



sartorius stedim
biotech

Sartobind[®] Product Guide on Improved Versions



turning science **into solutions**

Product change list

Arrow indicates change recommendation from existing to optimized capsule

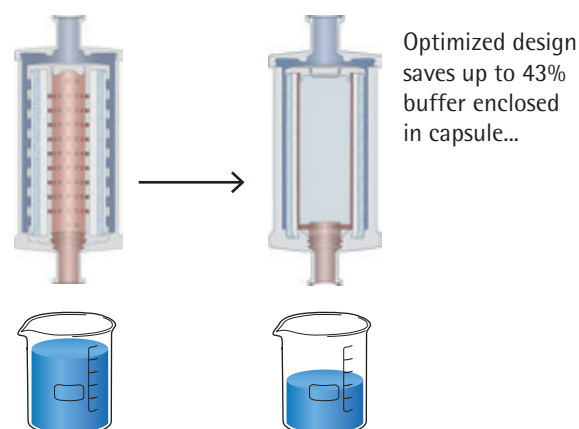
Size	Available	Bed height (mm)	Membrane volume (mL)	Optimized design	Connector type	Quantity per pack	Q (Quaternary amine)
Existing version to be discontinued in Oct. 2015 and July 2017							
New improved version							
96-well plate	Yes	0.8	0.019	-	-	2	99IEXQ42GC-----V
96-well plate	Yes	0.8	0.019	-	-	10	99IEXQ42GC-----D
Pico	Yes	4	0.08	-	Luer female	10	92IEXQ42DD-11--D
Nano 1 mL	until July 2017	4	1	-	Luer female	1	92IEXQ42DN-11
Nano 1 mL	Yes	4	1	Yes	Luer female	1	96IEXQ42DN-11
Nano 1 mL	until July 2017	4	1	-	Luer female	4	92IEXQ42DN-11--A
Nano 1 mL	Yes	4	1	Yes	Luer female	4	96IEXQ42DN-11--A
Mini	Yes	4	7	-	Hose barb	4	92IEXQ42D4-00--A
Mini	Yes	4	7	-	Sanitary 3/4"	4	92IEXQ42D4-SS--A
5"	until Oct. 2015	4	70	-	Hose barb	4	92IEXQ42D9-00--A
75 mL	Yes	4	75	Yes	Hose barb	4	96IEXQ42D9M00-A
5"	until Oct. 2015	4	70	-	Sanitary 1 1/2"	4	92IEXQ42D9-SS--A
75 mL	Yes	4	75	Yes	Sanitary 3/4"	4	96IEXQ42D9MFF--A
10"	until July 2017	4	180	-	Sanitary 1 1/2"	1	92IEXQ42D1-SS
200 mL	Yes	4	200	Yes	Sanitary 1 1/2"	1	96IEXQ42D1GSS
20"	until July 2017	4	360	-	Sanitary 1 1/2"	1	92IEXQ42D2-SS
400 mL	Yes	4	400	Yes	Sanitary 1 1/2"	1	96IEXQ42D2HSS
30"	until July 2017	4	540	-	Sanitary 1 1/2"	1	92IEXQ42D3-SS
600 mL	Yes	4	600	Yes	Sanitary 1 1/2"	1	96IEXQ42D3KSS
Mega 1.6 L	until July 2017	4	1620	-	Sanitary 1 1/2"	1	92IEXQ42DC3SS
Jumbo 2.5 L	Yes	4	2500	Yes	Sanitary 1 1/2"	1	96IEXQ42D3NSS
Nano 3 mL	Yes	8	3	Yes	Luer female	4	96IEXQ42EUC11--A
150 mL	Yes	8	150	Yes	Sanitary 3/4"	1	96IEXQ42E9BFF
400 mL	Yes	8	400	Yes	Sanitary 1 1/2"	1	96IEXQ42E1HSS
800 mL	Yes	8	800	Yes	Sanitary 1 1/2"	1	96IEXQ42E2LSS
1.2 L	Yes	8	1200	Yes	Sanitary 1 1/2"	1	96IEXQ42E3FSS
Jumbo 5 L	Yes	8	5000	Yes	Sanitary 1 1/2"	1	96IEXQ42E3ESS

* 5" STIC PA is available until July 2017

Optimized design of 4 mm (order number starting with 96) **reduces void volume 43% and improves dynamic binding capacity by 17%** compared to existing 4 mm design (order number starting with 92). The existing 8 mm products are already optimized. Compared to a standard filter design they display even up to 48% more dynamic binding capacity.

The membranes, method of membrane manufacturing, plastic materials, and bed height (4 mm) remain unchanged along with the recommended flow rate of 10–30 membrane volumes per minute.

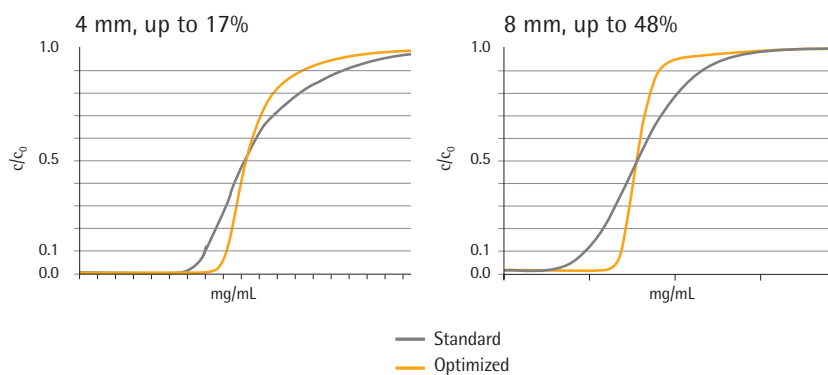
Sartobind® 96-well plates, pico and mini formats are kept unchanged.



S (Sulfonic acid)	STIC PA (Primary amine)	HIC (Phenyl)
99IEXS42GC-----V	99STPA42GC-----V	99HICP42GC-----V
99IEXS42GC-----D	99STPA42GC-----D	99HICP42GC-----D
92IEXS42DD-11--D	92STPA42DD-11--D	92HICP42DD-11--D
92IEXS42DN-11	-	-
96IEXS42DN-11	-	-
92IEXS42DN-11--A	92STPA42DN-11--A	-
96IEXS42DN-11--A	96STPA42DN-11--A	-
92IEXS42D4-00--A	-	-
92IEXS42D4-SS--A	-	-
92IEXS42D9-00--A	-	-
96IEXS42D9M00--A	-	-
92IEXS42D9-SS--A	92STPA42D9-FF--A*	-
96IEXS42D9MFF--A	96STPA42D9MFF--A	-
92IEXS42D1-SS	92STPA42D1-SS	-
96IEXS42D1GSS	96STPA42D1GSS	-
-	92STPA42D2-SS	-
96IEXS42D2HSS	96STPA42D2HSS	-
92IEXS42D3-SS	92STPA42D3-SS	-
96IEXS42D3KSS	96STPA42D3KSS	-
-	92STPA42DC3SS	-
-	96STPA42D3NSS	-
96IEXS42EUC11--A	-	96HICP42EUC11--A
96IEXS42E9BFF	-	96HICP42E9BFF
96IEXS42E1HSS	-	96HICP42E1HSS
96IEXS42E2LSS	-	96HICP42E2LSS
96IEXS42E3FSS	-	96HICP42E3FSS
96IEXS42E3ESS	-	96HICP42E3ESS



...and more dynamic binding capacity at 10% breakthrough



From

Conventional filter design

Q	Q	Q	Q	Bed height 8 mm
S	S	S	S	
HIC	HIC	HIC	HIC	
3 mL	150 mL	1,200 mL	5,000 mL	



1 mL	70 mL	180 mL	360 mL	540 mL	1,620 mL	Bed height 4 mm
Q	Q	Q	Q	Q	Q	
S	S	S	-	S	-	
STIC	STIC	STIC	STIC	STIC	STIC	

+>40% buffer savings | +>15% more binding capacity



Optimized design (48% more binding capacity)

Capsules with 400 and 800 mL membrane volumes added

Merged to one product line with optimized design

To

Q	Q	Q	Q	Q	Q	Bed height 8mm	Capture
S	S	S	S	S	S		
HIC	HIC	HIC	HIC	HIC	HIC		
3 mL	150 mL	400 mL	800 mL	1,200 mL	5,000 mL		



- Adenovirus
- Lentivirus
- Influenza A and B
- Vaccinia
- Virus Like Particles
- Conjugated Vaccines
- Adenoassociated Virus
- Baculovirus
- DNA plasmids
- Von Willebrand Factor vWF
- Factor VIII
- Aggregates
- Process related impurities
- Leached Ligands
- Endotoxins
- DNA
- Virus clearance
- Host cell proteins

1 mL	75 mL	200 mL	400 mL	600 mL	2,500 mL	4mm	Polishing
Q	Q	Q	Q	Q	Q	Bed height	
S	S	S	S	S	-		
STIC PA	STIC PA	STIC PA	STIC PA	STIC PA	STIC PA		

Chromatography as easy as filtration



◀ Sartobind App
www.sartorius-stedim.com/apps