample with Vivaspin® Turbo; now available in both Regenerated Cellulose (RC) and Polyethersulfone (PES) ultrafiltration membranes.

With ultrafiltration membranes – One size doesn't fit all.

Providing the fastest concentration times on the market with the only 4-15 mL device available with both key ultrafiltration membranes.





Two easy ways to get your free samples:

1. Fill in the fields on this card, take a picture with good light and send it to: **Leads@sartorius.com**
2. Or just scan this QR code and follow the steps.



Sample Request Form

Customer Name

Company | Organization

Address for Sample Delivery

City

State

Email

Phone Number

Volume of Sample for Concentration

Molecular Weight/Size of Your Molecule

Molecular Weight Cut off Currently in Use

General Application

This offer is valid while stocks last. Only one redemption per Lab.
Sample packs contain two devices of the same MWCO