

# Products

# Order Numbers

# SARTORIUS

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	-
mini	10 mL	4	■	Luer female	4	-	96IEXQ42D4R11--A	96IEXS42D4R11--A	96STPA42D4R11--A	-
mini	10 mL	4	■	Hose barb	4	-	96IEXQ42D4ROO--A	96IEXS42D4ROO--A	96STPA42D4ROO--A	-
mini	10 mL	4	■	Sanitary 3/4"	4	-	96IEXQ42D4RFF--A	96IEXS42D4RFF--A	96STPA42D4RFF--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
mini	20 mL	8	■	Luer female	4	-	96IEXQ42E4J11--A	96IEXS42E4J11--A	-	96HICP42E4J11--A
mini	20 mL	8	■	Hose barb	4	-	96IEXQ42E4JOO--A	96IEXS42E4JOO--A	-	96HICP42E4JOO--A
mini	20 mL	8	■	Sanitary 3/4"	4	-	96IEXQ42E4JFF--A	96IEXS42E4JFF--A	-	96HICP42E4JFF--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

# Sartobind® Selection Guide

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



## Germany

Sartorius Stedim Biotech GmbH  
August-Spindler-Strasse 11  
37079 Goettingen  
Phone +49 551 308 0

For further contacts, visit  
[www.sartorius.com](http://www.sartorius.com)

Specifications subject to change without notice.  
© 2021 Sartorius Stedim Biotech GmbH, August-Spindler-Strasse 11, 37079 Goettingen, Germany

Publication No.: SL-1099-e | Order No.: 85032-544-11 | Status: 06 | 17 | 2021



