

Microbiological Control  
Pharma QC Applications  
(Sales Training eBook)

Simplifying Progress

**SARTORIUS**

## Intelligent Tools for Microbiological Quality Control

Working in pharmaceutical manufacturing quality control is a big responsibility, customers are concerned for the patient's wellbeing. We understand that ensuring safety and compliance for microbiological processes is critical. Our microbiological QC portfolio is designed to give the best tools, ensuring results meet the required levels of safety and compliance. We offer intelligent solutions for unmet needs.

Reliability

### Microbial Enumeration

**USP <61>, <62> and <1231>**

Count the number of microbes in your sample using touch-free membrane transfer to agar.



### Sterility Testing

**USP <71> | EP 2.6.1 | Ph. Eur. 2.6.1**

Determine whether your sample is sterile. Use aseptic sampling for future quality assurance.



Compliance

### Rapid Microbial Detection

**USP <63>, | Ph. Eur. 2.6.7 | USP <1071>**

Detect mycoplasmas, as well as total bacteria and fungi with 3 hours (instead of 14 or 28 days) by qPCR.



### Microbial Air Monitoring

**USP <1116> EU GMP Annex 1**

Employ non-stop clean room monitoring for up to 8 hours.



Safety

Efficiency





Microbiological QC  
Solutions (all solutions  
for Pharma QC)

English

German

French



Short Introduction to  
Quality Control in the  
Biopharmaceutical  
Industry  
Video (slide show)

English



Guide to  
Pharmaceutical QC  
for Microbiology

English

EU GMP Guide, Annex 1   Room classification	
Classification	airflow at work surface (m/s)
A	0.3
B	0.1
C	0.1
D	0.1

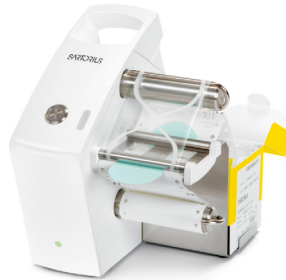
Short Explanation on  
Controlled  
Environments and  
Clean Room Classes  
Video (slide show)

English





Microsart® @filter and @media Plates



Microsart® e.motion Dispenser



Microsart® filtration Manifolds



# Microsart® @filter & @media

Our Microsart® @filter combined with the Microsart® @media is a unique system which effortlessly positions the membrane on the agar plate. With only a turn, it's locked and ready to incubate, reducing the risk of contamination.

## Key Benefits & Features

- + Convenience
- + Safety





Data Sheet Microsart®  
@filter

English

German

French

Spanish

Italian



Validation Guide  
Microsart® @filter  
Microsart® @media

English



Data Sheet Microsart®  
@media

English

German

French

Spanish

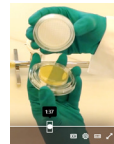
Italian

Russian



Microsart® @media  
Product Overview  
Training Video

English



Microsart® @filter and  
@media Full Workflow  
Training Video

English


Back to @filter @media

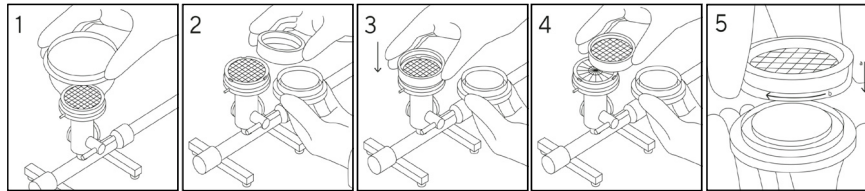


# Microsa

Our Microsart® @media unique system which plate. With only a tu risk of contamination

## Key Benefits & Feat

-  Convenience
-  Safety



1 Remove the funnel

2 Pick-up the plate lid

3 Apply the lid on filter

4 Membrane is transferred

5 Transfer filter on @media

## Touch-Free Membrane Transfer

Microsart® @media's innovative, adhesive lid transfers the filter and positions it precisely onto the agar plate to allow touch-free transfer of the membranes.



## Easy Colony Enumeration

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



# Microsart® @filter & @media

Our Microsart® @filter combines a unique system which effortlessly locks the plate. With only a turn, it's locked and the risk of contamination.

## Key Benefits & Features

- + Convenience
- + Safety



0,2 µm Minisart®  
for sterile venting

### Minimized Risk of Cross-Contamination

The integrated venting function of the new Microsart® Manifold minimizes the risk of cross-contamination during filtration.



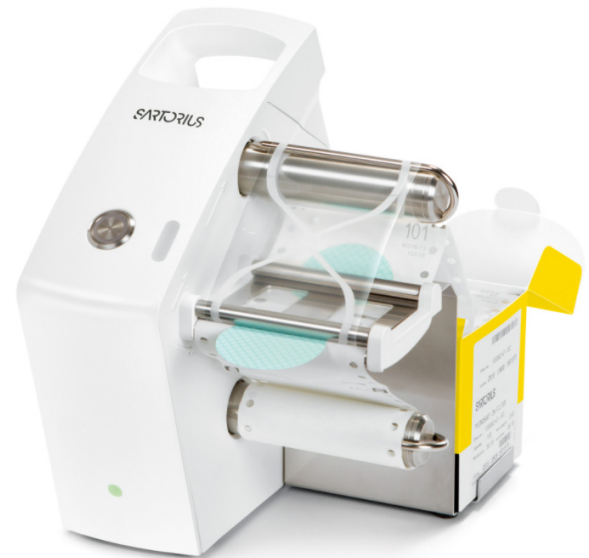


# Microsart® e.motion Dispenser

Thanks to its small footprint the Microsart® e.motion dispenser saves precious space under the laminar air flow which can be used for samples, media and funnels instead. The membrane filters are released fully automatically, at the touch of a button or hands-free or by foot switch!

## Key Benefits & Features

- + Convenience
- + Performance





Back to Microsart®  
e.motion

English

German

French

Italian



Microsart® e.motion  
Dispenser Manual

English

German

French

Spanish

Italian

Russian

Japanese

Chinese

Turkish

Polish



Data Sheet Microsart®  
e.motion Membrane  
Filters

English

German

French

Spanish

Italian



Back to Microsart® e.motion

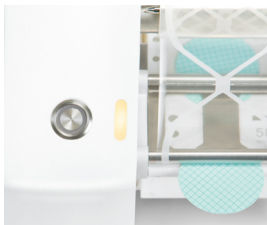


# Microsart®

Thanks to its small footprint saves precious space under the hood for samples, media and and for released fully automatically or by foot switch!

## Key Benefits & Features

-  Convenience
-  Performance



## Battery Driven

Microsart® e.motion dispenser can be purchased with battery or main power connection. Having no cable inside the laminar flow facilitates handling, cleaning and disinfection.

Color code indicates battery state.



## Automatic Dispensing

A sensor detects movements in front of the dispenser when approaching the forceps. The membrane is then released automatically facilitating handling.

## Remaining Membranes Info

Users can see any time how many membranes are left in the dispenser, helping streamlining their workflows with no stress.



# Microsart® e.motion Dispenser

Thanks to its small footprint it saves precious space under the hood for samples, media and funnels. Membrane filters are released fully automatically, or by foot switch!

## Key Benefits & Features

- + Convenience
- + Performance

# 38%

## Time Saving

Membrane filters are available in bands of 250 units, it will help users to save up to 38% of their time compared to single individually packed membrane filters.



# Microsart<sup>®</sup> Manifolds

The user can choose between different sizes of manifolds, from single branch to multi-branch. The different manifold sizes are connectable with each other by quick connection to enable the manifold to adapt to daily workload. Depending on the application and criticality of product to be tested, users can choose between different funnel and filtration units, from complete reusable to single sterile solutions.

## Key Benefits & Features

- + Convenience
- + Performance
- + Ergonomics
- + Adaptability





Microcart® Manifolds Datasheet

- English
- German
- French
- Italian



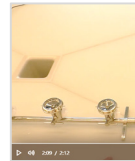
Microcart® Manifolds Video

- English



Microcart® Manifolds Manual

- English
- German
- French
- Spanish
- Italian
- Chinese



Microcart® Manifolds Video

- English





[Back to Microcart® Manifolds](#)






## Microsart® Manifolds

The user can choose between single-branch to multi-branch. They can be used together with each other by quick connectors to daily workload. Depending on the product to be tested, users can choose different filtration units, from complex

### Key Benefits & Features

-  Convenience
-  Performance
-  Ergonomics
-  Adaptability

**Assembly and disassembly without tools** facilitates handling and cleaning and sterilization of the manifolds.

**Quick connectors** on both sides of the manifolds improve usability and allow positioning the pump on left or right side.

**Silicone feet** prevent the manifold to slide on the working bench and facilitate overall handling.



# Microsart<sup>®</sup> Manifolds

The user can choose between different sizes of manifolds, from single branch to multi-branch. They can be used together with each other by quick connection to daily workload. Depending on the product to be tested, users can use different filtration units, from complete

## Key Benefits & Features

- + Convenience
- + Performance
- + Ergonomics
- + Adaptability

## Robustness

The Manifold is made of 100% stainless steel. The high quality of the stainless-steel material (Stainless Steel 316L - 1.4404) and the consequence in using one material type ensures a long working life of the manifolds.





# Microsart® Manifolds

The user can choose between different sizes of manifolds, from single branch to multi-branch. The difference lies in the number of ports with each other by quick connection to daily workload. Depending on the product to be tested, users can choose between different filtration units, from complete retention to low retention.

## Key Benefits & Features

- + Convenience
- + Performance
- + Ergonomics
- + Adaptability



**Low height** ensures efficient and ergonomic operation under the laminar flow, especially when heavy bottles must be repeatedly lifted over the funnel by the technician.

**Easy to operate valves** minimize the risks of strain injuries when numerous samples must be analyzed in a short period of time.



Marketing Resources

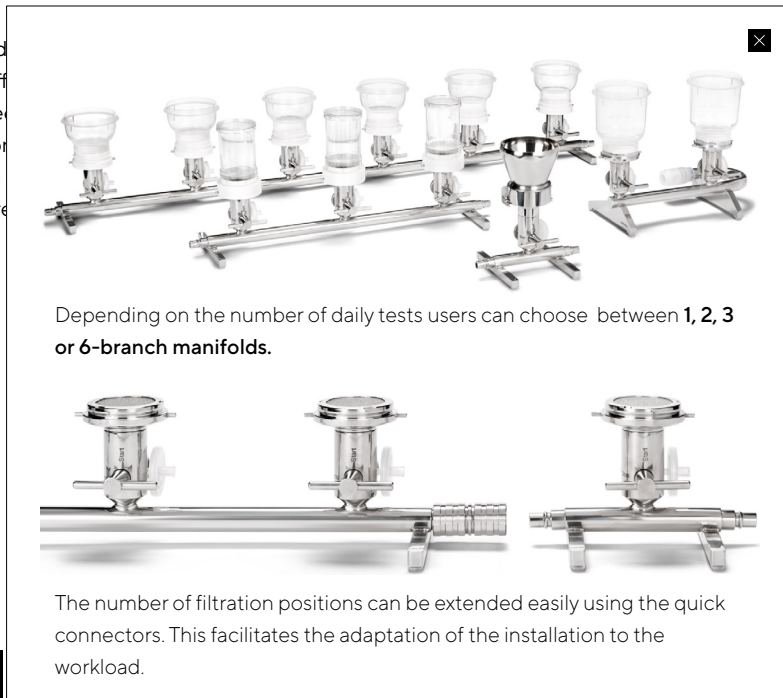


## Microsart® Manifolds

The user can choose between different configurations from single-branch to multi-branch. The different configurations are designed to work with each other by quick connectors to daily workload. Depending on the application and product to be tested, users can choose different filtration units, from complete replacement

### Key Benefits & Features

- + Convenience
- + Performance
- + Ergonomics
- + Adaptability



Marketing Resources





Sterisart® Pump



Sterisart® Disposable Canisters



# Sterisart® NF

State-of-the-art closed filtration system for lot-release sterility testing of sterile pharmaceuticals. The Sterisart® NF system is easy to use and ensures the maximum reliability of sterility test results. The portfolio includes several devices adapted to specific sterility testing needs and is fully compliant with all pharmacopeial and regulatory requirements.

## Key Benefits & Features

- + Convenience
- + Safety
- + Performance





Sterisart® NF Flyer

English



Video of Sterility Testing Workflow

English



Datasheet Sterisart® NF

English



Points to Consider When Validating Sterility Testing Canisters

English  
Portuguese



Sterisart® NF Brochure

English  
German



Why It Is Important to Have a Secondary Supplier

English  
Portuguese



Application Note Sterisart® Septum

English



Sterisart® Validation Guide

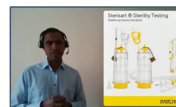
English



Sterisart® Sterility Testing eBook

English

Back to Sterisart® NF






Video Sterisart® Why Sartorius?

English

# Sterisart® NE

State-of-the-art closed filtration of sterile pharmaceuticals. The system is designed to ensure the maximum reliability and ensures the maximum reliability. The product portfolio includes several different models to meet various needs and is fully compliant with regulatory requirements.

## Key Benefits & Features

-  Convenience
-  Safety
-  Performance



The two different **color-coded clamps** on Sterisart® tubing eliminate the risk of confusion, making it easy for users to identify and fill the canisters with the correct media.

The extra-large, **grip-optimized filter caps** are directly attached to the vent filters. This is practical – and not only that. The caps connected by a band to the vent filters cannot fall off or touch any surfaces of the isolator, so they are safe from contamination.

Instead of being sealed in a separate, difficult-to-open package, the **easy-to-handle yellow plugs** for closing the canisters are right where users need them: in extra pockets directly inside the primary packaging. This facilitates the workflow and lets technicians stay focused on their core tasks.



Marketing Resources



# Sterisart® NE

State-of-the-art closed filtration of sterile pharmaceuticals. The device ensures the maximum recovery and ensures the maximum recovery. The portfolio includes several different models to meet your needs and is fully compliant with regulatory requirements.

## Key Benefits & Features

- + Convenience
- + Safety
- + Performance



A critical step in sterility testing is sampling from or supplementing the media in a canister. You usually do this by piercing a tube. This conventional method is complicated, involves the hazard of injury and increases the risk of false positive results.

The **septum technology** has been purpose-built to aseptically draw samples. The sample will remain safely protected, ensuring a sterile barrier against contamination. At the same time, the simple piercing mechanism minimizes risk of puncture injuries.

Our specially designed and patented **dual-needle**, with its **protective shield** and built-in venting filter, ensures safe and easy handling, even when working with cumbersome isolator gloves.



Marketing Resources



# Sterisart® N

State-of-the-art closed filtration of sterile pharmaceuticals. This unit and ensures the maximum recovery and our portfolio includes several designs to meet your needs and is fully compliant with regulatory requirements.

## Key Benefits & Features

- + Convenience
- + Safety
- + Performance



Our gamma-sterilized and **gas-impermeable** primary packaging enables users to omit the rinsing procedure and proceed right away with testing. This will save not only many steps, but also the additional costs incurred.

Oily substances and emulsions are difficult and take considerable time to filter. Sterisart® CA units with a cellulose acetate membrane deliver the **fastest flow rates** in customer tests.

The inlet port of the Sterisart® canisters has been designed with an **anti-foaming** dispensing channel. This facilitates rinsing and minimizes risks of venting filter clogging.

The venting filter is positioned with a **safety clearance** of 1 inch, or over 2½ cm, to effectively protect from contact with culture medium, even if foaming or the slightest vibration occurs.



Marketing Resources





# Sterisart® Universal Pump

The Sterisart® Universal Pump is available in two versions: as basic version 16419 and as an upgraded version 16420 with display and user software. The pump can be used in clean rooms, integrated into clean benches, or installed countersunk in the working surface of isolators. Its low, compact design has a space-saving footprint – a great benefit for most clean room benchtops and isolators.

## Key Benefits & Features

- + Convenience
- + Safety
- + Robustness





Sterisart® Universal Pump Brochure

English  
German



Sterisart® Universal Pump 16420 Manual

EN | DE  
French  
Russian



Sterisart® Easy Configuration Datasheet

English



Sterisart® Universal Pumps Equipment Qualification (IQ/OQ) Datasheet

English  
German



Sterisart® Universal Pump 16419 Manual

EN | DE



Sterisart® Universal Pumps Preventive Maintenance and Service Contracts

English  
German



Sterisart® Sterility Testing eBook

English



Video Sterisart® Why Sartorius?

English

Back to Sterisart® Pump



# Sterisart® Universal Pump

The Sterisart® Universal Pump is available in version 16419 and as an upgrade with the Sterisart® software. The pump can be used on clean benches, or installed count down benches. Its low, compact design has been designed for most clean room benches.

## Key Benefits & Features

- + Convenience
- + Safety
- + Robustness

Low profile



## Compact and User-friendly

As space is often limited in isolators and clean benches, the Sterisart® Universal pump has been designed with a compact footprint and a low profile. Thanks to its user-friendly construction, **it saves spaces**, ensuring excellent **ergonomic handling**, particularly in the areas in front of and above the pump. Despite the pump's compactness, it is easy to thread the tubing of the Sterisart® sterility test unit into the pump rotor. The special pump head design ensures excellent life of the tubing and, moreover, is easy to access for cleaning and disinfecting.

Marketing Resources



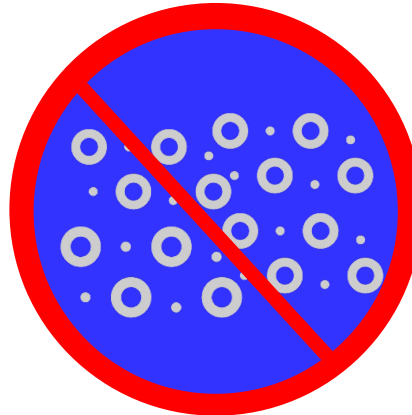
# Sterisart® Universal Pump

The Sterisart® Universal Pump is the latest version 16419 and as an upgrade to previous software. The pump can be used on benches, or installed count down units. Its low, compact design has made it suitable for most clean room benches.

## Key Benefits & Features

- + Convenience
- + Safety
- + Robustness

Zero particle emission



## Increased Safety and Reliability

A unique motor principle has resulted in a closed pump chassis design, eliminating the need for ventilation of the motor. Therefore, the pump will not release any particles or microbes into cleanroom or isolator environments or cause turbulence in laminar-flow areas. This results in increased safety and reliability.



Marketing Resources



# Sterisart® Universal Pump

The Sterisart® Universal Pump is an upgrade to version 16419 and as an upgrade to the software. The pump can be used on benchtop benches, or installed count down benches. Its low, compact design has made it a popular choice for most clean room benches.

## Key Benefits & Features

- + Convenience
- + Safety
- + Robustness



## Rugged and Maintenance-Free

As the pump chassis is made of stainless steel, it features excellent chemical compatibility with sterilants used in isolators. This, together with the unique type of motor, makes the pump highly rugged and virtually maintenance-free, considerably reducing service costs.



Marketing Resources





MD8 Airscan®



Gelatin Filters



# MD8 Airscan®

Monitors ambient air in cleanrooms and controlled areas in the (c) GMP environment for viable microorganisms. The MD8 Airscan® provides non-stop active air monitoring for at least 8 hours, using only one single gelatin membrane filter. Convenient, continuous monitoring no one else can offer.

## Key Benefits & Features

- + Convenience
- + Performance



Marketing Resources





MD8 Airscan®  
Datasheet

English  
German  
Russian



MD8 Airscan®  
Equipment Qualification  
(IQ/OQ) Datasheet

English  
German



MD8 Airscan®  
Sampling Head Manual

Multiple



MD8 Airscan®  
Equipment Qualification  
(IQ/OQ) Datasheet

English  
German



MD8 Airscan®  
Sampling Head Manual

English  
German  
French  
Spanish  
Italian  
Portuguese



MD8 Airscan®  
Product Overview  
Video

English



Dr. Eric Clement/Grahl  
Sartorius Manager

GMP Annex1 Air-  
Monitoring Webinar  
2021

English

Back to MD8 Airscan®







## MD8 Airscan

Monitors ambient air in cleanroom or GMP environment for viable particles. Provides non-stop active air monitoring with one single gelatin membrane filter. Monitoring no one else can.

### Key Benefits & Features

-  Convenience
-  Performance



The MD8 Airscan® Command Unit weighs **only 1.2 kg** (2.6 lbs). It includes 2 LEDs for displaying operating status and errors.

The specially designed bayonet connector allows the unit to function as a “stand-alone benchtop version”.

Quick connectors allow the command unit to be disconnected from one sampling head and connected to another one.

The MD8 Airscan® Command Unit can be connected with a thermal transfer printer.





Marketing Resources

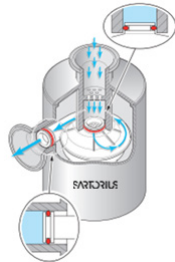


## MD8 Airscan

Monitors ambient air in cleanroom or GMP environment for viable particles. Provides non-stop active air monitoring with one single gelatin membrane filter. Monitoring no one else can.

### Key Benefits & Features

-  Convenience
-  Performance



Stores up to 300 test results from different sampling heads and displays operating status and errors.

Stainless-steel tri-clamp connectors facilitate installation of the MD8 Airscan® as an integrated unit in lab's air-line.

Flow path can be disinfected or sterilized inline with vaporized hydrogen peroxide (VHP).



Marketing Resources



# Gelatin Filters

Gelatin filter disposables are individually packaged, presterilized and ready-to-connect units, each consisting of a gelatin membrane filter and a holder. Gelatin filters in conjunction with the MD8 Airscan® Air Sampler are used for collection of airborne microbes and viruses.

## Key Benefits & Features

- + Convenience
- + Performance
- + Compliance





Gelatin Membrane Filters in Biosafe® Bags

English  
German



Continuous Air Monitoring Application Note

English  
German  
Japanese



Gelatin Filter Disposables Datasheet

English



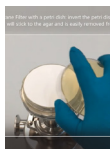
Air Monitoring in Cleanroom Environments Cording to EN 17141 and ISO 14698

English



Gelatin Filter Disposables Manual

Multiple



Gelatin Membrane Filters Product Features and Transfer to Agar Plate Video

English



Biosafe® Bags and Port Product Demonstration Video

English

Back to Gelatin Filters



Dr. Eric Christmann  
Senior Manager  
Global Air Monitoring




GMP Annex1 Air-Monitoring Webinar 2021

English

## Gelatin Filter

Gelatin filter disposables are ready-to-connect units, each consisting of a filter and a holder. Gelatin filters and Samplers are used for collection

### Key Benefits & Features

-  Convenience
-  Performance
-  Compliance



### Biosafe® Transfer Bags

Gelatin membrane filters (GMF) in Biosafe® Bags are aseptically transferred without any risk of contamination through the Biosafe® Port into the desired critical area, such as an isolator. The solution facilitates collapsing and streamlining of customers processes. (see video in marketing resources page).






Marketing Resources



# Gelatin Filters

Gelatin filter disposables are individual ready-to-connect units, each consisting of a filter and a holder. Gelatin filters in conjunction with the MD8 Airscan Sampler are used for collection of

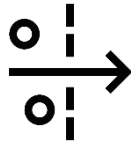
## Key Benefits & Features

-  Convenience
-  Performance
-  Compliance



### Long Term Monitoring

The MD8 Airscan® provides non-stop active air monitoring for at least 8 hours, using only one single gelatin membrane filter. Published experimental results have been proving the extraordinary performance of Gelatin Membrane Filters (see app. note).



### Absolute Retention

of organisms since the filters capture all bacteria laden particles in the sample regardless of particle size.



### Residual Moisture

in the filters prevents desiccation of the retained organisms, improving recovery rates.






Marketing Resources



## Gelatin Filter

Gelatin filter disposables are available in ready-to-connect units, each with a cap and a holder. Gelatin filters and Samplers are used for collecting

### Key Benefits & Features

-  Convenience
-  Performance
-  Compliance



### Regulatory Compliance

Biocontamination control strategies are a key requirement of recent standards and guidelines, such as the EN 17141 and the revision of the EU GMP Annex 1. The EN 17141, published in 2020, has been adopted in the European Union and United Kingdom, replacing the ISO 14698. The Annex E chapter of EN 17141 details guidance on culture-based microbiological monitoring methods and sampler verification. We have demonstrated that microbial air monitoring by Gelatin Membrane Filtration meets the requirements of the EN 17141 and the ISO 14698, which remains the standard outside the EU and the United Kingdom. (see application note).





Microsart® Mycoplasma



Microsart® ATMP Sterile Release





# Microsart<sup>®</sup> Mycoplasma

Viable but non-culturable (VBNC) Mycoplasma can go undetected when using the traditional method of mycoplasma culturing, risking false-negative results. Our Microsart<sup>®</sup> qPCR kits offer a method for early detection of mycoplasma contamination. These kits are fast, highly specific, sensitive, and compliant with international guidelines.

## Key Benefits & Features

- + Convenience
- + Performance
- + Compliance





Microsart® AMP  
Mycoplasma Datasheet

English



Correlation Between  
Colony Forming Units  
and Genome Copies of  
9 Mycoplasma Species -  
App. Note

English



Mycoplasma Assays for  
qPCR - The New Gold  
Standard - Flyer

English



Minimize the Risk of  
Mycoplasma  
Contamination in Your  
Cell Cultures - App. Note

English



Battle Card Microsart®  
Mycoplasma PCR Kits

English



Benchmarking of Two  
Mycoplasma Real-time  
PCR Kits According  
EP 2.6.7

English



Digital Tool - ATMP  
Sterile Release and  
Mycoplasma Testing

English




Back to Microsart® Mycoplasma



# Microsart®

Viable but non-culturable when using the traditional false-negative results. Our early detection of mycoplasma highly specific, sensitive, and

## Key Benefits & Features

-  Convenience
-  Performance
-  Compliance



## Reduced Pipetting Steps

with Our Master Mix The ready-to-use master mix has the DNA for internal control already included. It mitigates the risks of false-negative results and minimizes the number of reactions to run.

## Testing Right Away

Microsart® Mycoplasma qPCR kits can be used with any qPCR cyclers able to detect the fluorescent dyes FAM® and ROX®.




Our clear, **color-coded reagents** guide the user through the short and simple protocol.

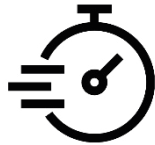


# Microsart® Mycoplasma

Viable but non-culturable (VBN) when using the traditional methods, leading to false-negative results. Our Microsart® enables early detection of mycoplasma with a highly specific, sensitive, and cost-effective assay.

## Key Benefits & Features

-  Convenience
-  Performance
-  Compliance



### Reliable Results Within 3 Hours

Users will avoid lengthy 28-day cultivation protocols when choosing Microsart® Mycoplasma qPCR Kits.



TaqMan® probes ensure the highest qPCR **specificity** without the need for a melting-curve analysis.



Lyophilized reagents ensure consistent quality **without the need for freezer** storage.



Validation standards for required mycoplasma species in Microsart® Mycoplasma qPCR kits are lyophilized to render them **non-infectious**.



Marketing Resources

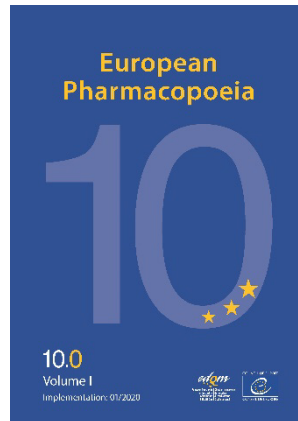


# Microsart® Mycoplasma

Viable but non-culturable (VBN) when using the traditional methods, leading to false-negative results. Our Microsart® Mycoplasma Kits enable early detection of mycoplasma contamination with highly specific, sensitive, and cost-effective results.

## Key Benefits & Features

- + Convenience
- + Performance
- + Compliance



## Validated Test Results in Compliance With International Standards

Microsart® AMP and ATMP Mycoplasma Kits are validated according to EP 2.6.7 for sensitivity, specificity, and robustness. They also fulfill the criteria of international authorities, such as USP and JP.



Marketing Resources



# Microsart® ATMP Sterile Release

Detection of bacterial and fungal contamination within 3 hours using the proven real-time PCR with our Microsart® ATMP Sterile Release Kit, a combination of Microsart® ATMP Bacteria and of Microsart® ATMP Fungi.

## Key Benefits & Features

- + Convenience
- + Performance
- + Compliance





Microsart® ATMP  
Bacteria | Fungi | Sterile  
Release  
Microsart® Research  
Bacteria | Fungi

English



Detection of Microbial  
Contaminations in High  
Cell Density Jurkat-,  
HPBMC and CHO-  
Cultures - App. Note

English



Short Introduction of  
Rapid Sterility and ATMP  
Video (Slide Show)

English



Equivalency of PCR-Based  
Rapid Sterility Testing and  
the Compendial Culture  
Method - App Note

English



Microsart® ATMP  
Bacteria Tutorial  
Video

English



Digital Tool - ATMP  
Sterile Release and  
Mycoplasma Testing

English




[Back to Microsart® ATMP Sterile Release](#)



# Microsart® ATMP qPCR Kit

Detection of bacterial and the proven real-time PCR Kit, a combination of Micro ATMP Fungi.

## Key Benefits & Features

-  Convenience
-  Performance
-  Compliance



### Reduced Pipetting Steps

with Our Master Mix

The ready-to-use master mix has the DNA for internal control already included. It mitigates the risks of false-negative results and minimizes the number of reactions to run.

### Testing Right Away

Microsart® Mycoplasma qPCR kits can be used with any qPCR cycle able to detect the fluorescent dyes FAM® and ROX®.

Our clear, **color-coded reagents** guide the user through the short and simple protocol.



Marketing Resources






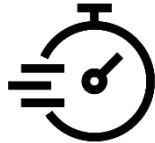


# Microsart® A

Detection of bacterial and fungal contamination using the proven real-time PCR with Microsart® ATMP Fungi Kit, a combination of Microsart® ATMP Bacteria and ATMP Fungi.

## Key Benefits & Features

-  Convenience
-  Performance
-  Compliance



### Reliable Results Within 3 Hours

Users will avoid lengthy 28-day cultivation protocols when choosing Microsart® Mycoplasma qPCR Kits.



TaqMan® probes ensure the highest qPCR **specificity** without the need for a melting-curve analysis.



Lyophilized reagents ensure consistent quality **without the need for freezer** storage.






Validation standards for required bacterial and fungal species in Microsart® ATMP Sterile Release are lyophilized to render them **non-infectious**.

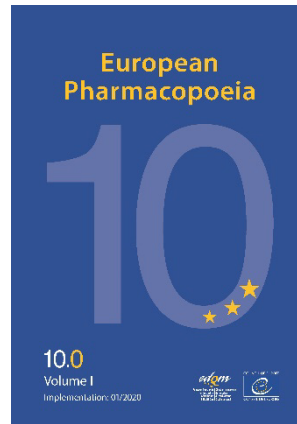


# Microsart® ATMP

Detection of bacterial and fungal contamination using the proven real-time PCR with Microsart ATMP Fungi Kit, a combination of Microsart ATMP Bacteria and ATMP Fungi.

## Key Benefits & Features

-  Convenience
-  Performance
-  Compliance



## Validated Test Results in Compliance With International Standards

Microsart® ATMP Sterile Release kits were developed in close cooperation with authorities and are validated according to EP 5.1.6 and USP <1223> for sensitivity, specificity, and robustness.

This test is not a replacement for the classic sterility test according to USP <71> or EP 2.6.1.



Marketing Resources



**Sartorius Lab Instruments GmbH & Co. KG**

Otto-Brenner-Strasse 20

37079 Goettingen

Phone +49 551 308 0


**Sartorius Corporation**

565 Johnson Avenue

Bohemia, NY 11716

Phone +1 631 254 4249

Toll-free +1 800 635 2906

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