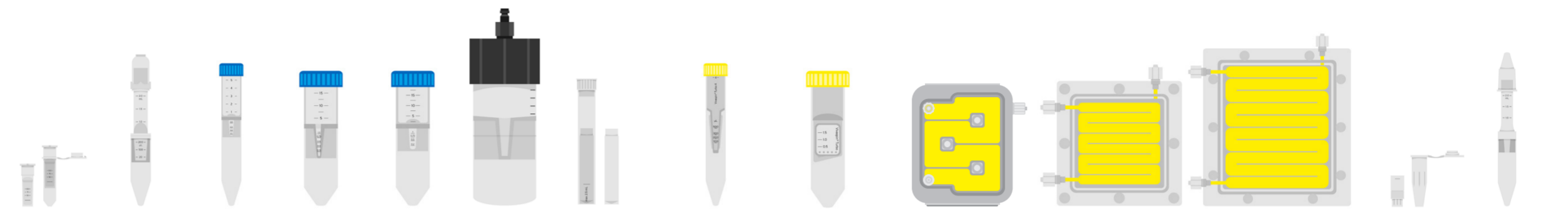


Laboratory Ultrafilters: Quick Reference

This guide summarizes the membrane materials, MWCOs and typical technical specifications for Sartorius laboratory ultrafilters. For further detail, please refer to the datasheets and instructions for use.

Membranes and MWCOs



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

¹Including Vivaspin® Endotest (20 kDa CTA).

²PCR grade (EO treated) options available for 30 kDa (Vivacon® 500), 100 kDa (Vivacon® 500 | 2) and 0.2 µm (Vivacon® 2).

³Cellulose acetate (CA) membrane.

Technical Specifications

	500	2	Vivaspin®				Filtrate	Vivaspin® Turbo		SU	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	98 13 116	100 24 100	126 38 138	45 12.4	125 16
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
Min. retentate volume (µL)	5	8	30	30	50	350	100	40-60	60-100	10,000	10,000	20,000	5	55
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 ⁶				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 ⁷	2,500 ⁷
Pressure														
Capacity (mL)					15	98								
Maximum pressure (bar)					5	5								
Pressure-Fuge max. RCF (g)					3,000 ⁴									
Pressure-Shake max. shake (rpm)						300								
TFF (Crossflow)														
Capacity, single cassette (mL)										500	500	2,500		
Capacity, two cassettes (mL)										1,000	1,000	5,000		
Max. retentate pressure (bar)										2.5	3.5	3.5		

⁴Max. RCF may differ for ultrafilters with ≥ 100 kDa MWCO.

⁵Max. RCF may differ for ultrafilters with ≥ 100 kDa MWCO and for different membrane materials.

⁶Capacity 11 mL for Vivaspin® Turbo 15 with PES membranes.

⁷Max. 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 retentate collection.