



Sterisart®
Sterility Testing
Redefining Standards

Simplifying Progress

SARTORIUS

Your benefits are all in the details:
Small innovations with great effects



1 Double safety for you and your sample: Sterisart® Septum for aseptic sampling and supplementation

A critical step in sterility testing is sampling from or supplementing the media in a canister. You usually do this by piercing a tube. This conventional method is complicated, involves the hazard of injury and increases the risk of false-positive results. This is also a destructive method that precludes the possibility of re-incubating the canisters. There is a better and safer way: septum closures.

The septum technology of our Sterisart® units has been purpose-built to draw samples for rapid sterility testing and has proven its effectiveness for more than ten years as the method of choice for aseptic transfer of liquids. Your sample will remain safely protected against contamination, as the septum ensures a sterile barrier. At the same time, the simple piercing mechanism minimizes your risk of puncture injuries.

2 Experience Easy Handling and User Safety

The quality of the needle plays a decisive role in determining the ease and safety of piercing closures on product containers and rinsing liquid bags, as well as of transferring culture media into the canisters.

Our specially designed and patented dual-needle, with its protective shield and built-in venting filter, ensures safe and easy handling, even when working with cumbersome isolator gloves.



🌐 **Learn more about our septum canisters at:**
<https://bit.ly/3bGkf26>



Start your test directly, without any need for rinsing: Gamma-sterilized and gas-impermeable primary packaging

When you use standard, gas-permeable primary packaging, you first have to go to the trouble of flushing the units with a rinsing liquid inside the isolator to remove all traces of residual gas.

Not so with our gamma-sterilized and gas-impermeable primary packaging. It enables you to omit the rinsing procedure and proceed right away with testing. This will save you not only many steps, but also the additional costs incurred.

3 No fear of foaming: Anti-foaming inlet and venting filter with safety clearance

An air filter is essential for sterile venting. But if it comes in contact with the culture medium, it can easily become clogged. At the same time, it can pose a risk to your sample as a result of microbial growth.

That's why we designed it with a safety clearance of 1 inch, or over 2½ cm, to effectively protect the filter from coming in contact with culture medium, even if foaming or the slightest vibration occurs.

4 Get a better grip: Large, pre-installed filter caps

Working with isolator gloves is a challenge because they considerably affect your feel for grip.

The extra-large, grip-optimized filter caps are directly attached to the vent filters. You'll find this practical – and not only that. The caps connected by a band to the vent filters cannot fall off or touch any surfaces of the isolator so they are safe from contamination.



5 Filter emulsions and other viscous substances the fastest way: Cellulose acetate specialty membranes

Oily substances and emulsions are difficult and take considerable time to filter. Besides that, several rinsing steps are normally needed to remove residual product.

Now there's a better, fast way: Sterisart® CA units with a cellulose acetate membrane deliver the fastest flow rates in

customer tests – convincing performance. As the pioneers of membrane filtration, we successfully developed a special membrane structure that it is less hydrophilic than are conventional cellulose acetate membranes. As a result, you'll save time during each filtration run and require fewer rinsing steps.



Optimize your workflow thanks to fewer manual steps: Just one primary package for all components

Working in an isolator calls for the highest dexterity and concentration – every additional manual step can be quite stressful.

Instead of being sealed in a separate, difficult-to-open package, the easy-to-handle yellow plugs for closing the canisters are right where you need them: in extra pockets directly inside the primary packaging. This facilitates your workflow and lets you stay focused on your core tasks.

The Sterisart® packaging is also colour- and bar-coded for easy identification and traceability.

6 Work with peace of mind knowing you can't go wrong: Two different color-coded tube clamps for clear identification

Making responsible decisions in final quality control for release of safe products can be highly stressful because every step has consequences – especially when you cannot tell one canister apart from the other. You can readily confuse these in busy-day routines.

The two different color-coded clamps on Sterisart® tubing eliminate the risk of confusion, making it easy for you to identify and fill the canisters with the correct media. Now you can work confidently without worrying about such risks.

Sterisart® sterility testing: Discover further benefits when you make the switch!

Choose the one that meets your requirements from over 20 Sterisart® units.

We offer all the adapters to fit your needs:

- Sterisart® units are compatible with any peristaltic pump on the market for sterility testing.
- Our Universal brand peristaltic pump is compatible with all commercially available sterility test units.



Be sure you stay within the limits: Use the peristaltic pump with the lowest particulate emissions

International regulations permit no more than 3,520 particles per m³ in ISO 5 class cleanrooms for sterility testing.

In a head-to-head comparison test at the highest speed, the particulate emissions of our Sterisart® Universal pump were far below this limit and thus even 60 times lower than that of the peristaltic pump of Competitor A.

With the Sterisart® Universal, you can be sure that you will be working in compliance with regulatory standards.

On-Site Demonstration

We would be happy to demonstrate the advantages of our Sterisart® sterility testing products in person. Just contact your regional Sartorius sales specialist.




<https://www.sartorius.com/en/applications/quality-control-testing/microbiological-quality-control/sterility-testing>

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
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 information, visit
www.sartorius.com