

# For Efficient Workflows in Microbiological Quality Control: Facilitated Microbial Enumeration Testing for Hundreds of Samples per Day



## How Do You Ensure Efficient Workflows while Dealing with Hundreds of Samples per Day?


A high number of samples during the daily routine analyses may result in:

- Increased manual handling efforts
- Crowded working space (filtration equipment, nutrient media, funnels, membrane filters & samples)
- Interruptions of the workflow (reloading the dispenser & the LAF)


Still, you need to complete your highly repetitive microbiological release testing fast, efficiently and reliably...




# For More Efficient Workflows and Less Interruptions in Your Routine Microbial Quality Control Testing

Need more Space in your working area for samples, funnels & media? 


Want less manual dispenser Reloading? 

Need Convenience in your working area? 

Want Full Shift working without interruptions? 

Want Flexibility of operation? 

Need Traceability? 

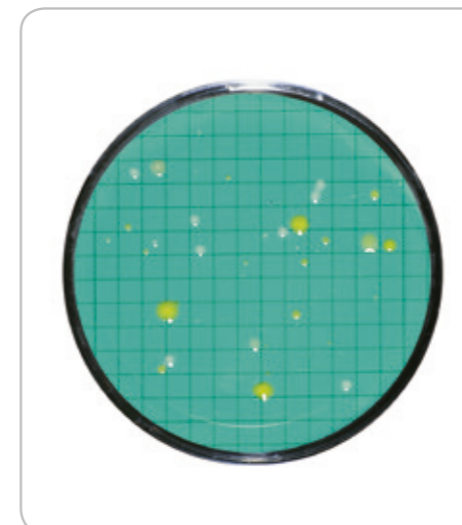
Interested in working Guideline Conform? 



## Benefit from Increased Efficiency and Smooth Workflows

The Microsart® e.motion Dispenser & Big Pack Filters will improve your highly repetitive and heavy workload by streamlining your workflow and simplifying your testing procedure.

- Continuous full shift working without interruptions for reloading the dispenser
- More space in the LAF for your samples due to the dispenser's small footprint and the increased capacity of the Big Pack Membrane Filter boxes and therefore ...
- ... saving you precious time as less reloading of the LAF is needed
- Guideline conform testing with the certified Microsart® e.motion membrane filters



## Microsart® e.motion Dispenser & Big Pack Filters: The Perfect Combination for Smooth and Most Efficient Workflows in a QC Lab with High Routine Sample Throughput.



With the compact design of the dispenser and the increased capacity of the Big Pack the new Microsart® e.motion product range facilitates your highly repetitive processes and heavy workloads day by day:

- Reduces the time spend for reloading the dispenser and the LAF
- Saves valuable space which is required for samples and filtration equipment
- Improves and simplifies your testing procedure by automatic, touch-free membrane filter dispensing

## Additional Information

Benefit from Sartorius' long-standing expertise in microbiological quality control and find out more about our comprehensive selection of membrane filters, dispensers and other dedicated products for microbial enumeration.



See attached file  
Data\_Microsart\_Emotion\_SM-2005-e



See attached file  
Data\_Microsart\_Emotion\_Membranes\_SM-2006-e



See attached file  
Broch\_Microsart\_Emotion\_Dispenser\_SM-1506-e



See attached file  
Broch\_Quality\_Control\_Guidelines\_SM-1505-e



# Sales and Service Contacts

For further contacts, visit [www.sartorius.com](http://www.sartorius.com)

## Europe

### Germany

Sartorius Lab Instruments  
GmbH & Co. KG  
Otto-Brenner-Strasse 20  
37079 Goettingen  
Phone +49.551.308.0

### France & Suisse Romande

Sartorius France  
2, rue Antoine Laurent de Lavoisier  
ZA de la Gaudrée  
91410 Dourdan  
Phone +33.1.70.62.50.00

### Austria

Sartorius Austria GmbH  
Modcenterstrasse 22  
1030 Vienna  
Phone +43.1.7965760.0

### Belgium

Sartorius Belgium N.V.  
Rue Colonel Bourg 105  
1030 Bruxelles  
Phone +32.2.756.06.90

### Finland & Baltics

Sartorius Biohit Liquid Handling Oy  
Laippatie 1  
00880 Helsinki  
Phone +358.9.755.951

### Hungary

Sartorius Hungária Kft.  
Kagyló u. 5.  
2092 Budakeszi  
Phone +3623.457.227

### Ireland

Sartorius Ireland Ltd.  
Unit 41, The Business Centre  
Stadium Business Park  
Ballycoolin Road  
Dublin 11  
Phone +353.1.8089050

### Italy

Sartorius Italy S.r.l.  
Via Torino 3/5  
20814 Varedo (MB)  
Phone +39.0362.5557.11

### Netherlands

Sartorius Netherlands B.V.  
Phone +31.30.60.53.001  
[info.netherlands@sartorius.com](mailto:info.netherlands@sartorius.com)

### Poland

Sartorius Poland sp.z o.o.  
ul. Wrzesinska 70  
62-025 Kostrzyn  
Phone +48.61.6473830

### Russian Federation

LLC "Sartorius RUS"  
Vasilyevsky Island  
5<sup>th</sup> line 70, Lit. A  
199178 St. Petersburg  
Phone +7.812.327.53.27

### Spain & Portugal

Sartorius Spain, S.A.  
Avda. de la Industria, 32  
Edificio PAYMA  
28108 Alcobendas (Madrid)  
Phone Spain +34.913.586.095  
Phone Portugal +351.800.855.800

### Switzerland

Sartorius Mechatronics Switzerland AG  
Ringstrasse 24a  
8317 Tagelswangen (ZH)  
Phone +41.44.746.50.00

### U.K.

Sartorius UK Ltd.  
Longmead Business Centre  
Blenheim Road, Epsom  
Surrey KT19 9QQ  
Phone +44.1372.737159

### Ukraine

LLS "Sartorius RUS"  
Post Box 440 "B"  
01001 Kiev, Ukraine  
Phone +380.44.411.4918

## Americas

### USA

Sartorius Corporation  
5 Orville Drive, Suite 200  
Bohemia, NY 11716  
Phone +1.631.254.4249  
Toll-free +1.800.635.2906

### Argentina

Sartorius Argentina S.A.  
Int. A. Ávalos 4251  
B1605ECS Munro  
Buenos Aires  
Phone +54.11.4721.0505

### Brazil

Sartorius do Brasil Ltda  
Avenida Senador Vergueiro 2962  
São Bernardo do Campo  
CEP 09600-000 - SP- Brasil  
Phone +55.11.4362.8900

### Canada

Sartorius Canada Inc  
1173 North Service Road West, D4  
Oakville, ON L6M 2V9  
Phone +1.905.569.7977  
Toll-Free +1.800.668.4234

### Mexico

Sartorius de México, S.A. de C.V.  
Libramiento Norte de Tepotzotlan s/n,  
Colonia Barrio Tlacateco,  
Municipio de Tepotzotlan,  
Estado de México,  
C.P. 54605  
Phone +52.55.5562.1102  
[leadsmex@sartorius.com](mailto:leadsmex@sartorius.com)

### Peru

Sartorius Peru S.A.C.  
Avenue Alberto del Campo 411  
Floor 12 - The Office  
15076 - San Isidro, Lima  
Phone +51.1.441 0158

## Asia | Pacific

### Australia

Sartorius Australia Pty. Ltd.  
Unit 5, 7-11 Rodeo Drive  
Dandenong South Vic 3175  
Phone +61.3.8762.1800

### China

Sartorius (Shanghai) Trading Co., Ltd.  
3rd Floor, North Wing, Tower 1  
No. 4560 Jinke Road  
Zhangjiang Hi-Tech Park  
Pudong District  
Shanghai 201210, P.R. China  
Phone +86.21.6878.2300

### Hong Kong

Sartorius Hong Kong Ltd.  
Unit 1012, Lu Plaza  
2 Wing Yip Street  
Kwun Tong  
Kowloon, Hong Kong  
Phone +852.2774.2678

### India

Sartorius Weighing India Pvt. Ltd.  
#69/2-69/3, NH 48, Jakkasandra,  
Nelamangala Tq  
562 123 Bangalore, India  
Phone +91.80.4350.5250

### Japan

Sartorius Japan K.K.  
4th Fl., Daiwa Shinagawa North Bldg.  
8-11, Kita-Shinagawa 1-chome  
Shinagawa-ku, Tokyo, 140-0001 Japan  
Phone +81.3.3740.5408

### Malaysia

Sartorius Malaysia Sdn. Bhd  
Lot L3-E-3B, Enterprise 4  
Technology Park Malaysia  
Bukit Jalil  
57000 Kuala Lumpur, Malaysia  
Phone +60.3.8996.0622

### Singapore

Sartorius Singapore Pte. Ltd  
10 Science Park Rd  
The Alpha #02-13/14  
Singapore Science Park II  
Singapore 117684  
Phone +65.6872.3966

### South Korea

Sartorius Korea Ltd.  
8th Floor, Solid Space B/D,  
PanGyoYeok-Ro 220, Bundang-Gu  
SeongNam-Si, GyeongGi-Do, 463-400  
Phone +82.31.622.5700

### Thailand

Sartorius (Thailand) Co. Ltd.  
129 Rama 9 Road,  
Huaykwang  
Bangkok 10310  
Phone +66.2643.8361-6



◀ [www.sartorius.com](http://www.sartorius.com)

## For More Efficient Workflows and Less Interruptions in Your Routine Microbial Quality Control Testing

Need more Space in your working area for samples, funnels & media?

Small Footprint by compact dispenser design: save up to 35% space in your working area compared to comparable dispensers.

In addition the new Big Pack 250 filter boxes will help you to reduce storing space.

Save 50 to 67% space compared to boxes with 100 or 150 filters and use the space for samples, funnels or media instead!

Want less manual dispenser Reloading?

Want Full Shift working without interruptions?


Need Traceability?

Interested in working Guideline Conform?






## For More Efficient Workflows and Less Interruptions in Your Routine Microbial Quality Control Testing




Need more Space in your working area for samples, funnels & media? 

Want less manual dispenser Reloading? 

Want Full Shift working without interruptions? 

Need Convenience in your working area?

The cableless & portable dispenser featuring:

-  Battery operation
-  Low weight and
-  Handle


makes you independent of power outlets and allows effortless lift up, removal or transport.


Need Traceability? 


Interested in working Guideline Conform? 




## For More Efficient Workflows and Less Interruptions in Your Routine Microbial Quality Control Testing

Need more Space in your working area for samples, funnels & media? 




Need Convenience in your working area? 

Want less manual dispenser Reloading? 

Want Full Shift working without interruptions? 

Want Flexibility of operation? 

The Microsart<sup>®</sup> e.motion dispenser offers 3 operating modes for flexible and error-free operation:


-  No-touch operation by sensor mode
-  Push-button or
-  Footswitch operation


Need Traceability? 


Interested in working Guideline Conform? 



## For More Efficient Workflows and Less Interruptions in Your Routine Microbial Quality Control Testing

Need more Space in your working area for samples, funnels & media? 

Want less manual dispenser Reloading? 

Need Convenience in your working area? 

Want Flexibility of operation? 

Want Full Shift working without 

### Need Traceability?

Number of remaining filters & all relevant information visible!


Printed on the paper of each membrane cell & visible through the dispensers' filter output are:

- Membrane filter type
- Diameter
- Lot number and
- Number of remaining filters

In addition, data, such as sterilization and expiration date, pore size and bar code on each box label make the usability of these membrane filters as convenient as possible.

## For More Efficient Workflows and Less Interruptions in Your Routine Microbial Quality Control Testing

Need more Space in your area for samples, funnels

Need Convenience in your working area? 

Want Flexibility of operation?

Want less manual dispenser Reloading?

The unique 250 filter package reduces the need to change the package every 100 or 150 filters, helping to increase your efficiency.





Packaging Type	Filters per box	Number of reloads per 1000 tests	Space required for 1000 tests
Standard Packaging	100 to 150	7 to 10	408 to 612 cm <sup>2</sup> (= 6 to 9 boxes)
Microsart <sup>®</sup> e.motion Big Pack	250	4	204 cm <sup>2</sup> (= 3 boxes)


without


g Guideline


## For More Efficient Workflows and Less Interruptions in Your Routine Microbial Quality Control Testing

Need more Space in your working area for samples, funnels & media? 

Need Convenience in your working area? 

Want Flexibility of operation? 

Want less manual dispenser Reloading? 

Want Full Shift working without interruptions? 


No workflow interruptions necessary with the big packs of 250 filters!


With a consumption of 250 filters per day, full shift working will be possible without reloading the dispenser.

Each membrane filter box allows 1.7 to 2.5 times more filtrations than the current products on the market before reloading.




# For More Efficient Workflows and Less Interruptions in Your Routine Microbial Quality Control Testing

Need more Space in your working area for samples, funnels & media? 

Want less manual dispenser Reloading? 

Want Full Shift working without interruptions? 

Need Convenience in your working area? 

Want Flexibility of operation? 

Interested in working Guideline Conform? 

Work guideline conform with our 0.45 µm membrane filters:

Certified quality according to ISO 7704. Find out more in the attached Guideline Landscape brochure.



See attached file  
Broch\_Quality\_Control\_Guidelines\_SM-1505-e

