



Microsart® @filter 100 | 250

Sterile single-use filter units featuring a click-fit closure

User Benefits

- Fully single-use design and sterile packaging help prevent secondary contamination
- Reliable results as no liquid remains after filtration
- Click-fit closure for fast routine analysis without the risk of leakage
- Transparent funnel for easy monitoring of the filtration progress
- Economical as the unit fits on multibranch manifolds – no added investment needed



Product Information

The key to manufacturing competitive products and maintaining conformity is effective quality assurance. 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. All components of the filtration system used in this method must comply with international guidelines, such as USP, EP or ISO standards.

Description

Ready-to-use, sterile Microsart® @filter units combine a funnel and a gridded membrane filter all in one unit. They have been specifically developed for the detection and enumeration of microorganisms in pharmaceuticals, cosmetics, food, beverages, water and other liquids. For sample volumes of up to 100 mL, Microsart® @filter 100 is the ideal device; larger sample sizes can be easily filtered by Microsart® @filter 250. Graduated marks on these units allow accurate sample volumes. In addition, their optimized design permits thorough rinsing of the system subsequent to filtration. No liquid is retained in the filter funnel. Optimal sealing to the 47 mm Microsart® Base is guaranteed by the bayonett closure. Microsart® @filter units, together with prefilled Microsart® @media agar dishes, constitute an innovative single-use system for transfer of membranes to these media without using tweezers.

Microsart® @vance

The Microsart® @filter units and Microsart® @media prefilled agar dishes are part of the Microsart® @vance product family that stands for cutting-edge products for microbiological analyses based on microbial enumeration. This product line excels in reducing workloads while eliminating the risk of contamination as only a few steps are needed from sampling to incubation.

Microsart® @filter packaging types

Microsart® @filter units are available in a choice of packages: either the filtration units are individually sterile packed in bags, packaged on trays for use in clean benches or in bags for use with the convenient Microsart® Funnel Dispenser. Both single sterile packed as well as the tray version are coming with one lid each.

Applications

Microbial enumeration, particle testing and microscopy

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	Fast for routine analysis; eliminates the risk of leakage
-------------------	---

Economy

Fits on Combisart®	Provides flexibility; no additional investment required
Transparent funnel material	Progress of an entire filtration run easy to track

Specifications

Materials	Funnel: Polypropylene Base: ABS (acrylonitrile butadiene styrene) Membrane filter: cellulose nitrate (cellulose ester); regenerated cellulose; choice of various colors and grids
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, prefilter 40 mm (particle testing only)
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 100, sterile single-use filter units with lid, 47 mm, 100 mL, packaged on trays, ideal for use in clean benches, 24 units

Pore size	Membrane filter* color grid color	Order no.
0.2	CN white black	16D01--10-07--TG
0.45, High Flow	CN white black	16D01--10-H6--TG
0.45, High Flow	CN gray white**	16D03--10-H6--TG
0.45	CN green dark green	16D02--10-06--TG
0.45	RC white (w/o grid)	16D05--10-06--TG***

Microsart® @filter 250, sterile single-use filter units with lid, 47 mm, 250 mL, packaged on trays, ideal for use in clean benches, 16 units

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

Microsart® @filter 100, sterile single-use filter units, 47 mm, 100 mL, stacked and packaged in bags, ideal for use with Microsart® Funnel Dispenser, 60 units

Pore size	Membrane filter* color grid color	Order no.
0.2	CN white black	16D01--10-07--BL
0.45, High Flow	CN white black	16D01--10-H6--BL
0.45, High Flow	CN gray white**	16D03--10-H6--BL
0.45	CN green dark green	16D02--10-06--BL
0.45	RC white (w/o grid)	16D05--10-06--BL***

Microsart® @filter 250, sterile single-use filter units, 47 mm, 250 mL, stacked and packaged in bags, ideal for use with Microsart® Funnel Dispenser, 48 units

Pore size	Membrane filter* color grid color	Order no.
0.2	CN white black	16D01--25-07--BK
0.45, High Flow	CN white black	16D01--25-H6--BK
0.45, High Flow	CN gray white**	16D03--25-H6--BK
0.45	CN green dark green	16D02--25-06--BK
0.65	CN gray white**	16D03--25-05--BK

Microsart® @filter 100, individually sterile packed, 100 ml capacity, with covers, 27 pieces

Pore size	Membrane filter* color grid color	Order no.
0.45 High Flow	CN white black	16D01--10-H6-ACG
0.45 High Flow	CN black white	16D03--10-H6-ACG
0.2	CN white black	16D01--10-07-ACG

Microsart® @filter 250, individually sterile packed, 250 ml capacity, with cover, 18 pieces

Pore size	Membrane filter* color grid color	Order no.
0.45 High Flow	CN white black	16D01--25-H6-ACF
0.45 High Flow	CN black white	16D03--25-H6-ACF
0.2	CN white black	16D01--25-07-ACF

Order no.

Microsart® Funnel Dispenser

Microsart® Funnel Dispenser for contamination-free removal of individual, sterile Microsart® @filter units	16A08
--	-------

Microsart® Base

	Order no.
Base, 47 mm, with frit, stainless steel, for attaching Microsart® @filter units to Combisart® and Microsart® Combi.jet manifolds	1ZU---0002

* CN = Cellulose Nitrate (Cellulose Ester), RC = Regenerated Cellulose

** 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

Europe

Germany

Sartorius Lab Instruments
GmbH & Co. KG
Weender Landstrasse 94-108
37075 Goettingen

Phone +49.551.308.0
Fax +49.551.308.3289

France & Suisse Romande

Sartorius France
2, rue Antoine Laurent de Lavoisier
ZA de la Gaudrée
91410 Dourdan

Phone +33.1.70.62.50.00
Fax +33.1.64.59.76.39

Austria

Sartorius Austria GmbH
Franzosengraben 12
1030 Vienna

Phone +43.1.7965760.0
Fax +43.1.7965760.24

Belgium

Sartorius Belgium N.V.
Leuvensesteenweg, 248/B
1800 Vilvoorde

Phone +32.2.756.06.71
Fax +32.2.253.45.95

Finland & Baltics

Sartorius Biohit Liquid Handling Oy
Laippatie 1
00880 Helsinki

Phone +358.9.755.95.1
Fax +358.9.755.95.200

Hungary

Sartorius Hungária Kft.
Kagyló u. 5.
2092 Budakeszi

Phone +3623.457.227
Fax +3623.457.147

Ireland

Sartorius Ireland Ltd.
Unit 41, The Business Centre
Stadium Business Park
Ballycoolin Road
Dublin 11

Phone +353.1.8089050
Fax +353.1.8089388

Italy

Sartorius Italy S.r.l.
Viale A. Casati, 4
20835 Muggiò (MB)

Phone +39.039.4659.1
Fax +39.039.4659.88

Netherlands

Sartorius Netherlands B.V.

Phone +31.30.60.53.001
Fax +31.30.60.52.917

info.netherlands@sartorius.com

Poland

Sartorius Poland sp.z o.o.
ul. Wrzesinska 70
62-025 Kostrzyn

Phone +48.61.6473830
Fax +48.61.6473839

Russian Federation

LLC "Sartorius ICR" and LLC "Biohit"
Uralskaya str. 4, Lit. B
199155, Saint-Petersburg

Phone +7.812.327.5.327
Fax +7.812.327.5.323

Spain & Portugal

Sartorius Spain, S.A.
Avda. de la Industria, 32
Edificio PAYMA
28108 Alcobendas (Madrid)

Phone Spain +34.902.123.367
Phone Portugal +351.800.855.800
Fax Spain +34.91.358.96.23
Fax Portugal +351.800.855.799

Switzerland

Sartorius Mechatronics Switzerland AG
Ringstrasse 24a
8317 Tagelswangen (ZH)

Phone +41.44.746.50.00
Fax +41.44.746.50.50

U.K.

Sartorius UK Ltd.
Longmead Business Centre
Blenheim Road, Epsom
Surrey KT19 9QQ

Phone +44.1372.737159
Fax +44.1372.726171

Ukraine

LLC "Biohit"
Post Box 440 "B"
01001 Kiev, Ukraine

Phone +380.44.411.4918
Fax +380.50.623.3162

America

USA

Sartorius Corporation
5 Orville Drive, Suite 200
Bohemia, NY 11716

Phone +1.631.254.4249
Toll-free +1.800.635.2906
Fax +1.631.254.4253

Argentina

Sartorius Argentina S.A.
Int. A. Avalos 4251
B1605ECS Munro
Buenos Aires

Phone +54.11.4721.0505
Fax +54.11.4762.2333

Brazil

Sartorius do Brasil Ltda
Avenida Senador Vergueiro 2962
São Bernardo do Campo
CEP 09600-000 - SP- Brasil

Phone +55.11.4362.8900
Fax +55.11.4362.8901

Canada

Sartorius Canada Inc.
2179 Dunwin Drive #4
Mississauga, ON L5L 1X2

Phone +1.905.569.7977
Toll-Free +1.800.668.4234
Fax +1.905.569.7021

Mexico

Sartorius de México S.A. de C.V.
Circuito Circunvalación Poniente
No. 149
Ciudad Satélite
53100, Estado de México
México

Phone +52.5555.62.1102
Fax +52.5555.62.2942

Asia | Pacific

Australia

Sartorius Australia Pty. Ltd.
Unit 5, 7-11 Rodeo Drive
Dandenong South Vic 3175

Phone +61.3.8762.1800
Fax +61.3.8762.1828

China

Sartorius Scientific
Instruments (Beijing) Co., Ltd.
33 Yu An Road, Airport Industrial Park
Zone B,
Shunyi District, Beijing 101300,
P.R.China

Phone +86.10.8042.6300
Fax +86.10.8042.6486

Hong Kong

Sartorius Hong Kong Ltd.
Unit 1012, Lu Plaza
2 Wing Yip Street
Kwun Tong
Kowloon, Hong Kong

Phone +852.2774.2678
Fax +852.2766.3526

India

Sartorius Weighing India Pvt. Ltd.
#69/2-69/3, NH 48, Jakkasandra,
Nelamangala Tq
562 123 Bangalore, India

Phone +91.80.4350.5250
Fax +91.80.4350.5253

Japan

Sartorius Japan K.K.
4th Fl., Daiwa Shinagawa North Bldg.
8-11, Kita-Shinagawa 1-chome
Shinagawa-ku, Tokyo, 140-0001 Japan

Phone +81.3.3740.5408
Fax +81.3.3740.5406

Malaysia

Sartorius Malaysia Sdn. Bhd
Lot L3-E-3B, Enterprise 4
Technology Park Malaysia
Bukit Jalil
57000 Kuala Lumpur, Malaysia

Phone +60.3.8996.0622
Fax +60.3.8996.0755

Singapore

Sartorius Singapore Pte. Ltd
1 Science Park Road,
The Capricorn, #05-08A,
Singapore Science Park II
Singapore 117528

Phone +65.6872.3966
Fax +65.6778.2494

South Korea

Sartorius Korea Ltd.
8th Floor, Solid Space B/D,
PanGyoYeok-Ro 220, Bundang-Gu
SeongNam-Si, GyeongGi-Do, 463-400

Phone +82.31.622.5700
Fax +82.31.622.5799

Thailand

Sartorius (Thailand) Co. Ltd.
129 Rama 9 Road,
Huaykwang
Bangkok 10310

Phone +66.2643.8361-6
Fax +66.2643.8367



◀ www.sartorius.com