SARTURIUS

Product Datasheet

Sartolab[®] Single-Use Polystyrene (PS) Storage Bottles



Product Information

Sartorius single-use polystyrene (PS) storage bottles are designed for safe, secure storage of tissue culture media and sera, buffers, and other aqueous solutions. The bottles are supplied as presterilized and ready to use units in a wide variety of capacities.

- Single-use polystyrene bottles for storage of media, buffers, and other aqueous solutions
- Bottles featuring a low-profile design for easy gripping with one hand
- Ergonomically designed cap to provide a rigid grip and one-hand handling of each bottle

- The bottles with an indented bottom are stackable to save space in the refrigerator and in the freezer
- Caps (45 mm) with airtight seal to minimize the risk of contamination
- The bottles can be used with all Sartolab® RF | BT vacuum filtration units with a 45 mm neck size
- Graduation marks on the bottles ensure accuracy and easy reading
- Designated writing field for clear labeling of samples
- The materials used in the production of the Sartolab® bottles are medical-grade for the highest quality, without any animal origin
- Sterile and endotoxin-free

Specifications

Choice of capacity	150 mL, 250 mL, 500 mL, 1,000 mL
Bottle neck size	45 mm
Color	Clear transparent bottle with bright yellow cap
Compatibility	All Sartolab® RF BT vacuum filtration units with a 45 mm neck
Autoclavable	No
Sterilization method	E-Beam (beta) irradiation
Packaging	Individually packaged, presterilized
Storage temperature*	-20° C to +60 °C
Operating temperature	0° C to +70° C
Operating pressure	-350 to -750 mbar

*Storage at temperatures of -20°C to 0°C is dependent on the liquid type and the storage conditions. We recommend conducting a storage test for the solution before use

Materials

1 Tatorialo					
Bottle	Polystyrene (PS)				
Screw cap	Polyethylene (HDPE)				
Product packaging	PET PE and PE PA multilayer films	٨			
Box for transportation	Corrugated cardboard	۵			

Suitable Liquids

Liquid types

Aqueous solutions, e.g., buffer or cell culture media

- It is not advisable to use the following fluids: Acids, aldehydes, aliphatic amines, aromatic amines, alkalis, esters, hydrocarbons, and ketones
- Fluids and temperatures may impair the properties of the product, e.g., mechanical strength, color, appearance, and dimensional stability

Alcohols may be suitable for some applications. We recommend testing the product with the solution.

Product Symbols

Symbol	Description			
Qty:	Quantity			
REF	Order number			
LOT	Batch number			
Ω	Use by			
STERILE R	Sterilized by irradiation			
②	Do not re-use			

Ordering Information

Order No.	Name	Nominal capacity (mL)	Height (mm)	Diameter (mm)	Weight with cap (g)	Quantity per pkg.
180-22E	Sartolab® bottle, 150 mL	150	77	87	49	12
180-23E	Sartolab® bottle, 250 mL	250	99	87	57	12
180-24E	Sartolab® bottle, 500 mL	500	114	112	87	12
180-25E	Sartolab® bottle 1,000 mL	1000	169	116	133	12

Germany

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



 ⊕ For further information, visit www.sartorius.com