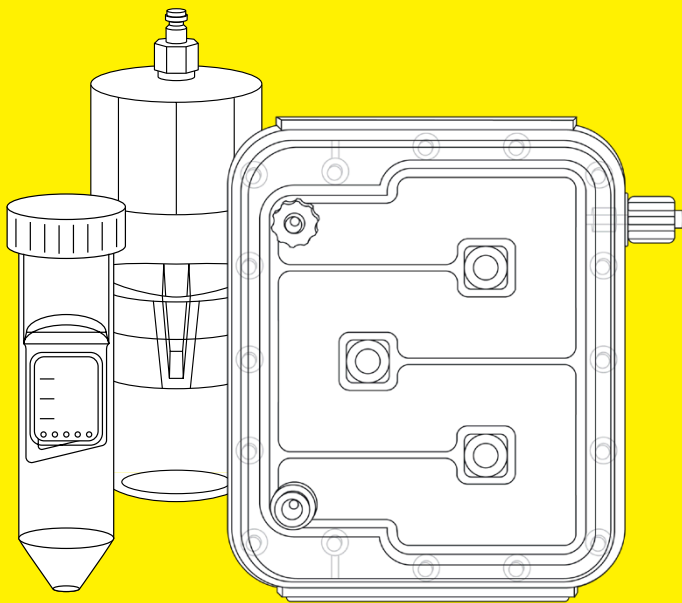


3 Steps to Choose Your Ultimate Ultrafilter



1



Confirm the **capacity** and minimum retentate volume will meet your required feed and final volumes.

2



Determine a suitable operating **method** for your sample and the equipment available in your laboratory.

3




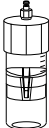

Select a **membrane** MWCO and material to support optimal process speed and molecule recovery.

SARTORIUS

Find Your Ultrafilter:



3 Steps to Choose Your Ultimate Ultrafilter

Capacity		Method	Membrane										
	Vivaspin® and Vivaspin® Turbo	Spin using a centrifuge	PES:	3	5	10	30	50	100	300	1,000 kDa	0.2 µm	
	Concentrate 0.1 - 90 mL feeds up to 100X or more		RC:	2	5	10	30	50	100		kDa		
	Vivaspin®	Pressure using compressed gas	PES:	3	5	10	30	50	100	300	1,000 kDa	0.2 µm	
	Concentrate 5 - 100 mL feeds up to 100X and more												
	Vivaflow®	TFF using a peristaltic pump	PES:	3	5	10	30	50	100	300	1,000 kDa	0.2 µm	
	Concentrate 100 - 2,500 mL feeds up to 100 - 250X		RC:	2	5	10		50	100	300	kDa		

Note: Not all methods and membranes are shown, and not all membranes are available in all products. Please refer to the laboratory ultrafiltration brochure and relevant datasheets for more information.

SARTORIUS

Find Your Ultrafilter:

