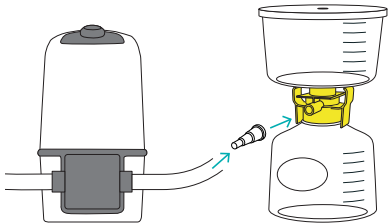
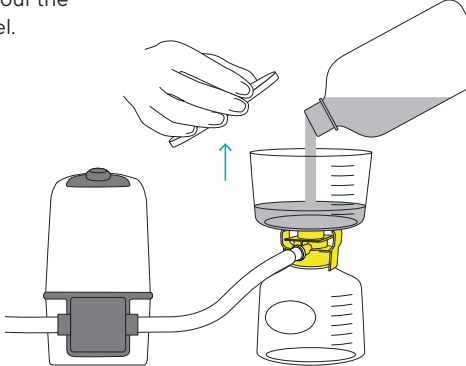
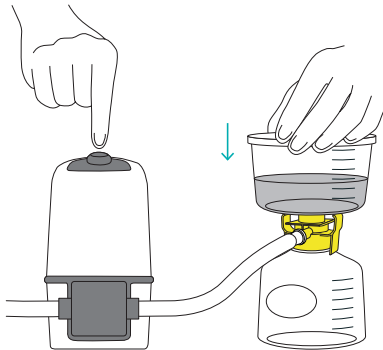
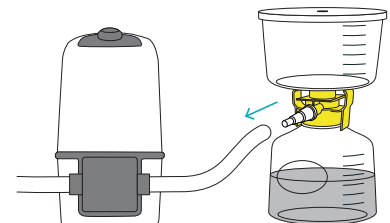
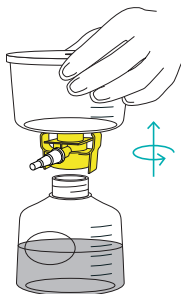
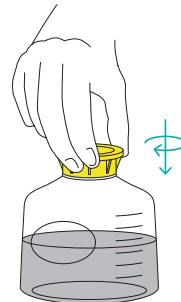


Sartolab® RF 150 – 1000 Vacuum Filtration Units Sartolab® BT 150 – 1000 Bottle Top Filters

Stand-Alone Filtration

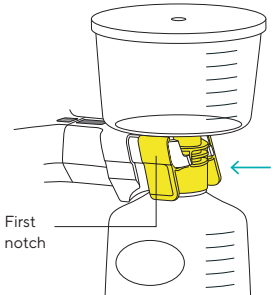
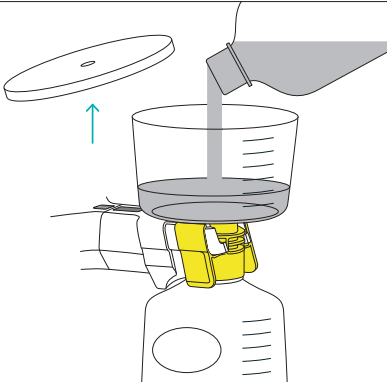
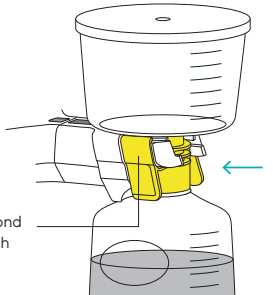
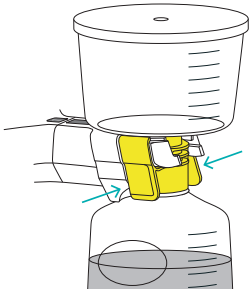
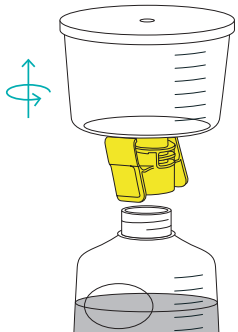
<p>1. Unpack the filtration device. For bottle top filters (Sartolab® BT), screw the funnel onto a bottle with a 45 mm neck.</p> <p>Unpack the tube connector and attach it to the yellow vacuum adapter. Connect the tube connector to the vacuum tube of the vacuum source.</p> 	<p>2. Remove the lid and pour the sample into the funnel.</p> 	<p>3. Place the lid back on. Apply vacuum to start filtration.</p> 
<p>4. To end filtration, turn off the vacuum source and detach the tubing from the connector.</p> 	<p>5. To detach the bottle, grasp the funnel at the top and unscrew it, along with the yellow adapter, then lift it off the bottle.</p> <p>Discard the funnel and the lid in accordance with your local legal regulations for waste disposal.</p> 	<p>6. For vacuum filtration units (Sartolab® RF), unpack the sterile screw cap and securely close the bottle for storage.</p> <p>For bottle top filters (Sartolab® BT), close the bottle with an appropriate screw cap.</p> 



Sartolab® RF 150 – 1000 Vacuum Filtration Units

Sartolab® BT 150 – 1000 Bottle Top Filters

Multiple Filtration Runs Using the Sartolab® MultiStation

<p>1. Unpack the filtration device. For bottle top filters (Sartolab® BT), screw the funnel onto a bottle with a 45 mm neck.</p> <p>Insert the vacuum filtration unit horizontally into the bracket (to first notch) of the MultiStation.</p>  <p>First notch</p>	<p>2. Remove the lid and pour the sample to be filtered into the funnel.</p> 	<p>3. Place the lid back on and push the filtration unit to the second notch.</p> <p>This automatically starts filtration when vacuum is applied to the MultiStation.</p>  <p>Second notch</p>
<p>4. To end filtration, gently squeeze the release tabs to move the vacuum filtration unit from the second notch up to the first notch and disengage the vacuum source from the device.</p> <p>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.</p> 	<p>5. To detach the bottle, grasp the funnel and unscrew it, along with the yellow adapter, then lift it off the bottle.</p> <p>Discard the funnel and the lid in accordance with your local legal regulations for waste disposal.</p> 	<p>6. For vacuum filtration units (Sartolab® RF), unpack the sterile screw cap and securely close the bottle for storage.</p> <p>For bottle top filters (Sartolab® BT), close the bottle with an appropriate screw cap.</p> 