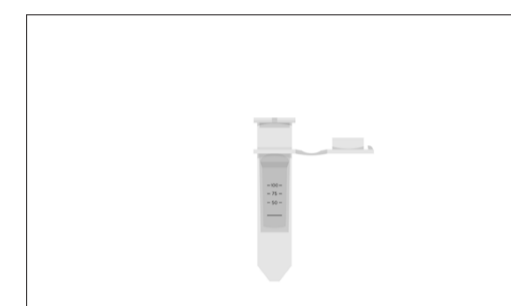


## Laboratory Ultrafilters

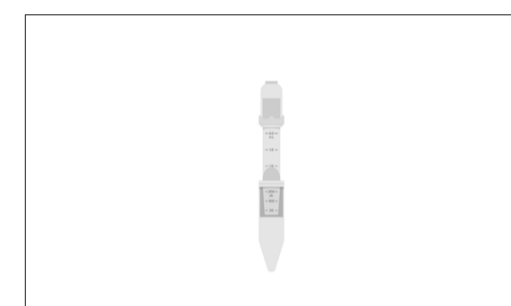
From Centrifugation to Crossflow: Unmatched Versatility and Performance for Any UF and DF Application



Vivaspin® 500

Feed volume: 0.1 – 0.5 mL

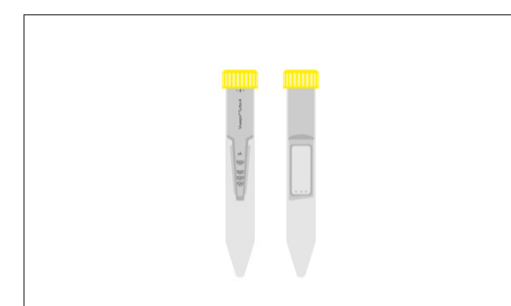
PES	25pc	100pc
3 kDa	V50191	V50192
5 kDa	V50111	V50112
10 kDa	V50101	V50102
30 kDa	V50121	V50122
50 kDa	V50131	V50132
100 kDa	V50141	V50142
300 kDa	V50151	V50152
1,000 kDa	V50161	V50162
0.2 µm	-	V50172



Vivaspin® 2

Feed volume: 0.4 – 3 mL

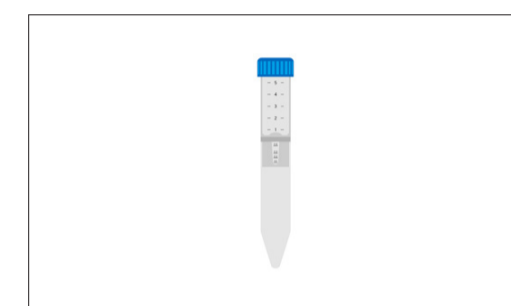
PES	25pc	100pc
3 kDa	V50291	V50292
5 kDa	V50211	V50212
10 kDa	V50201	V50202
30 kDa	V50221	V50222
50 kDa	V50231	V50232
100 kDa	V50241	V50242
<b>RC</b>		
2 kDa	V502H91	V502H92
5 kDa	V502H11	V502H12
10 kDa	V502H01	V502H02
30 kDa	V502H21	V502H22
<b>CTA</b>		
10 kDa	-	V502V2
20 kDa	-	V502X2



Vivaspin® Turbo 4

Feed volume: 1 – 4 mL

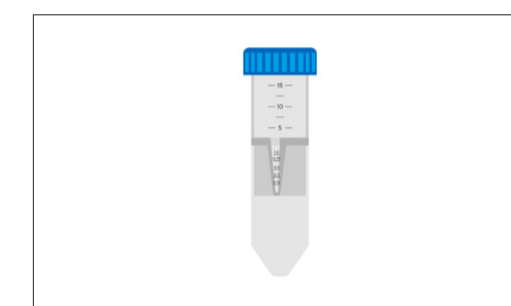
PES	25pc	100pc
3 kDa	V504T91	V504T92
5 kDa	V504T11	V504T12
10 kDa	V504T01	V504T02
30 kDa	V504T21	V504T22
50 kDa	V504T31	V504T32
100 kDa	V504T41	V504T42



Vivaspin® 6

Feed volume: 2 – 6 mL

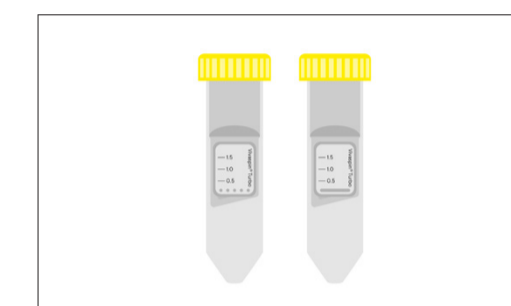
PES	25pc	100pc
3 kDa	V50691	V50692
5 kDa	V50611	V50612
10 kDa	V50601	V50602
30 kDa	V50621	V50622
50 kDa	V50631	V50632
100 kDa	V50641	V50642
300 kDa	V50651	V50652
1,000 kDa	V50661	V50662
0.2 µm	-	V50672



Vivaspin® 15R

Feed volume: 2 – 15 mL

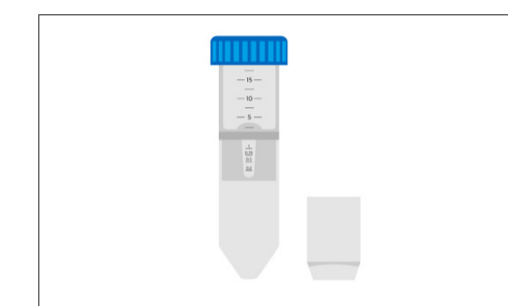
RC	12pc	48pc
2 kDa	V515RH91	V515RH92
5 kDa	V515RH11	V515RH12
10 kDa	V515RH01	V515RH02
30 kDa	V515RH21	V515RH22



Vivaspin® Turbo 15

Feed volume: 4 – 15 mL

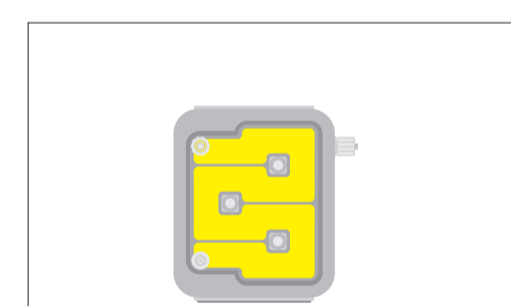
PES	12pc	48pc
3 kDa	V515T91	V515T92
5 kDa	V515T11	V515T12
10 kDa	V515T01	V515T02
30 kDa	V515T21	V515T22
50 kDa	V515T31	V515T32
100 kDa	V515T41	V515T42
<b>RC</b>		
5 kDa	V515TR11	V515TR12
10 kDa	V515TR01	V515TR02
30 kDa	V515TR21	V515TR22
50 kDa	V515TR31	V515TR32
100 kDa	V515TR41	V515TR42



Vivaspin® 20

Feed volume: 5 – 20 mL

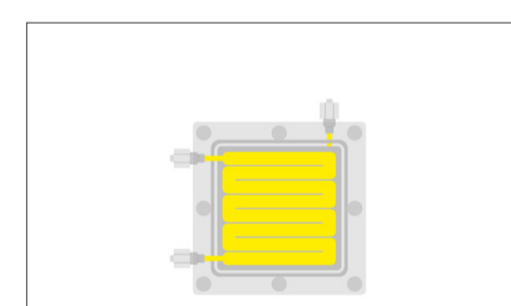
PES	12pc	48pc
3 kDa	V52091	V52092
5 kDa	V52011	V52012
10 kDa	V52001	V52002
30 kDa	V52021	V52022
50 kDa	V52031	V52032
100 kDa	V52041	V52042
300 kDa	V52051	V52052
1,000 kDa	V52061	V52062
0.2 µm	-	V52072



Vivaflow® SU

Feed volume: 0.1 - 1 L

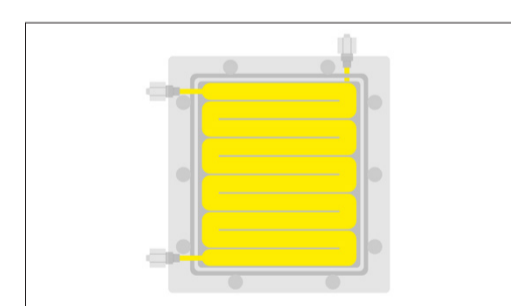
PES	2pc <sup>1</sup>
5 kDa	VF-S050P0005-IV
10 kDa	VF-S050P0010-IV
30 kDa	VF-S050P0030-IV
50 kDa	VF-S050P0050-IV
100 kDa	VF-S050P0100-IV
300 kDa	VF-S050P0300-IV
1,000 kDa	VF-S050P1000-IV
0.2 µm	VF-S050P2000-IV
<b>RC</b>	
2 kDa	VF-S050H0002-IV
5 kDa	VF-S050H0005-IV
10 kDa	VF-S050H0010-IV
30 kDa	VF-S050H0030-IV
100 kDa	VF-S050H0100-IV
300 kDa	VF-S050H0300-IV



Vivaflow® 50R

Feed volume: 0.1 - 1 L

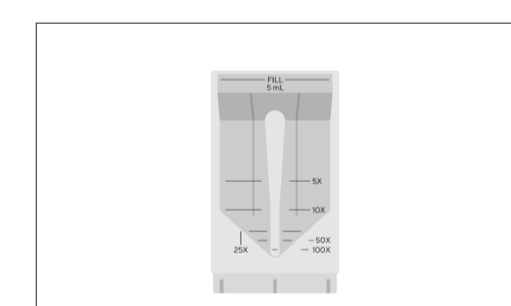
RC	1pc
5 kDa	VF05H1
10 kDa	VF05H0
30 kDa	VF05H2
100 kDa	VF05H4



Vivaflow® 200

Feed volume: 0.5 - 5 L

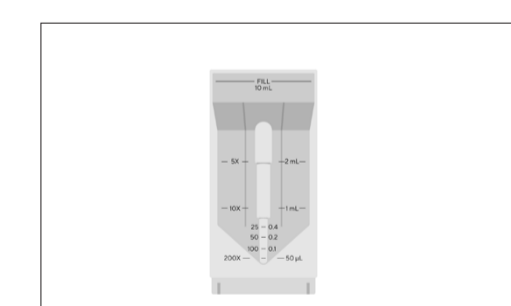
PES	1pc
3 kDa	VF20P9
5 kDa	VF20P1
10 kDa	VF20P0
30 kDa	VF20P2
50 kDa	VF20P3
100 kDa	VF20P4
0.2 µm	VF20P7
<b>RC</b>	
2 kDa	VF20H9
5 kDa	VF20H1
10 kDa	VF20H0
30 kDa	VF20H2
100 kDa	VF20H4



Vivapore® 5

Feed volume: 1- 5 mL

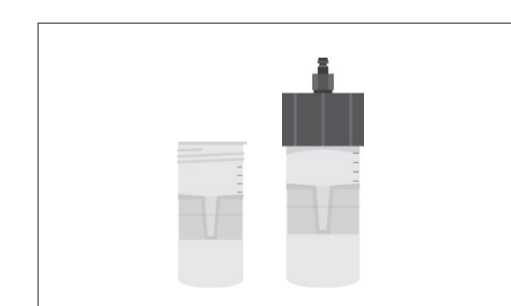
PES	30pc
7.5 kDa	VP-S005P0008-3



Vivapore® 10

Feed volume<sup>2</sup>: 2 - 20 mL

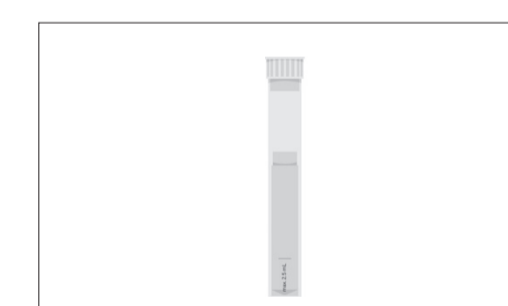
PES	30pc
7.5 kDa	VP-S010P0008-3



Vivaspin® 100

Feed volume: 20 – 98 mL

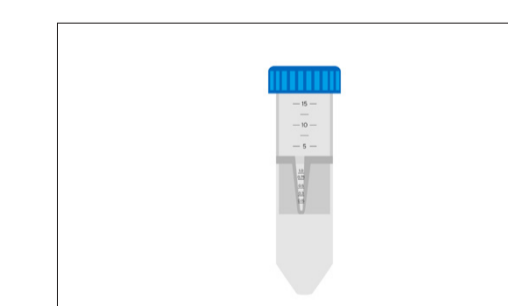
PES	2pc	10pc
5 kDa	VC1011	VC1012
10 kDa	VC1001	VC1002
30 kDa	VC1021	VC1022
50 kDa	VC1031	VC1032
100 kDa	VC1041	VC1042
300 kDa	VC1051	VC1052
1,000 kDa	VC1061	VC1062
0.2 µm	-	VC1072



Vivaspin® Filtrate

Feed volume: 0.5 - 2.5 mL

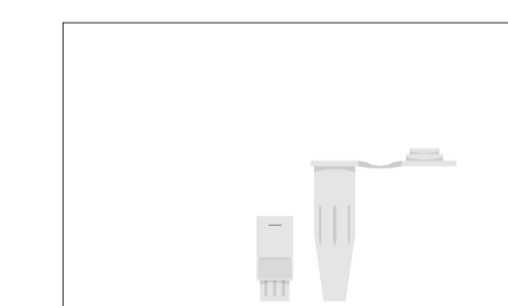
CTA	12pc
5 kDa	13229-E
10 kDa	13239-E
20 kDa	13249-E
<b>PES</b>	
100 kDa	13269-GE
300 kDa	13279-E



Vivaspin® Endotest

Feed volume: 2 – 15 mL

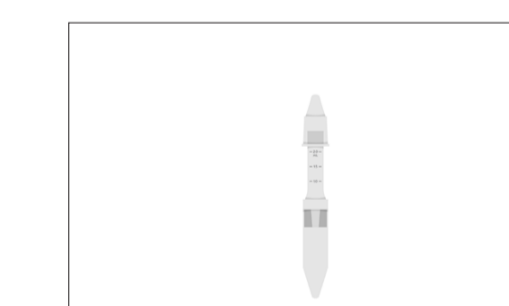
CTA	12pc
20 kDa	V515RXETO



Vivacon® 500

Feed volume: 0.1 – 0.5 mL

RC	25pc	100pc
2 kDa	VN01H91	VN01H92
10 kDa	VN01H01	VN01H02
30 kDa	VN01H21	VN01H22 <sup>3</sup>
50 kDa	VN01H31	VN01H32
100 kDa	VN01H41	VN01H42 <sup>3</sup>
<b>CA</b>		
125 kDa	-	VN01H82



Vivacon® 2

Feed volume: 0.4 – 2 mL

RC	100pc
2 kDa	VN02H92
10 kDa	VN02H02
30 kDa	VN02H22
50 kDa	VN02H32
100 kDa	VN02H42 <sup>3</sup>
<b>CA</b>	
125 kDa	VN02H82
0.2 µm	VN02C75ALE <sup>4</sup>

CTA = cellulose (tri)acetate  
 PES = polyethersulfone  
 RC = regenerated cellulose (including Hydrosart® RC)

<sup>1</sup> For feed volumes >0.5 to ≤1 L, replace -IV with -5V at the end of the order number.

<sup>2</sup> For feed volumes >10 to ≤20 mL, use an expansion reservoir (order no. VP-ARV0010-D).

<sup>3</sup> For PCR grade (EO treated) options, add ETO at the end of the order number.

<sup>4</sup> PCR grade (EO treated) product, 500pc with CA membrane and 2X permeate tubes, caps and lids per ultrafilter.