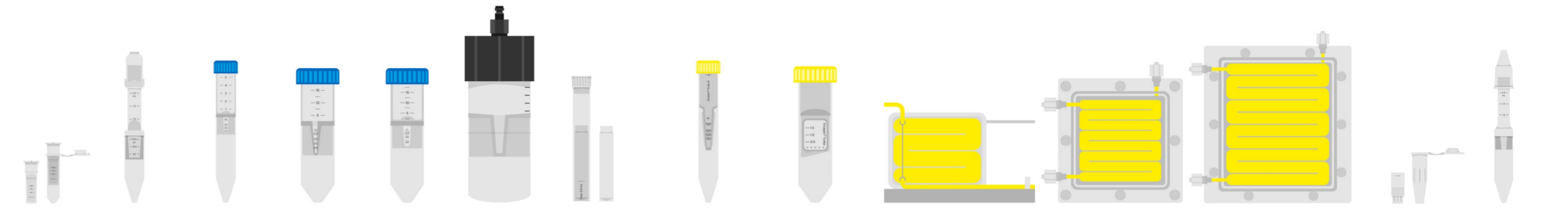


Laboratory Ultrafiltration Quick Reference Guide

This guide serves as a quick reference to the Sartorius lab ultrafiltration range. It summarizes the membrane materials, MWCOs and typical technical specifications for each product. For further detail, please refer to the brochure, data sheets and instructions for use.

Membranes and MWCOs



	500	2	Vivaspin®				Filtrate	Vivaspin® Turbo		50	Vivaflow®		Vivacon®	
			6	15R	20	100		4	15		50R	200	500	2
Polyethersulfone (PES)														
3 kDa	■	■	■		■			■	■	■		■		
5 kDa	■	■	■		■	■		■	■	■		■		
7.5 kDa														
10 kDa	■	■	■		■	■		■	■	■		■		
30 kDa	■	■	■		■	■		■	■	■		■		
50 kDa	■	■	■		■	■		■	■	■		■		
100 kDa	■	■	■		■	■	■	■	■	■		■		
300 kDa	■		■		■	■	■							
1,000 kDa	■		■		■	■				■				
0.2 µm	■		■		■	■				■		■		
Regenerated cellulose (HY RC)														
2 kDa		■		■									■	■
5 kDa		■		■					■		■	■		
10 kDa		■		■					■		■	■	■	■
30 kDa				■					■		■	■	■ ³	■ ³
50 kDa		■							■				■	■
100 kDa									■		■	■	■ ³	■ ³
Cellulose Triacetate (CTA)														
5 kDa		■												
10 kDa		■												
20 kDa		■		■ ¹										
125 kDa ¹													■ ^{2,3}	■ ^{2,3}

¹Vivaspin® Endotest
²Cellulose Acetate (CA)
³PCR grade versions also available

Technical Specifications

	500	2	Vivaspin®				Filtrate	Vivaspin® Turbo		50	Vivaflow®		Vivacon®	
			6	15R	20	100		4	15		50R	200	500	2
Dimensions														
L Ø or L W H (mm)	50 11	126 17	122 17	116 30	116 30	123 62	93 14	122.5 17	118 27	25 107 84	24 100 100	38 126 138	45 12.4	125 16
Active membrane area (cm ²)	0.5	1.2	2.5	3.9	6	23.5	0.79	3.2	7.2 or 8.6	50	50	200	0.32	0.95
Dead-stop volume (µL)	5	8	30	30	50	350	100	40-60	60-100	-	-	-	5	55
Hold-up volume (µL)	< 5	< 10	< 10	< 20	< 20	< 250	< 5	< 10	< 10	< 500	< 500	< 1,000	< 5	10
Centrifuge (Swing-Out)														
Capacity (mL)		3	6	15	20	90	2.5	4	15					
Maximum RCF (g)		4,000	4,000	3,000	4,000 ⁴	2,000	2,500	4,000 ⁴	4,000 ⁵					
Centrifuge (Fixed Angle)														
Capacity (mL)	0.5	2	6	12.5	14		2.5	4	9				0.5	2
Maximum RCF (g)	12,000	8,000	8,000 ⁴	6,000	6,000		2,000	7,500 ⁴	4,000 ⁵				2,500 ⁶	2,500 ⁶
Pressure														
Capacity (mL)					15	98								
Maximum pressure (bar)					5	5								
Pressure-Fuge max. RCF (g)					3,000 ⁴	n/a								
Pressure-Shake max. shake (rpm)					n/a	300								
Crossflow														
Capacity, single cassette (mL)										500	500	2,500		
Capacity, multiple cassettes (mL)										3,000	1,000	5,000		
Maximum pressure (bar)										3	4	4		

⁴Max. RCF may differ for devices with ≥ 100 kDa MWCO
⁵Max. RCF may differ for devices with ≥ 100 kDa MWCO and different membrane materials
⁶Max. RCF may differ dependent on MWCO, or different target molecule types | sizes. RCF indicated in this table is also the maximum value for reverse spin recovery of the retentate