

# Combisart® Kits

Parallel Vacuum Filtration  
With the Combisart®  
Manifold and 3 Glass or  
Polycarbonate Funnels



## Benefits

- Three-way valve allow the vacuum for each filter holder to be sterilely vented
- Low height of the manifold ports is particularly advantageous for working on a clean bench
- Easy-to-use setup
- Ideally suited for larger volume filtration with up to 0.75 L

## Product Information

Made of high-grade stainless steel (B.S. 304S31 | AISI 304); the three-branch manifold accommodates any type of vacuum funnel. Stainless steel three-way valves allow the vacuum for each filter holder to be sterilely vented.

The top and bottom part of the glass filter holders are easily and securely fastened together using the metal clamp. The centering rim on the filter support ensures correct positioning of the membrane filter. The reusable and practical filter holder is made of autoclavable plastic and, thus, ideal for microbiological and analytical testing outside the laboratory.

## Glass and Polycarbonate Material

The Glass Filter Holders are available in either 30 mL or 250 mL with two different filter supports per volume. These are a glass frit support and a PTFE-coated screen support. Both are equipped with a glass funnel, a glass base, a small clamp to tighten upper and lower part and a silicone O-ring in between. The polycarbonate filter holders can be screwed into the manifold and have a funnel capacity of 250 mL. They can be connected to a flushing tube via the hose nipple inlet on top of the holders.

## Applications

Typical applications for the glass filter holders are based in the field of ultracleaning or clarifying of up to liter volumes of solvents for analytical determinations (HPLC). The polycarbonate filter holders are mainly used for sterility testing, when the membrane filter is to be removed from the holder for incubation in culture media.



## Technical Specifications

	Specifications	Description
<b>Combisart®</b>	Dimension (L×H×D)	435 mm × 103 mm × 120 mm
	Max. Operating Pressure	Vacuum or max. 2 bar (29 psi) pressure
	Inlets	TR 20 × 2 female threads
	Outlets	Hose Nipple DN 10
	Sterilization	By autoclaving (134 °C max.) By dry heat (180 °C max.)
<b>Adapters for Glass Filter Holders</b>	Material	Base - AISI 304 stainless steel Stopper - Silicone
	Max. Operating Pressure	Vacuum only
	Outlets	TR 20 × 2 mm male thread
	Sterilization	By autoclaving (134 °C max.) By dry heat (180 °C max.)
<b>Glass Filter Holder 30 mL</b>	Base outlet	12 mm diameter
	Parts and Materials	Borosilicate glass funnel and base, PTFE   glass filter support (16306) or PTFE   stainless steel filter support, coated with PTFE (16315), Silicone O-ring 45 × 3 mm, Anodized aluminum clamp
	Chemical Compatibility	As for glass, PTFE and silicone, Silicone O-ring can be replaced by a fluorelastomer O-ring (00118)
	Funnel Capacity	30 mL
	Max Operating Pressure	Only for vacuum
	Filtration Area	3 cm <sup>2</sup>
	Suitable membrane filter diameter	25 mm
	Sterilization	By autoclaving (134 °C max.) By dry heat (180 °C max.)

	Specifications	Description
<b>Polycarbonate Filter Holder 250 mL</b>	Funnel Capacity	250 mL
	Max Operating Pressure	Vacuum or 2 bar (29 psi) pressure max.
	Filtration Area	12.5 cm <sup>2</sup>
	Suitable membrane filter diameter	47 mm
	Sterilization	By autoclaving (121 °C max.)
<b>Glass Filter Holder 250 mL</b>	Base outlet	115 mm diameter
	Parts and Materials	Borosilicate glass funnel and base, Silicone rubber lid, PTFE   glass filter support (16307) or PTFE   stainless steel filter support, coated with PTFE (16316), Silicone O-ring 45 × 3 mm, Anodized aluminum clamp
	Chemical Compatibility	As for glass, PTFE and silicone, Silicone O-ring can be replaced by a fluorelastomer O-ring (00124)
	Funnel Capacity	250 mL
	Max Operating Pressure	Only for vacuum
	Filtration Area	12.5 cm <sup>2</sup>
	Suitable membrane filter diameter	47 mm   50 mm
	Sterilization	By autoclaving (134 °C max.) By dry heat (180 °C max.)

## Ordering Information

Order No.	Description
16842-KIT-01	Combisart® Kit with three glass filter holders, 30 mL
16842-KIT-02	Combisart® Kit with three glass filter holders, 30 mL and PTFE coated filter support
16842-KIT-03	Combisart® Kit with three glass filter holders, 250 mL
16842-KIT-04	Combisart® Kit with three glass filter holders, 250 mL and PTFE coated filter support
16842-KIT-05	Combisart® Kit with three polycarbonate filter holders, 250 mL



### 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 more information, visit**  
[www.sartorius.com](http://www.sartorius.com)

Specifications subject to change without notice.

© 2024 Sartorius Stedim Biotech GmbH, August-Spindler-Strasse 11, 37079 Goettingen, Germany

Status: 02 | 2024