

Microsart® e.motion Dispenser

Sets Things in Motion

Benefits

- **Capacity of up to 250 membrane filters**
Efficient workflow, fewer interruptions for reloading
- **3 operating modes**
Highest flexibility of operation
- **Cableless operation and portability**
Easy to transport for convenient handling



Product Information

The ability to accurately detect and quantify microorganisms in liquid samples is of prime importance for any quality control laboratory. This is why membrane filtration is the established method of choice for liquid testing as it provides reliable and reproducible results. Facilitate your microbiological testing procedures with the new Microsart® e.motion dispenser: The membrane filters are released from their sterile packaging fully automatically at the touch of a button, by hands-free motion detection or even by using a foot switch: the choice is yours!

Applications

Microbiological quality control of liquid samples:

- Beverages (beer, wine, soft drinks, bottled water)
- Cosmetics
- Pharmaceutical analysis (WFI, purified water, microbial limit testing of non-sterile products, bioburden testing)
- Environmental analysis (water monitoring)

The Microsart® e.motion dispenser ensures efficient and ergonomic workflows for reliable and fast membrane filter dispensing. Supplied along with the new Big Packs of 250 membrane filters, this convenient dispenser eliminates the need for frequent reloading of filters. The clean, compact design of the dispenser allows quick and easy cleaning using standard disinfectants. In addition, the Microsart® e.motion dispenser offers the easiest and most flexible membrane filter dispensing: choose among three different operating modes. Whichever mode you opt for, 100% performance is always guaranteed.

Choose the mode most convenient for you:

- Push button
- Foot switch or
- Sensor mode

The dispenser's low weight and convenient handle make it easy to remove or transport. Moreover, the battery-operable Microsart® e.motion dispenser model operates without a cable, providing higher flexibility for work areas without a power outlet and eliminates cable clutter.

A window in the dispenser maintains the visibility of the information on the membrane filter band, such as lot number, membrane filter type and the quantity of filters. This ensures better scheduling of tests and easy traceability.

Membrane Filters

Microsart® e.motion membrane filters give you the highest flexibility: they can be used with most commercially available membrane filter dispensers.

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

The new Big Packs allow enable you to work a complete shift without the need to stop and reload membrane filters in the Microsart® e.motion dispenser. 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 Specifications

Dimensions (L x H x W)	233 mm x 230 mm x 172 mm 9.2" x 9" x 6.8"
Weight	
Model 16713----PS	2.4 kg 5.3 lbs.
Model 16713----BO	2.6 kg 5.7 lbs.
Operating voltage	100 V 240 V
Frequency	50 – 60 Hz
Max. power consumption	23 W
Certificates	CE mark EMC Directive EN 61326-1 EN 61010

Ordering Information

Order No.	Pore Size	Membrane Filter Color Grid Color	Diameter	Qty. per Pkg.
Big Packs - Membrane Filters, Cellulose Mixed Ester (Cellulose Nitrate)				
114H6Z-47----SJR	0.45 µm High-Flow	White Black	47mm	4 x 250
130H6Z-47----SJR	0.45 µm High-Flow	Gray* White	47mm	4 x 250
Membrane Filters, Cellulose Mixed Ester (Cellulose Nitrate)				
11407Z-47----SCM	0.2 µm	White Black	47 mm	3 x 100
11407Z-50----SCM	0.2 µm	White Black	50 mm	3 x 100
114H6Z-47----SCM	0.45 µm High-Flow	White Black	47 mm	3 x 100
114H6Z-50----SCM	0.45 µm High-Flow	White Black	50 mm	3 x 100
11406Z-47----SCM	0.45 µm	White Black	47 mm	3 x 100
11406Z-50----SCM	0.45 µm	White Black	50 mm	3 x 100
11404Z-47----SCM	0.8 µm	White Black	47 mm	3 x 100
11404Z-50----SCM	0.8 µm	White Black	50 mm	3 x 100
11403Z-47----SCM	1.2 µm	White Black	47 mm	3 x 100
11403Z-50----SCM	1.2 µm	White Black	50 mm	3 x 100
11402Z-47----SCM	3 µm	White Black	47 mm	3 x 100
139H6Z-47----SCM	0.45 µm High-Flow	White Green	47 mm	3 x 100
13906Z-47----SCM	0.45 µm	White Green	47 mm	3 x 100
13906Z-50----SCM	0.45 µm	White Green	50 mm	3 x 100
13806Z-47----SCM	0.45 µm	Green Dark green	47 mm	3 x 100
13806Z-50----SCM	0.45 µm	Green Dark green	50 mm	3 x 100
130H6Z-47----SCM	0.45 µm High-Flow	Gray* White	47 mm	3 x 100
130H6Z-50----SCM	0.45 µm High-Flow	Gray* White	50 mm	3 x 100
13006Z-47----SCM	0.45 µm	Gray* White	47 mm	3 x 100
13006Z-50----SCM	0.45 µm	Gray* White	50 mm	3 x 100
13005Z-47----SCM	0.65 µm	Gray* White	47 mm	3 x 100
13005Z-50----SCM	0.65 µm	Gray* White	50 mm	3 x 100
13004Z-47----SCM	0.8 µm	Gray* White	47 mm	3 x 100
13004Z-50----SCM	0.8 µm	Gray* White	50 mm	3 x 100

Membrane Filters, Other Materials

15406Z-47----SCM	0.45 µm Polyethersulfone	White	47 mm	3 x 100
18406Z-50----SCM	0.45 µm Regenerated Cellulose	White	50 mm	3 x 100
28006Z-47----SCM	0.45 µm PVDF	White	47 mm	3 x 100

* Gray membrane filters turn black when wet

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

Accessories	
Order No.	Description
1ZE----0028	Pedal switch for Microsart® e.motion dispenser. Continuous adjustment, single-pole, 6.0 (2.5) A 250 V, with built-in strip terminal and VDE* seal of approval. Pedal plate and housing made of thermoplastic material; antiskid thanks to 4 rubber feet. Protection rating: IP 22 according to DIN 40050
SB-12-01-0277	Holding clamp
16625	Stainless steel tweezers with blunt-edged tips


* VDE = Association of German Engineers

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 further contacts, visit
www.sartorius.com