

Sartobind® Rapid A

Simplify mAb Capture
With Sartobind® Rapid A



Product Information

Sartobind® Rapid A is a membrane chromatography platform designed for modern monoclonal antibody purification processes. Independent of the chromatography system, it offers a scalable solution for capturing Fc-region containing biomolecules, ideal for process intensification in single- or multi-column setups. Its agarose-based membrane combines convective and diffusive mass transport for higher flow rates at reduced back pressure, shortening process times without sacrificing binding capacity. Benefits include reducing set-up cost by up to 80%, ensuring up to 200 cycles for a full membrane lifespan, and high binding capacities of 40 g/L.

Benefits

- Ready-for-use consumables, quick and easy setup
- Scalable, cost-efficient portfolio
- Enables process intensification, shorter timelines
- Membrane designed for shorter residence time and diverse mAb sizes
- Designed for full utilization of the membrane's lifespan
- Expert, tailored consultancy services

Technical Specifications







Attribute	Sartobind® Rapid A 96-Well Plate	Sartobind® Rapid A Nano	Sartobind® Rapid A Mini	Sartobind® Rapid A 75 mL	Sartobind® Rapid A 200 mL	Sartobind® Rapid A Cassette
Membrane volume	19 µL/well	1.2 mL	10 mL	75 mL	200 mL	800 mL/cassette
Bed height [mm]	0.8	4	4	4	4	4
Dimensions [mm]	128×85.5×25	37×31	70×54.5	190×77	350×100	634×387×49
Connector	N/A	Female luer	Female luer; sanitary 3/4", 25 mm	Sanitary 3/4", 25 mm	Sanitary 1 1/2", 50.5 mm	Via manifold accessory: sanitary 1 1/2", 50.5 mm
Operating pressure limit	350 mbar	4 bar	4 bar	4 bar	4 bar	2 bar
Shelf life	2 Y	2 Y	2 Y	2 Y	2 Y	1 Y
DBC 10% [mg/mL]	38	38	38	38	38	38
Approximate weight	N/A	10 g	65 g	400 g	760 g	4.9 kg (6.0 kg wet)



Ordering Information

Product Name	Quantity	Description Adapter Inlet and Outlet	Order Number
Sartobind® Rapid A, 96-well plate	2	-	99R-PA19GC----V
Sartobind® Rapid A, Nano 1.2 mL, caps. 1 pc.	1	Female luer	96R-PA19C-A11ZX
Sartobind® Rapid A, Nano 1.2 mL, caps. 2 pcs.	2	Female luer	96R-PA19D-T11--V
Sartobind® Rapid A, Nano 1.2 mL, caps. 4 pcs.	4	Female luer	96R-PA19D-T11--A
Sartobind® Rapid A, Mini 10 mL	1	Female luer	96R-PA19D4R11
Sartobind® Rapid A, Mini 10 mL	1	Sanitary 3/4", 25 mm	96R-PA19D4RFF
Sartobind® Rapid A, 75 mL	1	Sanitary 3/4", 25 mm	96R-PA19D9MFF
Sartobind® Rapid A, 200 mL	1	Sanitary 1 1/2", 50.5 mm	96R-PA19D1GSS
Sartobind® Rapid A, 0.8 L cassette	1	Via manifold accessory: Sanitary 1 1/2", 50.5 mm	98R-PA19D-L

Peripherals and Accessories

Product Name	Order Number
Manifold set, 1 1/2" sanitary clamp, inlet and outlet adapter plate 	29Z-S00001
Pilot filter holder 	2ZGL--0005
Process filter holder 	2ZGL--0006
Double process filter holder 	2ZGL--0007
Drip pan for pilot filter holder	2ZGL--0008
Drip pan for process and double process filter holder	2ZGL--0015
Separation plate for serial or parallel fluid distribution	2ZGL--0013
Pressure safety device, to prevent excessive pressure during compressed air draining of the cassettes	2ZGL--0014
Adapter luer male to UNF 10 - 32 female, PEEK, 1 pc.	1ZA---0004
Adapter UNF 10 - 32 female to sanitary 3/4", 25 mm, polyoxymethylene, 2 pcs.	1ZAOGV0003

Germany

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

USA

Sartorius Stedim North America Inc.
565 Johnson Avenue
Bohemia, NY 11716
Toll-Free +1 800 368 7178

 **For more information, visit**
www.sartorius.com