



Microsart® e.motion Membrane Filters

Benefits

Reliable colony count

Various filter colors and distinctly visible gridlines provide the best contrast with the colonies

Increased efficiency

New Big Packs allow full shift working without reloading

Visible traceability

All relevant information on each filter box and each membrane filter cell



Product Information

Sartorius membrane filters not only feature the widest selection, three different filter colors, two different diameters and a variety of pore sizes. The various filter colors and distinctly visible gridlines provide the best contrast with the colonies for easy and reliable quantification and identification. The membrane filter band specially designed for the Microsart® e.motion dispenser can be conveniently inserted and easily changed, as needed.

Sartorius' experience in the design and continuous enhancement of membrane filters are also reflected in the results they deliver:

- High recovery rates of microorganisms
- The morphology and color of microorganisms be clearly identified
- Extremely fast filtration rate
- Unique grid lines make counting easy
- Compliance with the currently valid international standards, regulations and guidelines

For complete traceability, the membrane filter type, diameter, lot number and a sequential number (from 250 to 1) are printed on the paper of every membrane cell. The membrane filters' quality and compliance with current standards and regulations, such as ISO 7704 for all 0.45 µm membrane filters, are confirmed by the certificate supplied inside each box of membrane filters.

High Throughput

Your time is limited when you have to handle highly repetitive processes and heavy workloads day by day. In this case an efficient and ergonomic workflow is needed. The new Big Packs with a capacity of 250 membrane filters allow full shift working without the necessity of reloading the Microsart® e.motion dispenser. For efficient workflows and less interruptions!

Benefit from Sartorius' long-standing expertise in microbiological quality control and choose from a wide range of dedicated products for microbial enumeration – and more. Our comprehensive selection of membrane filters, filtration units, filtration systems and culture media will help you reach the highest levels of safety with reproducible, reliable results for in-process and final release testing – while streamlining your workflow and simplifying your testing procedures.

Technical Data

Microsart® e.motion Membrane Filter

Design

The membrane filters have a diameter of 47 or 50 mm, and are white, green or gray, and gridded. The grid divides the filter area into 130 squares; each measures 3.1 × 3.1 mm.

Sterilization

Gamma irradiation

Growth Test acc. to ISO 7704

Bacteria recovery tests performed with Sartorius membrane filters have shown that the growth and development of sensitive bacteria are not adversely affected or inhibited by the grid lines, the sterilization process or any extractable substances in the membrane.

Sterility Test

No growth

Thickness

acc. to DIN 53105, 115–145 µm

Thermal Resistance

130°C max.

Chemical Compatibility

Aqueous solutions (pH 4–8), hydrocarbons and various other organic solvents.

In addition, the certificate included in each package of Sartorius membrane filters confirms that they have been manufactured by applying the latest GMP standards. The development, production and distribution of these filters are subject to our stringent quality management system that has been certified for compliance with DIN|ISO 9001.

Typical Performance Data of Various Membrane Filter Types

Pore size		0.2 µm ¹⁾	0.45 µm ²⁾	0.45 µm ²⁾ High Flow	0.65 µm
Water flow rate per cm ² at 1 bar acc. to DIN 58355	in ml/min	20	70	100	130
Retention of coliforms	in %	100	100	100	n.a.
Recovery rate acc. to ISO 7704 as a batch release criterion	in %	≥ 90	≥ 90	≥ 90	≥ 90

Ordering Information

Order No.	Pore Size	Membrane Filter Color Grid Color	Diameter	Qty. per Pkg.
Big Packs				
114H6Z-47----SJR	0.45 µm High-Flow	White Black	47mm	4 × 250
130H6Z-47----SJR	0.45 µm High-Flow	Gray* White	47 mm	4 × 250
Membrane Filters, Cellulose Mixed Ester (Cellulose Nitrate)				
11407Z-47----SCM	0.2 µm	White Black	47 mm	3 × 100
11407Z-50----SCM	0.2 µm	White Black	50 mm	3 × 100
114H6Z-47----SCM	0.45 µm High-Flow	White Black	47 mm	3 × 100
114H6Z-50----SCM	0.45 µm High-Flow	White Black	50 mm	3 × 100
11406Z-47----SCM	0.45 µm	White Black	47 mm	3 × 100
11406Z-50----SCM	0.45 µm	White Black	50 mm	3 × 100
11404Z-47----SCM	0.8 µm	White Black	47 mm	3 × 100
11404Z-50----SCM	0.8 µm	White Black	50 mm	3 × 100
11403Z-47----SCM	1.2 µm	White Black	47 mm	3 × 100
11403Z-50----SCM	1.2 µm	White Black	50 mm	3 × 100
11402Z-47----SCM	3 µm	White Black	47 mm	3 × 100
139H6Z-47----SCM	0.45 µm High-Flow	White Green	47 mm	3 × 100
13906Z-47----SCM	0.45 µm	White Green	47 mm	3 × 100
13906Z-50----SCM	0.45 µm	White Green	50 mm	3 × 100
13806Z-47----SCM	0.45 µm	Green Dark green	47 mm	3 × 100
13806Z-50----SCM	0.45 µm	Green Dark green	50 mm	3 × 100
130H6Z-47----SCM	0.45 µm High-Flow	Gray* White	47 mm	3 × 100
130H6Z-50----SCM	0.45 µm High-Flow	Gray* White	50 mm	3 × 100
13006Z-47----SCM	0.45 µm	Gray* White	47 mm	3 × 100
13006Z-50----SCM	0.45 µm	Gray* White	50 mm	3 × 100
13005Z-47----SCM	0.65 µm	Gray* White	47 mm	3 × 100
13005Z-50----SCM	0.65 µm	Gray* White	50 mm	3 × 100
13004Z-47----SCM	0.8 µm	Gray* White	47 mm	3 × 100
13004Z-50----SCM	0.8 µm	Gray* White	50 mm	3 × 100
130H6Z-47----SCM	0.45 µm High-Flow	Gray* White	47 mm	3 × 100
Membrane Filters, Other Materials				
15406Z-47----SCM	0.45 µm Polyethersulfone	White	47 mm	3 × 100
18406Z-50----SCM	0.45 µm Regenerated Cellulose	White	50 mm	3 × 100
Dispenser				
Order No.	Description			
16713----BO	Microsart® e.motion dispenser with integrated battery and power supply			
16713----PS	Microsart® e.motion dispenser with power supply only			

* Gray membrane filters turn black when wet

Sales and Service Contacts

For further contacts, visit www.sartorius.com

Europe

Germany

Sartorius Lab Instruments
GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 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
Modecenterstrasse 22
1030 Vienna

Phone +43.1.7965760.0
Fax +43.1.7965760.24

Belgium

Sartorius Belgium N.V.
Rue Colonel Bourg 105
1030 Bruxelles

Phone +32.2.756.06.90
Fax +32.2.481.84.11

Finland & Baltics

Sartorius Biohit Liquid Handling Oy
Laippatie 1
00880 Helsinki

Phone +358.9.755.951
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.
Via Torino 3/5
20814 Varedo (MB)

Phone +39.0362.5557.11
Fax +39.0362.5557.99

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 RUS"
Vasilyevsky Island
5th line 70, Lit. A
199178 St. Petersburg

Phone +7.812.327.53.27
Fax +7.812.327.53.23

Spain & Portugal

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

Phone Spain +34.913.586.095
Phone Portugal +351.800.855.800
Fax Spain +34.913.589.623
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

LLS "Sartorius RUS"
Post Box 440 "B"
01001 Kiev, Ukraine

Phone +380.44.411.4918
Fax +380.50.623.3162

Americas

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
1173 North Service Road West, D4
Oakville, ON L6M 2V9

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.
Libramiento Norte de Tepetzotlan s/n,
Colonia Barrio Tlacateco,
Municipio de Tepetzotlan,
Estado de México,
C.P. 54605

Phone +52.55.5562.1102
Fax +52.55.5562.2942
leadsmex@sartorius.com

Peru

Sartorius Peru S.A.C.
Avenue Alberto del Campo 411
Floor 12 - The Office
15076 - San Isidro, Lima

Phone +51.1.441 0158
Fax +51.1.422 6100

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 (Shanghai) Trading Co., Ltd.
3rd Floor, North Wing, Tower 1
No. 4560 Jinke Road
Zhangjiang Hi-Tech Park
Pudong District
Shanghai 201210, P.R. China

Phone +86.21.6878.2300
Fax +86.21.6878.2882

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
10 Science Park Rd
The Alpha #02-13/14
Singapore Science Park II
Singapore 117684

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