SARTURIUS

Sterisart® Family

Sterility Testing Made Easy and Reliable



Benefits

- Compact, ergonomic design to cover all standard sterility testing applications
- Specially engineered needles for safe and easy piercing of stoppers and containers
- Anti-foaming inlet
- Raised sterile venting
- Tethered, large, grip-optimized filter caps
- Septum port that guarantees aseptic sampling and supplementation
- Recess in packaging that houses directly accessible outlet plugs

Product Information

The development and manufacturing of injectable, sterile medical products under GMP conditions is one of the most challenging and sensitive issues in the (bio-)pharmaceutical industry. Before a batch of the parenterals concerned can be released by the quality control department (QC), proof of sterility must be provided by performing the sterility test according to USP <71>, Ph. Eur. 2.6.1 and JP 4.06.

International Guidelines and Requirements

According to international pharmacopeias, parenterals (such as vaccines, mAbs, intravenous infusions, antibiotics, diabetes medication) that are injected into the human or animal body are subject to sterility testing. However, other products must be tested for sterility as well, for example:

- Ophthalmics
- Immunodiagnostics (urine | blood, etc.)
- Intermediates | APIs
- Cell culture media
- Cell banks
- Virus banks
- Medical instruments (e.g., scalpels)
- Ointments
- Creams

Various analytical methods are used depending on the volume to be analyzed; e.g., membrane filtration or direct inoculation.

Generally, the sterility of each batch of a sterile drug is verified by performing the sterility test at the latest on such pharmaceuticals packaged in their final containers. Given the role of the test method in the federal laws of some countries (e.g., U.S. CFR 610.12), this qualitative test for the presence or absence of microorganims, yeasts and fungi also has additional, legal significance.

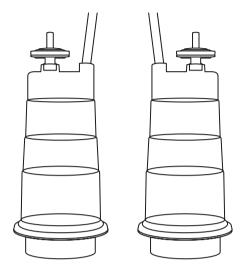
Quality Control

As Sartorius is a recognized supplier of products and services for the pharmaceutical industry, product safety and quality are its number one priorities. To meet the quality requirements according to ISO 9001, representative samples of each lot of Sterisart® systems are subjected to destructive testing. In-process and final quality control tests further ensure high product safety.

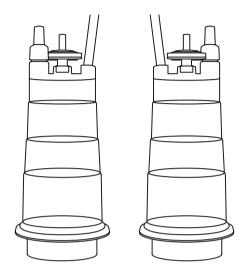
In addition to the bacteria challenge test and the Method Suitability test, the bubble point, flow rate, thickness and wetting time of the membrane are all checked.

Technical Specifications

Basic Structure of Sterisart® RC Units with a Regenerated Cellulose Membrane



Standard configuration, order no. ending in GBD



Configuration with septum, order no. ending in GSD

Sample Type	Product Container		Spike	Usage	Order Number
LVPs	Closed glass bottles with a septum				16466GBD
LVPs	Closed glass bottles with a septum	AS THE STATE OF TH		septum	16466GSD
LVPs, SVPs	Open containers e.g. glass ampoules, glass bottles, collapsible plastic bags				16467GBD

According to the United States Pharmacopeia:

^{*}LVPs: Large Volume Parenterals > 100 mL

^{*}SVPs: Small Volume Parenterals < 100 mL

Basic Structure of Sterisart® RC Units with a Regenerated Cellulose Membrane

Sample Type	Product Container	 Spike	Usage	Order Number
LVPs, SVPs	Open containers e.g. glass ampoules, glass bottles, col- lapsible plastic bags		septum	16467GSD
SVPs	Prefilled syringes with or without a spike			16469GBD
SVPs	Prefilled syringes with or without a spike		septum	16469GSD
Lyophilisates, soluble powders, liquid antibiotics	Closed glass bottles with a septum			16475GBD
Lyophilisates, soluble powders, liquid antibiotics	Closed glass bottles with a septum		septum	16475GSD
SVPs	Closed glass bottles with a septum			16476GBD
SVPs	Closed glass bottles with a septum		septum	16476GSD

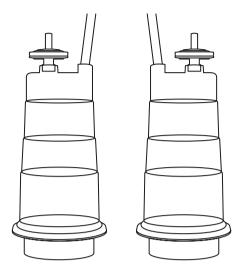
According to the United States Pharmacopeia: *LVPs: Large Volume Parenterals > 100 mL *SVPs: Small Volume Parenterals < 100 mL

Basic Structure of Sterisart® RC Units with a Regenerated Cellulose Membrane

Sample Type	Product Container		Spike	Usage	Order Number
LVPs, SVPs, eye drops	Closed containers; plastic containers with blow-fill seals; e.g. bottles, ampoules	And			16477GBD
LVPs	One-connector system for testing tube assemblies and bags. Fits product containers with male Luer lock or female Luer slip connectors				16468GBD
LVPs	Two-connector system for testing product containers with a male Luer lock; double-needle spike for simultaneous transfer of rinsing liquid				16478GBD

According to the United States Pharmacopeia: *LVPs: Large Volume Parenterals >100 mL *SVPs: Small Volume Parenterals <100 mL





Standard configuration, order no. ending in GBD

Basic Structure of Sterisart® CA Units with a Cellulose Acetate Membrane

Sample Type	Product Container	Spike	Usage	Order Number
LVPs	Closed glass bottles with septum		septum	16466CA-GSD
LVPs, SVPs	Open containers; e.g., glass ampoules, glass bottles, collaps- ible plastic bags		septum	16467CA-GSD

According to the United States Pharmacopeia: *LVPs: Large Volume Parenterals >100 mL *SVPs: Small Volume Parenterals <100 mL

Products for Direct Aseptic Transfer of Liquid Samples to Each Required Culture Medium

Sample Type	Product Container	Spike	Figure	Order Number
Liquids difficult to filter; medical materials	Closed glass bottles with a septum		+	16472GBD
Liquids difficult to filter; medical materials	Closed glass bottles with a septum open glass bottles			16471GBD

Sample Preparation

Sample Type	Product Container		Usage	Order Number
Powder with low solubility in closed glass bottles with a septum	Tube assembly with metal double needles of two different lengths			16470GBD

Optional Accessories

Sample Type	Product Container	Usage	Order Number
filled with rinsing solutions and	Spike with 0.2 µm sterilizing-grade filter, 4 cm, stainless steel, individually sterile-packaged, gamma- sterilized		16596HNK

Applications

Applications for Different Membranes and Systems

Sterisart® RC with a regenerated cellulose membrane is particularly suitable for testing aggressive, aqueous products and antibiotics.

Sterisart® CA with a cellulose acetate membrane has been specially designed to analyze the sterility of difficult-to-filter, viscous substances, such as emulsions, using membrane filtration.

Transfer sets: If the membrane filtration method cannot be used for sterility testing of liquid products, the Sartorius sterile transfer sets can be used for aseptic transfer of these products to liquid culture media for performing the direct inoculation method.

Overview of the Sterisart® Systems

Gamma-sterilized systems with regenerated cellulose (RC) membrane 16466-GBD, 16467-GBD, 16468-GBD, 16469-GBD, 16470-GBD, 16471-GBD, 16472-GBD, 16475-GBD, 16476-GBD, 16477-GBD, 16478-GBD,

Gamma-sterilized systems

with a septum and a regenerated cellulose (RC) membrane 16466-GSD, 16467-GSD, 16469-GSD, 16475-GSD, 16476-GSD

Gamma-sterilized systems with a cellulose acetate (CA) membrane 16466-CA-GSD; 16467-CA-GSD

Note:

The primary packaging of the systems as well as the primary packing of the transfer kits are gas-tight. This enables them to be used directly in isolators and prevents unnecessary rinsing steps.

EXPAND® Training

Available on request in your company or in our Training center.



Sterisart® Universal Pumps for Transfer of Liquids to the Sterisart® Units

16420 with a display and an integrated barcode scanner 16419 Basic version

Order number	Description		
1Z0004	Sterisart® Easy Configuration Software Drag & drop software, designed to increase process reliability for sterility testing		
1ZG0023	Drainage container cover for Sterisart® sterility test systems		
1ZE0033	Trolley for Sterisart® Universal Pump		
1ZGF0020	Tray for 10 Sterisart® units		
1ZE0039	Trolley for pump		
1ZA0002	Drain tubing		
1ZG0028	Drainage container and cover for Sterisart® Sterility units		
1ZGL0033	Bottle holder without wing nut		
1ZF0007	Wing nut		

Order number	Description
1ZGD0031	Stainless steel cover for the rotor
1EE0010	External barcode scanner
1ZE0050	Isolator installation kit
1ZG0024	Drainage container cover for competitor's sterility test consumables
1ZG0014	Adapter for Sterisart® systems in Equinox pumps; pkg. of 2
1ZA0028	Drainage container for Sterisart® pump 16419 16420
	16419 16420

Service

EXTEND® instrument services for Sterisart® Universal pumps

- Installation qualification and operational qualification (IQ | OQ)
- Preventive maintenance



General Technical System Details and Regulatory Requirements

The Sterisart® systems comply with GMP requirements with respect to sterility testing as well as to Regulation (EC) No. 1907/2006 (REACH) with particular reference to the plasticizer DEHP; ED/108/2014, ED/67/2008.

Sterilization

Gamma-sterilized at 25 kGy in compliance with DIN EN 552 and ISO 11137

Sterilization indicator on each carton; color change: from orange to red

Shelf Life

Gamma-sterilized units:

Sterile for up to 3 years after the date of manufacture; guaranteed for a minimum of 6 months after delivery





Dimensions and Weight	
Box (W×D×H)	28.1×27.3×24.6 cm
Weight of carton with 10 Sterisart® units	Approx. 2.2 kg
Gas-tight individual packaging (W× D× H)	5.5×13×27cm
Veight of Sterisart® unit vith packaging	Approx. 185 g
From bottom to sterile air filter ed cap attached (H×D)	Approx. 13.8 × 5.7 cm
Veight of the Sterisart® unit vithout primary packaging	Approx. 150 g, depending on the system version

Packaging				
Primary packaging material	OPA PE sealing foil			
Transparent film	A-PET/PE			
Transport packaging material	Polyamide (PA) and polyethylene (PE)			
Transparent film	PETG (polyethylene terephthalate)			





Sterisart® System

Method Suitability tests are performed on Sterisart® systems in compliance with the international pharmacopeias; USP <71> and Ph. Eur. 2.6.1.; an extractables profile has been created in compliance with the USP 23 and Ph. Eur. 3. The detection limits are below the requirements as specified for "water for injection".

Equal distribution of the product to the Sterisart® containers with a max. validated deviation of 10 %

Regenerated cellulose (RC) membrane				
Pore size	0.45 µm nominal; specified according to international pharmacopeias; USP <71>; Ph. Eur. 2.6.1			
Effective filter area	15.7 cm ²			
Integrity test	2.5 bar			
Membrane thickness	Approx. 150 – 170 μm			
Transparent film	PETG (polyethylene terephthalate)			
Cellulose acetate (CA) membrane				
Pore size	0.45 µm nominal; specified according to international pharmacopeias; USP <71>; Ph. Eur. 2.6.1			
Effective filter area	14.5 cm ²			
Integrity test	1.5 bar			
Membrane thickness	115 – 145 µm			
Transparent film	PETG (polyethylene terephthalate)			

Hydrophobic (sterile venting) filters	
Membrane	0.2 µm polytetrafluoroethylene (PTFE); validated according to HIMA for the retention of Brevundimonas diminuta
Burst pressure of the sterile air filter	At least 6 bar
Housing	Acrylic-based multipolymer
Water permeability pressure penetration pressure	>3 bar
Transparent film	PETG (polyethylene terephthalate)

Sterisart® container	
Upper part lower part	Styrene acrylonitrile (SAN)
Burst pressure of the housing	>5 bar
Max. operating pressure	3 bar at 20 °C
Max. operating temperature	50°C
Capacity	120 mL (50 mL, 75 mL and 100 mL graduated marks)
Red caps	Silicone
Integral membrane	Membrane incorporated in the system by a special clamping technology
Syringe holder	Styrene acrylonitrile (SAN) Sterisart® system, 16469
Female Luer lock and Luer slip connectors	Acrylonitrile butadiene styrene (ABS) Sterisart® system 16478



Accessory Materials Kit	
Wing nuts	Polyethylene (PE)
Tubing and Spikes	
Tubing material and length	PVC, 80 cm; Additionally in available as silicone tubing, order no. 16469; 60 cm
Sampling needle	Polycarbonate and stainless steel (Sterisart® system 16467)
Double-needle spike	Acrylonitrile butadiene styrene (ABS) and stainless steel
Needle for sterile venting	Polypropylene (PP) and stainless steel; PTFE membrane; Housing made of methyl acrylate- butadiene-styrene (MBS)

		6 mm
		26 mm
141 mm	135 mm	90 mm
		6 mm 13 mm



	Length	Outer Diameter
Needle for sterile venting of Sterisart® systems 16467, 16477 and 16468	40 mm	1.6 mm
Needle for Sterisart® system 16467	52 mm	1.5 mm
Needle for Sterisart® system 16468	60 mm	1.5 mm
Double-needle spike (long) for Sterisart® system 16466	41 mm	2.8 mm
Double-needle spike (short) for Sterisart® system 16476	21 mm	2.8 mm
Double-needle spike for Sterisart® systems 16469, 16471, 16475, 16478	35 mm	2.8 mm
Double-needle spike for Sterisart® system 16475	23 mm	2.8 mm
Double-needle spike (short) for Sterisart® system (one connector) 16472	22 mm	2.8 mm
Double-needle spike (short) for Sterisart® system (one connector) 16470	23 mm	2.8 mm
Double-needle spike (long) for Sterisart® system (one connector) 16470	37 mm	2.8 mm

Septum Material (only in systems ending in GSD)		
Protective cap	Polyethylene	
Septum material	Polyisoprene and acryloni trile butadiene styrene (ABS)	

Germany

Sartorius Lab Instruments GmbH & Co. KG Otto-Brenner-Strasse 20 37079 Goettingen Phone +49 551 308 0



Representation For For Further Information, visit www.sartorius.com

USA

Sartorius Corporation 565 Johnson Avenue Bohemia, NY 11716 Phone +1 631 254 4249 Toll-free +1 800 635 2906