

Touch-Free Membrane Transfer Using Microsart® @filter and @media

Quality control (QC) is a crucial component of drug manufacturing, ensuring the safety and quality of pharmaceutical products. A key component of QC is the detection and quantification of microbial contaminants. Membrane filtration is the regulatory-preferred method for microbiological QC in pharmaceutical and biotechnological applications.

Microsart® @filter are ready-to-use, sterile filtration units consisting of a funnel, a filter base and a membrane filter. They have been specifically developed for the detection and enumeration of microorganisms in pharmaceuticals, cosmetics, food, beverages, water and other liquids. Microsart® @media are pre-filled media dishes with different agar media types, sterile packaged and ready-to-use. When used in conjunction with Microsart® @filter, the Microsart® @media system offers a unique touch-free transfer mechanism, that effortlessly positions the membrane filter onto the agar plate, minimizing the risk of secondary contamination.



Instructions for Using Microsart® @media

1 Remove the Funnel



Do not close the vacuum until the funnel has been removed.



2 Pick Up the Microsart® @media Plate Adhesive Lid



3 Apply the Lid onto the Membrane Filter



To ensure perfect uptake of the membrane filter, the lid should be placed completely straight onto the base from above.



4 Lift the Lid, Including the Membrane Filter



5 Transfer the Membrane Filter onto the Microsart® @media Dish



Position the lid on the Microsart® @media dish at a slight angle before placing the membrane filter completely onto the agar. If something goes wrong, you can simply repeat the process as often as needed. Rotate the lid clockwise to close the locking mechanism.



Subsequently, the Microsart® @media plates can be placed in the incubator and incubated upside down at the required temperature.



[Click here to watch the full video](#)

Easy Colony Enumeration

Microsart® @media's convenient, liftable interior lid provides easy access to colonies after incubation.



[Click here to watch the full video](#)

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
3874 Research Park Dr.
Ann Arbor, MI 48108
Phone +1 734 769 1600



For further information, visit

www.sartorius.com