

# Microsart® Funnel 100 | 250

The sterile disposable  
filter funnel



## Benefits

- Reliable results – Use a new, sterile funnel for each test for certain avoidance of cross contamination!
- Time-saving – Just change the funnel, rather than spending time sanitizing it!
- Simpler handling – No more holding of hot funnels! And, you can see when filtration has been completed, particularly useful when using manifolds in routine testing.

In microbiological quality control, sterility of the equipment used for processing samples is a necessary basic requirement. The re-usable funnels made of stainless steel or other materials which are used for membrane filtration are usually sanitized between samples by flaming or with hot water. Both of these methods can be insufficiently reliable when not properly performed. Alternatively, the funnels could be sterilized by autoclaving, but this is too laborious for routine use. A disposable filter funnel is the ideal combination of reliability and time saving.

## Description

Microsart® Funnel 100 | 250 are sterile 100 mL and 250 mL respectively plastic funnels. They allow quick performance of the filtrations required in the routine testing of water, food and beverages, pharmaceutical and cosmetic products.

A Sartorius gridded membrane is placed on a stainless steel filter support. A Microsart® Funnel is simply and practically fitted on. The sample is filtered.

The funnel is made of polypropylene and thus is elastic enough for optimal sealing with a Click-Fit fastening. Graduations are marked as 20, 50 and 100 mL and 50, 100, 200 and 250 mL respectively for sufficiently exact sample volumes. The large inner diameter ensures a high flow rate. The optimized shape allows a thorough rinsing of the system subsequent to filtration. No liquid is retained in the filter funnel.

### Microsart® Funnel Dispenser

Funnel dispenser for secure removal of single, sterile Microsart® Funnel.

### Microsart® Stainless Steel Manifolds

Stainless steel manifolds, 1-, 2-, 3- and 6-branch. Made of high-grade stainless steel.

Stainless steel three-way valves (taps) allow the vacuum for each filter holder to be individually controlled and each holder to be sterilely vented.

## Order Information

### Microsart® Funnel and Accessories

#### Microsart® Funnel 100 | 250

Order No.	Order information	No.   pack
16A07--10-----N	Microsart® Funnels, sterile in 5 sealed bags	100
16A07--25-----N	Microsart® Funnels, sterile in 6 sealed bags	96

#### Accessories

Order No.	Description
16A08	Microsart® Funnel Dispenser
168M1-MS	Microsart® 1-branch stainless steel manifold
168M2-MS	Microsart® 2-branch stainless steel manifold
168M3-MS	Microsart® 3-branch stainless steel manifold
168M6-MS	Microsart® 6-branch stainless steel manifold

Microsart® Funnel 100 | 250 are a further Sartorius step towards quicker and simpler microbiological testing. They supplement products such as:

### Sartorius Stedim Biotech Nutrient Pad Sets

Nutrient Pads are 47 mm diameter absorbent pads which have been uniformly impregnated with a standard culture medium, then dried and sterilized in Petri dishes. A simple wetting with sterile water reactivates them for use. They are supplied as a set, complete with sterile, individually packaged membrane filters, of pore size and colour appropriate to the requirements of the individual medium, for optimum colony growth and convenience of colony counting.

A special packaging in aluminum bags protects the sensitive formula constituents of the nutrient pads during transport and storage from fluctuations in humidity and temperature. As a result, it guarantees the high quality of the NPS throughout their entire shelf life of 24 months.

#### The following are three examples of the 30 different types available

Ready to use units	No.   pack	Order No.
Endo Nutrient Pad Sets for E. coli and coliforms	100	14053--47-----N
Standard Nutrient Pad Sets for colony counts	100	14055--47-----N
Wort Nutrient Pad Sets for yeasts and moulds	100	14058--47-----N

## The Microsart® e.motion Dispenser

Sterile, individually packed membranes are the preferred type for microbiological testing. The Microsart® e.motion Dispenser releases the membrane filters from their sterile packaging fully automatically at the touch of a button or hands-free – a dispensing operation is triggered when the optical sensor detects approaching tweezers.

The membrane filters suitable for use in the Microsart® e.motion are sterile-sealed, without protective paper on top of each filter, in a specially designed individual package on a band. Various types of different membrane filters are available. Each box contains 100 membranes and it is designed so that it is easy to open and seal for storage.

## Order Information

### Microsart® e.motion Dispenser and Membrane Filters

Order No.	Description
16713----PS	Microsart® e.motion Dispenser, power supply
16713----BO	Microsart® e.motion Dispenser, Battery operated

Order No.	Pore size	Color Grid Diameter
11407Z-47----SCM*	0.2 µm	White   Black 47 mm
13006Z-47----SCM*	0.45 µm	Gray   White 47 mm
114H6Z-47----SCM*	0.45 µm	High Flow White   Black 47 mm
139H6Z-47----SCM*	0.45 µm	High Flow White   Green 47 mm
11406Z-47----SCM*	0.45 µm	White   Black 47 mm
13806Z-47----SCM*	0.45 µm	Green   Dark green 47 mm
13906Z-47----SCM*	0.45 µm	White   Green 47 mm
13005Z-47----SCM*	0.65 µm	Gray   White 47 mm
13004Z-47----SCM*	0.8 µm	Gray   White 47 mm
11403Z-47----SCM*	1.2 µm	White   Black 47 mm

Other types on request.

\*Materials supplied: Package with 3 boxes, each containing 100 membranes individually, sterile packaged


Details of these and other Sartorius Products for microbiological quality control are available on request.

**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 information, visit  
[www.sartorius.com](http://www.sartorius.com)