

Microsart® @media

Advanced System for Touch-free Membrane Transfer



Benefits

Innovative Membrane Transfer Concept

- Easy-to-use system
- Safe and reliable
- Time-saving

Product Information

The key to manufacturing competitive products and maintaining compliance is effective quality assurance and control in the highly regulated pharmaceutical industry. This is why microorganisms in liquids are quantified by the membrane filtration method. This method allows accurate quantification of bacteria, yeasts and molds when low counts in a high sample volume are anticipated. The method of transferring the membrane filter onto an agar medium is a critical step that can be a source of secondary contamination and lead to false-positive test results. Therefore, it is essential to reduce the risk of exogenous contamination entailed by the use of tweezers to transfer the membrane.

Description

The Microsart® @filter units and Microsart® @media introduce an innovative concept for membrane transfer to agar media. Microsart® @media are agar media dishes for Microbial Limit Testing. They are pre-filled with different agar media types, sterile-packaged and ready to use together with Microsart® @filter filtration units. Moreover, the Microsart® @media features an innovative patented lid. This active lid enables touch-free transfer of the membrane, without using any tweezers. In addition, the convenient, liftable interior lid provides easy access to select colonies after incubation for further analysis.

Applications

Microsart® @media enhance the safety of microbiological quality control in the pharmaceutical and biotech industries for

- Microbial Limit Testing according to the USP (Chapter <61>) and EP (Chapter 2.6.12), JP 4.05 part 1
- Bioburden Testing
- Water analyses as Purified Water or Water for Injection
- Efficient, advanced workflows in quality assurance laboratories

Innovative Membrane Transfer Concept

The active lid of Microsart® @media enables touch-free handling of the membrane and, therefore, reduces the risk of secondary contamination.

Easy-to-Use System

As a result of the combined development of the Microsart® @filter and Microsart® @media, the active lid of Microsart® @media fits perfectly onto the Microsart® @filter base and thus allows effortless and reliable transfer of the membrane filter onto the agar. Moreover, the innovative click-fit closure of the Microsart® @filter permits fast and easy removal of the funnel after filtration. This eliminates the risk of repetitive strain injury and the need for expensive additional devices to facilitate removal.

Safe and Reliable

Touch-free membrane transfer rules out membrane manipulation and handling, thus minimizing major sources of secondary contamination. This results in best growth conditions, outstanding recovery rates and reliable results.

Time-saving

The Microsart® filtration units are ready to use. In combination with the Microsart® @media membrane transfer concept, just a few quick steps are needed to proceed from sampling to incubation. As a result, this saves time and the cost of labor, while delivering more reliable results.

Summary

Microsart® @media stands for an innovative membrane transfer concept for microbiological analysis based on microbial enumeration. This product line excels in minimizing the risk of secondary contamination as just a few work-saving, touch-free steps are all it takes to accelerate your workflow. Microsart® @media meet the most stringent quality assurance standards, offering a convincing solution for reliable results and simple, time-saving handling.

Technical Data

Dimensions	Dish diameter	68.8 mm
	Dish height	14.9 mm
	Agar area	13.2 cm ²
Materials	Polypropylene	
	Agar media: R2A, TSA, Sabouraud, Plate Count Agar, Sabouraud+ (with Lecithin and Polysorbate 80), TSA+ (with Lecithin and Polysorbate 80)	
	Inhibitor-free glue	
Sterilization	Gamma irradiation within a range of 13.9 kGy to 25.0 kGy	
Lot certificate	Sterility, growth promotion, pH	
Shelf life	36 weeks at 2 °C to 8 °C	

Ordering Information

Microsart® @media prefilled agar media dishes, sterile double packaged and ready-to-use; quantity of 100 per box with 10 bags, each containing 10 media dishes

Media type	Target microorganisms	Order no.	Typical incubation time and temperature
Microsart® @media TSA (Tryptic Soy Agar)	Aerobic and anaerobic flora	14313--47----ACN	48 to 72 hrs. (USP) or 1 to 5 days (EP) at 30 °C to 35 °C
Microsart® @media SDA (Sabouraud Dextrose)	Yeasts and molds	14314--47----ACN	5 to 7 days at 20 °C to 25 °C
Microsart® @media R2A	Stressed heterotrophic bacteria in water	14322--47----ACN	5 to 7 days at 20 °C to 28 °C
Microsart® @media TSA (Tryptic Soy Agar) with Lecithin & Polysorbate	Heterotrophic microorganisms	14315--47----ACN	48 to 72 hrs. (USP) or 1 to 5 days (EP) at 30 °C to 35 °C
Microsart® @media SDA (Sabouraud Dextrose Agar) with Chloramphenicol	Yeasts and molds	14316--47----ACN	5 to 7 days at 20 °C to 25 °C
Microsart® @media PCA (Plate Count Agar)	Total count of aerobic bacteria in water	14317--47----ACN	48 to 72 hrs. at 32 °C to 35 °C (see also APHA water)

Accessories

	Order no.
Microsart® e.jet transfer pump	166MP-4
Microsart 2 branch Manifold	168M2-MS
Microsart 3 branch Manifold	168M3-MS
Microsart 6 branch Manifold	168M6-MS
Silicone pressure tubing, for pressure side; length in meters must be specified	1ZAS--0007
Venting filter 0.2 µm for Microsart Manifolds	S7575-----FXOSK

Microsart® @filter 100, sterile single-use filtration units with lid, 47 mm, 100 mL, packaged on trays, ideal for use in clean benches, 24 units

Pore size in µm	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

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

Pore size in µm	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

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

Pore size in µm	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

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

Pore size in µm	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-packaged filtration units, 100 ml capacity, with protective cover, 27 units

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

Microsart® @filter 250, individually sterile-packaged filtration units, 250 ml capacity, with protective cover, 18 units

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

* CN = Cellulose nitrate ** gray membranes; black after wetting

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