

Microsart[®] @Filter 100 | 250

Sterile Single-Use Filtration
Units Featuring
a Click-Fit Closure

Benefits

- Fully single-use design and sterile packaging help prevent secondary contamination
- Click-fit closure for effortless funnel removal
- Reliable results as no liquid remains after filtration
- Transparent funnel for easy monitoring of the filtration progress
- Filtration units with widest selection of integrated membrane filter types



Product Information

Microorganisms in liquids are quantified by the membrane filtration method as it allows accurate enumeration of bacteria, yeasts and molds when low counts in a high sample volume are anticipated. In addition, inhibitors, such as antibiotics or preservatives, can be removed by rinsing the membrane with buffer so that microbial growth is not suppressed. Microsart[®] @Filter filtration units have been specifically developed for the detection and enumeration of microorganisms in pharmaceuticals, cosmetics, beverages, water and other liquids.

Description

Microsart® @Filter units combine the benefits of membrane filtration with the advantages of a ready-to-use filtration unit. The sterile filtration unit consists of a funnel, a membrane filter and a base support. It is simply put on a stainless steel manifold in order to directly filter the sample. No decontamination of the stainless steel base or pump adaptor before usage necessary. After the filtration the funnel is exceptionally easy to remove, thanks to its click-fit closure.

Microsart® @Filter is available with either 100 or 250 mL volume funnels. Graduated marks on these units allow accurate sample volumes. In addition, their optimized design and transparency permits thorough rinsing subsequent to filtration so no liquid is retained in the funnel.

Microsart® @Filter – membrane selection

For each application the matching membrane filter type:

- Cellulose Nitrate (= Cellulose Mixed Ester) (CN) for aqueous solutions (pH 4–8)
- Polyvinylidene fluoride (PVDF) for aqueous solutions (pH 1–12), hydrocarbons and several other organic solvents and alcoholic solvents
- Polycarbonate (PC)
- Regenerated Cellulose (RC)

Microsart® @Filter packaging types

- individually packaged (with lid)
- packaged on trays (with lid)
- packaged in bags (without lid)

Microsart® @Media – Forceps-free, Effortless Membrane Transfer

Microsart® @Filter units*, together with the prefilled Microsart® @Media agar dishes, constitute an innovative single-use system for transfer of membranes to the media without needing forceps or forces.

The active lid of the media dish fits perfectly onto the filter unit base. The membrane filter is effortlessly fixed by the integrated adhesive ring in the lid and automatically positioned correctly on the agar plate. Just one turn is all it takes to lock the filter unit in place. Then just incubate and count as usual. Perform sampling on up to incubation in just a few easy steps while minimizing the risk of secondary contamination.

Applications

Microbial enumeration, bioburden testing, microbial limit testing and other microbial test of liquid samples.

A small sampling of the many advantages you will benefit from when using Microsart® @Filter units:

Reliable results

Sterile-packaged	No sterilization at the point of use required
Fully single-use base and funnel	Preparation- and sterilization-free procedure reduces the risk of secondary contamination
Optimized design and materials	No liquid remains after filtration; eliminates the need for rinsing

Easy use

Click-fit closure	Allows effortless and quick funnel removal
-------------------	--

Traceability

Transparent funnel material	Progress of an entire filtration run easy to track
-----------------------------	--

Specifications

Materials	Funnel: Polypropylene Base: ABS (acrylonitrile butadiene styrene) Filter support (PVDF version only): PET nonwoven (polyester) Membrane filter types: <ul style="list-style-type: none">▪ Cellulose Mixed Ester (CN)▪ Polycarbonate (PC)▪ Polyvinylidene fluoride (PVDF)▪ Regenerated cellulose (RC) Choice of various colors, grids and pore sizes
Capacity	100 mL (graduated marks at 20, 50 and 100 mL) 250 mL (graduated marks at 50, 100, 200 and 250 mL)
Filter diameter	47 mm
Filtration area	13.2 cm ²
Max. operating pressure	Vacuum only
Sterilization	Gamma irradiation
Lot certificate	Recovery rate, sterility and performance test

Ordering Information

Microsart® @Filter With CN Filter

Microsart® @Filter 100, with lid, packaged on trays, ideal for use in clean benches

Pore size	Membrane filter* color grid color	Funnel volume	Qty./pkg.	Order no.
0.2 µm	CN white black	100 mL	24	16D01--10-07--TG
0.45 µm, High Flow	CN white black	100 mL	24	16D01--10-H6--TG
0.45 µm, High Flow	CN gray white**	100 mL	24	16D03--10-H6--TG
0.45 µm	CN green dark green	100 mL	24	16D02--10-06--TG

Microsart® @Filter 250, with lid, packaged on trays, ideal for use in clean benches

Pore size	Membrane filter* color grid color	Funnel volume	Qty./pkg.	Order no.
0.2 µm	CN white black	250 mL	16	16D01--25-07--TF
0.45 µm, High Flow	CN white black	250 mL	16	16D01--25-H6--TF
0.45 µm, High Flow	CN gray white**	250 mL	16	16D03--25-H6--TF
0.45 µm	CN green dark green	250 mL	16	16D02--25-06--TF
0.65 µm	CN gray white**	250 mL	16	16D03--25-05--TF

Microsart® @Filter 100, stacked and packaged in bags, ideal for use with Microsart® Funnel Dispenser

Pore size	Membrane filter* color grid color	Funnel volume	Qty./pkg.	Order no.
0.2 µm	CN white black	100 mL	60	16D01--10-07--BL
0.45 µm, High Flow	CN white black	100 mL	60	16D01--10-H6--BL
0.45 µm, High Flow	CN gray white**	100 mL	60	16D03--10-H6--BL
0.45 µm	CN green dark green	100 mL	60	16D02--10-06--BL

Microsart® @Filter 250, stacked and packaged in bags, ideal for use with Microsart® Funnel Dispenser

Pore size	Membrane filter* color grid color	Funnel volume	Qty./pkg.	Order no.
0.2 µm	CN white black	250 mL	48	16D01--25-07--BK
0.45 µm, High Flow	CN white black	250 mL	48	16D01--25-H6--BK
0.45 µm, High Flow	CN gray white**	250 mL	48	16D03--25-H6--BK
0.45 µm	CN green dark green	250 mL	48	16D02--25-06--BK
0.65 µm	CN gray white**	250 mL	48	16D03--25-05--BK

Microsart® @Filter 100, individually sterile packaged, with lid

Pore size	Membrane filter* color grid color	Funnel volume	Qty./pkg.	Order no.
0.45 µm, High Flow	CN white black	100 mL	27	16D01--10-H6-ACG
0.45 µm, High Flow	CN black white	100 mL	27	16D03--10-H6-ACG
0.2 µm	CN white black	100 mL	27	16D01--10-07-ACG

Microsart® @Filter 250, individually sterile packaged, with lid

Pore size	Membrane filter* color grid color	Funnel volume	Qty./pkg.	Order no.
0.45 µm, High Flow	CN white black	250 mL	18	16D01--25-H6-ACF
0.45 µm, High Flow	CN black white	250 mL	18	16D03--25-H6-ACF
0.2 µm	CN white black	250 mL	18	16D01--25-07-ACF

Microsart® @Filter With PVDF Filter

Microsart® @Filter 100, stacked and packaged in bags, ideal for use with Microsart® Funnel Dispenser

Pore size	Membrane filter* color grid color	Funnel volume	Qty./pkg.	Order no.
0.45 µm	PVDF white no grid	100 mL	60	16D04--10-06--BL***

Microsart® @Filter 100, individually sterile packaged, with lid

Pore size	Membrane filter* color grid color	Funnel volume	Qty./pkg.	Order no.
0.45 µm	PVDF white no grid	100 mL	27	16D04--10-06-ACG***

Microsart® @Filter 100, with lid, packaged on trays, ideal for use in clean benches

Pore size	Membrane filter* color grid color	Funnel volume	Qty./pkg.	Order no.
0.45 µm	PVDF white no grid	100 mL	24	16D04--10-06--TG***

Microsart® @Filter With Polycarbonate Filter

Microsart® @Filter 100, stacked and packaged in bags, ideal for use with Microsart® Funnel Dispenser

Pore size	Membrane filter* color grid color	Funnel volume	Qty./pkg.	Order no.
0.45 µm	PC white no grid	100 mL	60	16D06--10-06--BL***

Microsart® @Filter With Regenerated Cellulose Filter

Microsart® @Filter 100, with lid, packaged on trays, ideal for use in clean benches

Pore size	Membrane filter* color grid color	Funnel volume	Qty./pkg.	Order no.
0.45 µm	RC white (w/o grid)	100 mL	24	16D05--10-06--TG***

Microsart® @Filter 100, stacked and packaged in bags, ideal for use with Microsart Funnel Dispenser

Pore size	Membrane filter* color grid color	Funnel volume	Qty./pkg.	Order no.
0.45 µm	RC white (w/o grid)	100 mL	60	16D05--10-06--BL***

Accessories

Microsart® Funnel Dispenser	Order no.
Microsart® Funnel Dispenser for contamination-free removal of individual, sterile Microsart® @Filter units	16A08

Equipment

Manifolds for use with Microsart® @Filter units	Order no.
1 place Microsart Manifold	168M1-MS
2 place Microsart Manifold	168M2-MS
3 place Microsart Manifold	168M3-MS
6 place Microsart Manifold	168M6-MS

* Microsart @Media can be used only with the CN membrane filter version of the Microsart @Filter

** Gray membranes, black after wetting

*** This Microsart® @Filter type cannot be used with Microsart® @Media.

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