SVISCISVS

Product Datasheet

Sartobind® Lab IEX 15, 75 and 100

Laboratory Scale Ion Exchange Membrane Adsorbers



Product Information

The Sartobind[®] Lab units with ion exchange (IEX) membranes are ready-to-use for laboratory experiments. Due to their large pore size of >3 µm the membranes are ideal for capture of large proteins and viruses, without observing size exclusion effects. The Lab IEX units can be operated manually by a syringe or on a FPLC system.

These units combine the advantages of Sartobind[®] membrane adsorbers with a syringe filter design, allowing for rapid and easy sample purification.

Benefits

- Ready-to-use units
- Reusable*
- High flow rate: up to 30 bed volumes per minute
- High dynamic binding capacity for large proteins and viruses
- Fast and convenient operation by syringe
- Compatible with peristaltic pumps or liquid chromatography systems
- Connect in series for increased binding capacity

For process development studies intended for later scale-up, we recommend to use Sartobind® Nano 1 or 3 mL capsules**.

Applications

Capture

Proteins

Viruses

- Oligonucleotides
- Plasmids

Contaminant removal

- DNA
- Viruses
- Host cell proteins
- Leached protein A
- Endotoxins

Technical Data

Specifications

Materials of construction		Membrane ligands		
Housing	Polypropylene		Strong basic anion exchanger	Quaternary ammonium (Q) R-CH₂-N⁺(CH₃)₃
Membrane matrix	Stabilized reinforced cellulose, nominal pore size >3 μm		Strong acidic cation exchanger	Sulfonic acid (S) R-CH ₂ -SO ₃ ⁻
			Weak basic anion exchanger	Diethylamine (D) R-CH2-N(C2H5)2
		SARTOBIND® Q15 Let no 80	StarroBIND* Q75 Latro B-	SARTOBIND® Q100 Lot no. 81543425 ↓
		Lab IEX 15	Lab IEX 75	Lab IEX 100
Membrane Area [cm²]		15	75	100
Number of layers		3	15	5
Bed height [mm]		0.8	4.0	1.4
Membrane volume [mL]		0.41	2.1	2.8
Membrane diameter [mm]		25	25	50
Typical dynamic binding capacity 10% for Q [mg/unit]		12	60	80
Typical dynamic binding capacity 10% for S [mg/unit]		10.5	52.5	70
Typical dynamic binding capacity 10% for D [mg/unit]		-	60	-
Recommended flow rate [mL/min]		2 - 12	10 - 63	14 - 84
Recommended flow rate [membrane volume/min]		5 - 30	5 - 30	5 - 30
Ligand density [µeq/cm²]		2 - 5 (Q, S), 2 -	4 (D) 2 - 5 (Q, S), 2 - 4 (D)	2 - 5 (Q, S), 2 - 4 (D)
Maximum pressure at 20°C		6 bar 0.6 MPa	87 psi 6 bar 0.6 MPa 87 p	osi 6 bar 0.6 MPa 87 psi
pH stability		Q	S	D
Short term		2-14	3 - 14	2 - 14
Long term		2-12	4 - 13	4 - 13

^{**} Sartorius Data sheet: Sartobind® Q and S, 4 and 8 mm bed height, Order No.: 85034-535-76

Dimensions and Connections

	Lab IEX 15	Lab IEX 75	Lab IEX 100
Dimensions (height × diameter) [mm]	25 × 36	28 × 36	31 × 66
Frontal surface area [cm²]	5	5	20
Connector inlet	Luer female	Luer female	Luer female
Connector outlet	Luer male	Luer male	Luer male
Approximate weight [g]	9	10	35

Ordering Information

Sartobind[®] Lab IEX units (reusable)

order number Description		Quantity	
93IEXQ42GB-12A	Sartobind [®] Lab Q15, Luer female and male connectors	4	
93IEXS42GB-12A	Sartobind [®] Lab S15 Luer female and male connectors	4	
93IEXQ42DB-12V	Sartobind [®] Lab Q75, Luer female and male connectors	2	
93IEXS42DB-12V	Sartobind [®] Lab S75, Luer female and male connectors	2	
93IEXD42DB-12V	Sartobind [®] Lab D75, Luer female and male connectors	2	
93IEXQ42BC-12	Sartobind [®] Lab Q100, Luer female and male connectors	1	
93IEXS42BC-12	Sartobind [®] Lab S100, Luer female and male connectors	1	
93IEXS42BC-12	Sartobind $^{\odot}$ Lab S100, Luer female and male connectors		

Each package of Lab IEX 15, 75 and 100 contains 1 × Luer male and 1 × Luer female adapters to UNF 10-32, PEEK, for operation via an FPLC system.

Accessories

Order number	Description	Quantity
1ZA0004	Luer male adapter to UNF 10-32 female, PEEK	1
1ZA0005	Luer female adapter to UNF 10-32 female, PEEK	1
17002140	$2 \times M6$ female to female Luer and $2 \times M6$ female to male Luer adapters, black Tefzel	1

Related Products

96-Well Plates and Accessories

Order number	Description	Quantity 2 (24 × 8-strips)	
99IEXQ42GCV	Sartobind® Q 96-well plate, 2 units		
99IEXQ42GCD	Sartobind® Q 96-well plate, 10 units	10 (120 × 8-strips)	
99STPA42GCV	Sartobind STIC® PA 96-well plate, 2 units	2 (24 × 8-strips)	
99STPA42GCD	Sartobind STIC® PA 96-well plate, 10 units	10 (120 × 8-strips)	
99IEXS42GCV	Sartobind® S 96-well plate, 2 units	2 (24 × 8-strips)	
99IEXS42GCD	Sartobind® S 96-well plate, 10 units	10 (120 × 8-strips)	
99HICP42GCV	Sartobind® Phenyl 96-well plate, 2 units	2 (24 × 8-strips)	
99HICP42GCD	Sartobind® Phenyl 96-well plate, 10 units	10 (120 × 8-strips)	
VW96VAC01	Vac96 vacuum manifold	1	
VW96VAA02	Vac96 liquid trap and reservoir	1	
VW96VAA04	96 deep well collection plate 2 mL (square wells)	25	
VW96VAA05	Replacement seal for Vac96 vacuum manifold	25	
16612	Vacuum pump, 98%, 220 V, 50 Hz	1	
16615	Vacuum pump, 98%, 110 V, 60 Hz	1	

Germany

USA

August-Spindler-Strasse 11 37079 Goettingen Phone +49 551 308 0

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

Learn more on

www.sartorius.com/chromatography-membranes