



sartorius

Sterisart[®] sterility testing
redefines standards

Minimize all remaining risks
in routine testing



We're sure about our test results – are you?!

Sterisart[®] sterility test units:
Make the switch for convincing proof.



Double safety for you and your sample:

Sterisart® Septum for aseptic sampling and injection

A critical step in sterility testing is sampling from or injecting into a canister. You usually do this by piercing a tube. This conventional method is not only complicated and involves the hazard of injury. It also increases the risk of false-positive results as this compromises the integrity of the closed system. There is a better and safer way: septum closures.

The septum technology of our Sterisart® units has been proving 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.

Your benefits are all in t
Small innovations with g

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.



Start your test directly, without any need for rinsing:

Gamma-sterilized and gas-impermeable primary packaging

When you use standard, gas-impermeable 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 your budget for rinsing liquid.



the details:
great effects

No fear of foaming:

Air 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 cm, or over 2½ inches, to effectively protect the filter from coming in contact with culture medium, even if foaming or the slightest vibration occurs.

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.

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 single extra manual step can add up to stress.

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.



Free Validation Service*

Save time and money in new or revalidation of your products with our Sterisart® units – we'll take the workload off your hands:

Option a)

We will perform validation of your products directly through Sartorius.

Option b)

You can validate your products on your own, and we'll match your expenses by supplying you with the equivalent in free products.

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

Customized Configuration

Get exactly what you need in your sterility test units – you name it, we'll configure it for you:

- Choose your favorite just right for your requirements from over 20 standard units.
- Any special requests? Let's customize your individual test units together with you.

Independent Flexibility

Be independent in the choice of your sterility testing equipment – 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.

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.



www.sartorius.com/sterisart



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

*This is a limited-time offer if Sartorius is selected as a primary supplier of sterility test units for an agreed term per each product validation. All details of our offer are available at www.sartorius.com/sterisart

Specifications subject to change without notice. Copyright Sartorius Lab Instruments GmbH & Co. KG. Printed in the EU on paper bleached without chlorine. Publication No.: SL-1126-e161101 · Order No.: 85037-556-13 · Ver. 11 | 2016