

Why It Is Important to Have a Secondary Supplier for USP <71> | EP 2.6.1 | JP 4.06 Sterility Testing

Supply chains are not insulated from today's risks and can leave you prone to operational disruptions. Your supplier's risk is your risk!

Delays in product release can lead to a build-up of inventory, impact the time to market, the shelf-life of your product, and its accessibility to patients. These challenges can cause significant financial burden and divert valuable resources to deal with such complications. Finding and validating a second supplier during a crisis is by no means a trivial process.

In the vital and highly regulated pharmaceutical industry, a secondary supplier for sterility testing can help you mitigate such uncertainty and avoid an absolute dependence on a single source. Let us help you manage your risks by being prepared!



Sartorius Sterisart® Benefits

Unique Septum Version for Aseptic Sampling

Sample extraction for sub-culturing and identification, or supplementation (inactivation of antibiotics).

Sharp Stainless Steel Needles

Convenient and safe piercing of septum on bottles, bags and vials. The dual-needle houses an in-built sterile venting filter.

Pre-Installed Color-Coded Tube Clamps

For easy and reliable identification.

Product- | Lot Number Identification

Traceability of the product for documentation | security.

Grip Optimized Filter Caps

Extra-large, grip optimized filter caps are tethered to the sterile venting filters to improve speed and convenience.

More Than 20 Canister Types

Gamma sterilized units in gas-impermeable packaging, for testing of ampoules, vials, bags, pre-filled syringes, bottles, lyophilized powders and medical devices.



Sales and Service Contacts

For further contacts, visit
www.sartorius.com

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