

# Sartobind<sup>®</sup> Rapid A

Plug. Play. Change the Game.  
Simplify mAb Capture  
With Sartobind<sup>®</sup> Rapid A



## Product Information

Resin-based packed bed chromatography for mAb capture hampers downstream processing with low productivity, laborious cleaning requirements, and high costs. Now, the introduction of Sartobind<sup>®</sup> Rapid A makes disposable “plug and play” a reality. A unique one-batch, one-membrane approach reduces bioburden risk and lowers CoGs compared to resin-based processes. Plus, rapid cycling translates to >10× higher productivity and linear scalability from PD to commercial scale.

## Benefits

- Eliminate column packing, cleaning validation and storage
- Reduce high cost of bulk resin and underutilization
- Accelerate speed to clinic with >10× productivity vs. resin-based batch operations
- Adopt a fully proven scalable solution from PD to commercial scale
- Cut risk of bioburden linked to columns

## Relevant Applications

- mAb capture
- Purification

## Relevant Process Steps

- Chromatography | capture and polish

# Technical Specifications







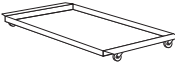
Attribute	Sartobind® Rapid A 96- Well Plate	Sartobind® Rapid A Pico	Sartobind® Rapid A Nano	Sartobind® Rapid A Mini	Sartobind® Rapid A 75 mL	Sartobind® Rapid A 200 mL	Sartobind® Rapid A Cassette
Membrane Volume	19 µL/well	0.08 mL	1.2 mL	10 mL	75 mL	200 mL	800 mL
Bed Height	0.8 mm	4 mm	4 mm	4 mm	4 mm	4 mm	4 mm
Dimensions [mm] h × d			37 × 31	70 × 54.5	190 × 77	350 × 100	634 × 387 × 49 w × h × d
Connector		Female luer	Female luer	Female luer; sanitary ¾", 25 mm	Sanitary ¾", 25 mm	Sanitary 1½", 50.5 mm	Via manifold accessory: sanitary 1½", 50.5 mm
Approximate Weight			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, Pico 0.08 mL	10	-	92R-PA19DD-11--D
Sartobind® Rapid A, Nano 1.2 mL	2	Female luer	96R-PA19D-T11--V
Sartobind® Rapid A, Nano 1.2 mL	4	Female luer	96R-PA19D-T11--A
Sartobind® Rapid A, Mini 10 mL	1	Female luer	96R-PA19D4R11
Sartobind® Rapid A, Mini 10mL	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
Sartocheck® 5 Plus Filter Integrity Tester	26787---FT---P
Adapter Luer male to UNF 10 - 32 female, PEEK, 1 piece	1ZA---0004
Adapter UNF 10 - 32 female to sanitary 3/4", 25 mm, polyoxymethylene, 2 pieces	1ZA0GV0003

**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](http://www.sartorius.com)