

Sartobind[®] Lab Protein A

Laboratory Scale Affinity
Membrane Adsorber



Benefits

- Purification as simple as filtration
- 10x higher flow rates than conventional columns
- High dynamic binding capacity (up to 90 mg/hr, 6 cycles)
- Ultimate flexibility for handling by syringe, pump or FPLC system
- Compatible with conventional buffer systems (e.g. PBS for equilibration | loading | washing, and 0.1 M glycine/HCl pH 2.8 – 3.5 for elution)

Product Information

Sartobind[®] Lab Protein A affinity membrane adsorbers are ideal for the rapid purification of antibodies, including IgGs and mAbs. The large pore size allows 10-fold faster flow rates than conventional resin-based columns, resulting in cycle times of only 10 minutes. This ensures a higher throughput than with any other Protein A matrix. Sartobind[®] Lab Protein A devices can be operated by syringe or, with the included UNF 10-32 adapters, on any peristaltic pump or liquid chromatography system.

Technical Specifications

Construction

Height × Diameter	27 × 36 mm
Housing	Polypropylene (PP)
Membrane	Stabilized Regenerated Cellulose
Ligand	Recombinant Protein A
Inlet Connector	Female Luer Lock
Outlet Connector	Male Luer Lock

Membrane Specifications

Frontal Surface Area	5 cm ²
Nominal Adsorption Area	100 cm ²
Bed Height	4 mm
Membrane Volume	2 mL
Ligand Density	1.5 mg/mL
Binding Capacity*	5 – 7.5 mg/mL 10 – 15 mg/unit

Operating Conditions

Recommended Flow Rate	5 – 10 mL/min
Maximum Pressure	0.6 MPa 6 bar 87 psi
pH Stability	3 – 9 (long term) 2 – 10 (short term)
Chemical Stability**	Stable in all common chromatography buffers, except oxidizing or reactive reagents

* Binding capacity for 1 mg/mL polyclonal human IgG sample at a flow rate of 10 mL/min with 10% breakthrough.

** A pH <3 is sometimes required to elute all strongly bound IgG. Under these conditions, proteins may hydrolyze.

Applications

Ultra rapid purification of antibodies (e.g. IgG, mAbs).

Ordering Information

Description	4 pc
Sartobind® Lab Protein A	93PRAP06HB-12--A


Each package contains 1x Luer male and 1x Luer female adapters to UNF 10–32, PEEK, for operation via an FPLC system.

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Straße 20
37079 Göttingen
Phone +49 551 308 0

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906

 For further information, visit
www.sartorius.com