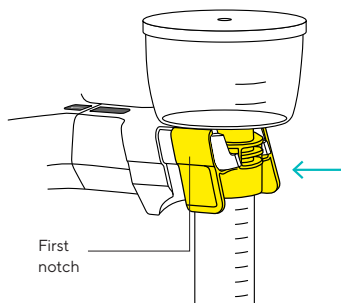


Sartolab® RF 50 Vacuum Filtration Unit (180E01-----E8 | 180F01-----E8)

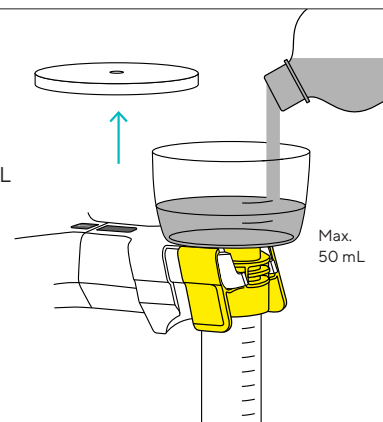
1. Check that the filtrate tube is correctly tightened. Hand-tighten it, if necessary.

Insert the vacuum filtration unit horizontally into the bracket (to first notch) of the MultiStation.

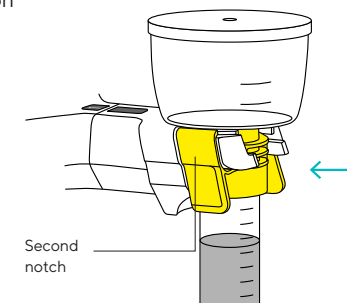


2. Remove the lid and pour the sample to be filtered into the funnel.

Caution
Do not pour more than 50 mL liquid into the funnel as this is the maximum capacity of the filtrate tube; otherwise, excess liquid might leak into the vacuum tubing.

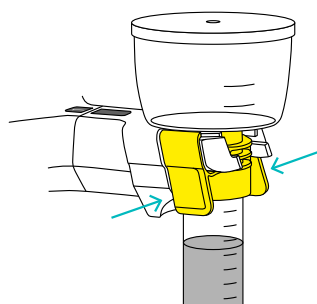


3. Place the lid back on and push the filtration unit to the second notch. This automatically starts filtration when vacuum is applied to the MultiStation.



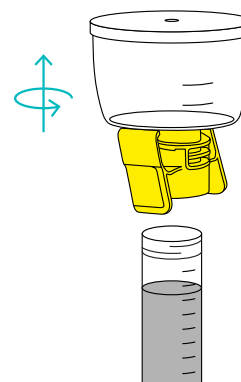
4. To end filtration, gently squeeze the release tabs to remove the vacuum filtration unit from the second notch to the first notch and disengage the vacuum from the device.

To remove the vacuum filtration unit from the MultiStation, squeeze the release tabs and pull the vacuum filtration unit away from the first notch, sliding it off the bracket.

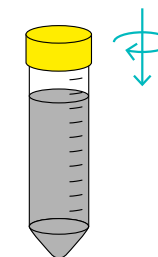


5. To detach the conical tube from the unit, grasp the funnel and unscrew it, along with the yellow adapter, then lift it off the tube.

Discard the upper part in accordance with your local legal regulations for waste disposal.



6. Use the screw cap to seal the filtrate tube for storage of the filtered sample.



Germany

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



Discover the complete
Sartolab® RF | BT product line:
www.sartorius.com/SartolabRFBT

For further information: www.sartorius.com/SartolabRFBTmanuals

Specifications subject to change without notice.
Copyright Sartorius Lab Instruments GmbH & Co. KG.
Material No.: 1000088974
Status: 04 | 2021