

Sartobind®

Ion Exchange and Affinity Chromatography
Solutions from Lab to Process



Research

Sartobind® Lab and Vivapure® units
and kits for:

- Scouting Purification
- Preparative Purification
- Proof of Concept

Learn more about membrane
chromatography for Research:

<https://sar.to/9ad9j>

Development

Sartobind® 96 Well Plates
and Nano, for:

- Molecule Development
- Buffer Screening
- Scalable Process Development

Learn more about membrane
chromatography for Development
and Production:

<https://sar.to/g4rwi>

Production

Sartobind® Capsules
and Cassettes for:

- Pilot Scale Production
- Process Scale Production

Ordering Information

Sartobind® Rapid A

	Size	MV	Inlet Connector	Qty	Order Number
	Lab	0.5 mL	Female Luer	1	SBLRA025EL-1
				4	SBLRA025EL-A
	96-Well Plate	0.019 mL	-	2	99R-PA19GC-----V
	Nano	1.2 mL	Female Luer	1	96R-PA19C-A11ZX
				2	96R-PA19D-T11--V
				4	96R-PA19D-T11--A
	Mini	10 mL	Female Luer	1	96R-PA19D4R11
			Sanitary 3/4", 25 mm	1	96R-PA19D4RFF
	5"	75 mL	Sanitary 3/4", 25 mm	1	96R-PA19D9MFF
	10"	200 mL	Sanitary 1 1/2", 50.5 mm	1	96R-PA19D1GSS
	Cassette	800 mL	Sanitary 1 1/2", 50.5 mm ¹	1	98R-PA19D-L

Sartobind® IEX

	Size	MV	Inlet Connector	Qty	Order Number Q	S
	Lab	0.41 mL	Female Luer	4	93IEXQ42GB-12--A	93IEXS42GB-12--A
		2.1 mL		2	93IEXQ42DB-12--V	93IEXS42DB-12--V
		2.8 mL		1	93IEXQ42BC-12	93IEXS42BC-12
	96-Well Plate	0.019 mL	-	2	99IEXQ42GC-----V	99IEXS42GC-----V
				10	99IEXQ42GC-----D	99IEXS42GC-----D
	Nano	3 mL	Female Luer	4	96IEXQ42EUC11--A ²	96IEXS42EUC11--A ²
	Mini	20 mL	Female Luer	4	96IEXQ42E4J11--A ²	96IEXS42E4J11--A ²
			Hose Barb	4	96IEXQ42E4JOO--A ²	96IEXS42E4JOO--A ²
			Sanitary 3/4", 25 mm	4	96IEXQ42E4JFF--A ²	96IEXS42E4JFF--A ²
	5"	150 mL	Sanitary 3/4", 25 mm	1	96IEXQ42E9BFF ²	96IEXS42E9BFF ²
	10"	400 mL	Sanitary 1 1/2", 50.5 mm	1	96IEXQ42E1HSS ²	96IEXS42E1HSS ²
	20"	800 mL	Sanitary 1 1/2", 50.5 mm	1	96IEXQ42E2LSS ²	96IEXS42E2LSS ²
	30"	1,200 mL	Sanitary 1 1/2", 50.5 mm	1	96IEXQ42E3FSS ²	96IEXS42E3FSS ²
	Jumbo	5,000 mL	Sanitary 1 1/2", 50.5 mm	1	96IEXQ42E3ESS ²	96IEXS42E3ESS ²
	Cassette	1,600 mL	Sanitary 1 1/2", 50.5 mm ¹	1	98IEXQ42E-P ²	98IEXS42E-P ²

¹ Via manifold.

² 8 mm bed height. For 4 mm bed height, replace 42E with 42D when placing your order.

Your Contact for Research:

Your Contact for Development and Production: