

Instructions for Use

Sartobind[®] Q and S

Void Volume Optimized Capsules and Cassettes
With 4 | 8 mm Bed Height



1000032237



SARTORIUS

Table of Contents

1	About these Instructions	6
1.1	Validity	6
1.2	Related Documents	8
1.3	Target Groups	8
1.4	Symbols Used	9
2	Safety	10
2.1	Intended Use	10
2.2	Modifications to the Product	12
2.3	Qualifications of Personnel	13
2.4	Personal Protective Equipment	13
2.5	Leaking Liquids from the Product	13
2.6	Components Under Pressure	14
2.7	Product weight	14
3	Product Overview	15
3.1	Operating Principle	15
4	Installation	21

5	Operation	27
5.1	Venting the Capsule	27
5.2	Venting the Nano Capsule	28
5.3	Cleaning and Equilibration	30
5.4	Autoclaving	31
5.5	Gamma Irradiation	32
5.6	Recommended Flow Rates and Equilibration Buffer Volumes	33
5.7	Buffer Conditions	36
5.8	Selection of pH Conditions	36
5.9	Contaminant Removal From Therapeutic Proteins And Other Sources in Flow-Through Mode	37
5.10	Sample Preparation	38
5.11	Washing	39
5.12	Elution	39
5.13	Draining	39
5.14	Regeneration and Storage	40
5.15	Operation of the Sartobind® Nano With Peristaltic Pumps or Liquid Chromatography (LC) Systems	40
5.16	Scaling Up	41

6	Testing of Membrane Samples Enclosed With 5 L Capsules	45
6.1	Testing Sartobind® Q 30 mm Discs	45
6.2	Testing Sartobind® S 30 mm Discs	50
7	Integrity Test by Diffusion	55
7.1	Installation	55
7.2	Operation Procedure	56
8	Troubleshooting	61
9	Technical Data	66
9.1	Bed Height 4 mm	66
9.2	Bed Height 8 mm	68
9.3	Materials	70
9.4	Binding Capacity	71
9.5	Chemical Stability	72
9.6	Storage Conditions	72
10	Quality Assurance	73

11	Ordering Information	74
11.1	Products 4 mm Bed Height.....	74
11.2	Products 4 mm Bed Height, Gamma Irradiatable.....	78
11.3	Products 4 mm Bed Height, Gamma Irradiated	80
11.4	Products 8 mm Bed Height.....	80
11.5	Products 8 mm Bed Height, Gamma Irradiatable.....	83
11.6	Accessories	86
12	Dimensions and Connections	88

1 About these Instructions

1.1 Validity

These instructions are part of the product; they must be read in full and retained. These instructions apply to the following versions of the product:



4 mm: Nano 1 mL

Mini 10 mL

75 mL

200 mL

400 mL

8 mm: Nano 3 mL

Mini 20 mL

150 mL

400 mL

800 mL



600 mL
1.2 L



Jumbo 2.5 L
Jumbo 5 L



Cassette 0.8 L
Cassette 1.6 L

1.2 Related Documents

In addition to these instructions, please observe the following documents:

- Operating instructions of the device in which the product is used
- Pilot Filter Holder Instructions for Use
- Process | Double Process Filter Holder Instructions for Use
- Validation Guide of the respective product

1.3 Target Groups

These instructions are addressed to the following target groups. The target groups must possess the knowledge specified below.

Target Group	Knowledge and Qualifications
Operator	The operator is familiar with the product and the associated work processes. The operator understands the hazards which may arise when working with the product, and knows how to prevent them.

1.4 Symbols Used

1.4.1 Warnings in Operation Descriptions

CAUTION

Denotes a hazard that may result in moderate or minor injury if it is **not** avoided.


NOTICE

Denotes a hazard that may result in property damage if it is **not** avoided.

1.4.2 Other Symbols

- ▶ Required action: Describes activities that must be carried out. The actions in the sequence must be carried out in succession.
- ▷ Result: Describes the result of the activities carried out.

2 Safety

 Use of the products in applications not specified or not described in this manual, may result in improper function, personal injury, or damage of the product or material. The products are supplied as non-sterile unless otherwise explicitly described. The membrane is dried from glycerol.

2.1 Intended Use

The product is intended exclusively for use in accordance with these instructions. Any other use is considered improper.

The membrane chromatography products also described as membrane adsorbers are intended and validated for single use to avoid carryover as well as tedious and costly cleaning validation procedure. However, it is technically possible to reuse after cleaning in place depending on application, character of sample and process. Additional cleaning and validation steps will be needed to assure constant binding capacity and flow rate after each cycle.

The **4 mm** product line has been developed for the highest possible flow rate and is typically applied for flow-through polishing applications in which binding capacity is typically no limitation.

The **8 mm** product line is used when bind & elute applications or the highest possible binding capacity is needed. If you are not sure what to use, the 8 mm should be used as this bed height provides the best ratio of void volume to membrane volume and the maximum achievable dynamic binding capacity.

For **scale-up** we recommend to stay within one bed height size as this is the least complex procedure. Nevertheless you may change to a different bed height later by keeping residence time constant (see chapter “5.16 Scaling Up”, page 41).

Sartobind® Nano capsules (1 and 3 mL) have been developed for working with small sample volumes. They are perfect for small scale applications, and also for screening purposes and laboratory-scale bind & elute and flow-through purifications.

Sartobind® Mini capsules (10 and 20 mL) have been developed for first scale up trials and preclinical production. These product sizes close the gap between the Nano and the 75 | 150 mL size.

Sartobind® capsules (75 and 150 mL) have been developed for intermediate and pilot scale.

Sartobind® capsules (200 mL up to Jumbo 5 L) have been developed for production purposes in the biopharmaceutical industry.

Sartobind® cassettes (0.8 or 1.6 L) are used in the Pilot Scale Filter Holder of up to 20.8 L membrane volume for the biopharmaceutical production.

2.2 Modifications to the Product

If the product is modified: Personnel may be put at risk. Product-specific documents and product approvals may lose their validity. If you have any queries regarding modifications to the product, contact Sartorius.

2.3 Qualifications of Personnel

Personnel who do **not** possess adequate knowledge about how to use the product safely may injure themselves and other personnel.

2.4 Personal Protective Equipment

Personal protective equipment protects against risks arising from the product. If the personal protective equipment is missing or is unsuitable for the work processes on the product: Personnel may be injured. The following personal protective equipment must be worn:

- Protective work clothing
- Safety gloves
- Safety glasses

2.5 Leaking Liquids from the Product

If the product is damaged or incorrectly installed:
Liquids can leak from the product.

- ▶ Do **not** exceed the maximum pressure (see chapter “9 Technical Data”, page 66).
- ▶ Perform a visual inspection before use.
- ▶ Ensure correct installation.

2.6 Components Under Pressure

Operating above the maximum operating pressure can lead to the bursting of pipes and plastic components. This can result in leakage of media and injuries due to bursting components.

Escaping medium can lead to infections.

- ▶ Wear appropriate personal protective equipment.
- ▶ Do **not** exceed the maximum operating pressure (see chapter “9 Technical Data”, page 66).

2.7 Product weight

The product may be heavy. There is a risk of injury when lifting and transporting the product, e.g. from the product being dropped.

- ▶ Wear appropriate personal protective equipment.
- ▶ If required get help from others when lifting and transporting the product.

3 Product Overview

3.1 Operating Principle

The capsules and cassettes with 4 and 8 mm bed height are ion exchange chromatography products based on macroporous membranes. They can be used for chromatographic separation in downstream processing of viruses and proteins. The ion exchange ligands are coupled to a membrane which is fitted into a plastic housing ready to use. The products are constructed with optimized fluid channels.

The capsules contain a central core and the cassettes a spacer element to minimize void volume. To set up and operate the Sartobind® Jumbo we recommend the trolley (see chapter “11.6 Accessories”, page 86). Capsules and cassettes with strong basic and strong acidic ion exchange Membrane Adsorber (MA) are available. These products are intended for single use to avoid carryover as well as tedious and costly cleaning validation procedures. However, it is technically possible to reuse after cleaning in place (see also section “9.5 Chemical Stability”, page 72).

Their major application is for single use as they are validated for contaminant removal from proteins in flow-through mode (negative chromatography) to bind DNA, residual protein, host cell proteins (HCP), endotoxins, viruses and aggregates. The products can also be used to capture large proteins, viruses such as adeno-, lentiviruses and virus like particles (VLP).

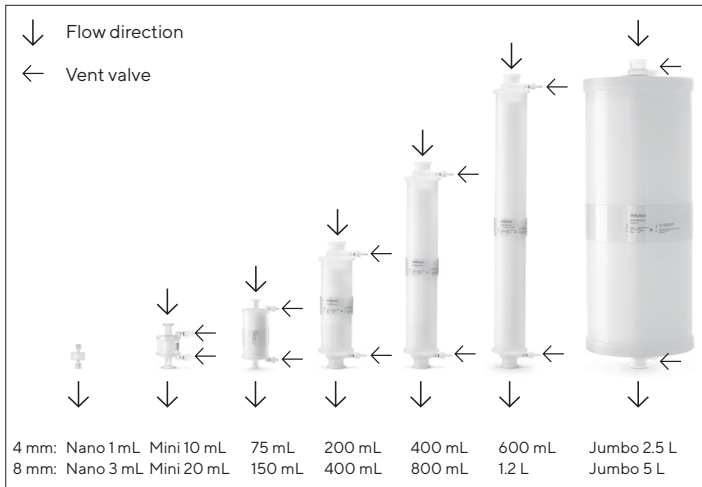


Fig. 1: Flow direction and position of vent valves of 4 and 8 mm capsules



Fig. 2: Flow direction and position of vent valve connection of 4 and 8 mm cassettes

For Mini and 75 | 150 mL products the central core is made from a solid polypropylene cylinder.

For the larger capsules it is made from a self-contained air filled polypropylene cylinder.

The interior of the core is inaccessible for gases and fluids. The 2 flat membrane stacks of the cassettes are separated by a central spacer element.

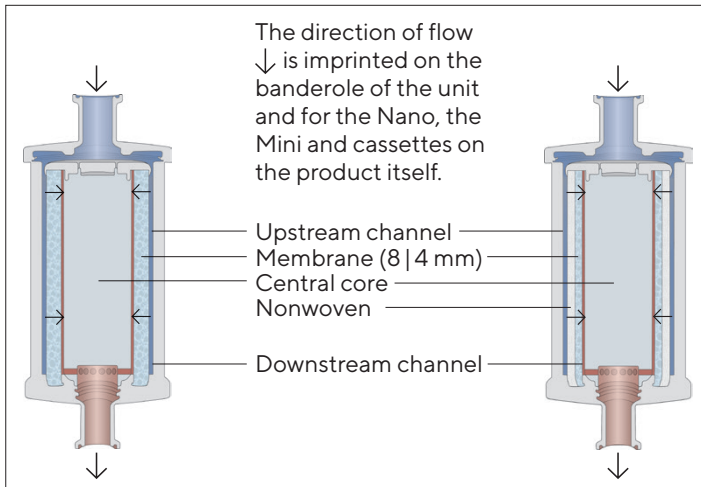


Fig. 3: Construction and flow path inside the 8 mm (left) and 4 mm (right) capsules.

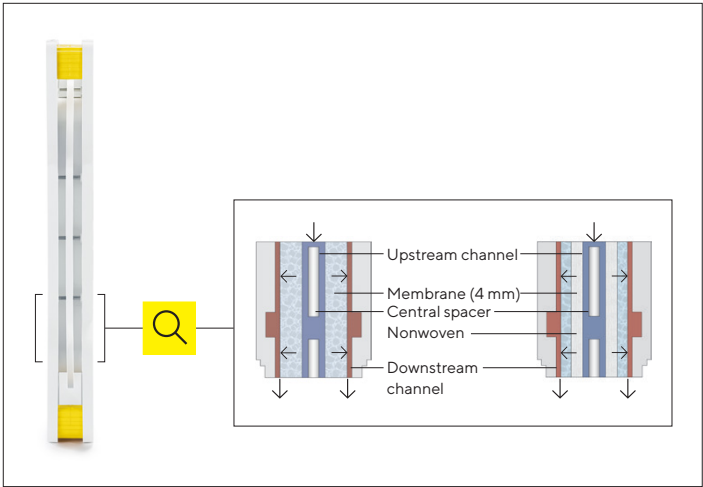


Fig. 4: Side view cassette; highlighted section see Fig. 5

Fig. 5: Construction and flow path inside the 8 mm (left) and 4 mm (right) cassettes

4 Installation

NOTICE

Malfunctions due to damaged products. A damaged product can cause malfunctions.

- ▶ Visually inspect the product before use.
 - ▶ Close vent valve before use by screwing the valve clockwise.
 - ▶ For the cassettes close the clamps at the manifold set.
 - ▶ Do **not** keep or place the products with the connector directly on the floor.
 - ▶ In case of damage: Replace the product.
-

Sartobind® Q and S Jumbo

- ▶ Take the capsule including the styrene foam end protectors, out of the box and place it upright on the end protectors.
- ▶ Move the Jumbo trolley (accessory) in place.
- ▶ Remove upper foam protection and transparent bag.
- ▶ Lift the Jumbo directly onto the trolley (inlet is up and the arrow imprinted on the banderole is pointing down).
- ▷ It is recommended to connect the Jumbo with the trolley with the 3 screws delivered with the trolley.

- ▷ The protective caps on inlet and outlet should stay until you use the unit.
- ▶ Store the caps when you plan to autoclave (see chapter “5.4 Autoclaving”, page 31).
- ▷ The Jumbo 2.5 and 5 L carry protective caps on vent valves as well.
- ▶ Remove before venting.

Sartobind® Q and S Capsules | Cassettes

- ▶ Protect inlet and outlet connectors when unpacking.
- ▶ Install the product in an upright position according to the process flow.
- ▷ The inlet is up.
- ▷ The flow is guided to the upstream channel (i.e. the solution enters the product) passing through the membrane layers to the downstream channel and to the outlet of the product (see Fig. 3).
- ▶ Install the capsule and cassette(s) in-line with a prefilter (0.2 µm or 0.45 µm) in front of the product to prevent blockage or pressure build-up.

- ▶ For Sartobind® cassettes use an appropriate cassette holder and one Manifold Set (see chapter “11.6 Accessories”, page 86).
- ▶ If you plan to use a different filter holder from other manufacturers, please contact Sartorius for technical advice.

Installing Sartobind® Phenyl Cassettes

- ▶ Unpack the Manifold Set containing one inlet and one outlet plate.
- ▶ Place the “INLET” marked plate at one end of the holder.
- ▷ “THIS SIDE UP” mark on the manifold should be readable on the top.
- ▶ Place the manifold marked with “OUTLET” at the other end of the holder, so that “THIS SIDE UP” is readable from the top.
- ▷ The fluid channels of both plates are oriented to the same direction.

- ▶ Place the cassettes (4, see Fig. 6) in the lowest possible position in the holder.
- ▶ The cassettes (4) used for chromatographic separation must originate from the same lot.
- ▶ Put the desired number of Sartobind® cassettes (4) between the manifolds (see Fig. 6).
- ▷ “THIS SIDE UP” is readable on top.
- ▶ Adjust the clamping force for cassettes in Pilot and Process holders to a minimum of 25 kN (optimal range: 25 – 30 kN).
- ▷ In the Pilot holder up to 13 cassettes and a manifold set can be installed.
- ▶ Close all DRAIN and VENT valves (1) of the manifold plates manually using the pinch clamp.

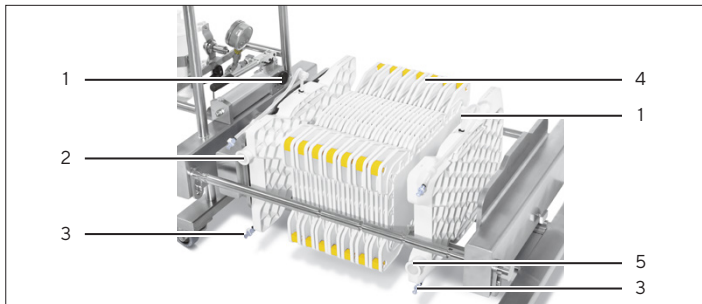


Fig. 6: Insert cassette(s) between the manifold inlet and outlet plates on the Pilot Filter Holder.

Pos.	Name
1	Vent valve
2	Inlet manifold
3	Drain valve
4	Cassettes
5	Outlet manifold

- ▶ Connect the inlet and outlet plates with 1½ inch tri-clamp to the process solution.
- ▷ Maximum pressure for the set-up of 1 to 13 cassette(s) is 2 bar (0.2 MPa, 29 psig).
- ▷ During flushing the clamping force may reduce.
- ▶ In order to avoid dripping during operation, it is recommended to re-adjust the clamping force to a minimum of 25 kN before you continue with equilibration.
- ▶ Make sure that the pump peak pressure caused by pulsation stays below this limit.

5 Operation

5.1 Venting the Capsule

All capsules except Nano have vent valves (see Fig. 1).

Procedure

- ▶ Remove the entire air from the unit before use.
- ▷ The vent valves are equipped with hose barb connectors for the fluid spilled out during venting.
- ▶ Check vent valve position.
- ▷ When turning anticlockwise, the valve is open, when turning clockwise, the valve is closed.
- ▶ Before opening the vent valve, connect the valves with flexible tubing (inner diameter 6 mm) to waste.
- ▶ **NOTICE** Insufficient closing of the valve due to excessive pressure! During venting of capsules do **not** exceed 0.05 MPa (0.5 bar | 7.3 psig).
- ▶ Open the vent valve screw $\frac{1}{8}$ turn to left until all air is replaced by fluid.
- ▶ For venting the cassettes, tubes with quick connectors are attached to the inlet and outlet manifolds and closed with a pinch clamp.

5.2 Venting the Nano Capsule

- ▶ Fill a 10 – 20 mL Luer syringe with equilibration buffer and connect to the capsule. Hold capsule upright (outlet is up) and expel air as shown in Fig. 7.
- ▶ If you still detect any air in the filled unit, close the outlet, hold the syringe up and move the plunger slightly up and down that air bubbles can ascend into the syringe.
- ▶ Alternatively, connect a second empty syringe to the top of the Nano and expel air and buffer into that syringe, disconnect the upper syringe to push out air and reconnect to the Nano, turn it and purge the solvent back and forth.
- ▷ Very small air bubbles observed directly below the inlet of the Nano do **not** disturb performance.
- ▷ The capsule function will **not** be influenced as long as the small air bubbles remain outside of the membrane bed.

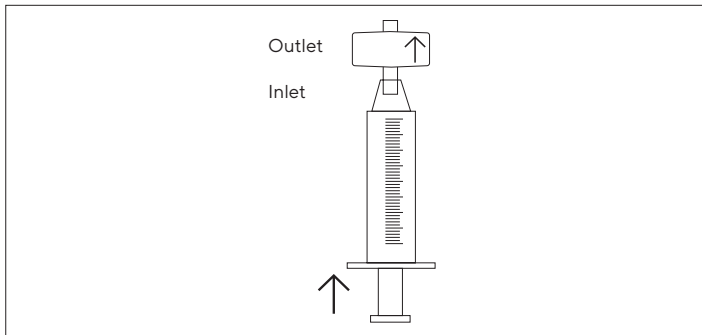


Fig. 7: Filling the Sartobind® Nano with a Luer syringe for air removal

5.3 Cleaning and Equilibration

Non-gamma-irradiated Sartobind® products have to be sanitized in place directly before use.

For **gamma-irradiated Sartobind®** Q capsules (see chapter “11 Ordering Information”, page 74), sanitization is optional.

- ▶ For sanitization use 30 membrane volumes (MV) of 1 N NaOH solution at the flow rate of 1 MV/min at 20°C. Preferentially, work at room temperature as low temperature increases viscosity of solvents. Also cold NaOH can cause swelling of the cellulose matrix significantly which will result in pressure increase.
- ▶ Flush with 10 MV of 1 N NaCl at 5 MV/min.
- ▶ Flush with 10 MV of equilibration buffer at 5 MV/min.

5.4 Autoclaving

NOTICE

Risk of damage to the product from improper handling!

- ▶ Do **not** perform in-line steam sterilization for the product.
 - ▶ Do **not** autoclave the cassettes. The cassette material is not compatible to autoclaving.
-

The capsules can be autoclaved once at 121°C for 30 minutes at 1 bar (0.1 MPa| 14.5 psig). Pre-wet the capsule with equilibration buffer. **Do not use pure water.** The protective caps enclosed in the Jumbo delivery must be reinstalled on inlet and outlet connectors of the Jumbo. Close valves immediately after sterilisation. For autoclaving of Sartobind® Jumbo refer to separate autoclaving instructions enclosed in delivery.

5.5 Gamma Irradiation

NOTICE

Damage to the product due to repeated gamma irradiation!

- ▶ Gamma irradiation may only be performed once.
-

Sartobind® Q with the product number starting 97 (see chapter 11.2 on page 78 and 11.5 on page 83) and Sartobind® Q Cassettes can also be gamma irradiated. The recommended dose for gamma irradiation is 25kGy and the maximum dose is 50kGy.

Products that are gamma irradiated can be identified by a colored indicator:

red indicator: gamma irradiated

orange indicator: **not** gamma irradiated

Once Sartobind® Q was irradiated, further autoclaving steps are prohibited, and vice versa.









5.6 Recommended Flow Rates and Equilibration Buffer Volumes

Membrane adsorbers can be run at much higher flow rate per volume than resin columns. As a rule of thumb, flow rates of 5 membrane volumes per minute are recommended for 8 mm bed height and 10 – 30 membrane volumes for 4 mm bed height. This recommendation is only a guideline as buffers and samples have different compositions and viscosities. Please test your respective flow rates with a small scale product to ensure that the flow rate fits with your pump capacities and the pressure limits.

Lower flow rates than the recommended ones can also be used but will typically **not** improve binding capacity or overall performance. Cold room temperature increases buffer viscosity and possibly back pressure.

The equilibration buffer volume is typically 10 membrane volumes depending on the type of buffer.

For the cassettes, flow rate and equilibration volumes have to be multiplied with the number of cassettes in use.

								
Membrane volume (MV)	1 mL	3 mL	10 mL	20 mL	75 mL	150 mL	200 mL	400 mL
Bed height (mm)	4	8	4	8	4	8	4	8
Rec. flow rate (L/min)	0.02	0.015	0.2	0.1	1.5	0.75	4	2
Rec. equilibration volume* (L)	0.01	0.03	0.1	0.2	0.75	1.5	2	4

* Refer to "5.3 Cleaning and Equilibration", page 30

** Multiply with number of used cassettes



400 mL



800 mL



600 mL



1.2 L



2.5 L



5 L



0.8 L



1.6 L

4

8

4

8

4

8

4

8

8

4

12

6

50

25

16**

8**

4

8

6

12

25

50

8**

16**

5.7 Buffer Conditions

In the majority of applications an equilibration buffer concentration of 10 mM provides sufficient buffering capacity and prevents the protein of interest from precipitation. The ionic strength should be kept as low as possible to avoid reduction of binding capacity.

The buffer should have a pKa within 0.5 pH units of the pH used. It should be filtered with 0.2 μm or 0.45 μm filters before use and the quality of water and chemicals should be of high purity.

NOTICE

Risk of reversible swelling and decreased permeability of the membrane due to the use of pure water!

▶ Do not apply pure water.

5.8 Selection of pH Conditions

In ion exchange chromatography a charged molecule is bound to oppositely charged groups attached to the insoluble matrix. This binding is reversible and induced by an increase of the salt concentration in the elution buffer.

The pH value at which a biomolecule has no net charge is the isoelectric point: pI . If the pH of the buffer is below the isoelectric point (rule of the thumb at least 1 pH unit), a protein has a positive net charge and will bind to a cation exchanger (e.g. Sartobind® S). If the pH of the buffer is above its isoelectric point (at least 1 pH unit), it will bind to an anion exchanger (e.g. Sartobind® Q).

5.9 Contaminant Removal From Therapeutic Proteins And Other Sources in Flow-Through Mode

For contaminant removal from products such as monoclonal antibodies, pH conditions in the range of pH 6 to 8 are used in order to bind highly negatively charged DNA, endotoxins, contaminating proteins, some host cell proteins and viruses with an anion exchanger. The product of interest, the monoclonal antibody with isoelectric points (pI) of 8–9.5 for example, will **not** bind and pass through Sartobind® Q. The influence of the flow rate on the performance is very low.

To remove contaminating proteins and aggregates with Sartobind® S in flow-through mode, process impurities have to be charged positively to bind to S while the target protein stays

negative. When the pH of the buffer is above the pI, the protein product flows through without binding.

5.10 Sample Preparation

Procedure

- ▶ Adjust the sample to the starting buffer conditions and prefilter it through a 0.2 µm membrane filter e.g. Sartopore® 2 XLG capsule.
- ▶ For small volumes in the mL range, use a 0.2 µm Minisart® filter with Luer outlet (order number 16532-K for polyethersulfone or 16534-K for cellulose acetate membrane).

NOTICE

Unfiltered feed will block the Membrane Adsorber and lead to capacity loss and increased back pressure!

- ▶ We recommend inline filtering during operation.
 - ▶ When the pressure increases replace the prefilter.
-

5.11 Washing

Procedure

- ▶ When using capsules in bind & elute mode, wash with equilibration buffer after sample loading.

5.12 Elution

Procedure

- ▶ To elute target protein or the virus, virus like particle (VLP) use buffers with appropriate salt concentration.

5.13 Draining

A dual air regulator system is recommended to prevent over-pressure of the Sartobind® products. The first regulator should reduce line air pressure to 2 bar.

The second regulator, positioned immediately upstream of the Sartobind®, should reduce the 2 bar regulated supply pressure to the <1 bar (14.5 psig) for a capsule and 0.5 bar (7.3 psig) for 1 to 13 cassettes draining pressure.

Procedure

- ▶ You may drain the capsule or cassette by application of air or nitrogen pressure (<1 bar | 14.5 psig) to the inlet of the product.

5.14 Regeneration and Storage

After elution wash with equilibration buffer. If necessary, use 1 N NaOH, 1 N HCl or 70% ethanol for 1 hour for regeneration and store in 20% ethanol in equilibration buffer.

5.15 Operation of the Sartobind® Nano With Peristaltic Pumps or Liquid Chromatography (LC) Systems

Procedure

- ▶ After the unit is filled completely with equilibration buffer, close the outlet of the Sartobind® Nano and remove the syringe.
- ▶ Start your LC system or peristaltic pump at a low flow rate.
- ▶ When fluid emerges, stop the pump, connect the tubing to the inlet of the Sartobind® Nano.

- ▶ Make sure that no air is introduced.
- ▶ Remove the cap from outlet.
- ▶ Run the pump until fluid emerges from the outlet of the unit and stop it.
- ▶ Connect the outlet of the unit via Luer adapter to the LC detector and proceed with loading.
- ▶ If your system pressure is too high, refer to your LC system manual to remove any flow restrictor after the UV cell, as the system may generate a pressure above the allowed maximum pressure.
- ▷ As membrane adsorbers are typically run at much higher flow rates than columns, there is no risk of bubble formation in the UV cell when removing the flow restrictor.

5.16 Scaling Up

Run break through experiments for the target compound (contaminants) to be bound on the membrane matrix. After optimization of the binding conditions, the purification step can be scaled up to a larger capsule.

Recommendations:

Maintain

- Bed height (stay within the same bed height when scaling up)
- Linear flow (when using capsules with same bed height, the flow rate will scale up linear when keeping MV/min constant)
- Sample concentration

Increase (see scaling factors in the following table)

- Sample load volume
- Volumetric flow rate
- Membrane volume

Scale up calculations are done preferably by keeping the bed height constant and adjusting the membrane volume. This will make the calculation simple. Other methods for scale up via residence time will lead to same results. Residence time is calculated by the membrane volume divided by the flow rate. You may change from 4 mm to 8 mm by keeping the residence time constant which is equal to the flow rate expressed in membrane volumes per minute but not above 5 MV/min.

When using Sartobind® Nano 1 or 3 mL, the scale up factor for flow rate and binding capacity is equal to the multiplication factor of membrane volumes for the listed scale up products:

Size	Membrane Volume (mL)	Factor to increase* (from Nano)	Bed height (mm)	Membrane Volume (mL)	Factor to increase* (from Nano)	Bed height (mm)
Nano	1	-	4	3	-	8
Mini	10	10	4	20	6.7	8
5"	75	75	4	150	50	8
10"	200	200	4	400	133	8
20"	400	400	4	800	266	8
30"	600	600	4	1200	400	8
Jumbo	2500	2500	4	5000	1667	8
Cassette	800	800	4	1600	533	8
Cassettes**	10400	10400	4	20800	6933	8

* Flow rate and binding capacity; ** 13 Cassettes as example

Example: After breakthrough experiments with the Nano 3 mL, you realize that a 1500-fold higher binding capacity is needed for a large scale run. Then you choose the 5 liter Jumbo capsule. To determine the running conditions of the Jumbo and to keep consistent upscaling, adjust the flow rate by a factor of ~1670. To assure the scale up, additional experiments with the 150 mL capsule (increase by a factor of 50) support this scale up calculation.

In the example above one and 13 cassettes are used, but any number of cassettes between one or 13 which is the maximum in the pilot filter holder can be applied. Then the factor to increase is adapted proportionally. For a larger scale above 20 L, a process and a double process filter holder is available as accessory (see chapter “11.6 Accessories”, page 86). To assure the scale, intermediate sizes are recommended to assure the design.

NOTICE

Keep sample concentration constant in lab and production scale.

- ▶ Adjustments might be required due to additional volumes from tubing and the system.
-

6 Testing of Membrane Samples Enclosed With 5 L Capsules

Membrane quality is always tested as described in chapter “8 Troubleshooting”, page 61. Additionally the 5 L Jumbo is delivered with membrane samples of same membrane lot that was used in the capsule. This allows again for testing of membrane with user specific tests before use of capsule. Static Binding capacity tests can be performed as follows:

6.1 Testing Sartobind® Q 30 mm Discs

Using the appropriate conditions, bovine serum albumin can be selectively bound to membrane adsorber 30 mm discs. By increasing the salt concentration of the buffer solution, the target substance can be recovered for subsequent quantification. BSA in excess is loaded at a batch experiment on Sartobind® Q membrane disc. The total capacity is obtained after eluting the bound BSA by increasing salt concentration. The amount of BSA is determined by spectrophotometer.

6.1.1 Materials for Test of Q Membrane Discs

Membrane sample:

Sartobind® Q membrane, 30 mm diameter
(membrane area: 7.1 cm²)

Test protein: bovine serum albumin (BSA)
Cat.No. 11930.03 SERVA

Buffers:

Equilibration buffer: 20 mM Tris|HCl,
(pH 7.4, conductivity 1.80 mS/cm)

Test solution: 0.2% BSA in equilibration buffer (2 mg/mL)

Elution buffer: 1 M NaCl in equilibration buffer
(Temperature of buffers: 20 – 25°C)

Equipments:

Glass or plastic dish

pH meter

Conductivity meter (temperature coefficient 2.00%/°C
at 25°C reference temperature)

Spectrophotometer set at 280 nm

Orbital or reciprocal shaker set at approx. 80 rpm

Operation temperature: 20 – 25°C

6.1.2 Method for Test of Q Membrane Discs

Equilibration

Apply 2 mL/cm² equilibration buffer (= 14.2 mL per 30 mm membrane disc) into the glass dish.

Place a membrane sample into the equilibration buffer (make sure that the membrane is wetted completely) and shake three times for 5 min each at 80 rpm.

Remove the equilibration buffer after each step by vacuum pump.

Loading

Apply 2 mL/cm² test solution into the dish. Close the dish and shake for 12 – 18 h at 80 rpm.

Remove excess test solution by vacuum pump as much as possible.

Washing

Apply 2 mL/cm² of equilibration buffer and shake for 15 min.

Remove equilibration buffer by vacuum pump.

Repeat this step once.

Elution

Place each membrane disc in a separate Petri dish. Apply 10 mL elution buffer onto each membrane disc with a pipet. Shake for 1 h at 80 rpm (= elution solution). The volume of the elution solution = V_{elute} (mL).

Extinction measurements

Zero the instrument with equilibration buffer.

Measure the extinction of the test solution = E_{test} .

Zero the instrument with elution buffer.

Measure extinction of the elution solution = E_{elute} .

6.1.3 Calculation of Membrane Binding Capacity

Binding capacity (mg/cm²)

$$= \frac{E_{\text{elute}} + C_{\text{test}} + V_{\text{elute}}}{E_{\text{test}} + MA}$$

$$= \frac{E_{\text{elute}} + 2 + 10}{1.25 + 7.1}$$

$$= E_{\text{elute}} + 2.25$$

E_{elute} : extinction 280 nm of the elution solution

C_{test} : concentration of the test solution (2 mg/mL)

V_{elute} : volume of the elution solution (20 mL)

E_{test} : extinction 280 nm of the test solution (= 1.25),
d (cell width) = 1 cm

MA: membrane area (= 7.1 cm²)

6.2 Testing Sartobind® S 30 mm Discs

Using the appropriate conditions, lysozyme can be selectively bound to membrane adsorber 30 mm discs. By increasing the salt concentration of the buffer solution, the target substance can be recovered for subsequent quantification. Lysozyme in excess is loaded at a batch experiment on Sartobind® S membrane disc. The total capacity is obtained by eluting the bound protein by increasing salt concentration. The amount of lysozyme is determined by spectrophotometer.

6.2.1 Materials for Test of S Membrane Discs

Membrane sample:

Sartobind® S membrane, 30 mm diameter

(membrane area: 7.1 cm²)

Test protein: lysozyme Cat.No. L-6776 Sigma

Buffers:

Equilibration buffer: 10 mM potassium phosphate,
(pH 7.0, conductivity 1.75 mS/cm)

Test solution: 0.2% lysozyme in equilibration buffer
(2 mg/mL)

Elution buffer: 1 M NaCl in equilibration buffer
(Temperature of buffers: 20 – 25°C)

Equipments:

Plastic or glass dish

pH meter

Conductivity meter (temperature coefficient 2.00%/°C
at 25°C reference temperature)

Spectrophotometer set at 280 nm

Orbital or reciprocal shaker set at approx. 80 rpm

Operation temperature: 20 – 25°C

6.2.2 Method for Test of S Membrane Discs

Equilibration

Apply 2 mL/cm² equilibration buffer into the dish. Place a membrane sample into the equilibration buffer (make sure that the membrane is wetted completely) and shake three times for 5 min each at 80 rpm. Remove the equilibration buffer after each step by vacuum pump.

Loading

Apply 2 mL/cm² test solution into the dish. Close the dish and shake for 12 – 18 h at 80 rpm.

Remove excess test solution by vacuum pump as much as possible.

Washing

Apply 2 mL/cm² of equilibration buffer and shake for 15 min.

Remove equilibration buffer by vacuum pump.

Repeat this step once.

Elution

Place each membrane disc in a separate Petri dish. Apply 10 mL elution buffer onto each membrane disc with a pipet. Shake for 1 h at 80 rpm (= elution solution). The volume of the elution solution = V_{elute} (mL).

Extinction measurements

Zero the instrument with equilibration buffer. Measure the extinction of the test solution = E_{test} .

Zero the instrument with elution buffer. Measure extinction of the elution solution = E_{elute} .

6.2.3 Calculation of Membrane Binding Capacity

Binding capacity (mg/cm^2)

$$= \frac{E_{\text{elute}} + C_{\text{test}} + V_{\text{elute}}}{E_{\text{test}} + \text{MA}}$$

$$= \frac{E_{\text{elute}} + 2 + 10}{5.0 + 7.1}$$

$$= E_{\text{elute}} + 0.56$$

E_{elute} : extinction 280 nm of the elution solution

C_{test} : concentration of the test solution (2 mg/mL)

V_{elute} : volume of the elution solution (20 mL)

E_{test} : extinction 280 nm of the test solution (= 5.0)

MA: membrane area (= 7.1 cm^2)

7 Integrity Test by Diffusion

The integrity of a membrane adsorber can be tested by a diffusion test.

The testing procedure describes the diffusion test for pre and post use. The test is intended to discriminate between defective and intact products and to detect major bypasses, large holes and faulty assembly.

7.1 Installation

Procedure

- ▶ Install adsorber as shown in Fig. 8.

The test procedure has been developed and checked with the Sartochek[®] instrument family e.g. Sartochek[®] 4 Plus (26288) or 4 (16288). The use of Sartochek[®] instruments older than Sartochek[®] 4 will generate faulty data.

- ▶ Note that the test procedure with other vendor's integrity testers can require a different set up.

7.2 Operation Procedure

7.2.1 Pre-washing of Device

Prior to integrity testing, the capsule needs to be prewashed, to remove any glycerol. The washing solution should have room temperature.

NOTICE

Risk of reversible swelling and decreased permeability of the membrane due to the use of pure water!

- ▶ Do **not** apply pure water.

Procedure

- ▶ Keep the unit in an upright position for proper venting and open the vent screw on top of the product until all air is replaced by the testing solvent.
- ▶ Pre-wash the product with 30 membrane volumes (MV) of buffer or 0.9% NaCl in water at recommended flow rate.

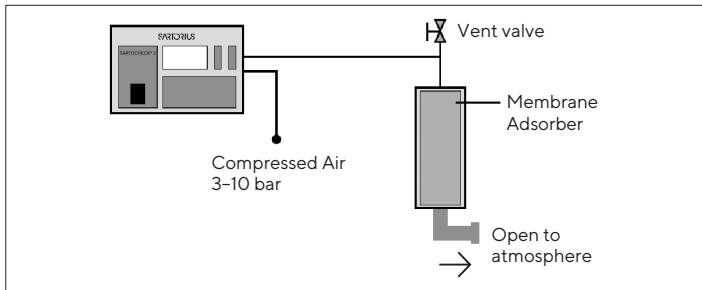


Fig. 8: Set up of diffusion test with Sartocheck®.

7.2.2 Diffusion Measurement With Sartocheck®

- Choose “Programming” in the main menu
- Choose “Diffusion Test”

Procedure

- ▶ Choose the test pressure, stabilization and testing time for your product from the table (next page).
- ▶ If you set the Net Volume to zero, Sartocheck® automatically measures the upstream void volume including tubing.

Test Parameters 4 mm

Size	Bed height (mm)	Membrane Volume (MV) (mL)	Test pressure mbar (psig)	Stabilization time (min)	Testing time (min)	Diffusion max. (mL/min)
Nano	4	1	200 (2.9)	2	1	15
Mini	4	10	200 (2.9)	2	1	15
5"	4	75	200 (2.9)	2	1	15
10"	4	200	200 (2.9)	3	1	15
20"	4	400	200 (2.9)	3	1	15
30"	4	600	200 (2.9)	3	1	15
Jumbo	4	2500	200 (2.9)	3	1	15
Cassette(s)*	4	800 – 10,400	200 (2.9)	5	1	15 – 195

* Diffusion max. per 4 or 8 mm cassette is 15 mL/min multiplied by number of cassettes

Test Parameters 8 mm

Size	Bed height (mm)	Membrane Volume (MV) (mL)	Test pressure mbar (psig)	Stabilization time (min)	Testing time (min)	Diffusion max. (mL/min)
Nano	8	3	200 (2.9)	2	1	15
Mini	8	20	200 (2.9)	2	1	15
5"	8	150	200 (2.9)	2	1	15
10"	8	400	200 (2.9)	3	1	15
20"	8	800	200 (2.9)	3	1	15
30"	8	1,200	200 (2.9)	3	1	15
Jumbo	8	5,000	200 (2.9)	5	1	15
Cassette(s)*	8	1,600–20,800	200 (2.9)	5	1	15–195

* Diffusion max. per 4 or 8 mm cassette is 15 mL/min multiplied by number of cassettes

7.2.3 Results and Evaluation for One Device

- Diffusion \leq Diffusion max.:
Test passed (diffusion value on the print out)
- Diffusion $>$ Diffusion max.:
Test failed (red text on the print out)

The maximum allowed diffusion values are per product.

If you set up for example 10 cassettes in a holder the values have to be added up and the max diffusion value will be 150 mL.

8 Troubleshooting

Problem	Possible cause	Action
Air bubbles can be seen	Incomplete air removal	Small air bubbles seen in the top of the unit do not interfere with the purification as long as they do not touch the membrane bed. If too much air is enclosed, repeat removal as described in chapter “5.1 Venting the Capsule”, page 27.
I installed the capsule upside down	Installation of capsule may be easier in the process flow	Validation has been done with a process flow from top to bottom. Thus it is clearly recommended to use capsules in the described flow direction (Feed enters capsule on top and leaves it on bottom).
I deviated from the CIP and flushing equilibration procedure		The capsules have been qualified and validated according the given procedure. If a deviation is necessary, the results may also deviate from the given validation data.
High back pressure	Material has not been filtered	Prefilter with 0.2 μm or 0.45 μm filter before processing through the unit (preferentially inline).

Problem	Possible cause	Action
High back pressure	Material has been filtered but was stored before purification	Proteins can form aggregates within hours or during operation. Thus we recommend to prefilter inline by attaching a 0.2 μm filter in front of the adsorber. When you observe again pressure built up, replace the filter.
	LC system generates high pressure	Remove restrictor after the UV cell.
	The adsorber is clogged membrane fouling	Replace unit. Perform a regeneration cycle. You may backflush within given flow and pressure limits.
	Viscosity swelling effects	Work at room temperature, avoid lower temperatures.
	Pure water leads to swelling of membrane	Add sodium chloride or use ionic buffers.
Target molecule is not bound	Conditions for binding are insufficient	Decrease conductivity, control other process parameters as type of buffer and pH.

Problem	Possible cause	Action
Binding capacity is not sufficient	Process conditions not optimized	Use larger adsorber product, or: connect two adsorbers (same size) in series (i.e connect outlet of first adsorber to inlet of second) to achieve higher binding capacity. As a rule of thumb the pressure doubles when the flow rate is kept constant and the number of membrane layers is doubled.
Reuse is needed	For economic or practical reasons	The major application of Sartobind® capsules is the single use and they are constructed in plastic housing for this. Also they are validated and certified only for one use. Technically they can be reused. The durability of the unit depends on the nature of sample and sample preparation, prefiltration as well as proper regeneration and application. Plastic materials and membranes allow CIP and long term storage if carefully treated. For reuse validation we assist you with our Validation Service. Please ask your local representative.

Problem	Possible cause	Action
Binding capacity decreases after several uses	Improper filtration	Prefilter with 0.2 μm filter before processing through the unit.
	Some molecule species binds tightly (e.g. DNA on Sartobind® Q) and cannot be removed with 1 N NaOH 1 h	Use capsule only once or use DNA endonucleases such as Denarase®.
	Protein or contaminants are still bound from last cycle	Run a 1 M NaCl buffer step to elute tightly bound proteins quantitatively. Then regenerate adsorber by loading with 1 N NaOH and keep it for 1 hour at room temperature (20°C).
	Wrong storage	Do not store in sodium hydroxide containing buffers. Store long term* in 20% ethanol buffer (e.g. equilibration buffer) solution and do not use oxidative chemicals in buffers.
A vertical line is seen on one capsule side when filled	Membrane edge visible	No action necessary. It can be visible the edge of the nonwoven touching the inner tube.

Problem	Possible cause	Action
I purged with air or nitrogen and lost flow and binding capability.	Air has entered into the pores	See troubleshooting “Applied bubble point instead of diffusion test” below.
Accidentally a bubble point test instead of diffusion test has been run	Operation error	The membrane has then to be purged extensively to remove all the air which has been pressed into the pores. If properly purged, the diffusion test can be run successfully and the product works as expected.
Cassettes system leaks or fails at integrity test	Wrong assembly	Position manifolds and cassettes at the lowest position in the holder otherwise seals are not perfectly aligned.

9 Technical Data

9.1 Bed Height 4 mm

Membrane volume (MV)	1 mL	10 mL	75 mL
Nominal membrane area	36.4 cm ²	364 cm ²	2,700 cm ²
Bed height	4 mm	4 mm	4 mm
Design	Cylindrical	Cylindrical	Cylindrical
Sartobind® Q typical 10% dynamic binding capacity*	29 mg	290 mg	2.16 g
Sartobind® S typical 10% dynamic binding capacity*	25 mg	250 mg	1.89 g
Maximum pressure bar (MPa, psig) at 20°C	4 (0.4, 58)	4 (0.4, 58)	4 (0.4, 58)
Maximum pressure during venting bar (MPa, psig) at 20°C	-	0.5 (0.05, 7)	0.5 (0.05, 7)
Nominal void volume (mL)	3.5	32	200
Nominal void volume (MV)	3.5	3.2	2.7
Approximate weight	10 g	65 g	400 g

1 mL membrane = 36.4 cm² membrane

Ion capacity per cm² of membranes: 2–5 µeq

Short term pH stability Q | S: 1-14 | 3-14 refers to cleaning in place and regeneration procedures during operation

200 mL	400 mL	600 mL	2.5 L	800 mL
7,300 cm ²	14,600 cm ²	22,000 cm ²	91,000 cm ²	29,000 cm ²
4 mm	4 mm	4 mm	4 mm	4 mm
Cylindrical	Cylindrical	Cylindrical	Cylindrical	Flat sheet
5.8 g	11.7 g	17.6 g	73 g	23.2 g
5.1 g	10.2 g	15.4 g	-	20.3 g
4 (0.4, 58)	4 (0.4, 58)	4 (0.4, 58)	3 (0.3, 43.5)	2 (0.2, 29)
0.5 (0.05, 7)	0.5 (0.05, 7)	0.5 (0.05, 7)	0.5 (0.05, 7)	0.5 (0.05, 7)
540	1,080	1,600	7,000	2,500
2.7	2.7	2.7	2.8	3.1
750 g	1.3 kg	1.9 kg	16 kg 20 kg wet 23 kg filled	4.9 kg 6.0 kg wet

Long term storage pH stability Q|S: 2-12|4-13 refers to overnight storage and longer. Preferably store units in 20% ethanol|buffer

* See section “9.4 Binding Capacity”, page 71

9.2 Bed Height 8 mm

Membrane volume (MV)	3 mL	20 mL	150 mL
Nominal membrane area	110 cm ²	728 cm ²	5,500 cm ²
Bed height	8 mm	8 mm	8 mm
Design	Cylindrical	Cylindrical	Cylindrical
Sartobind® Q typical 10% dynamic binding capacity*	88 mg	580 mg	4.4 g
Sartobind® S typical 10% dynamic binding capacity*	77 mg	500 mg	3.9 g
Maximum pressure bar (MPa, psig) at 20°C	4 (0.4, 58)	4 (0.4, 58)	4 (0.4, 58)
Maximum pressure during venting bar (Mpa, psig) at 20°C	-	0.5 (0.05, 7)	0.5 (0.05, 7)
Nominal void volume (mL)	4	32	200
Nominal void volume (MV)	1.3	1.6	1.3
Approximate weight	10 g	65 g	400 g

1 mL membrane = 36.4 cm² membrane

Ion capacity per cm² of membranes: 2–5 µeq

Short term pH stability Q | S: 1-14 | 3-14 refers to cleaning in place and regeneration procedures during operation

400 mL	800 mL	1.2 L	5 L	1.6 L
14,600 cm ²	29,000 cm ²	44,000 cm ²	182,000 cm ²	58,000 cm ²
8 mm	8 mm	8 mm	8 mm	8 mm
Cylindrical	Cylindrical	Cylindrical	Cylindrical	Flat sheet
11.7 g	23.3 g	35 g	145 g	46 g
10.2 g	20.4 g	31 g	127 g	41 g
4 (0.4, 58)	4 (0.4, 58)	4 (0.4, 58)	3 (0.3, 43.5)	2 (0.2, 29)
0.5 (0.05, 7)	0.5 (0.05, 7)	0.5 (0.05, 7)	0.5 (0.05, 7)	0.5 (0.05, 7)
540	1,080	1,600	7,000	2,900
1.4	1.4	1.3	1.4	1.8
760 g	1.3 kg	1.9 kg	16 kg 20 kg wet 23 kg filled	4.9 kg 6.5 kg wet

Long term storage pH stability Q | S: 2-12 | 4-13 refers to overnight storage and longer. Preferably store units in 20% ethanol | buffer

* See section “9.4 Binding Capacity”, page 71

9.3 Materials

Membrane materials

Matrix	Stabilized reinforced cellulose
Membrane thickness membrane volume = membrane area	275 μm 1 mL = 36.4 cm^2
Nominal pore size	> 3 μm
Ion exchanger ligand Q	Strong anion Q: trimethyl ammonium moiety ($-\text{N}^+(\text{CH}_3)_3$)
Ion exchanger ligand S	Strong cation S: sulfonic acid moiety ($-\text{SO}_3^-$)

Capsule materials

Outer cage, inner core, end caps, capsule housing	PP (polypropylene)
Nonwoven Non-gamma-irradiatable products	PP (polypropylene)
Gamma-irradiatable gamma-irradiated products	PET (polyethylene terephthalate)
O-ring in vent valve (except Nano)	EPDM (ethylene propylene diene monomer)

Cassette materials

Outer cage, seal, nonwoven

ABS (acrylonitrile butadiene styrene),
silicone, PET (polyethylene terephthalate), TPE (thermoplastic elastomers)

9.4 Binding Capacity

Data are based on dynamic binding capacity measurements 10% using 3 layers of 5 cm² membrane discs (15 cm² total area, membrane thickness of 275 μm) arranged in a holder and run at 10 mL/min.

	Typical dynamic binding capacity 10%	Reference protein and buffer
Q	0.8 mg/cm ² (29 mg/mL)	BSA (bovine serum albumin) in 20 mM Tris/HCl, pH 7.5
S	0.7 mg/cm ² (25 mg/mL)	Lysozyme in 10 mM potassium phosphate, pH 7.0

9.5 Chemical Stability

Stable for all commonly used buffers in chromatography

No oxidizing agents

9.6 Storage Conditions

Clean, dry place in closed bag and box at room temperature.

Protected from direct sunlight.

Store membrane samples enclosed in Jumbo shipment in a safe place as they **cannot** be ordered again.

10 Quality Assurance

The final Sartobind® products are tested for protein dynamic binding capacity and flow rate. Sartobind® membranes are tested for protein dynamic binding capacity, flow rate, thickness, and evenness.

Capsules, cassettes and membranes are manufactured in a controlled environment. The products meet all Sartorius standards for traceability, production and specifications as given here or exceeded them as certified in the quality assurance certificate enclosed. A validation and an extractables guide are available on request.

11 Ordering Information

11.1 Products 4 mm Bed Height

Order number	Description	Quantity
96IEXQ42DN-11	Sartobind® Q Nano 1 mL, 4 mm, Luer female connectors, 2 PEEK adapters Luer male to UNF 10-32 female	1
96IEXQ42DN-11--A	Sartobind® Q Nano 1 mL, 4 mm, Luer female connectors, 2 PEEK adapters Luer male to UNF 10-32 female	4
96IEXS42DN-11	Sartobind® S Nano 1 mL, 4 mm, Luer female connectors, 2 PEEK adapters Luer male to UNF 10-32 female	1
96IEXS42DN-11--A	Sartobind® S Nano 1 mL, 4 mm, Luer female connectors, 2 PEEK adapters Luer male to UNF 10-32 female	4

Order number	Description	Quantity
96IEXQ42D4R11--A	Sartobind® Q Mini 10 mL, 4 mm, Luer female connectors, 8 PEEK adapters Luer male to UNF 10 - 32 female	4
96IEXQ42D4RFF--A	Sartobind® Q Mini 10 mL, 4 mm, ¾" sanitary clamp	4
96IEXQ42D4ROO--A	Sartobind® Q Mini 10 mL, 4 mm, hose barb connectors	4
96IEXS42D4R11--A	Sartobind® S Mini 10 mL, 4 mm, Luer female connectors, 8 PEEK adapters Luer male to UNF 10 - 32 female	4
96IEXS42D4RFF--A	Sartobind® S Mini 10 mL, 4 mm, ¾" sanitary clamp	4
96IEXS42D4ROO--A	Sartobind® S Mini 10 mL, 4 mm, hose barb connectors	4
96IEXQ42D9MOO--A	Sartobind® Q 75 mL, 4 mm, hose barb connectors	4
96IEXQ42D9MFF--A	Sartobind® Q 75 mL, 4 mm, ¾" sanitary clamp	4

Order number	Description	Quantity
96IEXS42D9MOO--A	Sartobind® S 75 mL, 4 mm, hose barb connectors	4
96IEXS42D9MFF--A	Sartobind® S 75 mL, 4 mm, ¾" sanitary clamp	4
96IEXQ42D1GSS	Sartobind® Q 200 mL, 4 mm, 1½" sanitary clamp	1
96IEXS42D1GSS	Sartobind® S 200 mL, 4 mm, 1½" sanitary clamp	1
96IEXQ42D2HSS	Sartobind® Q 400 mL, 4 mm, 1½" sanitary clamp	1
96IEXS42D2HSS	Sartobind® S 400 mL, 4 mm, 1½" sanitary clamp	1
96IEXQ42D3KSS	Sartobind® Q 600 mL, 4 mm, 1½" sanitary clamp	1
96IEXS42D3KSS	Sartobind® S 600 mL, 4 mm, 1½" sanitary clamp	1
96IEXQ42D3NSS	Sartobind® Q Jumbo 2.5 l, 4 mm, 1½" sanitary clamp, 2 protective caps	1

Order number	Description	Quantity
98IEXQ42D-L	Sartobind® Q Cassette 0.8 L, 4 mm, 1½" sanitary clamp via manifold set (accessory)	1
98IEXS42D-L	Sartobind® S Cassette 0.8 L, 4 mm, 1½" sanitary clamp via manifold set (accessory)	1
98IEXQ42DGL	Sartobind® Q Cassette 0.8 L, 4 mm, gamma irradiated, 1½" sanitary clamp via manifold set (accessory)	1

11.2 Products 4 mm Bed Height, Gamma Irradiatable

Order number	Description	Quantity
971EXQ42D-U11	Sartobind® Q Nano 1 mL, 4 mm, Luer female connectors, 2 PEEK adapters Luer male to UNF 10-32 female	1
971EXQ42D-U11--A	Sartobind® Q Nano 1 mL, 4 mm, Luer female connectors, 2 PEEK adapters Luer male to UNF 10-32 female	4
971EXQ42D4R11	Sartobind® Q Mini 10 mL, 4 mm, Luer female connectors, 8 PEEK adapters Luer male to UNF 10-32 female	1
971EXQ42D4R11--A	Sartobind® Q Mini 10 mL, 4 mm, Luer female connectors, 8 PEEK adapters Luer male to UNF 10-32 female	4
971EXQ42D4RFF	Sartobind® Q Mini 10 mL, 4 mm, ¾" sanitary clamp	1
971EXQ42D4RFF--A	Sartobind® Q Mini 10 mL, 4 mm, ¾" sanitary clamp	4

Order number	Description	Quantity
97IEXQ42D4ROO	Sartobind® Q 75 mL, 4 mm, hose barb connectors	1
97IEXQ42D4ROO--A	Sartobind® Q 75 mL, 4 mm, hose barb connectors	4
97IEXQ42D9MFF	Sartobind® Q 75 mL, 4 mm, ¾" sanitary clamp	1
97IEXQ42D9MOO	Sartobind® Q 75 mL, 4 mm, hose barb connectors	1
97IEXQ42D1GSS	Sartobind® Q 200 mL, 4 mm, 1½" sanitary clamp	1
97IEXQ42D1GOO	Sartobind® Q 200 mL, 4 mm, hose barb connectors	1
97IEXQ42D2HSS	Sartobind® Q 400 mL, 4 mm, 1½" sanitary clamp	1
97IEXQ42D2HOO	Sartobind® Q 400 mL, 4 mm, hose barb connectors	1
97IEXQ42D3KSS	Sartobind® Q 600 mL, 4 mm, 1½" sanitary clamp	1
97IEXQ42D3KOO	Sartobind® Q 600 mL, 4 mm, hose barb connectors	1

11.3 Products 4 mm Bed Height, Gamma Irradiated

Order number	Description	Quantity
SBG500000	Sartobind® Q Nano 1 mL, 4 mm Luer female connectors, 2 PEEK adapters Luer male to UNF 10-32 female	4

11.4 Products 8 mm Bed Height

Order number	Description	Quantity
96IEXQ42EUC11--A	Sartobind® Q Nano 3 mL, 8 mm, Luer female connectors, 2 PEEK adapters Luer male to UNF 10-32 female	4
96IEXS42EUC11--A	Sartobind® S Nano 3 mL, 8 mm, Luer female connectors, 2 PEEK adapters Luer male to UNF 10-32 female,	4
96IEXQ42E4J11--A	Sartobind® Q Mini 20 mL, 8 mm, Luer female connectors, 8 PEEK adapters Luer male to UNF 10-32 female	4

Order number	Description	Quantity
96IEXQ42E4JFF--A	Sartobind® Q Mini 20 mL, 8 mm, ¾" sanitary clamp	4
96IEXQ42E4JOO--A	Sartobind® Q Mini 20 mL, 8 mm, hose barb connectors	4
96IEXS42E4J11--A	Sartobind® S Mini 20 mL, 8 mm, Luer female connectors, 2 PEEK adapters Luer male to UNF 10-32 female	4
96IEXS42E4JFF--A	Sartobind® S Mini 20 mL, 8 mm, ¾" sanitary clamp	4
96IEXS42E4JOO--A	Sartobind® S Mini 20 mL, 8 mm, hose barb connectors	4
96IEXQ42E9BFF	Sartobind® Q 150 mL, 8 mm, ¾" sanitary clamp	1
96IEXS42E9BFF	Sartobind® S 150 mL, 8 mm, ¾" sanitary clamp	1
96IEXQ42E1HSS	Sartobind® Q 400 mL, 8 mm, 1½" sanitary clamp	1
96IEXS42E1HSS	Sartobind® S 400 mL, 8 mm, 1½" sanitary clamp	1

Order number	Description	Quantity
96IEXQ42E2LSS	Sartobind® Q 800 mL, 8 mm, 1½" sanitary clamp	1
96IEXS42E2LSS	Sartobind® S 800 mL, 8 mm, 1½" sanitary clamp	1
96IEXQ42E3FSS	Sartobind® Q 1.2 L, 8 mm, 1½" sanitary clamp	1
96IEXS42E3FSS	Sartobind® S 1.2 L, 8 mm, 1½" sanitary clamp	1
96IEXQ42E3ESS	Sartobind® Q Jumbo 5 L, 8 mm, 1½" sanitary clamp, 2 protective caps, 15 membrane discs 30 mm	1
96IEXS42E3ESS	Sartobind® S Jumbo 5 L, 8 mm, 1½" sanitary clamp, 2 protective caps, 15 membrane discs 30 mm	1
98IEXQ42E-P	Sartobind® Q Cassette 1.6 L, 8 mm, 1½" sanitary clamp via manifold set (accessory)	1
98IEXS42E-P	Sartobind® S Cassette 1.6 L, 8 mm, 1½" sanitary clamp via manifold set (accessory)	1

Order number	Description	Quantity
98IEXQ42EGP	Sartobind® Q Cassette 1.6 L, 8 mm, gamma irradiated, 1½" sanitary clamp via manifold set (accessory)	1

11.5 Products 8 mm Bed Height, Gamma Irradiatable

Order number	Description	Quantity
97IEXQ42E-C11	Sartobind® Q Nano 3 mL, 8 mm, Luer female connectors, 2 PEEK adapters Luer male to UNF 10-32 female	1
97IEXQ42E-C11--A	Sartobind® Q Nano 3 mL, 8 mm, Luer female connectors, 2 PEEK adapters Luer male to UNF 10-32 female	4
97IEXQ42E4J11	Sartobind® Q Mini 20 mL, 8 mm, Luer female connectors, 2 PEEK adapters Luer male to UNF 10-32 female	1
97IEXQ42E4J11--A	Sartobind® Q Mini 20 mL, 8 mm, Luer female connectors, 8 PEEK adapters Luer male to UNF 10-32 female	4

Order number	Description	Quantity
97IEXQ42E4JFF	Sartobind® Q Mini 20 mL, 8 mm, ¾" sanitary clamp	1
97IEXQ42E4JFF--A	Sartobind® Q Mini 20 mL, 8 mm, ¾" sanitary clamp	4
97IEXQ42E4JOO	Sartobind® Q Mini 20 mL, 8 mm, hose barb connectors	1
97IEXQ42E4JOO--A	Sartobind® Q Mini 20 mL, 8 mm, hose barb connectors	4
97IEXQ42E9BFF	Sartobind® Q 150 mL, 8 mm, ¾" sanitary clamp	1
97IEXQ42E9BOO	Sartobind® Q 150 ml, 8 mm, hose barb connectors	1
97IEXQ42E1HSS	Sartobind® Q 400 mL, 8 mm, 1½" sanitary clamp	1
97IEXQ42E1HOO	Sartobind® Q 400 mL, 8 mm, hose barb connectors	1
97IEXQ42E2LSS	Sartobind® Q 800 mL, 8 mm, 1½" sanitary clamp	1




Order number	Description	Quantity
97IEXQ42E2LOO	Sartobind® Q 800 mL, 8 mm, hose barb connectors	1
97IEXQ42E3FSS	Sartobind® Q 1.2 L, 8 mm, 1½" sanitary clamp	1
97IEXQ42E3FOO	Sartobind® Q 1.2 L, 8 mm, hose barb connectors	1

11.6 Accessories






Order number	Description	Quantity
1ZA---0004	Adapter Luer male to UNF 10-32 female, PEEK	1
1ZAOGV0003	Adapter UNF 10-32 female to sanitary ¾", 25 mm, polyoxymethylene	2
5ZGI--0001	Holder for 1 × 200 to 1,200 mL (10-30") capsule, stainless steel, 3 legs	1
5ZALB-0002	Distribution adapter for 3 × 200 (10-30") to 1200 mL capsules, 1 × 2", 3 × 1½", sanitary, stainless steel	1
7ZAL-V0013	Reducing adapter 1½" (50.5 mm) to ¾" (25 mm), sanitary	1
7ZAL-V0010	Reducing adapter 2" (64 mm) to 1½" (50.5 mm), sanitary	1
9ZGL--0102	Trolley for Jumbo 2.5 or 5 L, stainless steel	1
26787---FT	Sartocheck® 5 Filter Tester	1
26787---FT---P	Sartocheck® 5 Plus Filter Tester	1
29Z-S00001	Manifold set for Sartoclear® Sartobind®, 1½" sanitary clamp	2

Order number	Description	Quantity
29Z-S00003	Manifold set for Sartobind®, gamma irradiated, 1½" sanitary clamp	2
2ZGL--0005	Pilot filter holder for Sartoclear® Sartobind®	1
2ZGL--0006	Process filter holder for Sartoclear® Sartobind®	1
2ZGL--0007	Double process filter holder for Sartoclear® Sartobind®	1
2ZGL--0008	Drip pan for Pilot Filter holder	1
2ZGL--0015	Drip pan for Process and double Process Filter Holder	1

12 Dimensions and Connections

			
Membrane volume			
4 mm	1 mL	10 mL	75 mL
8 mm	3 mL	20 mL	150 mL
Size	Nano	Mini	5"
Dimensions in mm	37×32 H×Ø	Luer: 70×55 Sanitary: 100×55 Hose barb: 110×55 H×Ø	Sanitary: 190×77 Hose barb: 204×77 H×Ø
Connectors	Luer female	- Luer female - Sanitary 3/4", 25 mm outer, 14 mm inner Ø - Hose barb 1/2", 12.7 mm*	- Sanitary 3/4" 25 mm outer, 14 mm inner Ø - Hose barb 1/2", 12.7 mm*
Gaskets	n.a.	3/4", inner Ø 16 mm	3/4", inner Ø 16 mm

n.a. = not available | * Recommended internal diameter of flexible tube: 1/2", 12.7 mm

				
200 mL 400 mL	400 mL 800 mL	600 mL 1.2 L	2.5 L 5 L	0.8 L 1.6 L
10"	20"	30"	Jumbo	Cassette
Sanitary: 347 × 100 Hose barb: 359 × 100 H × Ø	Sanitary: 579 × 100 Hose barb: 591 × 100 H × Ø	Sanitary: 807 × 100 Hose barb: 819 × 100 H × Ø	850 × 302 H × Ø	634 × 387 × 49 W × L × Ø
Sanitary 1½" 50.5 mm outer, 36 mm inner Ø	Sanitary 1½" 50.5 mm outer, 36 mm inner Ø	Sanitary 1½" 50.5 mm outer, 36 mm inner Ø	Sanitary 1½" 50.5 mm outer, 36 mm inner Ø	Via manifold: Sanitary 1½" 50.5 mm outer, 36 mm inner Ø
1½", inner Ø 35.8 mm	1½", inner Ø 35.8 mm	1½", inner Ø 35.8 mm	1½", inner Ø 35.8 mm	For manifold: 1½", inner Ø 35.8 mm

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen, Germany

Phone: +49 551 308 0
www.sartorius.com

© 2024
Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen, Germany

ML | Publication No.: SL-6210-e240501
DIR: 2624867-000-06

The information and figures contained in these instructions correspond to the version date specified below.

Sartorius reserves the right to make changes to the technology, features, specifications and design of the equipment without notice. Masculine or feminine forms are used to facilitate legibility in these instructions and always simultaneously denote all genders.

Copyright notice:

These instructions, including all components, are protected by copyright.

Any use beyond the limits of the copyright law is not permitted without our approval.

This applies in particular to reprinting, translation and editing irrespective of the type of media used.

Last updated:
05 | 2024

List of Sartorius Material Numbers Applying to EPA-FIFRA

96IEXQ42DN-11	96IEXS42D9MFF--A	97IEXQ42D4R11
96IEXQ42DN-11--A	96IEXQ42D1GSS	97IEXQ42D4R11--A
96IEXS42DN-11	96IEXS42D1GSS	97IEXQ42D4RFF
96IEXS42DN-11--A	96IEXQ42D2HSS	97IEXQ42D4RFF--A
96IEXQ42D4R11--A	96IEXS42D2HSS	97IEXQ42D4ROO
96IEXQ42D4RFF--A	96IEXQ42D3KSS	97IEXQ42D4ROO--A
96IEXQ42D4ROO--A	96IEXS42D3KSS	97IEXQ42D9MFF
96IEXS42D4R11--A	96IEXQ42D3NSS	97IEXQ42D9MOO
96IEXS42D4RFF--A	98IEXQ42D-L	97IEXQ42D1GSS
96IEXS42D4ROO--A	98IEXS42D-L	97IEXQ42D1GOO
96IEXQ42D9MOO--A	98IEXQ42DGL	97IEXQ42D2HSS
96IEXQ42D9MFF--A	97IEXQ42D-U11	97IEXQ42D2HOO
96IEXS42D9MOO--A	97IEXQ42D-U11--A	97IEXQ42D3KSS

97IEXQ42D3KOO	96IEXQ42E3FSS	97IEXQ42E9BOO
SBG500000	96IEXS42E3FSS	97IEXQ42E1HSS
96IEXQ42EUC11--A	96IEXQ42E3ESS	97IEXQ42E1HOO
96IEXS42EUC11--A	96IEXS42E3ESS	97IEXQ42E2LSS
96IEXQ42E4J11--A	98IEXQ42E-P	97IEXQ42E2LOO
96IEXQ42E4JFF--A	98IEXS42E-P	97IEXQ42E3FSS
96IEXQ42E4JOO--A	98IEXQ42EGP	97IEXQ42E3FOO
96IEXS42E4J11--A	97IEXQ42E-C11	
96IEXS42E4JFF--A	97IEXQ42E-C11--A	
96IEXS42E4JOO--A	97IEXQ42E4J11	
96IEXQ42E9BFF	97IEXQ42E4J11--A	
96IEXS42E9BFF	97IEXQ42E4JFF	
96IEXQ42E1HSS	97IEXQ42E4JFF--A	
96IEXS42E1HSS	97IEXQ42E4JOO	
96IEXQ42E2LSS	97IEXQ42E4JOO--A	
96IEXS42E2LSS	97IEXQ42E9BFF	