



sartorius stedim  
biotech

## Microporous PES Membranes High Efficiency Membranes for Filtration and Venting Applications



turning science **into solutions**

# Sartorius Stedim Biotech

## Strong Commitment to the Industry

Sartorius Stedim Biotech is a leading provider of cutting-edge membrane technologies for the industry around the world. Sartorius Stedim Biotech focuses on technologies to meet the rapidly changing requirements of the industry it serves. Strongly rooted in the scientific community and closely allied with customers and technology partners, the company is dedicated to its philosophy of "Turning science into solutions".

Headquartered in Aubagne, France, Sartorius Stedim Biotech is listed on the Eurolist of Euronext Paris. With its own manufacturing and R&D sites in Europe, North America and Asia and a global network of sales companies, Sartorius Stedim Biotech enjoys a worldwide presence. Its key membrane manufacturing and R&D site are located in Germany.

Sartorius Stedim Biotech enjoys a dedicated sales structure for OEM membranes to meet the customer needs. Whether located in Europe, America or Asia our sales representative can meet with you. Our experienced sales organization can help you to develop OEM solutions to your request and to define product specifications to your needs.





## Leading Membrane Technology

Sartorius Stedim Biotech produces a wide variety of microporous membranes that are especially designed, developed and manufactured to meet differing needs of biotechnology and medical device industries. Backed up by many decades of experience in membrane manufacturing and by utilizing the most state-of-the-art production equipment on the market, we guarantee excellent performance, consistent quality and a reliable supply of our membrane products.

All Sartorius Stedim Biotech membranes are manufactured in our premier production facilities in Göttingen, Germany. With our advanced casting technologies and strict environmental controls in place we ensure a stable and clean manufacturing process as well as a strong security of supply. Consistent process performance is secured by stringent qualification of all components and assemblies, manufacturing processes and personnel. Visits to our manufacturing facilities may be arranged on request.



## Quality Assurance from Sartorius

All Sartorius Stedim Biotech membranes are manufactured according to an ISO 9001 certified Quality Management System. Consistent high quality of Sartorius Stedim Biotech membranes is assured by careful selection of the raw materials and well-planned and validated production technologies, all of which results in high lot-to-lot reproducibility. The test procedures used are based both on external standard methods, such as the USP, EP and ASTM, and on in-house methods which are

the result of Sartorius' experience over the past 60 years. A lot is not released until all inprocess and final quality control data are available and fulfilled.

Sartorius Stedim Biotech's quality philosophy ensures that our customers receive products and services that meet the strictest quality requirements. The entire range of our membrane products is supplied with Certificates of Analysis.

## Broad Range of Applications

Microporous polyethersulfone (PES) membranes are available with a multitude of different pore sizes and performance characteristics making them ideally suited for a wide variety of applications including:

- Sterilizing Grade Filtration
- Mycoplasma Retentive Filtration
- Bioburden Reduction
- Prefiltration
- Clarification
- Venting & Gas Filtration

## Complete Traceability

All membrane products of Sartorius Stedim Biotech feature a comprehensive and easy to read out labeling which includes:

- Filter Type
- Order Code
- Lot Number
- Pore Size

The traceable lot number allows convenient retrieval of all data compiled on the materials used, production steps and QC tests.

# PES Membrane Portfolio – Quality Made in Germany

## Removal of Particles and Microorganisms from Liquids and Gases

Membrane processes are one of the most effective separation processes and they are steadily under development leading to new prospects of their applications. Sartorius Stedim Biotech polyethersulfone (PES) membranes are available in a wide variety of different pore sizes and structures as well as surface properties to serve nearly unlimited selectivity of separation. Their superior intra and inter lot consistency guarantees for reliable results. Further outstanding features like their excellent gamma compatibility and high mechanical and thermal resistance make Sartorius Stedim Biotech PES membranes the first choice for all major liquid and gas filtration applications, including medical devices.



The picture shows the wetting abilities of a hydrophilic, hydrophobic and oleophobic PES membrane (from right to left). The hydrophilic PES membrane wets out quickly and completely resulting in excellent flow rates and high throughputs. The hydrophobic PES membrane reliably repels water and other high surface tension fluids. The oleophobic PES membrane even resists wet-out with low surface tension fluids like oil or organic solvents assuring high performance in demanding venting applications.

## Hydrophilic PES Membranes

### Clarification | Sterile Filtration

Hydrophilic PES membranes are ideal for clarification and sterilization of aqueous liquids laden with particles, e.g. for preparation of pharmaceuticals or infusion solutions. All microorganisms and particles are reliably removed, without any effects on the ingredients, due to adsorption or decomposition. For optimal results, our hydrophilic PES membranes provide high flow rates and lowest adsorption characteristics. Different structures from symmetric to highly asymmetric allow you to select the membrane with the best combination of selectivity, flux and total throughput for your individual OEM device.

## Hydrophobic and Oleophobic PES Membranes

### Sterile Venting | Medical Use

These hydrophobic and oleophobic PES membranes have been developed to combine exceptionally high air flow rates with reliable liquid repelling properties. The membrane serves as a barrier against contamination from particles, aerosols, microorganisms, and other undesirable substances. In addition, venting membranes efficiently equalize pressure fluctuations that can occur during manufacturing processes or within a product during normal use, thus leading to increased process and product security. These high efficiency membranes are ideally suited for fast air filtration or sterile venting while preventing any liquid passage or condensate blockage.

## Applications

	Hydrophilic PES membranes	Hydrophobic and oleophobic PES membranes
Medical Filtration   Devices	IV therapy Ophthalmic Drug preparation Clinical reagents	Infusion Sets IV filter Transducer protectors Catheters Tubing systems Insufflation filters
Life Sciences	Buffer, media and sera filtration	Vent caps Liquid stopper devices
Pharmaceutical   Biotechnology Industry	Cell separation Water treatment Sterile filtration	Sterile vents for containers, bioreactors, fermenters
Food & Beverage Industry	Clarification Concentration Sterile filtration	Liquid bottling Food packaging
Electronics Industry		Sensors Electric motors

## Custom Made Membranes and Dimensions

OEM membranes often need custom dimensions in order to suit individual customer application requirements. From the master rolls, stored immediately after production, smaller slit rolls, sheets or disks of different format can be generated. Based on technical feasibility and quantity, a customized membrane development is possible. Please contact us at [unisart@sartorius-stedim.com](mailto:unisart@sartorius-stedim.com) to get more information.

## Do You Need Assistance to Choose the Best Suitable Membrane?

Please do not hesitate to contact us at [unisart@sartorius-stedim.com](mailto:unisart@sartorius-stedim.com). Our team of experts and scientists will assist you with the choice of the right membrane for your OEM application.



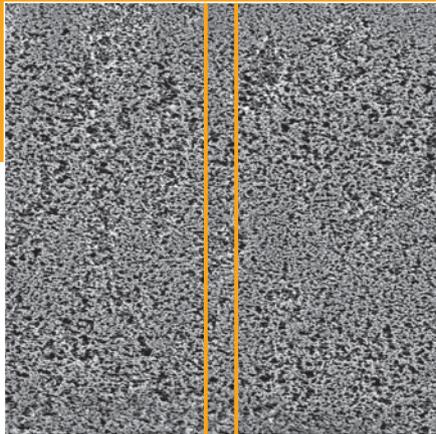
# Hydrophilic PES Membranes for Liquid Filtration

Hydrophilic PES membranes have a high internal porosity. They perform well at high flux with an excellent throughput of aqueous solutions over the entire pH range of 1–14. Their low level of extractables makes them suitable for environmental analysis. Thanks to their low non-specific protein binding the PES membranes are recommended for filtering biological and pharmaceutical solutions.

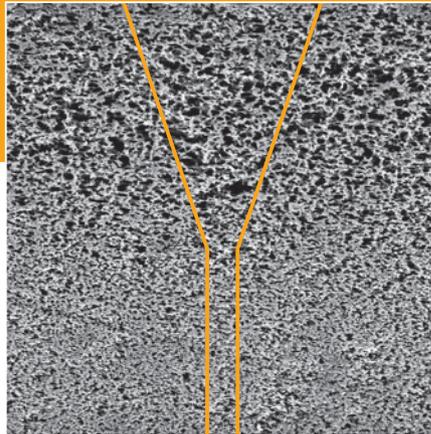
Hydrophilic PES membranes are available in different membrane structures. The symmetric membrane structure allows for high physical stress levels in the final product manufacturing, still ensuring the integrity of the membrane. The asymmetric membrane structure features higher flow rates in combination with relatively high physical stress levels in final device manufacturing. If total throughput is most important for the application, a highly asymmetric membrane is the membrane of choice.

Regardless of the membrane structure, 0.1–0.22  $\mu\text{m}$  rated PES membranes provide sterile filtration making the membranes well suited for biological sample preparation or sterile filtration of culture media and buffers.

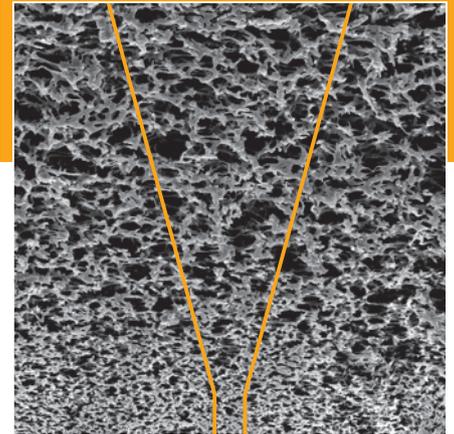




Symmetric PES membrane



Asymmetric PES membrane



Highly asymmetric PES membrane

## Typical Performance Characteristics

Material Type	Nominal Pore Size [μm]	Structure*	Water Flow Rate [ml/(cm <sup>2</sup> min bar)]	Bubble Point Sartochek® [bar]	Burst Pressure [bar]	Thickness [μm]	Retention (10 <sup>7</sup> /cm <sup>2</sup> Filter Area)
15442	5	AS	> 40**	0.3****	≥ 0.2	160	N/A
15402	3	AS	> 25**	0.5****	≥ 0.2	160	N/A
15413	1.2	AS	> 12**	0.8	≥ 0.3	160	N/A
15405	0.65	AS	> 80	1.5	≥ 0.4	160	N/A
15445	0.5	AS	90	2.3	≥ 0.4	160	N/A
15456	0.45	AS	> 40	2.4	≥ 0.5	160	N/A
15406	0.45	S	45	2.6	≥ 0.4	150	100% Serratia marcescens
15467	0.3	AS	> 23	> 2.9	≥ 0.4	150	N/A
15407	0.2	S	25	> 3.5	≥ 0.4	150	100% Brevundimonas diminuta
15407MI	0.2	AS	> 30	> 3.2	≥ 0.4	140	100% Brevundimonas diminuta
15427EP	0.22	HAS	> 34	> 3.55	≥ 0.4	160	100% Brevundimonas diminuta
15458	0.1	S	9	> 2.5****	≥ 0.5	150	100% Brevundimonas diminuta; LRV ≥ 7 Acholeplasma laidlawii

\* S = symmetric, AS = asymmetric, HAS = highly asymmetric

\*\* air flow rate [l/(m<sup>2</sup> s) at 200 Pa]

\*\*\* with isopropyl alcohol | water 60 vol% | 40 vol%

\*\*\*\* visual

### Structure

- Symmetric, asymmetric, highly asymmetric

### Adsorption, Non-specific

- ~ 10 μg/cm<sup>2</sup> for γ-globulin
- < 8 μg/cm<sup>2</sup> for BSA

### pH Stability

- Resistant to aqueous solutions pH 1–14

### Thermal Resistance

- 200 °C max.

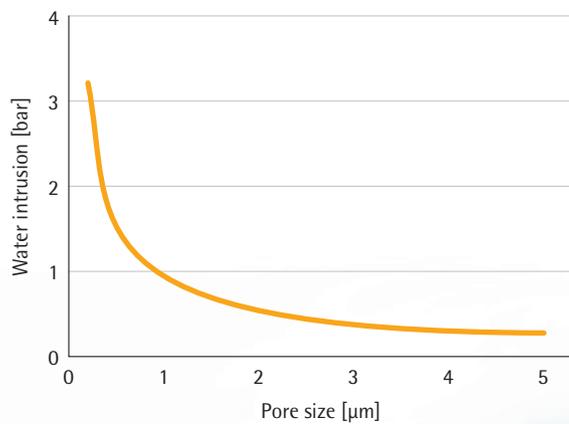
### Biocompatibility According to USP Standard

- Passes USP Biological Tests Plastic Class VI for biocompatibility and cytotoxicity
- Tests are partially performed on final devices

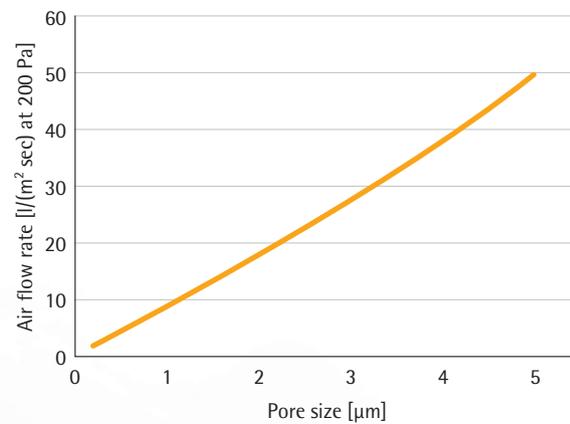
# Hydrophobic and Oleophobic PES Membranes for Venting & Gas Filtration

These PES membranes are treated to make their hydrophobic or oleophobic surfaces perfectly liquid repelling. Even under high humidity or high moisture, outstanding air permeability is guaranteed to assure rapid venting at low differential pressure. Due to their high mechanical stability, the membranes can easily be welded and integrated into a wide variety of devices. An additional specific feature of these membranes is their excellent gamma compatibility. In contrast to other common polymeric matrices for venting and gas filtration, such as PTFE, Sartorius Stedim Biotech PES membranes can be sterilized by gamma irradiation making them also best suited for the most demanding applications like medical devices.

### Water Intrusion Pressure



### Air Flow Rate





## Typical Performance Characteristics

Material Type	Surface Property	Nominal Pore Size* [μm]	Thickness [μm]	Air Flow Rate [l/(m <sup>2</sup> sec) at 200 Pa]	Water Intrusion Pressure		IPA** Wettability
					[bar]	[psi]	
15C77	hydrophobic	0.2	135	1.4	2.8	41	+
15D07MI	oleophobic	0.2	135	1.4	2.8	41	-
15D13	oleophobic	1.2	160	14	0.5	7.3	-
15D02	oleophobic	3.0	160	25	0.4	5.8	-
15D42	oleophobic	5.0	160	50	0.3	4.4	-

\* Additional pore sizes are available, please contact us at [unisart@sartorius-stedim.com](mailto:unisart@sartorius-stedim.com)

\*\* isopropyl alcohol

### Thermal Resistance

– 200 °C max.

### Biocompatibility According to USP Standard

– Passes USP Biological Tests Plastic Class VI for biocompatibility and cytotoxicity

# Technical Information

## Sterilization Methods

All Sartorius Stedim Biotech PES membrane products show excellent gamma compatibility and high thermal resistance giving you the freedom to choose the sterilization method that best meets the requirements of your device:

- Gamma irradiation ( $\leq 50$  kGy)
- Autoclaving at 121 °C or 134 °C
- EtO sterilization

## Membrane Sealing Methods

Due to their good tensile strength and high thermal resistance, all PES membrane products can easily be sealed onto plastic support or housing using most common welding methods like heat, radio frequency, ultrasonic and adhesives.



## Storage and Handling

By respecting the following recommendations you will ensure that your membrane products will always be at their optimum condition before use.

- The membranes are stored in their original package until use.
- Do not expose the membrane to direct sunlight or chemical vapors and keep it away from sources of heat.
- Constant temperatures between 15–25 °C and a relative humidity of max. 70% are best for storage and handling of the membrane. Very dry or humid storage conditions might alter the wettability and handling properties of the membrane. This change is only temporarily and can be overcome by conditioning the membrane at relative humidity between 40–70% for approximately 12 hours before processing the membrane.
- After unpacking the membrane, please avoid any direct contact of the membrane to materials that have the potential to release chemicals or additives. Adsorption or absorption of such substances could affect essential membrane properties like wettability. For the same reason, any direct contact with the membrane should be avoided.

## Would You Like to Use Membranes with Other Polymeric Matrices?

Microporous membranes can be composed of various polymers that differ from one another in their chemical and physical properties. Together with the characteristics of the filter pores these polymer properties govern the results in many filtration applications. The table below provides an overview which other polymeric matrices are available at Sartorius Stedim Biotech in addition to our PES portfolio.

Polymer	Features	Typical Applications
Cellulose acetate (CA)	high flow rates thermal stability very low non-specific adsorption	protein filtration, biological and clinical analysis, sterility tests
Surfactant free cellulose acetate (SFCA)	excellent wettability very low non-specific adsorption low content of extractables	removal of particles and microorganisms from aqueous solutions, sterile filtration
Cellulose nitrate (CN)	very high protein and DNA binding	cell retention, buffer filtration, microbiological testing
Regenerated cellulose (RC)	strong chemical resistance low protein binding	particle removal from organic and aqueous media, ultracleaning of solutions for HPLC
Polyamide (PA)	chemically resistant to alkaline solutions and organic solvents	particle removal in solutions for HPLC, filtration of protein-free culture media

Please contact us at [unisart@sartorius-stedim.com](mailto:unisart@sartorius-stedim.com) for a complete listing of our membrane products.

# Sales and Service Contacts

For further contacts, visit [www.sartorius-stedim.com](http://www.sartorius-stedim.com)

## Europe

**Germany**  
Sartorius Stedim Biotech GmbH  
August-Spindler-Strasse 11  
37079 Goettingen  
Phone +49.551.308.0

Sartorius Stedim Systems GmbH  
Robert-Bosch-Strasse 5-7  
34302 Guxhagen  
Phone +49.5665.407.0

**France**  
Sartorius Stedim FMT S.A.S.  
ZI des Paluds  
Avenue de Jouques - CS 91051  
13781 Aubagne Cedex  
Phone +33.442.845600

Sartorius Stedim France SAS  
ZI des Paluds  
Avenue de Jouques - CS 71058  
13781 Aubagne Cedex  
Phone +33.442.845600

**Austria**  
Sartorius Stedim Austria GmbH  
Modecenterstrasse 22  
1030 Vienna  
Phone +43.1.7965763.18

**Belgium**  
Sartorius Stedim Belgium N.V.  
Rue Colonel Bourg 105  
1030 Bruxelles  
Phone +32.2.756.06.80

**Hungary**  
Sartorius Stedim Hungária Kft.  
Kagyló u. 5  
2092 Budakeszi  
Phone +36.23.457.227

**Italy**  
Sartorius Stedim Italy S.r.l.  
Via dell'Antella, 76/A  
50012 Antella-Bagno a Ripoli (FI)  
Phone +39.055.63.40.41

**Netherlands**  
Sartorius Stedim Netherlands B.V.  
Phone +31.30.60.25.080  
[filtratie.nederland@sartorius-stedim.com](mailto:filtratie.nederland@sartorius-stedim.com)

**Poland**  
Sartorius Stedim Poland Sp. z o.o.  
ul. Wrzesinska 70  
62-025 Kostrzyn  
Phone +48.61.647.38.40

**Russian Federation**  
LLC "Sartorius Stedim RUS"  
Vasilyevsky Island  
5<sup>th</sup> line 70, Lit. A  
199178 St. Petersburg  
Phone +7.812.327.53.27

**Spain**  
Sartorius Stedim Spain, S.A.U.  
Avda. de la Industria, 32  
Edificio PAYMA  
28108 Alcobendas (Madrid)  
Phone +34.913.586.098

**Switzerland**  
Sartorius Stedim Switzerland AG  
Ringstrasse 24 a  
8317 Tagelswangen  
Phone +41.52.354.36.36

**U.K.**  
Sartorius Stedim UK Ltd.  
Longmead Business Centre  
Blenheim Road, Epsom  
Surrey KT19 9 QQ  
Phone +44.1372.737159

**Ukraine**  
LLC "Sartorius Stedim RUS"  
Post Box 440 "B"  
01001 Kiev, Ukraine  
Phone +380.44.411.4918

## Americas

**USA**  
Sartorius Stedim North America Inc.  
5 Orville Drive, Suite 200  
Bohemia, NY 11716  
Toll-Free +1.800.368.7178

**Argentina**  
Sartorius Argentina S.A.  
Int. A. Ávalos 4251  
B1605ECS Munro  
Buenos Aires  
Phone +54.11.4721.0505

**Brazil**  
Sartorius do Brasil Ltda  
Avenida Senador Vergueiro 2962  
São Bernardo do Campo  
CEP 09600-000 - SP- Brasil  
Phone +55.11.4362.8900

**Mexico**  
Sartorius de México, S.A. de C.V.  
Libramiento Norte de Tepetzotlan s/n,  
Colonia Barrio Tlacateco,  
Municipio de Tepetzotlan,  
Estado de México,  
C.P. 54605  
Phone +52.55.5562.1102  
[leadsmex@sartorius.com](mailto:leadsmex@sartorius.com)

**Peru**  
Sartorius Peru S.A.C.  
Avenue Alberto del Campo 411  
Floor 12 - The Office  
15076 - San Isidro, Lima  
Phone +51.1.441 0158

## Asia | Pacific

**Australia**  
Sartorius Stedim Australia Pty. Ltd.  
Unit 5, 7-11 Rodeo Drive  
Dandenong South Vic 3175  
Phone +61.3.8762.1800

**China**  
Sartorius Stedim Biotech (Beijing) Co. Ltd.  
No. 33 Yu'an Road  
Airport Industrial Park Zone B  
Shunyi District, Beijing 101300  
Phone +86.10.80426516

Sartorius Stedim (Shanghai)  
Trading Co., Ltd.  
3rd Floor, North Wing, Tower 1  
No. 4560 Jinke Road  
Zhangjiang Hi-Tech Park  
Pudong District  
Shanghai 201210, P.R. China  
Phone +86.21.6878.2300

Sartorius Stedim Biotech (Beijing) Co. Ltd.  
Guangzhou Representative Office  
Unit K, Building 23  
Huihua Commerce & Trade Building  
No. 80 Xianlie Middle Road  
Guangzhou 510070  
Phone +86.20.37618687 | 37618651

**India**  
Sartorius Stedim India Pvt. Ltd.  
#69/2-69/3, NH 48, Jakkasandra  
Nelamangala Tq  
562 123 Bangalore, India  
Phone +91.80.4350.5250

**Japan**  
Sartorius Stedim Japan K.K.  
4th Fl., Daiwa Shinagawa North Bldg.  
8-11, Kita-Shinagawa 1-chome  
Shinagawa-ku, Tokyo, 140-0001 Japan  
Phone +81.3.4331.4300

**Malaysia**  
Sartorius Stedim Malaysia Sdn. Bhd.  
Lot L3-E-3B, Enterprise 4  
Technology Park Malaysia  
Bukit Jalil  
57000 Kuala Lumpur, Malaysia  
Phone +60.3.8996.0622

**Singapore**  
Sartorius Stedim Singapore Pte. Ltd.  
10 Science Park Rd  
The Alpha #02-13/14  
Singapore Science Park II  
Singapore 117684  
Phone +65.6872.3966

**South Korea**  
Sartorius Korea Biotech Co., Ltd.  
8th Floor, Solid Space B/D,  
PanGyoYeok-Ro 220, Bundang-Gu  
SeongNam-Si, GyeongGi-Do, 463-400  
Phone +82.31.622.5700



▶ [www.sartorius-stedim.com](http://www.sartorius-stedim.com)