

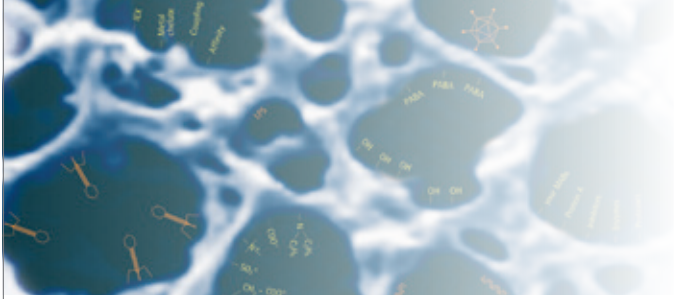


sartorius stedim
biotech

Sartobind[®] selection guide



turning science into solutions



Step 1: Select ligand

Single-use	Re-use	High salt condition	Ligand and binding capacity	For Polishing	For Capture
Anion exchange					
•	•	-	Q, quaternary ammonium 29 mg/mL (BSA)	DNA, host cell proteins, viruses, endotoxins	Large proteins (e.g. FVIII), viruses, VLP
Anion exchange					
•	-	•	STIC PA, primary amine 50 mg/mL (BSA at 17 mS/cm)	DNA, host cell proteins, viruses, endotoxins	-
Cation exchange					
•	•	-	S, sulfonic acid 26 mg/mL (lysozyme)	Host cell proteins, aggregates	Large proteins, viruses, VLP
Hydrophobic interaction					
•	•	•	Phenyl, hydrophobic interaction 15 mg/mL (IgG)	Aggregates, viruses	Large proteins, viruses, VLP

- Applicable
- Not applicable



Step 2: Screen buffer type, composition, pH and conductivity



96-well plate

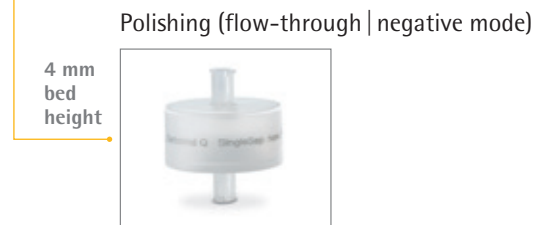


pico 0.08 mL

Step 3: Determine binding capacity



nano 3 mL

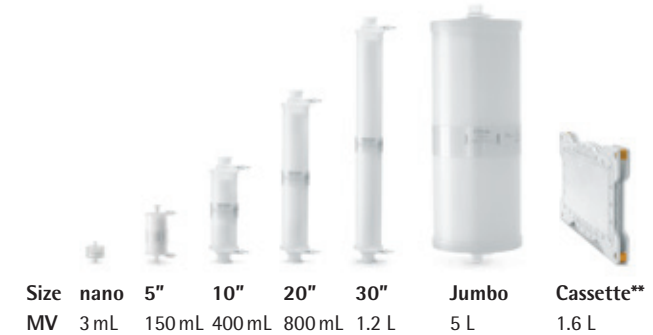


nano 1 mL

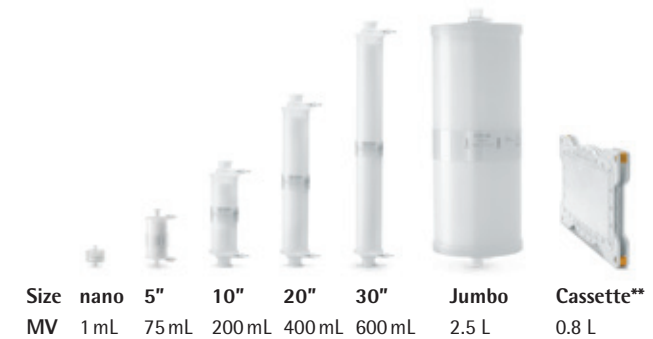


Step 4: Select product for capture or polishing

Bed height 8 mm Flow rate 5 MV*/min Void volume 1.3-1.8 MV



Bed height 4 mm Flow rate 10-30 MV*/min Void volume 2.7-3.5 MV



* Membrane volume

** option γ -irradiated

Each size is available in 4 or 8 mm bed height.
For capture choose 8 mm, for flow-through polishing 4 mm.

Products

Order numbers

Size	MV*	Bed height (mm)	Optimized design	Connector type	Quantity per pack	Q (Sterile validated)	Q (Quaternary ammonium)	S (Sulfonic acid)	STIC PA (Primary amine)	HIC (Phenyl)
96-well plate	0.019 mL	0.8	-	-	2	-	99IEXQ42GC -----V	99IEXS42GC -----V	99STPA42GC -----V	99HICP42GC -----V
96-well plate	0.019 mL	0.8	-	-	10	-	99IEXQ42GC -----D	99IEXS42GC -----D	99STPA42GC -----D	99HICP42GC -----D
Pico	0.08 mL	4	-	Luer female	10	-	92IEXQ42DD-11 --D	92IEXS42DD-11 --D	92STPA42DD-11 --D	92HICP42DD-11 --D
Nano	1 mL	4	•	Luer female	1	-	96IEXQ42DN-11	96IEXS42DN-11	-	-
Nano	1 mL	4	•	Luer female	4	-	96IEXQ42DN-11 --A	96IEXS42DN-11 --A	96STPA42DN-11 --A	-
5"	75 mL	4	•	Hose barb	4	-	96IEXQ42D9MOO --A	96IEXS42D9MOO --A	-	-
5"	75 mL	4	•	Sanitary 3/4"	4	-	96IEXQ42D9MFF --A	96IEXS42D9MFF --A	96STPA42D9MFF --A	-
10"	200 mL	4	•	Sanitary 1 1/2"	1	-	96IEXQ42D1GSS	96IEXS42D1GSS	96STPA42D1GSS	-
20"	400 mL	4	•	Sanitary 1 1/2"	1	-	96IEXQ42D2HSS	96IEXS42D2HSS	96STPA42D2HSS	-
30"	600 mL	4	•	Sanitary 1 1/2"	1	-	96IEXQ42D3KSS	96IEXS42D3KSS	96STPA42D3KSS	-
Jumbo	2.5 L	4	•	Sanitary 1 1/2"	1	-	96IEXQ42D3NSS	-	96STPA42D3NSS	-
Cassette	800 mL	4	•	Sanitary 1 1/2***	1	98IEXQ42DGL	98IEXQ42D-L	98IEXS42D-L	98STPA42D-L	-
Nano	3 mL	8	•	Luer female	4	-	96IEXQ42EUC11 --A	96IEXS42EUC11 --A	-	96HICP42EUC11 --A
5"	150 mL	8	•	Sanitary 3/4"	1	-	96IEXQ42E9BFF	96IEXS42E9BFF	-	96HICP42E9BFF
10"	400 mL	8	•	Sanitary 1 1/2"	1	-	96IEXQ42E1HSS	96IEXS42E1HSS	-	96HICP42E1HSS
20"	800 mL	8	•	Sanitary 1 1/2"	1	-	96IEXQ42E2LSS	96IEXS42E2LSS	-	96HICP42E2LSS
30"	1.2 L	8	•	Sanitary 1 1/2"	1	-	96IEXQ42E3FSS	96IEXS42E3FSS	-	96HICP42E3FSS
Jumbo	5.0 L	8	•	Sanitary 1 1/2"	1	-	96IEXQ42E3ESS	96IEXS42E3ESS	-	96HICP42E3ESS
Cassette	1.6 L	8	•	Sanitary 1 1/2***	1	98IEXQ42EGP	98IEXQ42E-P	98IEXS42E-P	-	98HICP42E-P

- Available
- Not available
- * Membrane Volume
- ** Via manifold



Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen, Germany

Phone +49.551.308.0
Fax +49.551.308.3289

www.sartorius-stedim.com