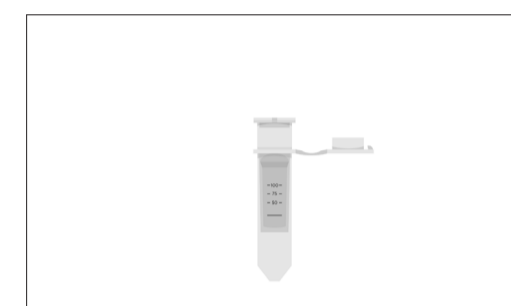


Laboratory Ultrafilters

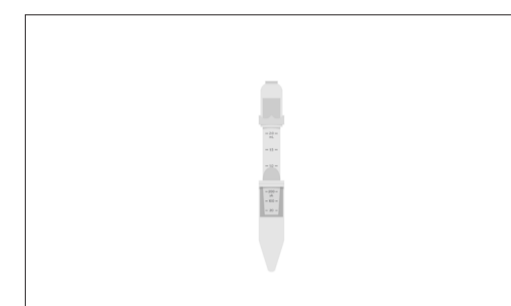
For Any Ultrafiltration and Diafiltration Application



Vivaspin® 500

Initial volume: 0.1 – 1 mL

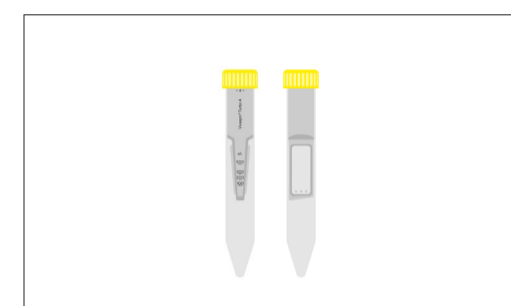
PES	25pc	100pc
3 kDa	VS0191	VS0192
5 kDa	VS0111	VS0112
10 kDa	VS0101	VS0102
30 kDa	VS0121	VS0122
50 kDa	VS0131	VS0132
100 kDa	VS0141	VS0142
300 kDa	VS0151	VS0152
1,000 kDa	VS0161	VS0162
0.2 µm	-	VS0172



Vivaspin® 2

Initial volume: 0.4 – 3 mL

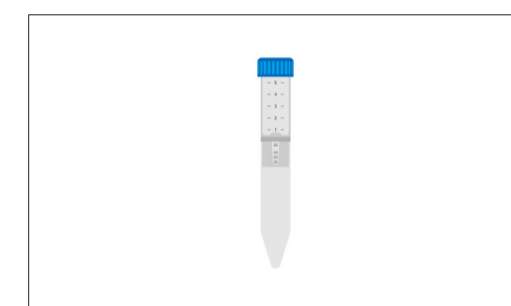
PES	25pc	100pc
3 kDa	VS0291	VS0292
5 kDa	VS0211	VS0212
10 kDa	VS0201	VS0202
30 kDa	VS0221	VS0222
50 kDa	VS0231	VS0232
100 kDa	VS0241	VS0242
HY		
2 kDa	VS02H91	VS02H92
5 kDa	VS02H11	VS02H12
10 kDa	VS02H01	VS02H02
30 kDa	VS02H21	VS02H22
CTA		
10 kDa	-	VS02Y2
20 kDa	-	VS02X2



Vivaspin® Turbo 4

Initial volume: 1 – 4 mL

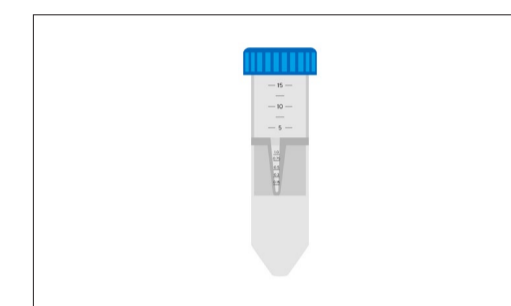
PES	25pc	100pc
3 kDa	VS04T91	VS04T92
5 kDa	VS04T11	VS04T12
10 kDa	VS04T01	VS04T02
30 kDa	VS04T21	VS04T22
50 kDa	VS04T31	VS04T32
100 kDa	VS04T41	VS04T42



Vivaspin® 6

Initial volume: 2 – 6 mL

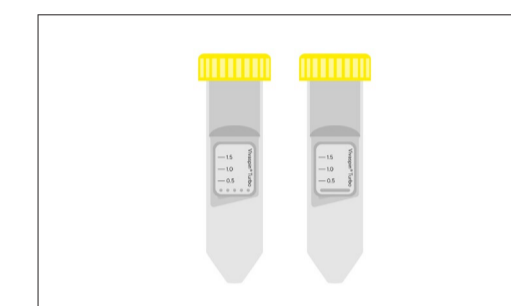
PES	25pc	100pc
3 kDa	VS0691	VS0692
5 kDa	VS0611	VS0612
10 kDa	VS0601	VS0602
30 kDa	VS0621	VS0622
50 kDa	VS0631	VS0632
100 kDa	VS0641	VS0642
300 kDa	VS0651	VS0652
1,000 kDa	VS0661	VS0662
0.2 µm	-	VS0672



Vivaspin® 15R

Initial volume: 2 – 15 mL

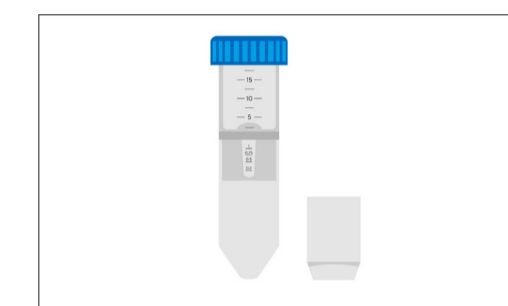
HY	12pc	48pc
2 kDa	VS15RH91	VS15RH92
5 kDa	VS15RH11	VS15RH12
10 kDa	VS15RH01	VS15RH02
30 kDa	VS15RH21	VS15RH22



Vivaspin® Turbo 15

Initial volume: 4 – 15 mL

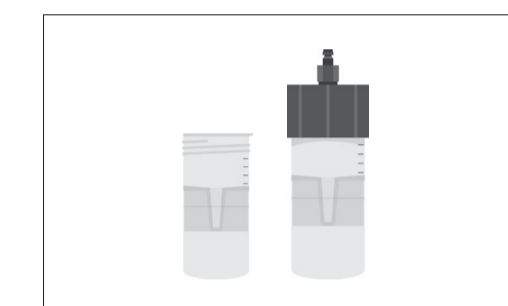
PES	12pc	48pc
3 kDa	VS15T91	VS15T92
5 kDa	VS15T11	VS15T12
10 kDa	VS15T01	VS15T02
30 kDa	VS15T21	VS15T22
50 kDa	VS15T31	VS15T32
100 kDa	VS15T41	VS15T42
RC		
5 kDa	VS15TR11	VS15TR12
10 kDa	VS15TR01	VS15TR02
30 kDa	VS15TR21	VS15TR22
50 kDa	VS15TR31	VS15TR32
100 kDa	VS15TR41	VS15TR42



Vivaspin® 20

Initial volume: 5 – 20 mL

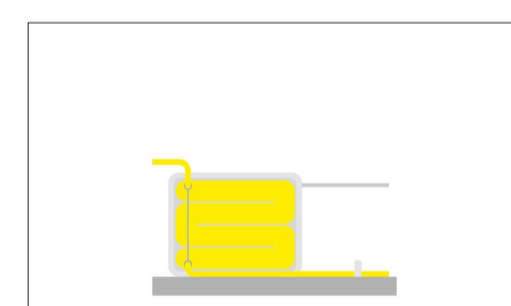
PES	25pc	100pc
3 kDa	VS2091	VS2092
5 kDa	VS2011	VS2012
10 kDa	VS2001	VS2002
30 kDa	VS2021	VS2022
50 kDa	VS2031	VS2032
100 kDa	VS2041	VS2042
300 kDa	VS2051	VS2052
1,000 kDa	VS2061	VS2062
0.2 µm	-	VS2072



Vivaspin® 100

Initial volume: 20 – 98 mL

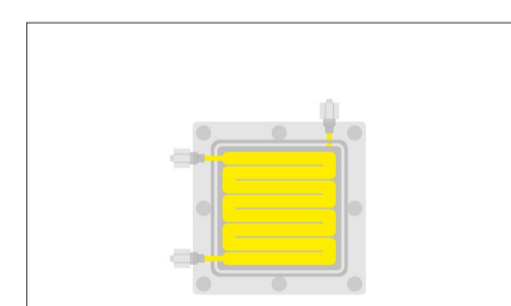
PES	2pc	10pc
5 kDa	VC1011	VC1012
10 kDa	VC1001	VC1002
30 kDa	VC1021	VC1022
50 kDa	VC1031	VC1032
100 kDa	VC1041	VC1042
300 kDa	VC1051	VC1052
1,000 kDa	VC1061	VC1062
0.2 µm	-	VC1072



Vivaflow® 50

Initial volume: 0.1 – 3 L

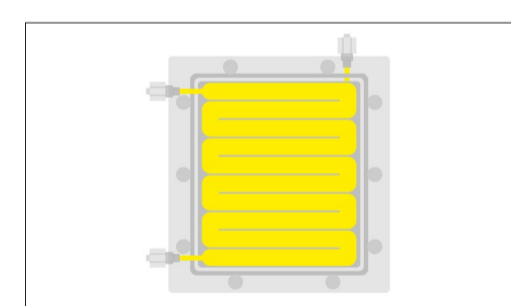
PES	2pc
3 kDa	VF05P9
5 kDa	VF05P1
10 kDa	VF05P0
30 kDa	VF05P2
50 kDa	VF05P3
100 kDa	VF05P4
1,000 kDa	VF05P6
0.2 µm	VF05P7



Vivaflow® 50R

Initial volume: 0.1 – 1 L

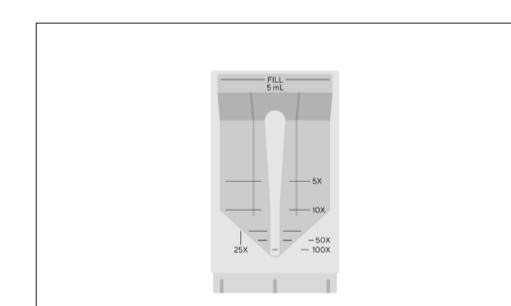
HY	2pc
5 kDa	VF05H1
10 kDa	VF05H0
30 kDa	VF05H2
100 kDa	VF05H4



Vivaflow® 200

Initial volume: 0.5 – 5 L

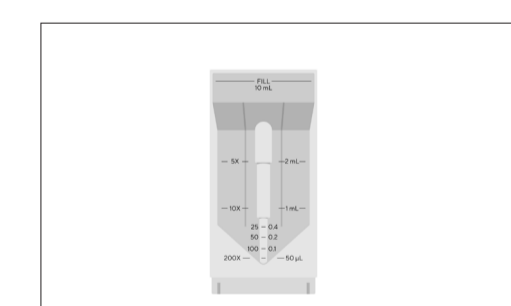
PES	2pc
3 kDa	VF20P9
5 kDa	VF20P1
10 kDa	VF20P0
30 kDa	VF20P2
50 kDa	VF20P3
100 kDa	VF20P4
0.2 µm	VF20P7
HY	
2 kDa	VF20H9
5 kDa	VF20H1
10 kDa	VF20H0
30 kDa	VF20H2
100 kDa	VF20H4



Vivapore® 5

Initial volume: 1 – 5 mL

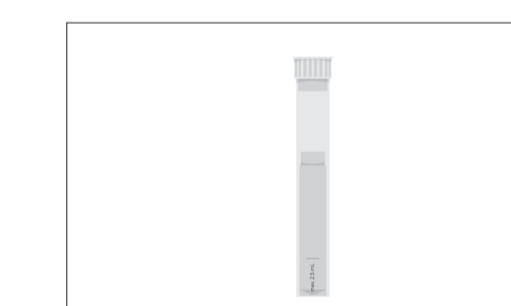
PES	30pc
7.5 kDa	VP-S005P0008-3



Vivapore® 10

Initial volume: 2 – 10 mL¹

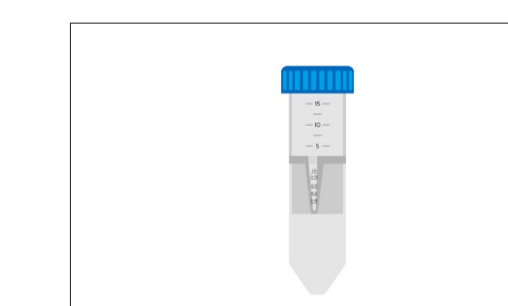
PES	30pc
7.5 kDa	VP-S010P0008-3



Vivaspin® Filtrate

Initial volume: 0.5 – 2.5 mL

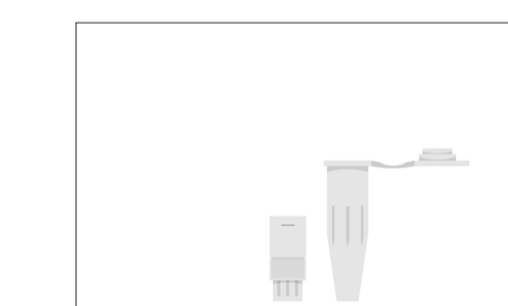
HY	12pc
5 kDa	13229-E
10 kDa	13239-E
20 kDa	13249-E
PES	
100 kDa	13269-GE
300 kDa	13279-E



Vivaspin® Endotest

Initial volume: 2 – 15 mL

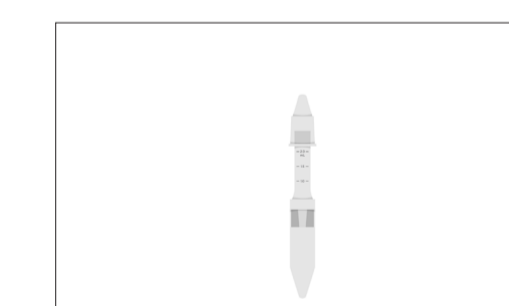
CTA	12pc
20 kDa	VS15RXETO



Vivacon® 500

Initial volume: 0.1 – 0.5 mL

HY	25pc	100pc
2 kDa	VN01H91	VN01H92
10 kDa	VN01H01	VN01H02
30 kDa	VN01H21	VN01H22 ^{2,3}
50 kDa	VN01H31	VN01H32
100 kDa	VN01H41	VN01H42 ²
CTA		
125 kDa	VN01H81	VN01H82 ^{2,3}



Vivacon® 2

Initial volume: 0.4 – 2 mL

HY	25pc	100pc
2 kDa	VN02H91	VN02H92
10 kDa	VN02H01	VN02H02
30 kDa	VN02H21	VN02H22 ²
50 kDa	VN02H31	VN02H32 ²
100 kDa	VN02H41	VN02H42 ^{2,3}
CTA		
125 kDa	VN02H81	VN02H82 ²

PES = Polyethersulfone
HY = Hydrosart® regenerated cellulose
RC = Regenerated cellulose
CTA = Cellulose triacetate

¹ Up to 20 mL sample capacity when used with an expansion reservoir
² PCR grade version available (100pc pack size)
³ PCR grade version available (500pc pack size)