

The new Microsart[®] e.motion Dispenser – Sets Things in Motion

Simplifying Progress



To start click here



Membrane filters on demand

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, hands-free or by foot switch: the choice is yours!



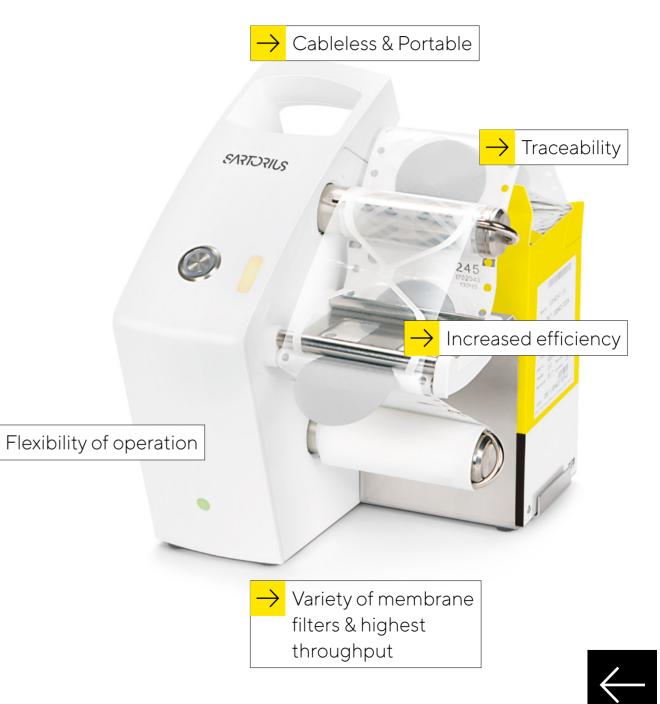




Speed and Convenience in Microbial Enumeration

In a high-throughput production environment, you need to complete your microbiological analyses fast, efficiently and reliably. Any contaminants in your liquid sample will have a direct impact on consumers' health and your product liability.

The fully automated Microsart[®] e.motion dispenser offers efficient and ergonomic workflows for reliable and fast membrane filter dispensing.





Flexibility of operation

The Microsart[®] e.motion dispenser offers the easiest and most flexible way of membrane filter dispensing.

You have the freedom of choice between 3 different operation modes! Easily switch between the operating modes just by pressing the button. The type of operation mode will be indicated by light.

Push-button

Simply push the button whenever you are ready for the next filtration to start.



- push-buttonsensor mode or
- foot-switch

Functional capability always 100% guaranteed!

Choose what is most convenient for you:





Sensormode

No-touch operation - one single movement over the sensor and a membrane filter is automatically dispensed.

Foot-switch

Hands-free operation with the foot-switch option.









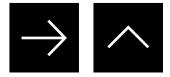
Sartorius offers the widest selection of membrane filters: three different filter colors, a varity of pore sizes and distinctly visible gridlines. Making guantification and identification of the colonies as easy and reliable as possible.

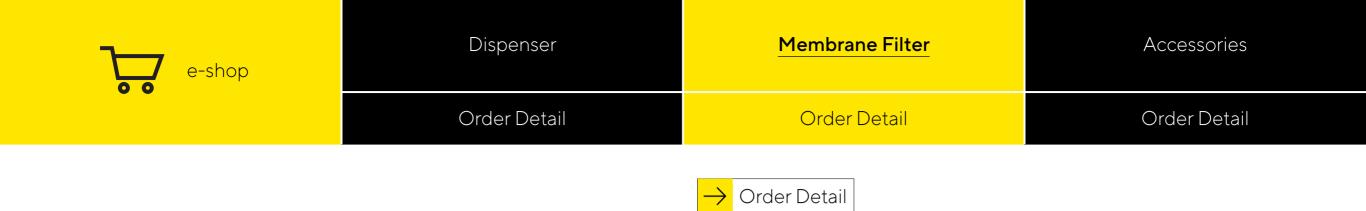
The membrane filter band specially designed for the Microsart[®] e.motion dispenser can be conveniently inserted, and changed easily as needed. In addition you have the highest flexibility: the Microsart® e.motion membranefilters are useable with most commercially available membrane filter dispensers!

High Throughput

The new Big Packs with a capacity of up to 250 membrane filters allow full shift working without the necessity of reloading the Microsart® e.motion dispenser. For efficient workflows and less interruptions!







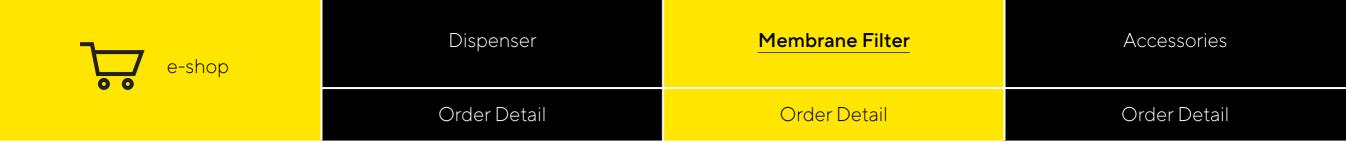
Increased efficiency

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. Save more than 35% of your time by using Microsart e.motion membranes.

The new Big Pack Microsart[®] e.motion membrane filters with a capacity of 250 filters allow full shift working without the necessity of reloading the Microsart[®] e.motion dispenser. For efficient workflows and less interruptions!







Traceability

For complete traceability, the membrane filter type, diameter, lot number and a sequential number are printed on the paper of each membrane cell. In addition, data, such as sterilization and expiration date, pore size and bar code, make handling of these membranes easy and flexible.

The membrane filters' 100% quality and compliance with current standards and regulations, such as ISO 7704 for all 0.45 µm membrane filters, are confirmed in a certificate coming with each box of membrane filters.









Membrane filter overview for Water testing

Correct choice according to current ISO standards

All gridded 0.45 μm Sartorius filters are ISO 7704 conform

Organism	Guideline	Filters per guideline	Sartorius dispenser-ready filters
Enterococci	ISO 7899 - 2:2000	0.45 µm	114H6Z-47SCM (CN, white, black grid)
E. Coli	ISO DIS 9308-1:2012	0.45 µm Cellulose-Esters	- 139H6Z-47SCM (CN, white, green grid) 114H6Z-47SJR (CN, white, black grid, big pack)
Legionella	ISO 11731:2017	direct placing on culture media: 0.2 or 0.45µm black Cellulose Nitrate Membranes For Concentration and Eluation: 0.2µm PC or PESU Membranes	130H6Z-47SCM (CN, 0.45μm, black, white grid) 130H6Z-47SJR (CN, 0.45μm, black, white grid, big pack)
Clostridium perfringens	ISO DIS 14189:2013	0.45 µm	139H6Z-47SCM (CN, white, green grid) 114H6Z-47SCM (CN, white, black grid) 114H6Z-47SJR (CN, white, black grid, big pack)
Pseudomonas aeruginosa	ISO 16266:2006	0.45 µm	114H6Z-47SCM (CN, white, black grid) 114H6Z-47SJR (CN, white, black grid, big pack)
Campylobacter	ISO 17995:2005	0.45 µm; Cellulose-Esters	114H6Z-47 SCM (CN, white, black grid) 114H6Z-47 SJR (CN, white, black grid, big pack)
Salmonella	ISO 19250:2010	0.45 µm	114H6Z-47SCM (CN, white, black grid) 114H6Z-47SJR (CN, white, black grid, big pack)
Anaerobes (Clostridia)	ISO 6461-2:1986	0.2 µm	11407Z-47SCM (CN, white, black grid)
Bacteriophages	ISO 10705 part 1-4	0.2 µm	11407Z-47SCM (CN, white, black grid)









Cableless & Portable

The dispenser's low weight and the convenient handle make it easy to remove or transport.

Moreover the Microsart[®] e.motion dispenser with battery operation can work cableless offering higher flexibility for working areas without power outlet. This allows convenient handling and eliminates cable clutter inside the laminar airflow cabinet.





Microsart[®] e.motion dispenser

Choose between the battery-operated Microsart[®] e.motion dispenser which includes a long-lasting lithium accu or the version with power supply operation only.

Description	Order Number
Microsart [®] e.motion dispenser, battery operation	16713 BO
Microsart [®] e.motion dispenser, power supply only	16713 PS

Dimensions & Weight

Dimensions: (L×W×H)mm: 233×172×230

Weight: 16713----BO: 2.4 kg 16713----PS: 2.6 kg







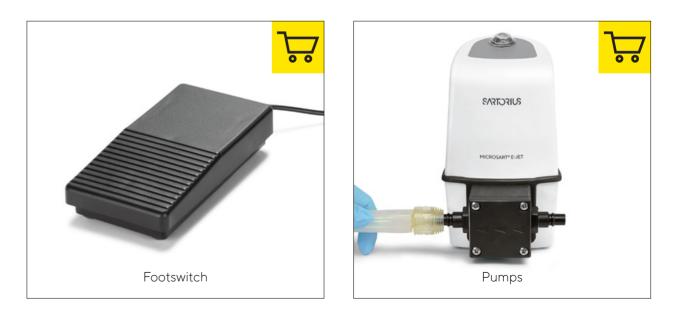
The right equipment for your application

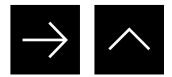
The Microsart[®] e.motion dispenser meets all requirements placed on advanced equipment for microbial enumeration. In addition you can find everything you need for your liquid filtration: from tweezers to vacuum pumps and manifolds.

Choose the products best suited for your application from our comprehensive range of membrane filters, filtration units and filtration systems to ensure:

- Total ease of use
- Quality results
- Excellent cost-effectiveness









Order Detail Accessories

Facilitate your microbiological testing procedures: Choose from our wide selection of manifolds and accessories as well as ready-to-use disposable units and funnels perfectly tailored to meet your specific requirements.

Please use the product configurator to find the equipment that meets your needs.

Description	Order Number
Foot switch for Microsart® e.motion dispenser	1ZE0028
Stainless-steel tweezers	16625
Microsart [®] e.jet Pump	166MP-4
Microsart [®] 3-branch manifold	168M3-MS
Microsart [®] 6-branch manifold	168M6-MS
Microsart [®] 100 ml Funnel	16A0710N
Microsart [®] 250 ml Funnels	16A0725N

