



Sterisart®

Redefining Industry Standards

Simplifying Progress

SARTORIUS

Our History- Sterility Testing

Sartorius, established in 1870, has a rich and diverse history. Founded as a company developing precision measuring instruments for academic research laboratories, it has transformed into a global enterprise supporting the biopharmaceutical industry through innovative technologies. These technologies enable our customers to develop therapies more efficiently, more economically, and ensures that these therapies remain safe for patient use.

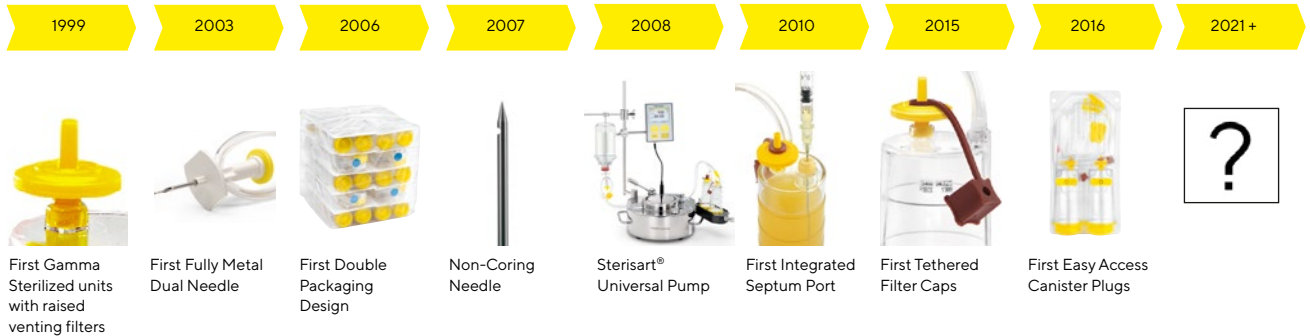
In 1927, Sartorius co-founded a company manufacturing membrane filters which for the first time made the sterilisation of heat-sensitive solutions possible. These membranes have been continually enhanced for a range of applications and remain at the core of our current offering.

In the late 1960's, identifying a major need to ensure that pharmaceuticals are free of microbial contamination, Sartorius partnered with leading scientific institutes and the German pharmaceutical industry to develop the appropriate methods and tools. Besides sterilising grade membranes, Sartorius developed membrane filters for use in the predecessor of the modern single-use sterility test system. The 'Schiller system', as it was known in Germany, was part of the Sartorius portfolio and was the first closed, reusable sterility test which was in widespread use in Europe in the early 70's. Its successor, the Sterisart® family, through small but judicious innovations, has been at the vanguard and continues to redefine the standards of sterility testing today.



Innovation Timeline

Two Decades of Innovation



We believe in continuous improvement and are committed to serving the needs of our customers. Through regular discussions with our customers we have continually adapted and bettered our range of sterility testing solutions. In doing so we have redefined industry standards.

Sterility Testing Canisters

Sterisart® NF System

For more information please click on the 

State-of-the-art closed filtration system for lot-release sterility testing of sterile pharmaceuticals.

The Sterisart® NF system is easy to use and ensures the maximum reliability of your sterility test results.

The portfolio includes a number of devices adapted to cater to your specific sterility testing needs and is fully compliant with all pharmacopeial and regulatory requirements.



See Brochure



See Flyer



See Datasheet



See Second Supplier Validation Guide



Key Features and Benefits:

- Covers all standard sterility testing applications
- Compact, closed, ergonomic design
- Specially constructed needles for safe and easy piercing
- Guaranteed sterile barrier for aseptic septum sampling
- Gas-impermeable, color- and bar-coded packaging
- Adapters for use of the Sterisart® NF system in all pumps
- Meticulously controlled quality testing
- Fully compliant with USP <71>, Ph. Eur. <2.6.1> and JP <4.06>

Sterility Testing Canisters

Sterisart® NF System

For more information

State-of-the-art
for lot-release of
pharmaceuticals.

The Sterisart® NF system is easy to use and ensures the maximum reliability of your sterility test results.

The portfolio includes a number of devices adapted to cater to your specific sterility testing needs and is fully compliant with all pharmacopeial and regulatory requirements.

Pre-installed, color coded clamps for intuitive handling.



Key Features and Benefits:

- Covers all standard sterility testing applications
- Compact, closed, ergonomic design
- Specially constructed needles for safe and easy piercing
- Guaranteed sterile barrier for aseptic septum sampling
- Gas-impermeable, color- and bar-coded packaging
- Adapters for use of the Sterisart® NF system in all pumps
- Meticulously controlled quality testing
- Fully compliant with USP <71>, Ph. Eur. <2.6.1> and JP <4.06>

Sterility Testing Canisters

Sterisart® NF System

For more information please click on the 

State-of-the-art closed filtration system

for lot-release of pharmaceuticals

The Sterisart® NF system ensures the maximum sterility test results

Raised sterile venting, with a safety clearance, to prevent accidental clogging and minimise secondary contamination.

The portfolio includes a number of devices adapted to cater to your specific sterility testing needs and is fully compliant with all pharmacopeial and regulatory requirements.




Key Features and Benefits:

- Covers all standard sterility testing applications
- Compact, closed, ergonomic design
- Specially constructed needles for safe and easy piercing
- Guaranteed sterile barrier for aseptic septum sampling
- Gas-impermeable, color- and bar-coded packaging
- Adapters for use of the Sterisart® NF system in all pumps
- Meticulously controlled quality testing
- Fully compliant with USP <71>, Ph. Eur. <2.6.1> and JP <4.06>

Sterility Testing Canisters

Sterisart® NF System

For more information please click on the 

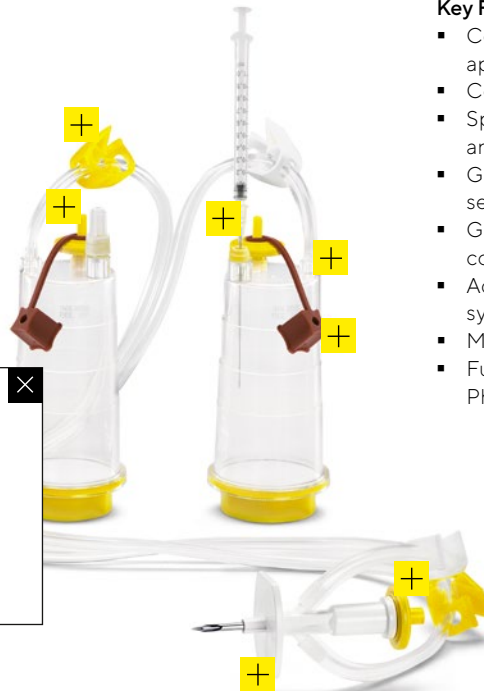
State-of-the-art closed filtration system for lot-release sterility testing of sterile pharmaceuticals.

The Sterisart® NF system is easy to use and ensures the maximum reliability of your sterility test results.

The portfolio includes a number of devices adapted for testing with all requirements.

Carefully chosen Sartochem® regenerated cellulose and cellulose acetate membranes to cover all sterility testing applications.

Robust membranes to meet low adsorption, chemical compatibility, flow rate and microbial retention requirements.



Key Features and Benefits:

- Covers all standard sterility testing applications
- Compact, closed, ergonomic design
- Specially constructed needles for safe and easy piercing
- Guaranteed sterile barrier for aseptic septum sampling
- Gas-impermeable, color- and bar-coded packaging
- Adapters for use of the Sterisart® NF system in all pumps
- Meticulously controlled quality testing
- Fully compliant with USP <71>, Ph. Eur. <2.6.1> and JP <4.06>

Sterility Testing Canisters

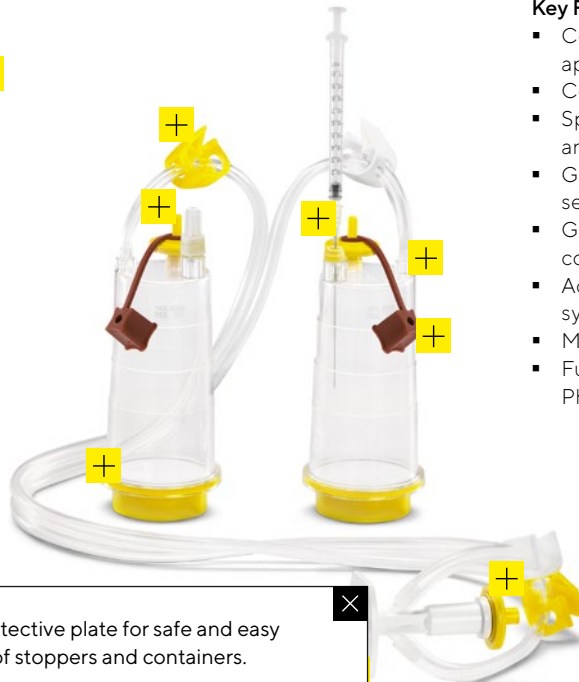
Sterisart® NF System


For more information please click on the 

State-of-the-art closed filtration system for lot-release sterility testing of sterile pharmaceuticals.

The Sterisart® NF system is easy to use and ensures the maximum reliability of your sterility test results.

The portfolio includes a number of devices adapted to cater to your specific sterility testing needs and is fully compliant with all pharmacopeial and regulatory requirements.



Large protective plate for safe and easy piercing of stoppers and containers. 

Key Features and Benefits:

- Covers all standard sterility testing applications
- Compact, closed, ergonomic design
- Specially constructed needles for safe and easy piercing
- Guaranteed sterile barrier for aseptic septum sampling
- Gas-impermeable, color- and bar-coded packaging
- Adapters for use of the Sterisart® NF system in all pumps
- Meticulously controlled quality testing
- Fully compliant with USP <71>, Ph. Eur. <2.6.1> and JP <4.06>

Sterility Testing Canisters

Sterisart® NF System

For more information please click on the 

State-of-the-art closed filtration system for lot-release sterility testing of sterile pharmaceuticals.

The Sterisart® NF system is easy to use and ensures the maximum reliability of your sterility test results.

The portfolio includes a number of devices adapted to cater to your specific sterility testing needs and is fully compliant with all pharmacopeial and regulatory requirements.



Specially constructed dual-needle with built-in sterile venting for closed sample containers.

Key Features and Benefits:

- Covers all standard sterility testing applications
- Compact, closed, ergonomic design
- Specially constructed needles for safe and easy piercing
- Guaranteed sterile barrier for aseptic septum sampling
- Gas-impermeable, color- and bar-coded packaging
- Adapters for use of the Sterisart® NF system in all pumps
- Meticulously controlled quality testing
- Fully compliant with USP <71>, Ph. Eur. <2.6.1> and JP <4.06>

Sterisart® NF System


Sterisart® NF System

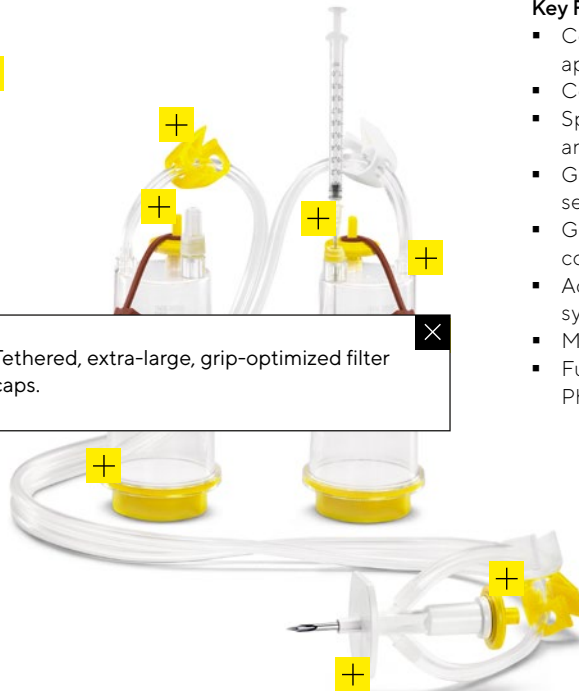
For more information please click on the 

State-of-the-art closed filtration system for lot-release sterility testing of sterile pharmaceuticals.

The Sterisart® NF system is easy to use and ensures the maximum reliability of your sterility test results.

The portfolio includes a number of devices adapted to cater to your specific sterility testing needs and is fully compliant with all pharmacopeial and regulatory requirements.

Tethered, extra-large, grip-optimized filter caps. 




Key Features and Benefits:

- Covers all standard sterility testing applications
- Compact, closed, ergonomic design
- Specially constructed needles for safe and easy piercing
- Guaranteed sterile barrier for aseptic septum sampling
- Gas-impermeable, color- and bar-coded packaging
- Adapters for use of the Sterisart® NF system in all pumps
- Meticulously controlled quality testing
- Fully compliant with USP <71>, Ph. Eur. <2.6.1> and JP <4.06>

Sterility Testing Canisters

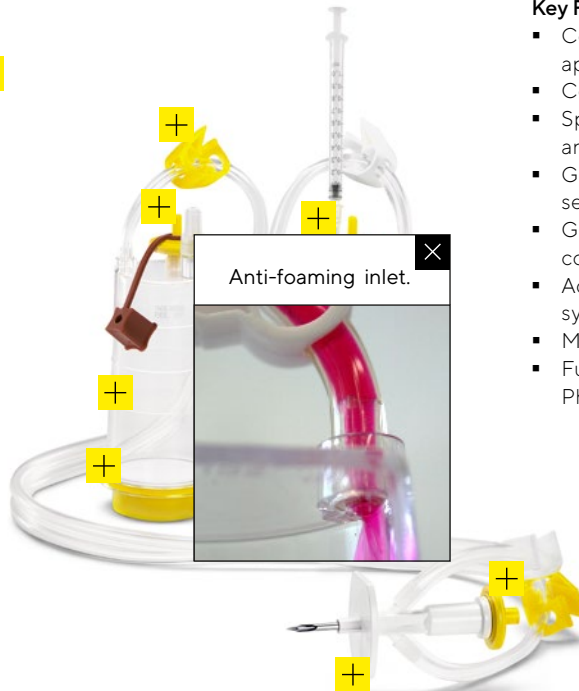
Sterisart® NF System

For more information please click on the 

State-of-the-art closed filtration system for lot-release sterility testing of sterile pharmaceuticals.

The Sterisart® NF system is easy to use and ensures the maximum reliability of your sterility test results.

The portfolio includes a number of devices adapted to cater to your specific sterility testing needs and is fully compliant with all pharmacopeial and regulatory requirements.



Key Features and Benefits:

- Covers all standard sterility testing applications
- Compact, closed, ergonomic design
- Specially constructed needles for safe and easy piercing
- Guaranteed sterile barrier for aseptic septum sampling
- Gas-impermeable, color- and bar-coded packaging
- Adapters for use of the Sterisart® NF system in all pumps
- Meticulously controlled quality testing
- Fully compliant with USP <71>, Ph. Eur. < 2.6.1> and JP <4.06>

Sterility Testing Canisters

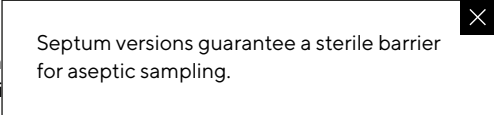
Sterisart® NF System

For more information please click on the 

State-of-the-art closed filtration system for lot-release sterility testing of sterile pharmaceuticals.

The Sterisart® NF system is easy to use and ensures the maximum reliability of sterility test results.

The portfolio includes a number of devices adapted to cater to your specific sterility testing needs and is fully compliant with all pharmacopeial and regulatory requirements.

 Septum versions guarantee a sterile barrier for aseptic sampling.



Key Features and Benefits:

- Covers all standard sterility testing applications
- Compact, closed, ergonomic design
- Specially constructed needles for safe and easy piercing
- Guaranteed sterile barrier for aseptic septum sampling
- Gas-impermeable, color- and bar-coded packaging
- Adapters for use of the Sterisart® NF system in all pumps
- Meticulously controlled quality testing
- Fully compliant with USP <71>, Ph. Eur. <2.6.1> and JP <4.06>

Canister Types



Yellow Base

Regenerated Cellulose



White Base

Cellulose Acetate

Material of construction

- Housing: Styrene acrylonitrile (SAN)
- Tubing: PVC (double lumen) & silicone
- Standard sampling needle: Polycarbonate & stainless steel
- Dual needle: Acrylonitrile butadiene styrene (ABS) & stainless steel
- Wing nuts: Polyethylene (PE)
- Filter plugs: Silicone
- Septum: Polyisoprene and acrylonitrile butadiene styrene (ABS)

Maximum operating pressure:

3 bar at 20 °C

Maximum operating temperature:

50 °C

Burst pressure of the housing:

> 5 bar

Capacity:

120 ml (graduations at 50 ml, 75 ml and 100 ml)

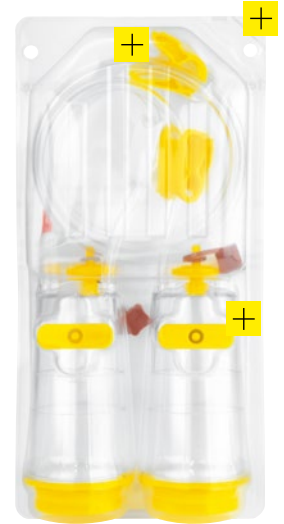
Venting filter

- Hydrophobic polytetrafluoroethylene (PTFE) membrane
- Penetration pressure > 3 bar

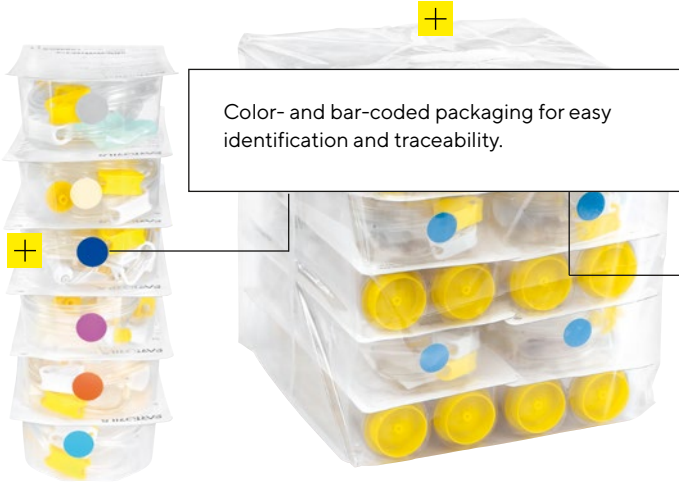
Sterilization:

Gamma irradiation

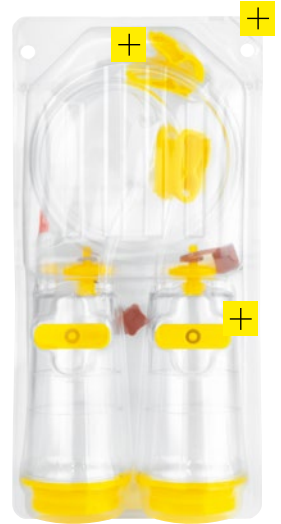
Packaging



Packaging



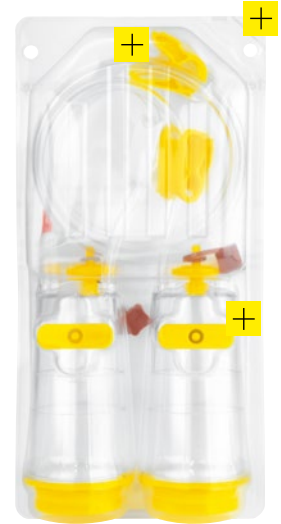
Color- and bar-coded packaging for easy identification and traceability.



Packaging



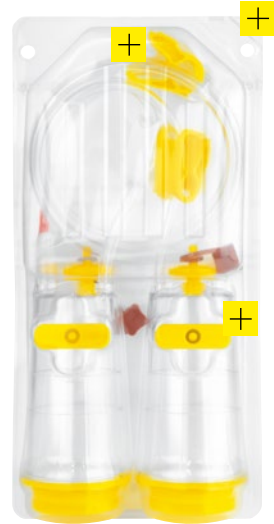
Double packed for easy and secure transfer into a clean-room, laminar flow hood or isolator.



Packaging



Gamma-sterilized to prevent secondary contamination and mitigate the risk of false positives.



Packaging



Gas-impermeable packaging to eliminate the risk of false-negatives, following VHP-decontamination in isolators.

Packaging



Corner reinforcement for easy-peeling and hanger holes.

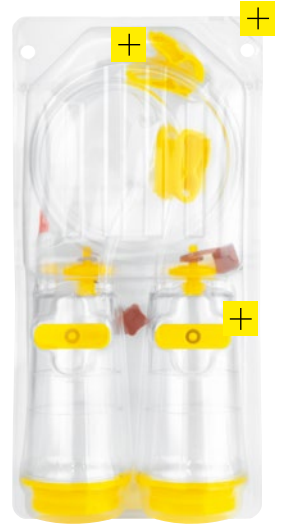


Packaging



Recess housing directly-accessible, large canister plugs.

Packaging



Sterisart® Septum

During growth incubation, sampling of the culture medium may be required for any one of the following reasons:

- The growth media is rendered turbid by microbial growth, following incubation, and necessitates the identification of the micro-organism as part of a root cause analysis.
- The product renders the growth medium turbid, prior to incubation, and requires sub-culturing | dilution.
- Samples are supplemented with agents to counteract anti-microbial components of the tested product.
- Samples are drawn to test for microbial contamination by rapid detection methods

Sampling via the tubing can compromise the integrity of the sterility test and preclude re-incubation of the sterility testing canisters. The aseptic sampling port not only eliminates the risk of introducing false positives through external contamination but also ensures operator safety and prevents accidental spillage during sampling.

Given the interest in rapid sterility testing solutions, Sartorius has partnered with Charles River to pair our Sterisart® Septum canisters with the Celsis® detection platform.



See Application Note

[Click here for a study on how the Sterisart® Septum facilitates the recurrent sterile extraction of samples](#)



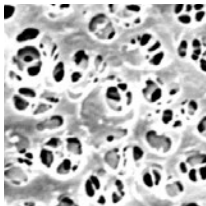
Learn More

[Learn more about our partnership with Charles River here:](#)

Membranes

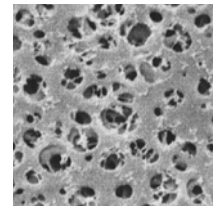
Regenerated Cellulose

Material	Regenerated cellulose, reinforced with non-woven cellulose
Properties	<ul style="list-style-type: none">▪ Combines excellent chemical resistance and thermal stability with very low adsorption characteristics▪ Hydrophilic
Chemical Compatibility	Aqueous solutions (pH 3-12) and organic solvents
Pore size	0.45 μm
Flow Rate for Water per cm^2 (DIN 58355:)	30 ml/min at $D_p = 1 \text{ bar}$ $\sim 15 \text{ psi}$
Thickness (DIN 53105)	150 – 170 μm
Wetting time	< 1 second, with deionised water




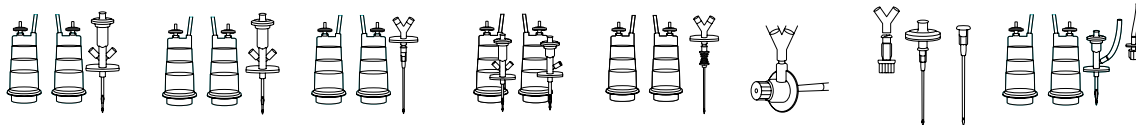
Cellulose Acetate

Material	Cellulose acetate
Properties	<ul style="list-style-type: none">▪ Combines high-flow rates and thermal stability with very low adsorption characteristics, making it ideal for difficult-to-filter, viscous substances▪ Hydrophilic
Chemical Compatibility	Aqueous solutions (pH 4-8), oils, alcohols and several organic solvents
Pore size	0.45 μm
Flow Rate for Water per cm^2 (DIN 58355:)	65 ml/min at $D_p = 1 \text{ bar}$ $\sim 15 \text{ psi}$
Thickness (DIN 53105)	115 – 145 μm
Wetting time	< 1 second, with deionised water



Ordering Information









For more information, please select each column's 



Regenerated Cellulose


Standard	16466-----GBD	16476-----GBD	16467-----GBD	16475-----GBD	16477-----GBD	16469-----GBD	16468-----GBD	16478-----GBD
Septum	16466-----GSD	16476-----GSD	16467-----GSD	16475-----GSD		16469-----GSD		

Cellulose Acetate

	16466-----CA-GSD		16467-----CA-GSD					
Sample Type								
Product Container	Closed	Closed	Open	Closed	Closed with bow-fill seals	Syringes	Male luer locks or female slip	Male luer locks



Ordering Information

For more information, please select each column's 

Sterility Testing Solutions for Liquid Transfer and Dilution



(for sample prep and dilution)

16470-----GBD



(for liquid transfer)

16471-----GBD



(for liquid transfer)

16472-----GBD

Sample Type



Closed container



Open Container



Open/Vented Container



Ordering Information

Regenerated Cellulose

Standard Canisters

16466-----GBD



Septum Canisters

16466-----GSD



Cellulose Acetate

Septum Canisters

16466-----CA-GSD



Connector: Long (41 mm) dual-needle with built-in venting filter

Designed for: Liquids in closed, large volume containers

Ordering Information

Regenerated Cellulose

Standard Canisters

16476-----GBD



Septum Canisters

16476-----GSD



Connector: Short (21 mm) dual-needle with built-in venting filter

Designed for: Liquids in closed, small volume containers

Ordering Information

Regenerated Cellulose

Standard Canisters

16467-----GBD



Septum Canisters

16467-----GSD



Cellulose Acetate

Septum Canisters

16467-----CA-GSD



Connector: Standard sampling needle (52 mm) Separate needle for sterile venting (40 mm)

Designed for: Liquids in open containers

Ordering Information

Regenerated Cellulose

Standard Canisters

16475-----GBD



Septum Canisters

16475-----GSD



Connector: Standard (35 mm) dual-needle with built-in venting filter & sampling dual-needle (23 mm)

Designed for: Soluble lyophilisates in closed containers

Ordering Information

Regenerated Cellulose

Standard Canisters

16477-----GBD



Non-coring needle

Connector: Non-coring needle (56 mm)

Separate needle for sterile venting (40 mm)

Designed for: Liquids in plastic containers with blow-fill seals

Ordering Information

Regenerated Cellulose

Standard Canisters

16469-----GBD



Septum Canisters

16469-----GSD



Connector: Standard (35 mm) dual-needle with built-in venting filter & syringe adaptor

Designed for: Liquids in pre-filled syringes

Ordering Information

Regenerated Cellulose

Standard Canisters

16468-----GBD (One-Connector System)



Connector: Standard sampling needle (60 mm) & detachable two-way female Luer adaptor. Separate needle for sterile venting (40 mm)

Designed for: Liquids in containers with a Luer connector

Ordering Information

Regenerated Cellulose

Standard Canisters

16478-----GBD (Two-Connector System)



Connector: Standard (35 mm) dual-needle with built-in venting filter & female Luer adaptor.

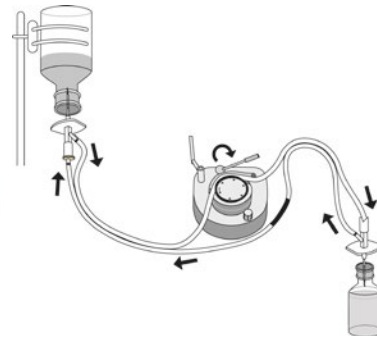
Designed for: Liquids in containers with a male Luer connector

Ordering Information

Sterility Testing Solutions for Liquid Transfer and Dilution

Sample Preparation

16470-----GBD



Connector: 37 mm and 23 mm dual-needle for sampling

Designed for: Sample preparation of poorly soluble lyophilized products

Ordering Information

Sterility Testing Solutions for Liquid Transfer and Dilution

Sample Preparation

16471-----GBD



Connector: 29 mm dual-needle for sampling

Designed for: Direct inoculation: Transfer of liquids that are difficult to filter
Sample pooling of small volumes prior to membrane filtration

Ordering Information

Sterility Testing Solutions for Liquid Transfer and Dilution

Sample Preparation

16472-----GSD



Connector: 17 mm dual-needle and 80 mm standard needle for sampling

Designed for: Direct inoculation: Transfer of liquids that are difficult to filter
Sample pooling of small volumes prior to membrane filtration

Quality Assurance

Routine testing of every lot includes:


- **Incoming Materials Testing**
Qualification of the membrane filters, plastic and tubing
- **Rigorous In-Process Controls**
Physical integrity test of the housing container, tubing and venting filter
- **Stringent Final Release Testing**
Includes a physical integrity test, bacterial challenge test and growth promotion test.

A quality assurance certificate is delivered with every unit for your records.

A comprehensive validation/ qualification guide is available on request.



Sterility Testing Pump

For more information please click on the 

The Sterisart® Universal Pump, can be used in isolators, laminar flow hoods, biosafety cabinets and cleanroom benches. Since space is often at a premium, the pump features a compact architecture with a space-saving footprint.

We have designed our pump focusing on your convenience. Combined with our Sterisart® sterility testing canisters, we offer you complete confidence in your sterility test results.



Key Features and Benefits:

- Versatile pump designed for bench-top use or integration into an isolator
- Compact, low-profile for comfortable use
- Pioneering closed pump chassis design with built-in passive cooling to prevent particulate emission
- Effortlessly detachable components to facilitate easy-access for cleaning
- Robust & durable: Rugged stainless-steel body that is compatible with all commonly used chemical sterilants and fully VHP-resistant
- Easy tube-placement with convenient, malfunction-proof, locking mechanism
- Open-design: compatible with other filtration devices on the market

Model No. 16420



See Brochure

Sterility Testing Pump

For more information please click on the [+](#)

The Sterisart® Universal Pump, can be used in isolators, laminar flow cabinets and clean rooms. The compact design and space saving footprint features a compact space-saving footprint.

Height adjustable bottle and bag holder.


We have designed our pump focusing on your convenience. Combined with our Sterisart® sterility testing canisters, we offer you complete confidence in your sterility test results.



Key Features and Benefits:

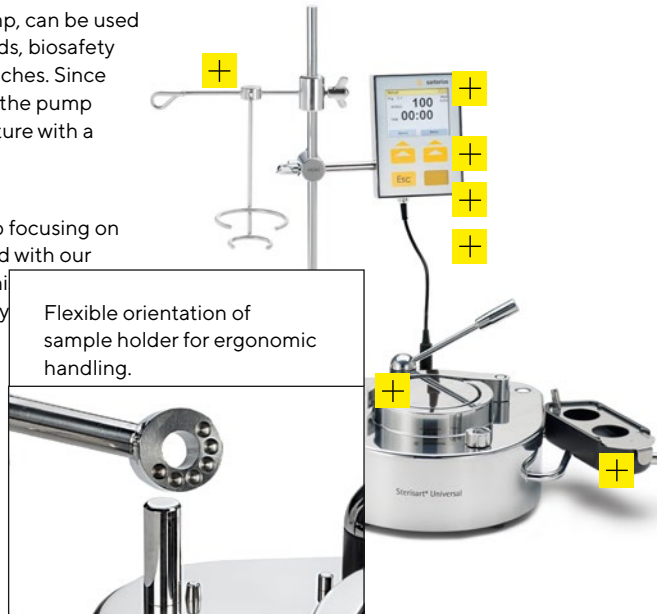
- Versatile pump designed for bench-top use or integration into an isolator
- Compact, low-profile for comfortable use
- Pioneering closed pump chassis design with built-in passive cooling to prevent particulate emission
- Effortlessly detachable components to facilitate easy-access for cleaning
- Robust & durable: Rugged stainless-steel body that is compatible with all commonly used chemical sterilants and fully VHP-resistant
- Easy tube-placement with convenient, malfunction-proof, locking mechanism
- Open-design: compatible with other filtration devices on the market

Sterility Testing Pump

For more information please click on the 

The Sterisart® Universal Pump, can be used in isolators, laminar flow hoods, biosafety cabinets and cleanroom benches. Since space is often at a premium, the pump features a compact architecture with a space-saving footprint.

We have designed our pump focusing on your convenience. Combined with our Sterisart® sterility testing canisters you complete confidence in your test results.




Flexible orientation of sample holder for ergonomic handling.

Key Features and Benefits:

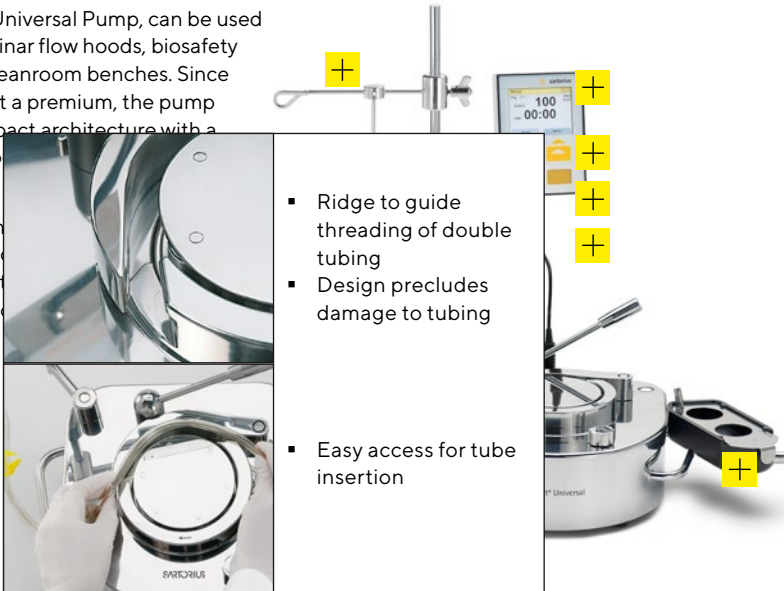
- Versatile pump designed for bench-top use or integration into an isolator
- Compact, low-profile for comfortable use
- Pioneering closed pump chassis design with built-in passive cooling to prevent particulate emission
- Effortlessly detachable components to facilitate easy-access for cleaning
- Robust & durable: Rugged stainless-steel body that is compatible with all commonly used chemical sterilants and fully VHP-resistant
- Easy tube-placement with convenient, malfunction-proof, locking mechanism
- Open-design: compatible with other filtration devices on the market

Sterility Testing Pump

For more information please click on the 

The Sterisart® Universal Pump, can be used in isolators, laminar flow hoods, biosafety cabinets and cleanroom benches. Since space is often at a premium, the pump features a compact architecture with a space-saving form factor.

We have designed your convenience. Sterisart® sterilization you complete confidence in test results.




- Ridge to guide threading of double tubing
- Design precludes damage to tubing

- Easy access for tube insertion

Key Features and Benefits:

- Versatile pump designed for bench-top use or integration into an isolator
- Compact, low-profile for comfortable use
- Pioneering closed pump chassis design with built-in passive cooling to prevent particulate emission
- Effortlessly detachable components to facilitate easy-access for cleaning
- Robust & durable: Rugged stainless-steel body that is compatible with all commonly used chemical sterilants and fully VHP-resistant
- Easy tube-placement with convenient, malfunction-proof, locking mechanism
- Open-design: compatible with other filtration devices on the market

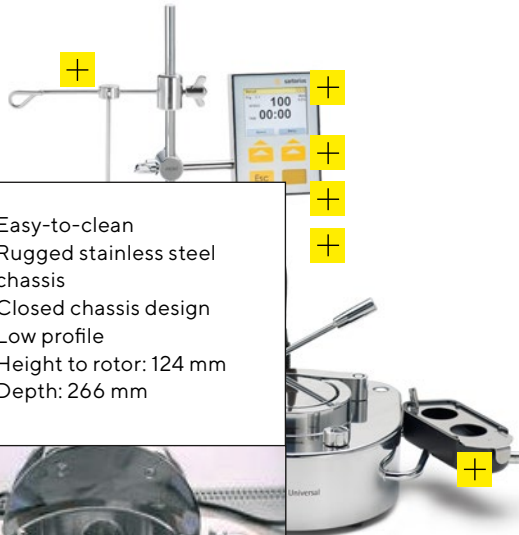
Sterility Testing Pump

For more information please click on the 

The Sterisart® Universal Pump, can be used in isolators, laminar flow hoods, biosafety cabinets and cleanroom benches. Since space is often at a premium, the pump features a compact architecture with a space-saving footprint.

We have designed our pump focusing on your convenience. Combined with our Sterisart® sterility testing canisters, we give you complete confidence in your sterility test results.


- Easy-to-clean
- Rugged stainless steel chassis
- Closed chassis design
- Low profile
Height to rotor: 124 mm
Depth: 266 mm



Key Features and Benefits:

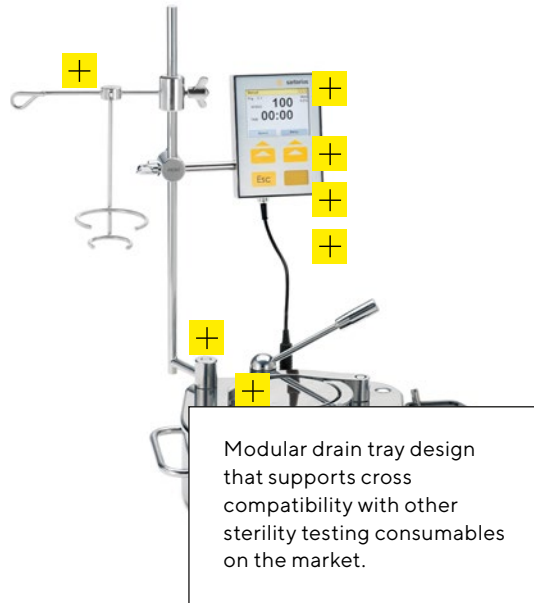
- Versatile pump designed for bench-top use or integration into an isolator
- Compact, low-profile for comfortable use
- Pioneering closed pump chassis design with built-in passive cooling to prevent particulate emission
- Effortlessly detachable components to facilitate easy-access for cleaning
- Robust & durable: Rugged stainless-steel body that is compatible with all commonly used chemical sterilants and fully VHP-resistant
- Easy tube-placement with convenient, malfunction-proof, locking mechanism
- Open-design: compatible with other filtration devices on the market

Sterility Testing Pump

For more information please click on the 

The Sterisart® Universal Pump, can be used in isolators, laminar flow hoods, biosafety cabinets and cleanroom benches. Since space is often at a premium, the pump features a compact architecture with a space-saving footprint.


We have designed our pump focusing on your convenience. Combined with our Sterisart® sterility testing canisters, we offer you complete confidence in your sterility test results.



Key Features and Benefits:

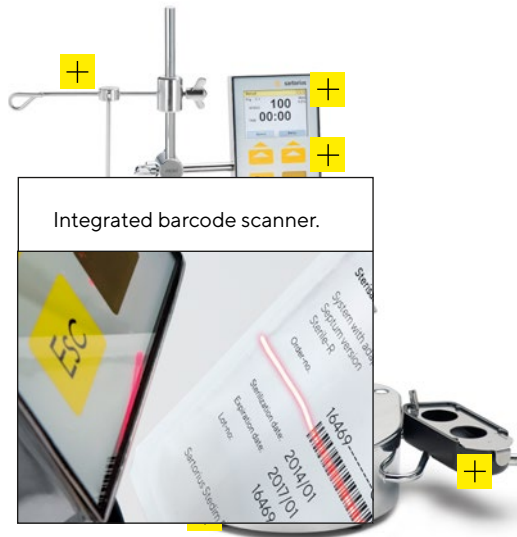
- Versatile pump designed for bench-top use or integration into an isolator
- Compact, low-profile for comfortable use
- Pioneering closed pump chassis design with built-in passive cooling to prevent particulate emission
- Effortlessly detachable components to facilitate easy-access for cleaning
- Robust & durable: Rugged stainless-steel body that is compatible with all commonly used chemical sterilants and fully VHP-resistant
- Easy tube-placement with convenient, malfunction-proof, locking mechanism
- Open-design: compatible with other filtration devices on the market

Sterility Testing Pump

For more information please click on the 

The Sterisart® Universal Pump, can be used in isolators, laminar flow hoods, biosafety cabinets and cleanroom benches. Since space is often at a premium, the pump features a compact architecture with a space-saving footprint.


We have designed our pump focusing on your convenience. Combined with our Sterisart® sterility testing canisters, we offer you complete confidence in your sterility test results.



Key Features and Benefits:

- Versatile pump designed for bench-top use or integration into an isolator
- Compact, low-profile for comfortable use
- Pioneering closed pump chassis design with built-in passive cooling to prevent particulate emission
- Effortlessly detachable components to facilitate easy-access for cleaning
- Robust & durable: Rugged stainless-steel body that is compatible with all commonly used chemical sterilants and fully VHP-resistant
- Easy tube-placement with convenient, malfunction-proof, locking mechanism
- Open-design: compatible with other filtration devices on the market

Sterility Testing Pump

For more information please click on the 

The Sterisart® Universal Pump, can be used in isolators, laminar flow hoods, biosafety cabinets and cleanroom benches. Since space is often at a premium, the pump features a compact architecture with a space-saving footprint.


We have designed our pump focusing on your convenience. Combined with our Sterisart® sterility testing canisters, we offer you complete confidence in your sterility test results.



Key Features and Benefits:

- Versatile pump designed for bench-top use or integration into an isolator
- Compact, low-profile for comfortable use
- Pioneering closed pump chassis design with built-in passive cooling to prevent particulate emission
- Effortlessly detachable components to facilitate easy-access for cleaning
- Robust & durable: Rugged stainless-steel body that is compatible with all commonly used chemical sterilants and fully VHP-resistant
- Easy tube-placement with convenient, malfunction-proof, locking mechanism
- Open-design: compatible with other filtration devices on the market

Sterility Testing Pump

For more information please click on the 

The Sterisart® Universal Pump, can be used in isolators, laminar flow hoods, biosafety cabinets and cleanroom benches. Since space is often at a premium, the pump features a compact architecture with a space-saving footprint.


We have designed our pump focusing on your convenience. Combined with our Sterisart® sterility testing canisters, we offer you complete confidence in your sterility test results.



Key Features and Benefits:

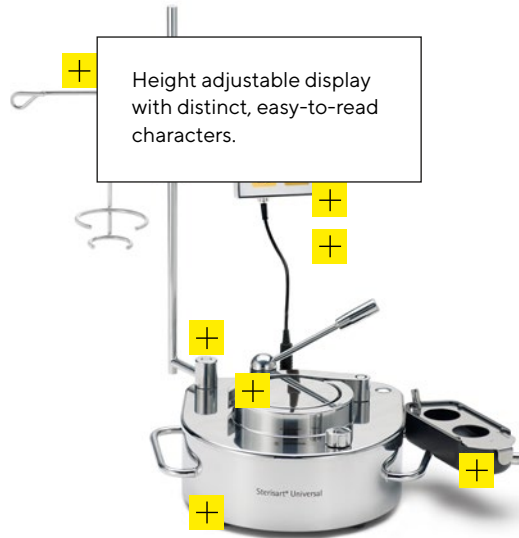
- Versatile pump designed for bench-top use or integration into an isolator
- Compact, low-profile for comfortable use
- Pioneering closed pump chassis design with built-in passive cooling to prevent particulate emission
- Effortlessly detachable components to facilitate easy-access for cleaning
- Robust & durable: Rugged stainless-steel body that is compatible with all commonly used chemical sterilants and fully VHP-resistant
- Easy tube-placement with convenient, malfunction-proof, locking mechanism
- Open-design: compatible with other filtration devices on the market

Sterility Testing Pump

For more information please click on the 

The Sterisart® Universal Pump, can be used in isolators, laminar flow hoods, biosafety cabinets and cleanroom benches. Since space is often at a premium, the pump features a compact architecture with a space-saving footprint.

We have designed our pump focusing on your convenience. Combined with our Sterisart® sterility testing canisters, we offer you complete confidence in your sterility test results.



Key Features and Benefits:

- Versatile pump designed for bench-top use or integration into an isolator
- Compact, low-profile for comfortable use
- Pioneering closed pump chassis design with built-in passive cooling to prevent particulate emission
- Effortlessly detachable components to facilitate easy-access for cleaning
- Robust & durable: Rugged stainless-steel body that is compatible with all commonly used chemical sterilants and fully VHP-resistant
- Easy tube-placement with convenient, malfunction-proof, locking mechanism
- Open-design: compatible with other filtration devices on the market

Accessories

Model No.	Description	
1ZE---0040	Communication Kit	+
1Z---0004	Sterisart® Easy Configuration Software	+
1ZW---0002	Ampoule breaker	+
1ZGD--0031	Protective stainless-steel rotor cover/shield	+
1ZA---0024	Syringe adaptor/ support for pre-filled syringes	+
1ZE---0033	Foot switch	
1EE---0009	Extension cable control unit/ display	
1ZG---0024	Stainless steel drain cover/ adaptor for Millipore sterility test units	+
1ZE---0050	Isolator installation kit	
1ZGF--0020	Transport tray (10-canister holder)	+



Accessories

Model

1ZE---0

1Z---0

1ZW---

1ZGD-

1ZA---

1ZE---0

1EE---

1ZG---0024 Stainless steel drain cover/ adaptor for Millipore sterility test units

1ZE---0050 Isolator installation kit

1ZGF--0020 Transport tray (10-canister holder)



1Z---0040 - Communication Kit

Contains:

- 1GB SD card
- USB male to female converter
- RS-232 to RS-232 converter
- RS-232 cable (approx. 2 m)
- Driver CD, USB to serial port
- Card reader, RS-232 to SD

Accessories

Model

1ZE---

1Z---

1ZW---

1ZGD-

1ZA---

1ZE---

1EE---

1ZG---

1ZE---

1ZGF---



1ZE---0004 - Sterisart® Easy Configuration Software

- Designed to increase the process reliability for sterility testing
- Intuitive drag-and-drop programming of sterility testing workflow
- Step-by-step guidance of the operator through each programmed sterility test
- User account management, audit trail and password protection
- Scanning, saving and printing of process-relevant material data
- FDA 21 CFR Part 11 compliant electronic records



See Datasheet

Accessories

Model No.
1ZE---0040
1Z---0004
1ZW---0002
1ZGD--0031
1ZA---0024
1ZE---0033
1EE---0009
1ZG---0024
1ZE---0050
1ZGF--0020

1ZW---0002 - Ampoule Breaker



Accessories

Model No.
1ZE---0040
1Z---0004
1ZW---0002
1ZGD--0031
1ZA---0024
1ZE---0033
1EE---0009
1ZG---0024
1ZE---0050
1ZGF--0020

1ZGD--0031 - Stainless Steel Rotor Cover



Accessories

Model

1ZE---0

1Z---0

1ZW---

1ZGD-

1ZA---

1ZE---0

1EE---

1ZG---

1ZE---0

1ZGF--



1ZA---0024- Syringe adaptor for pre-filled syringes

For liquids in pre-filled syringes

For use with 16469-----GBD or 16469-----GSD.

A foot-switch controls both the operation of the pump and the pinch-valve, which acts as a flow restrictor, through a communication hub. The valve regulates the flow of the rinsing fluid through the tubing and thereby eliminates the need to overturn the rinsing fluid bottle between sample injections. The Sterisart® syringe adaptor can also be attached to the main support rod of the pump to minimize the space used.

Accessories

Model

1ZE---0

1ZG---0024 Stainless steel drain cover/ adaptor for Millipore™ sterility test units

1Z---0

1ZW---

1ZGD-

1ZA---

1ZE---

1EE---

1ZG---

1ZE---

1ZGF---



Accessories

Model

1ZE---0

1Z---0

1ZW---

1ZGD-

1ZA---

1ZE---0

1EE---0

1ZG---

1ZE---0

1ZGF--

1ZGF--0020 - Transport tray (10-canister holder)



Service Support

Service life cycle management is a critical element when purchasing your equipment. This routinely involves:

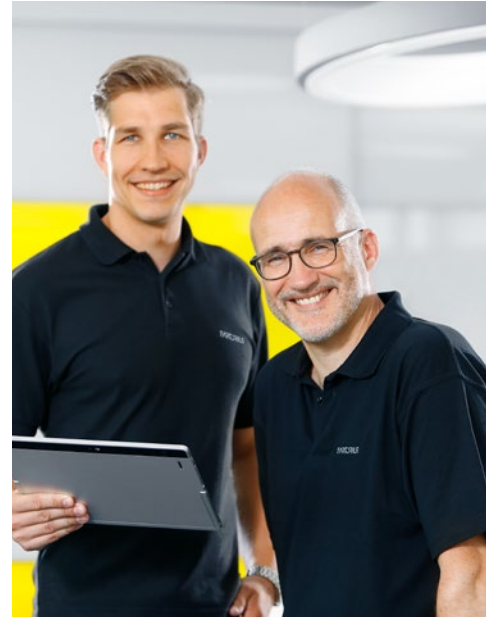
- Proper installation of the equipment
- Basic user training
- Routine preventative maintenance visits (IQ/OQ)

These services will increase the longevity of the unit, while reducing the downtime.

We serve customers around the globe – with service contracts and a full range of services to suit every need. If you have questions about our service offerings or are in need of technical support, we are here for you. Just provide us with detailed information in our contact form, and we will promptly get in touch with you.



[Click to Learn More](#)



Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Stedim North America Inc.
5 Orville Drive, Suite 200
Bohemia, NY 11716
Toll-Free +1 800 368 7178



For further information, visit
www.sartorius.com