



OEM-Products and Components

Weighing | Filtration | Final Filling | Pipetting | Connection and Disconnection |
Tubing | Aseptic Fluid Management | Fittings and Sensors

Simplifying Progress

SARTORIUS

<u>Introduction</u>	Weighing	Filtration	Final Filling	Tubing
Pipetting	Connects Disconnects	Aseptic Fluid Management	Contact	

OEM-Products

Customized Solutions for Your Specific Requirements

Discover the Sartorius portfolio of OEM products that offer you the flexibility to find the customized solution tailored to your individual requirements.

We offer an extensive portfolio of advanced filtration devices, membranes, weigh cells, and dispensing solutions, as well as connecting and disconnecting technologies.

Our experts are happy to support you in finding the most suitable solution to meet your specific needs.

By combining our technology leadership and manufacturing excellence with your unique requirements, we can help you find the performance and design characteristics ideally suited to your individual product requirement.

 **Find out more**

For more information, please visit
www.sartorius.com/en/products/oem

Introduction

Weighing

Filtration

Final Filling

Tubing

Pipetting

Connects | Disconnects

Aseptic Fluid Management

Contact

OEM Weigh Cells & High Resolution Load Cells

Usability Leading engineering and automation companies utilize our weigh cell components globally as a force or weighing sensor in production lines, industrial equipment or other types of machines



Benefits

- A monolithic weighing system in stainless steel housing
- Force compensation technology system with a resolution of 1 μg
- Easy to integrate, connect & configure
- Possibility for under floor weighing
- Fast warm-up by separation of weighing system and electronics
- Variety of dimensions for optimal fit where space is an issue
- Software customization configuration for specific applications
- Fully automated calibration and adjustment function for accurate results all the way
- IP protection
- Explosion protected weigh cell for use in hazardous areas
- Adjustable load receptor with overload protection
- RS232C interface for configuration and data transfer
- 10 years long term availability
- Global service and support

 Find out more

For more information, please visit

www.sartorius.com/en/products/oem/oem-weigh-cells

Introduction	<u>Weighing</u>	Filtration	Final Filling	Tubing
Pipetting	Connects Disconnects	Aseptic Fluid Management	Contact	

Basic OEM-Weigh Cell

The Basic series focus just on the essentials. Sartorius provides the mechanical and electronic solution. These weighing devices should be integrated by the system provider to overcome the influences of the ambient conditions.



Special solution OEM-Weigh Cells

The WZA25-NC has a capacity of 20 g with a readability of 10 µg combined with a small housing. A multiple weighing system can be arranged in linear fashion with a distance of 25 mm between the weighing pins. Head to Head configuration allow 12.5 mm track distance.



Explosion Protected IP44 OEM-Weigh Cells

Explosion protected weigh cell for use in hazardous areas

- Zone 1 and 2 with gasses of group IIC and temperature class T4
- Zone 21 and 22 for dust atmospheres ($T_{max} = 80^{\circ}C$)

High Precision OEM-Weigh Cells

High precision weigh cells with a readability of 1 µg for the weighing capacity of 20 g.

The weighing system has an internal calibration weight for easy checks in automated environments.



Introduction	Weighing	<u>Filtration</u>	Final Filling	Tubing
Pipetting	Connects Disconnects	Aseptic Fluid Management	Contact	

OEM Membranes

Microporous PES Membranes – High Efficiency Membranes for Filtration

Leading Membrane Technology

Sartorius 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 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.

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



Find out more

For more information, please visit
www.sartorius.com/en/products/oem/oem-membranes-and-devices

Introduction	Weighing	<u>Filtration</u>	Final Filling	Tubing
Pipetting	Connects Disconnects	Aseptic Fluid Management	Contact	

OEM Membranes

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 exceptional 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 PES membranes the first choice for all major liquid and gas filtration applications, including medical devices.

Hydrophilic PES Membranes

Custom Made Membranes and Dimensions

Hydrophobic and Oleophobic PES Membranes

Customized Membranes with Other Polymeric Matrices



Introduction	Weighing	<u>Filtration</u>	Final Filling	Tubing
Pipetting	Connects Disconnects	Aseptic Fluid Management	Contact	

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 low adsorption characteristics. Different structures from symmetric to highly asymmetric allow you to select the membrane with the combination of selectivity, flux, and total throughput ideally suited to your individual OEM device.



Introduction

Weighing

Filtration

Final Filling

Tubing

Pipetting

Connects | Disconnects

Aseptic Fluid Management

Contact

Hydrophobic and Oleophobic PES Membranes

Sterile Venting | Medical Use

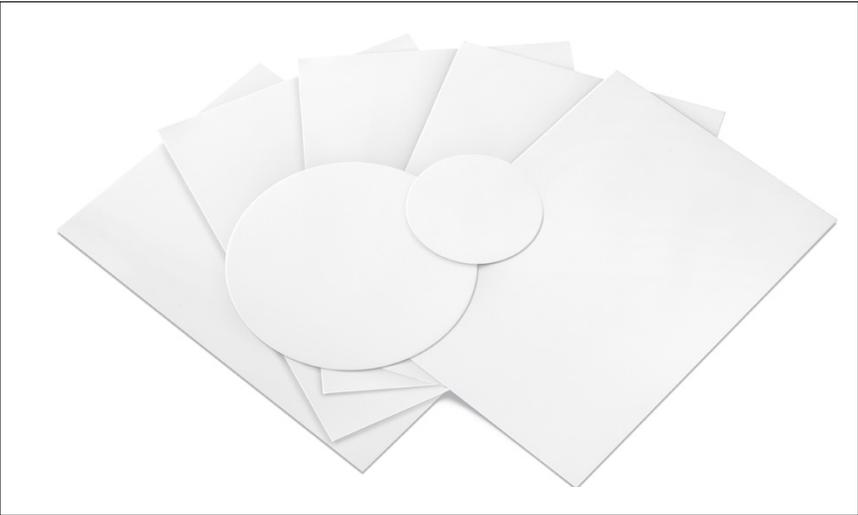
These hydrophobic and oleophobic PES membranes have been developed to combine exceptionally high air flow rates with reliable liquid repelling properties. The membrane serves as a barrier against contamination from particles, aerosols, microorganisms, and other undesirable substances. In addition, venting 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.



Introduction	Weighing	<u>Filtration</u>	Final Filling	Tubing
Pipetting	Connects Disconnects	Aseptic Fluid Management	Contact	

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.



Introduction	Weighing	<u>Filtration</u>	Final Filling	Tubing
Pipetting	Connects Disconnects	Aseptic Fluid Management	Contact	

Customized 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 provides an overview of different polymeric matrices that are available at Sartorius in addition to our PES portfolio.



Polymer	Features	Typical Applications
Cellulose acetate (CA)	<ul style="list-style-type: none"> high flow rates thermal stability very low non-specific adsorption 	<ul style="list-style-type: none"> protein filtration, biological and clinical analysis, sterility tests
Surfactant free cellulose acetate (SFCA)	<ul style="list-style-type: none"> excellent wettability very low non-specific adsorption low content of extractables 	<ul style="list-style-type: none"> removal of particles and microorganisms from aqueous solutions sterile filtration
Cellulose nitrate (CN)	<ul style="list-style-type: none"> very high protein and DNA binding cell retention 	<ul style="list-style-type: none"> buffer filtration, microbiological testing
Regenerated cellulose (RC)	<ul style="list-style-type: none"> strong chemical resistance low protein binding 	<ul style="list-style-type: none"> particle removal from organic and aqueous media ultracleaning of solutions for HPLC



Introduction

Weighing

Filtration

Final Filling

Tubing

Pipetting

Connects | Disconnects

Aseptic Fluid Management

Contact

Custom Made Syringe Filters

Sartorius Premium Filters for Removal of Particles and Microorganisms from Liquids and Gases

- Sartorius syringe filters made of a PP housing and membrane components featuring maximum chemical compatibility and minimum extractables to ensure excellent results.
- Sartorius filters made of acrylic housing provide high flow rates and low adsorption characteristics. This housing type could be color coded, e.g. for easy pore size identification.



Housing Material		γ -irradiable	EO-sterilizable
Acrylic-based Multipolymer	acrylic	yes	yes
Polypropylene	PP	no	yes
Polypropylene*	PP*	yes	yes

* γ -sterilizable PP is available on special request

Introduction	Weighing	<u>Filtration</u>	Final Filling	Tubing
Pipetting	Connects Disconnects	Aseptic Fluid Management	Contact	

Custom Made Syringe Filters – Membrane Material



Membrane Material		γ -irradiable	EO-sterilizable
Surfactant-free Cellulose Acetate	SFCA	yes	yes
Polyethersulfon, hydrophilic	PES	yes	yes
Polyethersulfon, hydrophobic	PES	yes	yes
Regenerated Cellulose	RC	yes	yes
Polyamide Nylon	PA	no	yes
Polytetrafluoroethylene	PTFE	no	yes
Glass Fiber	GF	yes	yes

Introduction	Weighing	<u>Filtration</u>	Final Filling	Tubing
Pipetting	Connects Disconnects	Aseptic Fluid Management	Contact	

OEM Filter Elements

Our filter elements with a multitude of different membrane and filter active fleece materials as well as various sterilization methods making them ideally suited for a broad range of applications include:

- Sterilizing Grade Filtration
- Bioburden Reduction
- Prefiltration and Clarification
- Air & Gas Filtration

Benefit from the Extensive Filtration Portfolio of Sartorius

Broad Range of Filter Materials

Flexible in Filter Designs & Connectors



Introduction	Weighing	<u>Filtration</u>	Final Filling	Tubing
Pipetting	Connects Disconnects	Aseptic Fluid Management	Contact	

Broad Range of Filter Material – Liquid Filtration

Polyethersulfone

Application:
 pH adjustment in Cell Culture |
 Fermentation Buffer Preparation |
 Form & Fill

Bioburden Reduction: 0.8 | 0.45 µm
 Sterilizing Grade: 0.45 | 0.2 µm

Filter Designs	Sterilization Methods	Delivered in sterile condition (autoclaved)
Mini Cartridges	In-line steamable or autoclavable	
Cartridges	In-line steamable or autoclavable	
T-Style Maxicaps®	Autoclavable or γ-irradiatable (hybrid version)	
Maxicaps®	Autoclavable	
	γ-irradiatable	
Midicaps®	Autoclavable	▪
	γ-irradiatable	
Capsule Size 4	Autoclavable	▪
	γ-irradiatable	
Capsule Size 5	Autoclavable	▪
	γ-irradiatable	



Introduction	Weighing	<u>Filtration</u>	Final Filling	Tubing
Pipetting	Connects Disconnects	Aseptic Fluid Management	Contact	

Broad Range of Filter Material – Liquid Filtration

Polypropylene

Application:

pH adjustment in Cell Culture |

Fermentation | Form & Fill | Solvents & Oily formulations

Retention Rate¹:

50 µm	3 µm
20 µm	1.2 µm
8 µm	0.65 µm
5 µm	0.45 µm

¹ Nominal

Filter Designs	Sterilization Methods
Mini Cartridges	In-line steamable or autoclavable
Cartridges	In-line steamable or autoclavable
T-Style Maxicaps®	Autoclavable
Midicaps®	Autoclavable
Capsule Size 4	Autoclavable
Capsule Size 5	Autoclavable



Introduction	Weighing	<u>Filtration</u>	Final Filling	Tubing
Pipetting	Connects Disconnects	Aseptic Fluid Management	Contact	

Broad Range of Filter Material – Liquid Filtration

Glass Fiber

Application:
Cell Removal | Clarification

Retention Rate¹:

- 1.2 µm
- 0.65 µm

¹ Nominal

Filter Designs	Sterilization Methods
Cartridges	In-line steamable or autoclavable
Maxicaps®	Autoclavable
Midicaps®	Autoclavable

PTFE

Application:
Sterile Filtration of Solvents

Sterilization Grade: 0.2 µm (LG)

Filter Designs	Sterilization Methods	Delivered in sterile condition (autoclaved)
Mini Cartridges	In-line steamable or autoclavable	
Junior	In-line steamable or autoclavable	
Cartridges	In-line steamable or autoclavable	
Maxicaps®	Autoclavable	▪
Midicaps®	Autoclavable	▪
Capsule Size 4	Autoclavable	▪
Capsule Size 5	Autoclavable	▪



Introduction	Weighing	<u>Filtration</u>	Final Filling	Tubing
Pipetting	Connects Disconnects	Aseptic Fluid Management	Contact	

Broad Range of Filter Material – Air & Gas Filtration

Hydrophobic Polyethersulfone

Application:

- Venting, inflating & leak testing of single-use bags | assemblies
- Venting of single-use Bioreactors | Fermentors
- Protection of integrity test devices

Sterilizing Grade: 0.2 µm

Filter Designs	Sterilization Methods
Maxicaps®	γ-irradiatable
Midicaps®	γ-irradiatable
Capsule Size 4	γ-irradiatable
Midisart®	γ-irradiatable or autoclavable

Hydrophobic Glass Fiber

Application:

Venting of tanks (non pressure resistant)

Bioburden Reduction | Particle Retention:
0.2 µm (LG)

Filter Designs	Sterilization Methods
Mini Cartridges	In-line steamable or autoclavable
Cartridges	In-line steamable or autoclavable



Introduction	Weighing	<u>Filtration</u>	Final Filling	Tubing
Pipetting	Connects Disconnects	Aseptic Fluid Management	Contact	

Broad Range of Filter Material – Air & Gas Filtration

PTFE

Application:

- Compressed Gases
- Venting of tanks (pressure resistant)
- Venting of production machinery
- Venting of stainless steel Bioreactors | Fermentors
- Protection of integrity test devices
- Venting of glass & plastic bottles

Sterilizing Grade: 0.2 µm

Filter Material (Membrane): PTFE

Filter Designs	Sterilization Methods	Delivered in sterile condition (autoclaved)
Mini Cartridges	In-line steamable or autoclavable	
Junior	In-line steamable or autoclavable	
Cartridges	In-line steamable or autoclavable	
Maxicaps®	Autoclavable	▪
Midicaps®	Autoclavable	▪
Capsule Size 4	Autoclavable	▪
Capsule Size 5	Autoclavable	▪
Midisart®	γ-irradiatable or autoclavable	

Application:

Venting of WFI loops (high temperature)

Sterilizing Grade: 0.2 µm

Filter Material (Membrane): PTFE

Filter Designs	Sterilization Methods
Cartridges	In-line steamable or autoclavable



Introduction	Weighing	<u>Filtration</u>	Final Filling	Tubing
Pipetting	Connects Disconnects	Aseptic Fluid Management	Contact	

Flexible in Filter Designs & Connectors

Cartridges



Adapter 25
2 Flange
Bayonet adapter with 226 double o-ring



Adapter 27
Bayonet adapter with 222 double o-ring



Adapter 28
3 Flange
Bayonet adapter with 222 double o-ring



Maxicaps®



Connector Type S
1½" Tri-Clamp
50.5 mm



Connector Type F
¾" Tri-Clamp
25 mm

Connector Type O
½" single stepped hose barb



T-Style Maxicaps®



Connector Type S
1½" Tri-Clamp
50.5 mm



Connector Type O
½" single stepped hose barb



Multi-use

Single-use



Introduction

Weighing

Filtration

Final Filling

Tubing

Pipetting

Connects | Disconnects

Aseptic Fluid Management

Contact

Flexible in Filter Designs & Connectors

Mini Cartridges



Adapter 15
Bayonet adapter
with o-ring



Adapter 18
Plug adapter
with double
o-ring



Midicaps®



**Connector
Type S**
1½" Tri-Clamp
50.5 mm



**Connector
Type F**
¾" Tri-Clamp
25 mm



**Connector
Type O**
½" single
stepped
hose barb



Capsules



**Connector
Type O**
¼" multiple
stepped hose
barb



**Connector
Type S**
¾" Tri-Clamp
25 mm



Multi-use

Single-use



Introduction

Weighing

Filtration

Final Filling

Tubing

Pipetting

Connects | Disconnects

Aseptic Fluid Management

Contact

Flexible in Filter Designs & Connectors

Midicaps®



**Standard Stepped
Hose Barb**
6-12 mm



**Small Stepped
Hose Barb**
4-12 mm



Thread
8" NPT



Tri-Clamp
3/4" Tri-Clamp
25 mm



Multi-use

Single-use



Introduction

Weighing

Filtration

Final Filling

Tubing

Pipetting

Connects | Disconnects

Aseptic Fluid Management

Contact

Midisart® 2000

Midisart® 2000 is a 20 cm² small-scale sterile air | gas PTFE disposable vent filter. Available in 0.2 µm or 0.45 µm, sterile or non-sterile.



Introduction

Weighing

Filtration

Final Filling

Tubing

Pipetting

Connects | Disconnects

Aseptic Fluid Management

Contact

Sartopore® Air Midisart®

The Sartopore® Air Midisart® with its revolutionary hydrophobic 0.2 µm PES membrane completely addresses the rigorous demands of small-scale, single-use, venting applications. Sartopore® Air Midisart is a Gamma-Irradiatable or Autoclavable Single-Use Venting Filter available in 20 cm².



Introduction

Weighing

Filtration

Final Filling

Tubing

Pipetting

Connects | Disconnects

Aseptic Fluid Management

Contact

Sartolab[®] P20

Sartorius Sartolab[®] P20 pressure sterile filtration devices represent an eco-friendly approach to aqueous filtration that offers a faster and more adaptable sterile filtration for 100 mL to 5 L sample volumes.

Benefit from the combination of a large filtration area (20 cm²) and the optimal design of the filter support guarantees high flow rates with high throughputs and automatic venting to ensure that the entire filter surface is used for effective filtration.



Housing Material		EO-sterilizable
Polycarbonate	PC	yes

Membrane Material		EO-sterilizable
Surfactant-free Cellulose Acetate	SFCA	yes
Polyethersulfon, hydrophilic	PES	yes

Introduction

Weighing

Filtration

Final Filling

Tubing

Pipetting

Connects | Disconnects

Aseptic Fluid Management

Contact

Minisart®

Minisart® syringe filters are ideal for sterilization and removal of particles from air and other gases, and are optimal for sterile venting of containers, bioreactors, fermenters and tubing systems in medical devices.



Introduction

Weighing

Filtration

Final Filling

Tubing

Pipetting

Connects | Disconnects

Aseptic Fluid Management

Contact

Final Filling

Biosafe® Solutions

The Biosafe® range of Aseptic Transfer Systems offers reliable and easy-to-use solutions for the secure transfer of components, fluids and powders while maintaining the integrity of the critical area in isolators, RABS and cleanrooms.

Octoplus FF® Single Use Bags

Octoplus FF® is a pre-sterilized, ready to use, single-use final filling set that can be configured with a large variety of components and connectors.



 Find out more

For more information, please visit
www.sartorius.com/en/products/fluid-management/final-filling

Aseptic Transfer Systems

The Biosafe® system consists in a hardware (alpha-port) combined with different types of consumables (beta-port) depending on the applications to be covered:

Application	Component transfer (stoppers, caps, plungers...)	Removal of components, waste or petri dishes	Fluid transfer
Solution	Biosafe® bags <ul style="list-style-type: none"> Gamma sterilizable or autoclavable Sent to component manufacturers or directly to the end-users 	Biosafe® bags gamma sterile	Rapid Aseptic Fluid Transfer bags (RAFT) <ul style="list-style-type: none"> Use the Biosafe® connection to transfer a single use tubing
			



Find out more

For more information, please visit

www.sartorius.com/en/products/fluid-management/final-filling



Introduction

Weighing

Filtration

Final Filling

Tubing

Pipetting

Connects | Disconnects

Aseptic Fluid Management

Contact

Final Filling Set

Octoplus FF[®] provides a completely single-use alternative to traditional stainless steel “break tank” or “buffer tank” systems in a wide range of applications. The systems are designed using 100% single-use components from the bulk vessel aseptic connection all the way to the final filling needle. Octoplus FF[®] can be designed to work with different types of dosing systems, and in any type of filling environment (traditional cleanrooms, RABS, isolators).

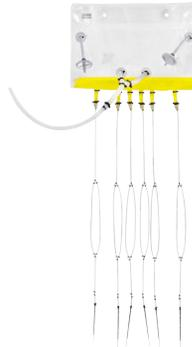
Application

Drug product repartition

Solution

Octoplus[®] Final Filling single use bag
Transferred to the filling line:

- through the Biosafe[®] Port
- manually prior VHP decontamination



Introduction	Weighing	Filtration	Final Filling	<u>Tubing</u>
Pipetting	Connects Disconnects	Aseptic Fluid Management	Contact	

Tuflux[®] Tubing

Benefit from the best of two worlds

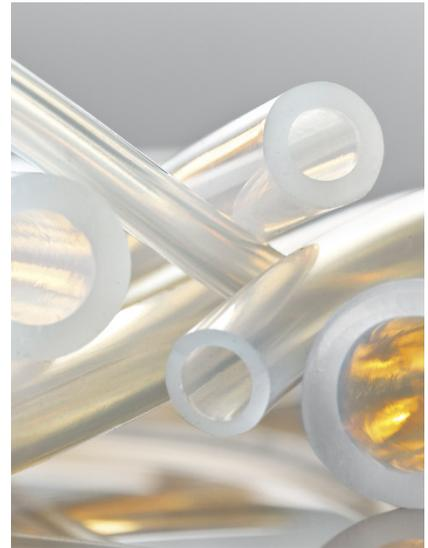
Liquid Transfer is an essential part of bio-manufacturing processes. For every transfer between different single-use, multi-use or hybrid systems, Tuflux[®] is the right solution. With Tuflux[®] Tubing, you benefit from extremely pure Silicone and TPE tubing for the transfer of your critical fluids.

Both tubing types have excellent extractable | profiles and guarantee that the cleanliness of the drug substances they carry is maintained across all process steps.

A broad range of standards

Tuflux[®] SIL & TPE are available in 7 dimensions from 1/8" (3.2 mm) to 3/4" (19.1 mm) for the internal dimensions and a wall thickness from 1.6 to 4.8 mm.

We ship from stock to ensure some of the most competitive lead times on the market. Our robust change control process provides reliable communication channels on any product changes.



Introduction	Weighing	Filtration	Final Filling	<u>Tubing</u>
Pipetting	Connects Disconnects	Aseptic Fluid Management	Contact	

Tuflux[®] SIL – Platinum Cured Silicone Tubing

Consistent Purity, Robustness and additional Ease of Use

Tuflux[®] SIL is designed to facilitate fluid transfer in the biopharmaceutical manufacturing processes. It is a highly resistant platinum cured silicone tubing. Tuflux[®] SIL is manufactured by Raumedic in ISO7 clean room and undergoes an extensive post-curing treatment after extrusion to obtain an extremely low level of siloxanes. With its 60 Shore A hardness Tuflux[®] SIL ensures solid compromise between pressure resistance, long pumping life time, high flow rates while still maintaining high flexibility of the tube without kinking.

Plasma coating (Low-Tack treatment) provides a less sticky outer surface of the Tuflux[®] SIL tubing in comparison with common non-coated silicone tubing products. This eases the unwinding of coils as well as handling with gloves in general. We ship from stock to ensure some of the most competitive lead times on the market. Our robust change control process provides reliable communication channels on any product changes.



- Excellent Pumping Performance
- Easy Handling
- Outstanding Purity



Introduction

Weighing

Filtration

Final Filling

Tubing

Pipetting

Connects | Disconnects

Aseptic Fluid Management

Contact

Tuflux[®] TPE Thermoplastic Elastomer Tubing

Comfortable Handling, Better Cell Growth and More Flexibility

Tuflux[®] TPE is the new tubing within the Sartorius Stedim Biotech portfolio made for liquid transfer in the biopharmaceutical industry.

It can be heat sealed and welded with Biosealer[®] TC & Biowelder[®] TC to maintain sterile conditions also with other TPE tubing formulations like C-Flex or Advata-Flex. Unique color-coding per Dimension, aligned with Biowelder[®] TC tube holders, minimizes the risk of operator errors.

Due to the absence of the Antioxidant Irgafos 168, there is no inhibition of Cell Growth even under worst case conditions.

- Ease of Use | Error Proving
- High flexibility for use with other materials
- Optimized for Cell Growth Results



Introduction

Weighing

Filtration

Final Filling

Tubing

Pipetting

Connects | Disconnects

Aseptic Fluid Management

Contact

Pipetting and Dispensing

Customizeable components and products for accurate, precise and reliable liquid handling tailored for your application.

Rline[®] Robotic Dispenser Module

Picus[®] Electronic Pipette

Roboline

Optifit and Safetyspace[™] Filter Tips

Mline[®] & Proline[®] Plus
Mechanical Pipettes

Automation Tips

Midi Plus[™] Pipetting Controller



Introduction

Weighing

Filtration

Final Filling

Tubing

Pipetting

Connects | Disconnects

Aseptic Fluid Management

Contact

Rline[®] Robotic Dispenser Module

The Sartorius Rline[®] Robotic Dispenser Module brings the proven technology of Sartorius electronic pipettes into the field of laboratory automation.

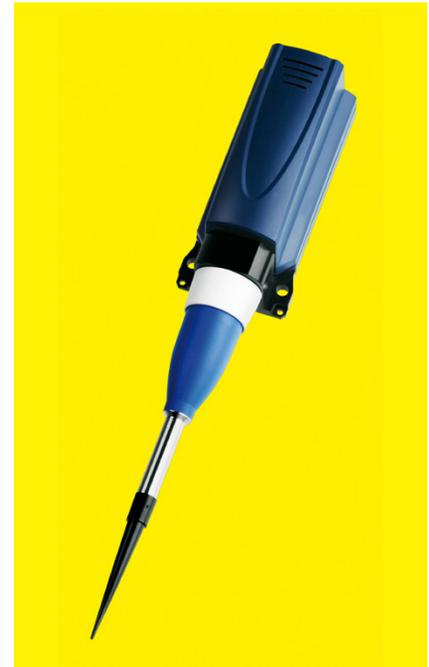
The unique and innovative features of the module ensure the accuracy and performance of sample processors and automated instrumentation:

- Outstanding accuracy and reliability with the enhanced DC-motor concept, electronic brake system and optical sensor
- Combination of Liquid Level Sensor with conductive robotic tips enables the detection of the liquid surface in the target vessel. This prevents erroneous attempts to aspirate liquid before the liquid surface has been reached.

Typical Applications

- Laboratory automation for research and clinical applications
- Laboratory robotics and automated workflows
- Industrial applications and production systems

The Rline[®] can be used as an off-the-shelf module or tailored to fit a specific application. To help you to get started with Rline[®], we also offer development kits!



Find out more

For more information, please visit
www.sartorius.com/en/products/oem/dispensing

Introduction

Weighing

Filtration

Final Filling

Tubing

Pipetting

Connects | Disconnects

Aseptic Fluid Management

Contact

Roboline

The Sartorius Roboline is the ideal companion to automate routine pipetting work. This small, quiet and compact liquid handling automate is easy to integrate as a part of a workflow.

Open programmability makes the Roboline easy to use and suitable for a wide range of applications. With this companion you can simply automate a single work stage or pipetting steps for a whole application.

The robotic Rline® dispensing module guarantees that your results are accurate and precise without human error.



Introduction

Weighing

Filtration

Final Filling

Tubing

Pipetting

Connects | Disconnects

Aseptic Fluid Management

Contact

Mline® & Proline® Plus Mechanical Pipettes

Mline®

Sartorius' mechanical pipette family – the Mline® – offers excellent ergonomics, performance and reliability in manual pipetting. It is designed to maintain high accuracy and precision in repetitive and long-lasting manual pipetting.

Excellent ergonomics and low pipetting forces protect laboratory workers from injury. Mline® has exceptionally light pipetting and tip ejection forces due to its patented spring mechanism.

Mline® covers the full volume range of 0.1 µl to 10 ml and is available in single- and multichannel models.

Proline® Plus

The Proline® plus mechanical pipette family is designed to offer comfort and quality for your everyday manual pipetting. It combines durable construction with ease and lightness of use, and is therefore the perfect choice for liquid-handling professionals and students alike. In addition, it has an exceptionally wide pipette range, including fixed volume pipettes for when volumes need to be ready-set to avoid errors.

Proline® plus has low pipetting forces, a comfortable handle and an ergonomic finger support for effortless pipetting. Good fit in hand minimizes the grip force needed to hold the pipette, thereby reducing the risk of strain injury.

Proline® plus covers the full volume range of 0.1 µl to 10 ml and is available in single- and multi-channel models, as well as fixed volume single channel models.



Find out more

For more information, please visit

www.sartorius.com/en/products/pipetting-and-dispensing



Introduction

Weighing

Filtration

Final Filling

Tubing

Pipetting

Connects | Disconnects

Aseptic Fluid Management

Contact

Midi Plus™ Pipetting Controller

The Midi Plus™ is a lightweight electronic cordless pipetting controller, which allows aspiration from bottles and tubes, without the arm and hand elevations required in the case of serological or volumetric pipettes.

It fits all commonly used 1 – 100 ml glass or plastic pipettes, but can also be used with Sartorius 5 ml and 10 ml disposable tips. The speed can be fine-tuned by applying varying finger pressure to the operating buttons. The Midi Plus™ is ideal, for example, in microbiological work: dispensing into a culture media dish can be performed carefully, drop by drop, without breaking the fine surface of the medium.

Midi Plus™ can be customized to fit to your brand including the pipette controller and packaging.



Find out more

For more information, please visit

www.sartorius.com/en/products/pipetting-and-dispensing



Introduction

Weighing

Filtration

Final Filling

Tubing

Pipetting

Connects | Disconnects

Aseptic Fluid Management

Contact

Picus[®] Electronic Pipette

Excellent Ergonomics

The uniquely lightweight and streamlined design of the Picus[®] pipette ensures an effortless pipetting experience. The pipette rests lightly in the user's hand due to the comfortable handle and finger hook, and minimal gripping force is needed. The conveniently located soft-touch operating buttons, and unique electronic tip ejection, help minimize muscle strain, further enhancing the ergonomics of the pipette.

Reliable Results

The electronic brake and piston control system guarantee accurate and precise pipetting results, independent of the user. Using the unique plate tracker increases reliability in microwell plate dispensing, by guiding the steps to pipette into the correct microplate wells.

Fast Execution of Pipetting Tasks

The unique adjustment wheel of Picus[®] pipettes offers extremely fast volume setting and menu navigation. The user can perform pipetting tasks quickly and easily with the extensive range of pipetting modes, from diluting and titrating to serial dispensing. The pipetting functions are quick to learn, using the intuitive user interface, available in a choice of language options: English, French, German, Russian and Chinese.

Picus[®] is available in single-channel models, covering a volume range of 0.2 – 10,000 μ l and in multi-channel models from 0.2 μ l to 1,200 μ l.

Picus software and visual details can be modified to fit your brand and application.



 Find out more

For more information, please visit
www.sartorius.com/en/products/pipetting-and-dispensing

Introduction

Weighing

Filtration

Final Filling

Tubing

Pipetting

Connects | Disconnects

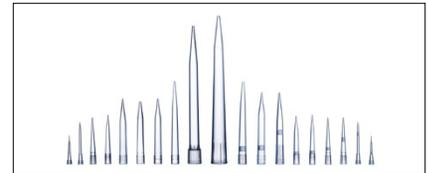
Aseptic Fluid Management

Contact

Optifit and Safetyspace™ Filter Tips

The pipette is a precision measuring instrument which creates a system with the pipette tip. Sartorius pipettes tips complete any OEM or private label pipetting solution with wide range of options to choose from. Safetyspace™ Filter tips bring peace of mind to any application and prevent cross-contamination and pipette contamination. Optifit tips are natural decision for most liquid handling applications. Low Retention tips enhance sample recovery when handling difficult solutions.

- Excellent to OEM and PL electronic and mechanical pipettes or automation dispensers from Sartorius
 - Picus®
 - Mline®, Proline Plus®, Proline®
 - Rline®
- Complies to stringent purity standards and has excellent chemical resistance
- Safetyspace™ filter tips are available from 10 to 1,200 µl volumes
- Optifit tips are available from 10 to 10,000 µl volumes
- Optifit tips are available in single tray, refill tower, refill pack and in Flexibulk format.



Find out more

For more information, please visit
www.sartorius.com/en/products/pipetting-and-dispensing



Introduction

Weighing

Filtration

Final Filling

Tubing

Pipetting

Connects | Disconnects

Aseptic Fluid Management

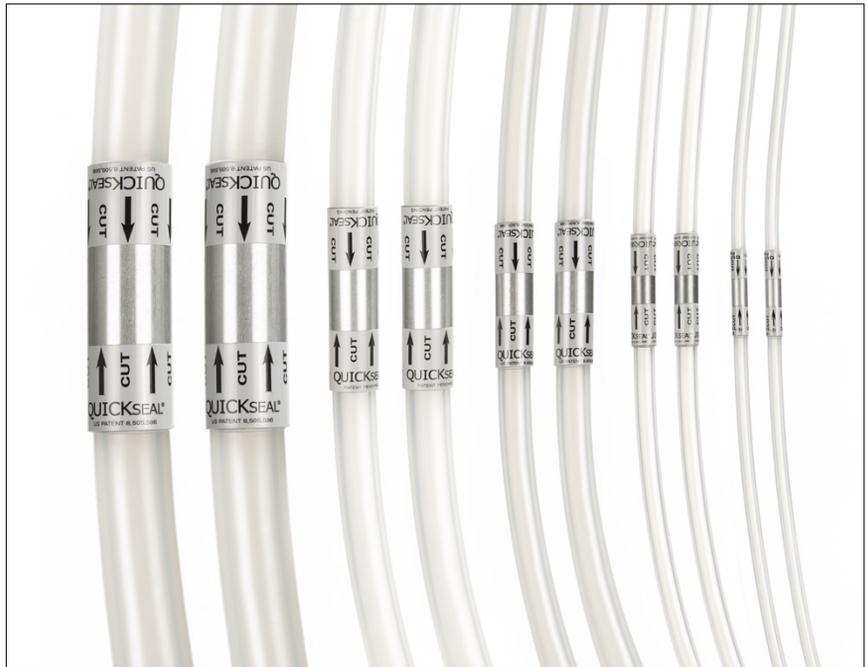
Contact

Quickseal®

Installed on an array of tube materials including platinum-cured silicone, C-Flex® and other thermoplastic elastomers, Quickseal® is an exceptionally quick aseptic tube separation technique. Extensively tested, Quickseal® is proven aseptic and stronger than the tubing it is part of.

Quickseal® can be included as a component on single-use tube, bag or bottle assemblies, sampling manifolds, single-use bioreactors and more.

Quickseal® reduces operating time and improves reliability for biopharmaceutical and other high-purity fluid handling industries.



Introduction

Weighing

Filtration

Final Filling

Tubing

Pipetting

Connects | Disconnects

Aseptic Fluid Management

Contact

Opta[®] SFT

The Opta[®] SFT Sterile Connector is a single use device that allows a sterile connection between two separate, pre-sterilized components in biopharmaceutical manufacturing processes. Opta[®] SFT Sterile Connectors are quick and easy to use, and are backed by extensive validation work and 100% in house integrity testing.



Introduction

Weighing

Filtration

Final Filling

Tubing

Pipetting

Connects | Disconnects

Aseptic Fluid Management

Contact

Clipster® Aseptic Disconnecter

The Clipster® aseptic disconnecter is a single use device that performs aseptic disconnections of tubing. The Clipster® aseptic disconnecter allows aseptic disconnection in non classified and classified environments while maintaining product sterility. It can be applied to multiple types and sizes of tubing. The Clipster® aseptic disconnecter may be sold as a stand alone product or preassembled on our Fluid Management bag assemblies.



Introduction

Weighing

Filtration

Final Filling

Tubing

Pipetting

Connects | Disconnects

Aseptic Fluid Management

Contact

Celsius® EVA Single Use Bags

Celsius® EVA chambers are specifically designed for controlled freezing, thawing and long term frozen storage of biopharmaceuticals such as process intermediates, bulk drug substances, product in clinical phases (pre, I, II +), vaccines.

The mechanical properties of the S71 film have been characterized between -80 °C and +45 °C. The film strength offers security during handling and thawing while flexibility ensures integrity during frozen storage and shipment. (Cleanrooms, RABS, Isolators).



Find out more

For more information, please visit
www.sartorius.com/en/products/fluid-management/freeze-thaw

Introduction

Weighing

Filtration

Final Filling

Tubing

Pipetting

Connects | Disconnects

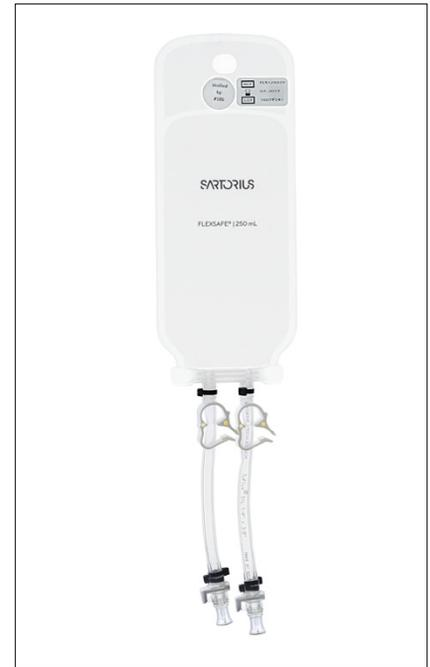
Aseptic Fluid Management

Contact

Chambers

Both the EVA Flexboy® Single Use Bags and the LDPE Flexsafe® Single Use Bags are available to the OEM market in sterile or non-sterile chambers. This can include tubing and connections of your preference or chambers with ports and no tubing.

Depending on your application or customer's needs, we have a variety of single use components and assemblies.



Introduction

Weighing

Filtration

Final Filling

Tubing

Pipetting

Connects | Disconnects

Aseptic Fluid Management

Contact

LDPE Flexsafe® Single Use Bags

The LDPE Flexsafe® Single Use Bags have a LLDPE Polyethylene contact layer which has excellent biocompatibility and ideal for media storage and cell culture. The versatile range of LDPE Flexsafe® Single Use 2D bags allows for storage, shipping and transfer of all biopharmaceutical fluids in all process steps.

The resins and additives used for LDPE Flexsafe® Single Use Bags are optimized to ensure excellent and reproducible growth. Available in sizes ranging from 5ml to 50L.



Introduction	Weighing	Filtration	Final Filling	Tubing
Pipetting	Connects Disconnects	<u>Aseptic Fluid Management</u>	Contact	

Flexboy® Single Use Bags

Standard Flexboy® bags are designed for the preparation, storage and transport of biopharmaceutical solutions, intermediates and final bulk products. They provide a single-use alternative to traditional glass, stainless steel and rigid plastic carboys in a large variety of applications.

The film used for the manufacture of all the Flexboy® bags combines multiple layers of different materials to provide a robust high barrier structure and a broad chemical compatibility. Flexboy® bags have been applied in a wide range of biopharmaceutical processes with a large variety of application fluids for more than decades. Available in sizes ranging from 5ml to 50L.



Introduction

Weighing

Filtration

Final Filling

Tubing

Pipetting

Connects | Disconnects

Aseptic Fluid Management

Contact

Mycap[®] Bottle Closures

Mycap[®] bottle closure systems make customized molded bottle closures available and affordable to everyone. The patent-pending Mycap[®] bottle closure systems fit almost any bottle, flask, tube or rigid container with a screw cap.

A platinum-cured silicone stopper with integral tubing is securely molded into the cap.



Find out more

For more information, please visit
www.sartorius.com/en/products/fluid-management/aseptic-bottle-closure

Introduction

Weighing

Filtration

Final Filling

Tubing

Pipetting

Connects | Disconnects

Aseptic Fluid Management

Contact

Takeone®

Takeone® aseptic sampling solutions are ideal to collect process fluids for analysis while protecting the product from adventitious agents. Takeone® collects perfectly representative samples removing the risk of false-positives for each process step.

Eliminate assembly operations where errors can occur or bioburden be introduced. Takeone® connects directly to process equipment to immediately take perfectly representative samples. Eliminate the risk of contamination from improper or ineffective cleaning and incorrect assembly of reusable parts.



Introduction

Weighing

Filtration

Final Filling

Tubing

Pipetting

Connects | Disconnects

Aseptic Fluid Management

Contact

Custom Manifolds

The Manifold system of single use bags or bottles are especially designed for fluid distribution from one process step to several volumes of bags with filtration, sampling and further connections capabilities.



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 contacts, visit
www.sartorius.com