



# Monitoring the Quality of Beverages, Water and Foods

Biosart® 100 Monitors  
and Biosart® 100 Media

Simplifying Progress

**SARTORIUS**

# Fast and Reliable: Biosart® 100 Monitors & Media

## Microbiological Quality Control With Biosart® 100 Monitors Increases Efficiency and Saves Time

The detection of microbial contamination in sample liquids such as final product, incoming inspection or during in-process testing plays a significant role in the quality assurance process. The requirements for a practical microbiological test method are that it permits quantitative and reproducible detection of trace contamination and that it can be performed efficiently and economically under routine conditions. These requirements are fulfilled optimally by the membrane filtration method. The use of ready-to-use disposable units simplifies the testing procedure and prevent crosscontamination of samples.

Biosart® 100 Monitors have been specifically designed for the detection and enumeration of microorganisms in beverages, water, pharmaceuticals, cosmetics, foods and other liquids. These sterile disposables with an incorporated membrane filter and cellulose pad are ready-to-use. After filtration, just remove the 100 mL funnel to convert the Monitor into a petri dish eliminating the need for membrane manipulation. Culture media for wetting the pad are available in individually sterilized, convenient plastic ampoules. Biosart® 100 Monitors are ready-to-use filter units designed to be placed onto the bases of a vacuum manifold, eliminating the cleaning and sterilization required of reusable funnels.

## High Flow Membranes

Biosart® 100 Monitors are also available with the new 0.45 µm High Flow membranes. The special pore structure allows shorter filtration times due to 30% higher flow rates.

## Compliance with International Standards

The membrane filtration method is worldwide accepted and the preferred method of choice for the analysis of microbial contamination in liquid samples. Biosart® 100 Monitors and Media are in compliance with the membrane filtration procedures referenced in the:

- European drinking water directive (Council Directive 98/83/EC on the quality of water)
- Standard Methods for the Examination of Water and Waste Water, 20th edition
- U.S. Environmental Protection Agency, 600/8-78-017.
- ISO Standard's microbiological methods, such as ISO 7704, ISO 9308-1, EN 12780, ISO 8199
- WHO Guidelines for Drinking Water Quality, 1997
- International Pharmacopoeia, such as the current editions of the USP and EP

The quality management system of Sartorius meets the requirements of the International Standard ISO 9001. For quality assurance all materials are selected carefully in accordance with current regulations and recommendations, such as the FDA CFR's and applicable current Good Manufacturing Practices.

## Biosart® 100 Monitors

### Specifications

|                         |  |
|-------------------------|--|
| Housing                 | Polystyrene  |
| Membrane filter         | Cellulose nitrate (cellulose ester): choice of white, green or grey, with grid; Regenerated cellulose: white; can be used as documentation |
| Plug and adapter        | Polyethylene   |
| Pad                     | Cellulose  |
| Capacity                | 100 mL, 10 mL graduations  |
| Pore size               | 0.2 µm, 0.45 µm or 0.8 µm  |
| Filter diameter         | 47 mm  |
| Filtration area         | 14.5 cm <sup>2</sup>   |
| Max. operating pressure | Vacuum only  |
| Outlet                  | 6.5 × 1.5 mm   |
| Lot certificates        | Recovery rate, sterility and specifications  |



Each lot is tested by Sartorius for accordance with established specifications before release, and each box includes a lot certificate.

## Easy Work Flow – Reliable Results

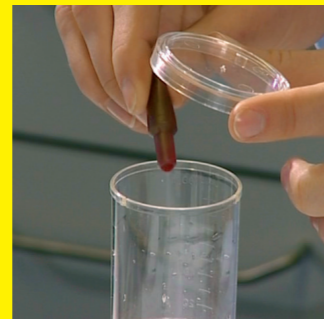
1 | Pour the Sample



2 | Apply Vacuum & Filter the Sample



3 | Add the Biosart® 100 Nutrient Media



4 | Close the Outlet



5 | Remove the Funnel



6 | Incubate the Petri Dish



# Microsart® Manifolds – The Filtration Stand That Adapts to Your Need

## Product Information

The ability to accurately detect and quantify microorganisms in liquid samples is of prime importance for any quality control laboratory. This is why membrane filtration is the established method of choice for liquid testing as it provides reliable and reproducible results. A suitable and thoughtful filtration equipment is at least as important as high quality membrane filters and reliable media plates. Facilitate your daily microbiological testing procedures with the new Microsart® Manifold: select the proper manifold for your individual microbiological testing procedure, choose between reusable and single use funnels or filtration units, and decide between different sizes of manifolds and enjoy the manifold that adapts to your needs.

## Applications

Microbiological quality control of liquid samples:

- Beverages (beer, wine, soft drinks, bottled water)
- Pharmaceutical analysis (WFI, purified water, Microbial Limit Testing of non-sterile products, bioburden testing)
- Environmental water testing (water monitoring)
- Cosmetics

### The Manifold That Adapts to Daily Needs

Depending on the number of daily tests you 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 your daily workload. Standardized quick connectors at both sides of the manifolds enable fast connections of either tubing, another manifold or end caps. Depending on your application and criticality of product to be tested, choose between different funnel and filtration units, from complete reusable to single sterile solutions. Simply connect the filter onto the manifold by using the correct adaptor. Simply push the filter adapter into the manifold to fix it and choose your favorite between 3 different working positions. This new and fast way to connect the respective filter adapter (base support) to the manifold eliminates the need of additional tools.

### Minimized Risk of Secondary Contamination by Design

The Manifold is made of 100% stainless steel. For sterilization of your working equipment, no disassembly is needed – simply put the complete manifold including the filter adapters into the autoclave for reliable sterilization. The easy to clean, compact design of the manifold allows quick and easy cleaning using standard disinfectants. The higher quality of the stainless steel material and the consequence in using one material type ensures a long working life of the manifold in your lab. During filtration, residual remaining liquid below the filter could potentially cross-contaminate the filter and therefore your results. The newly designed and integrated sterile venting step enables full drainage of liquid below the filter when using the Microsart® base support (Microsart® base). This integrated venting step eliminates the risk of secondary contamination and ensures a secure working procedure.

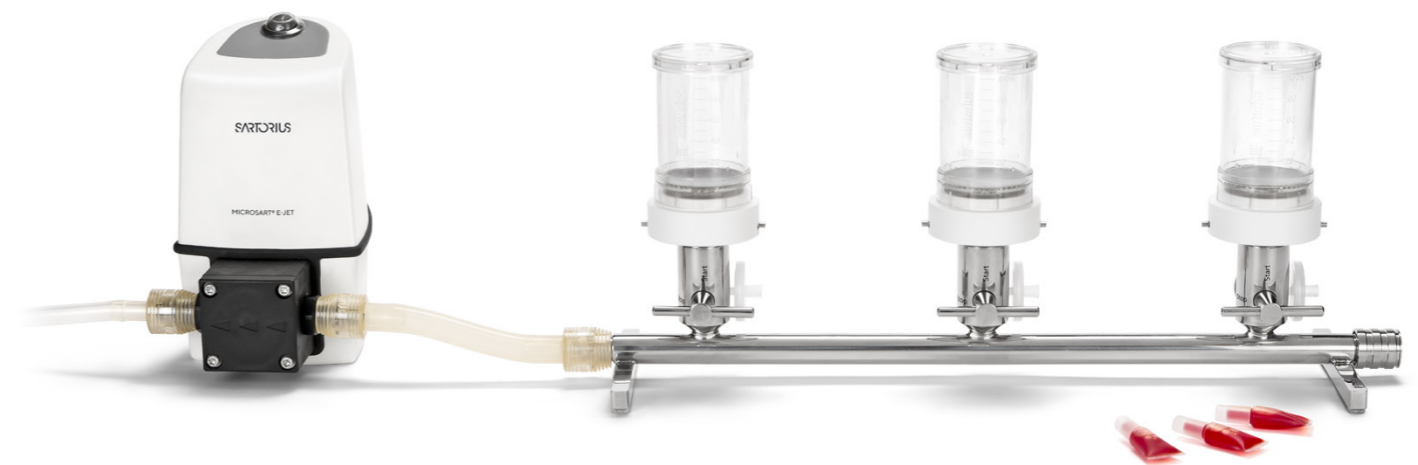
### Safe and Stable Ergonomic Working Procedures

The Manifolds rubber feet providing better stability at the workplace. The low height ensures efficient and ergonomic working under the laminar flow. Intuitive handling of the valves is given by clear start | stop feedback when turning the valves and indications on the manifold. Venting is automatically occurring when turning the valves to the stop position at the end of the filtration.

## Microsart® Manifolds

### Specifications

| Materials of construction                                 |                               |
|---|-------------------------------|
| Manifold and base supports                                | Stainless Steel 316L (1.4404) |
| Valves  | Monel                         |
| O-rings   | Silicon                       |
| Feet  | Silicon                       |
| Dimensions (L × H × W) in mm (without Funnels and Tubing) |                               |
| 1 branch  | 176 × 120 × 98                |
| 2 branch  | 246 × 130 × 118               |
| 3 branch  | 474 × 120 × 98                |
| 6 branch  | 924 × 120 × 98                |
| Weight in kg  |                               |
| 1 branch (for Microsart® products)                        | 0.275                         |
| 2 branch (for Microsart® products)                        | 0.675                         |
| 3 branch (for Microsart® products)                        | 0.725                         |
| 6 branch (for Microsart® products)                        | 1.400                         |
| Autoclaving conditions                                    |                               |
| 121 °C for 30 min   |                               |



# Ordering Information

## Biosart® 100 Monitors



### Superior Performance

- High flow rate
- High total throughput

### Safe and Reliable

- Sterile or individually, sterile packaged
- Consistently recovery
- Membranes meet ISO 7704
- Membranes available in various colors
- Without any hydrophobic adhesive areas

### Economical

- Ready to connect and easy to use
- Minimal amount of equipment needed

Biosart® 100 Monitors, 100 mL, 47 mm, individually packaged, sterile, 48 units

| Pore size | Membrane filter color   Grid color   | Order No.        |
|-----------|--------------------------------------|------------------|
| 0.2 µm    | Cellulose nitrate white   black      | 16401-47-07--ACK |
| 0.45 µm   | Cellulose nitrate white   black      | 16401-47-06--ACK |
| 0.45 µm   | Cellulose nitrate green   dark green | 16402-47-06--ACK |
| 0.45 µm   | Cellulose nitrate gray   white*      | 16403-47-06--ACK |

Biosart® 100 Monitors, 100 mL, 47 mm, packaged in trays, sterile, 48 units

|                      |                                      |                  |
|----------------------|--------------------------------------|------------------|
| 0.2 µm               | Cellulose nitrate white   black      | 16401-47-07----K |
| 0.45 µm<br>High Flow | Cellulose nitrate white   black      | 16401-47-H6----K |
| 0.45 µm              | Cellulose nitrate white   black      | 16401-47-06----K |
| 0.45 µm              | Cellulose nitrate green   dark green | 16402-47-06----K |
| 0.45 µm              | Cellulose nitrate gray   white*      | 16403-47-06----K |
| 0.8 µm               | Cellulose nitrate gray   white*      | 16403-47-04----K |
| 0.45 µm              | Regenerated cellulose white          | 16404-47-06----K |

Biosart® 100 Monitors, 100 mL, 47 mm, sterile, 48 units

|                      |                                 |                  |
|----------------------|---------------------------------|------------------|
| 0.45 µm<br>High Flow | Cellulose nitrate white   black | 16401-47-H6-V--K |
| 0.45 µm              | Cellulose nitrate white   black | 16401-47-06-V--K |
| 0.45 µm              | Cellulose nitrate gray   white* | 16403-47-06-V--K |
| 0.8 µm               | Cellulose nitrate gray   white* | 16403-47-04-V--K |

\* Gray membranes after wetting black

## Biosart® 100 Nutrient Media



### Safe and Reliable

- Presterilized media
- Certificate of quality for every batch
- In compliance with international standards
- Consistently recovery

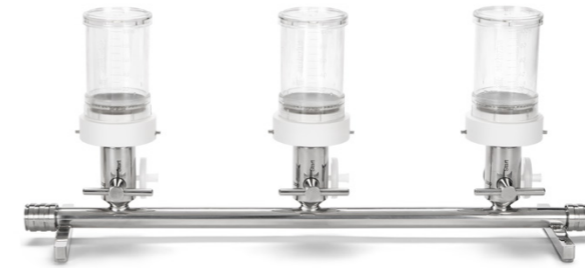
### Economical

- Ready-to-use
- Long shelf life

Biosart® 100 Nutrient Media, 2.5 mL, individually, sterile packaged in ampoules, 50 units

| Determination of                   | Media type                                  | Order No.        |
|------------------------------------|---|------------------|
| Total count                        | Caso (acc. USP)                             | 16400-02----CA-K |
|                                    | R2A (acc. EP)                               | 16400-02----RA-K |
|                                    | TGE   Total Count                           | 16400-02----TC-K |
|                                    | Total Count TTC                             | 16400-02----TZ-K |
| E. coli and coliforms              | m Endo                                      | 16400-02----EN-K |
|                                    | m FC  | 16400-02----MF-K |
|                                    | Lauryl Sulfate   Teepol                     | 16400-02----LS-K |
|                                    | Tergitol TTC                                | 16400-02----TT-K |
| Enterococci                        | KF Strep   Azide                            | 16400-02----KF-K |
| Pseudomonas aeruginosa             | Cetrimide                                   | 16400-02----CE-K |
| Yeasts and molds                   | Sabouraud (acc. USP)                        | 16400-02----SB-K |
|                                    | m Green yeast and mold   Schaufus Pottinger | 16400-02----MG-K |
|                                    | m Green yeast and mold selective            | 16400-02----GS-K |
|                                    | Wort  | 16400-02----WZ-K |
| Yeasts and molds and bacteria      | WL Nutrient   Wallerstein Nutrient          | 16400-02----WN-K |
| Bacteria in fermentation processes | WL Differential   Wallerstein Differential  | 16400-02----WL-K |
| Acid-tolerant microorganisms       | Orange Serum                                | 16400-02----OS-K |

## Microsart® Manifolds for Biosart® Monitors



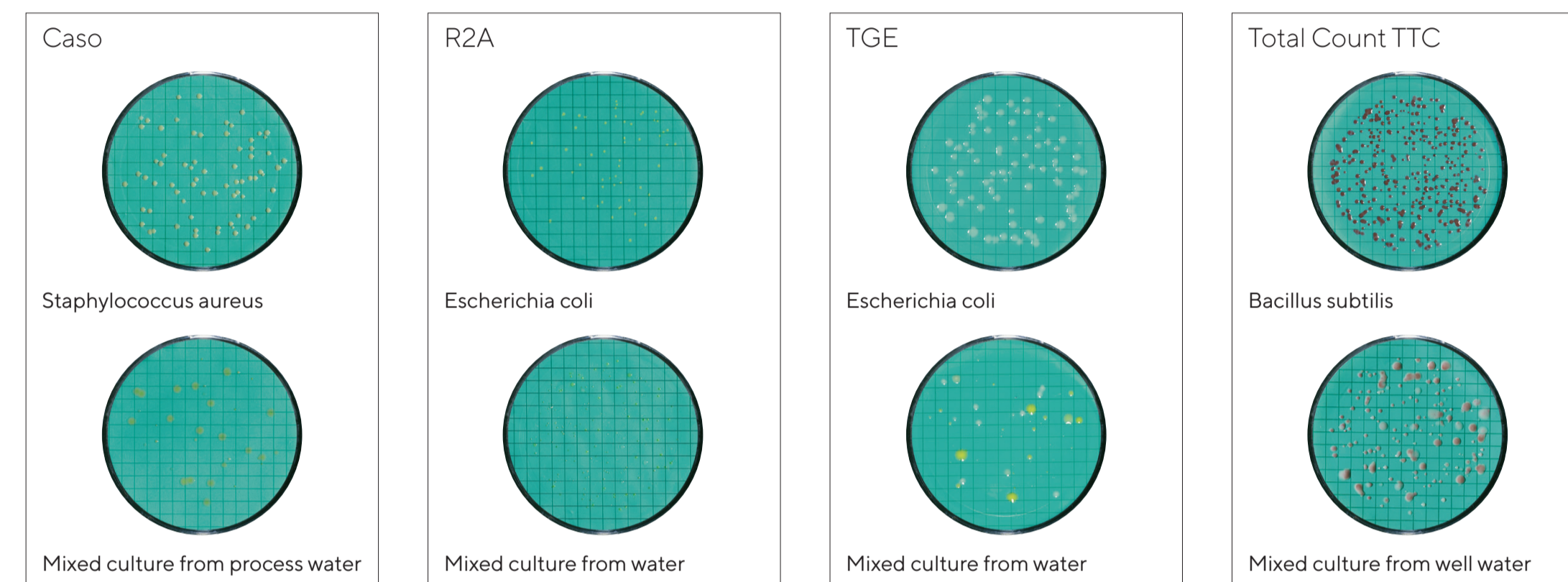
| Description  | Order No.   |
|--|-------------|
| Microsart® 1-branch stainless steel manifold with adapter for Biosart® Monitors  | 168M1-BS100 |
| Microsart® 2-branch stainless steel manifold with adapters for Biosart® Monitors | 168M2-BS100 |
| Microsart® 3-branch stainless steel manifold with adapters for Biosart® Monitors | 168M3-BS100 |
| Microsart® 6-branch stainless steel manifold with adapters for Biosart® Monitors | 168M6-BS100 |

## Biosart® 100 Monitor Adapter and Membrane Lifter

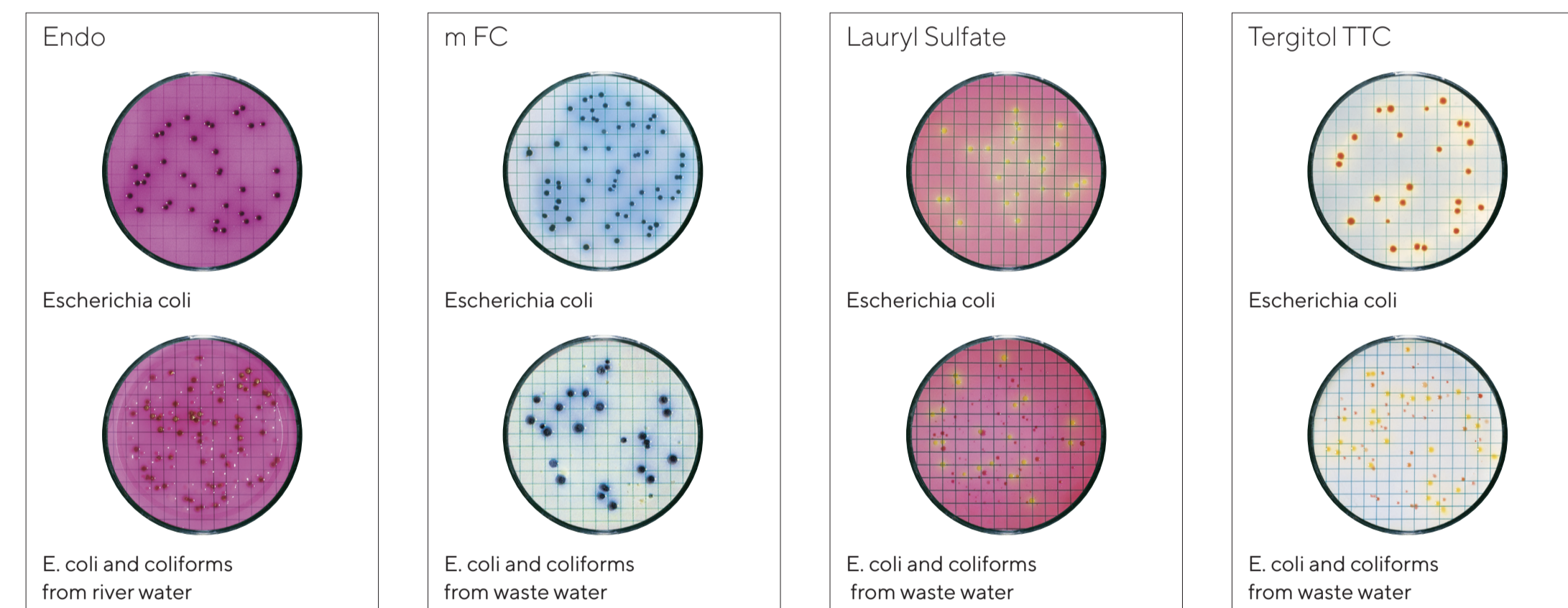
| Description                       | Adaptation  | Order No. |
|-----------------------------------|---|-----------|
| Biosart® 100 Adapter              | Adapter for use Biosart® 100 Monitors on Microsart® Manifolds | 16424     |
| Biosart® 100 Membrane Lifter, ABS | For easy transfer to agar                                     | 16417     |

## Biosart® 100 Nutrient Media

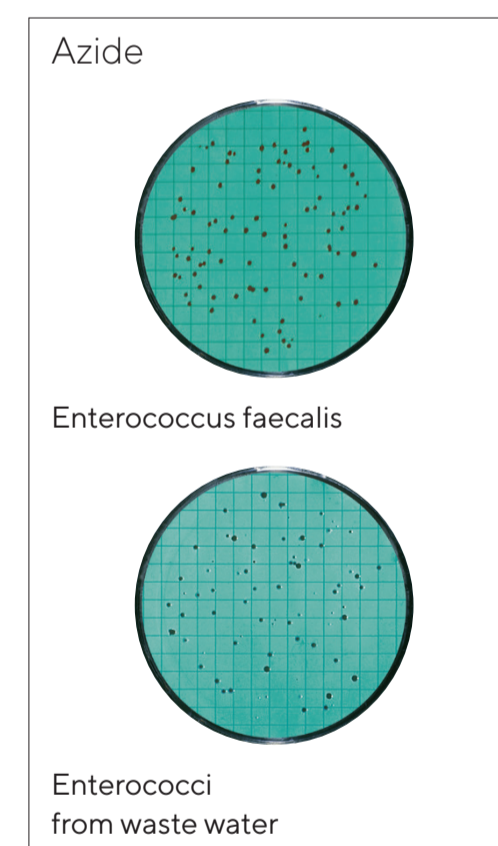
### Total colony count



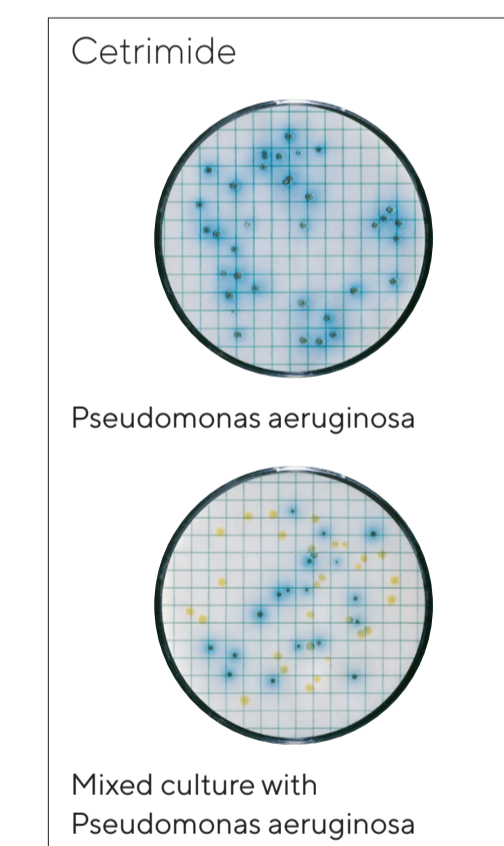
### E. coli and coliforms, Enterobacteria



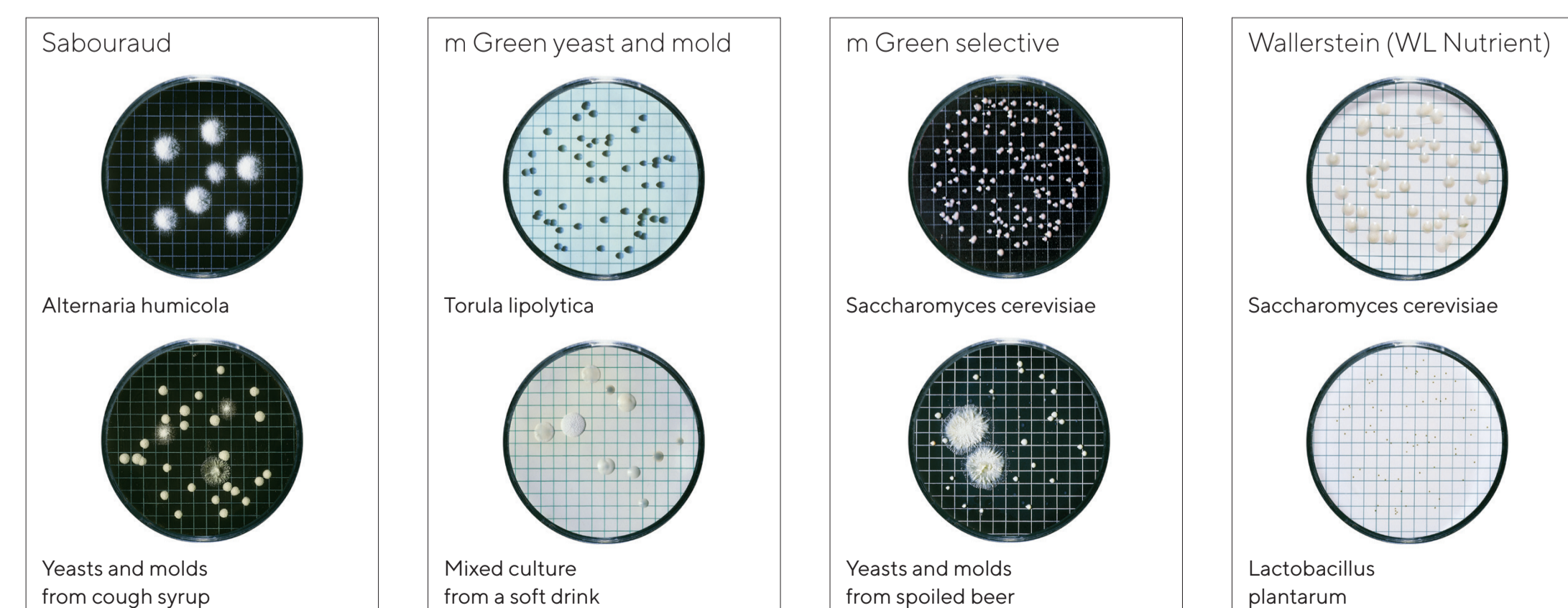
### Other faecal bacteria



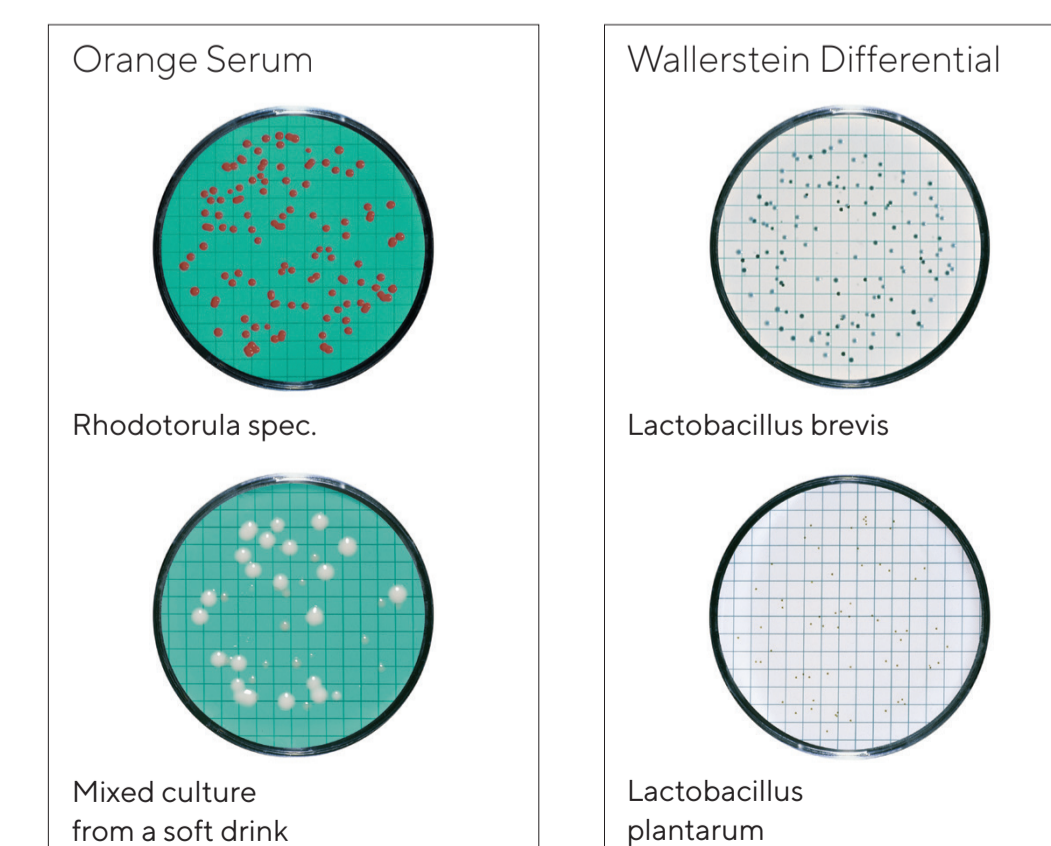
### Non-faecal, pathogenic bacteria



### Yeasts and molds



### Product-spoiling microorganisms



### Typical Application Examples

| Product  | Detection and enumeration of ...                          | Biosart® 100 Nutrient Media Type  |
|--|---|---|
| Beer   | Lactobacilli and Pediococci other beer spoiling organisms | Lactobacilli and Pediococci and   |
|  | Total colony count  | Total Count TTC   |
|  | Yeasts and molds  | Wallerstein Nutrient, Wort  |
| Foods  | Acid-tolerant microorganisms                              | Orange Serum  |
|  | Enterobacteria, E. coli and coliforms                     | Endo, m FC, Teepol   Lauryl Sulphate, Tergitol TTC  |
|  | Enterococci, Enterococcus faecalis                        | Azide   KF Strep  |
|  | Pseudomonas aeruginosa                                    | Cetrimide   |
|  | Total colony count  | Caso, TGE   Tryptone Glucose Extract  |
| Fruit juice  | Yeasts and molds  | Wort  |
|  | Enterobacteria, E. coli and coliforms                     | Endo, Tergitol TTC*   |
|  | Oenococcus and other product spoiling organisms           | Orange Serum, Wallerstein Differential  |
| Milk   | Yeasts and molds  | m Green yeast and mold   Schaufus Pottinger, Wallerstein Nutrient   |
|  | E. coli and coliforms                                     | Endo  |
| Pharmaceuticals, WFI, raw materials and cosmetics                  | Enterococci, Enterococcus faecalis                        | Azide   KF Strep  |
|  | Enterococci, Enterococcus faecalis                        | Azide   KF Strep  |
|  | Pseudomonas aeruginosa                                    | Cetrimide   |
|  | Total colony count  | Caso, R2A   |
|  | Yeasts and molds, Candida albicans                        | Sabouraud   |
| Soft drinks, concentrates  | Acid-tolerant microorganisms, Lactic-acid bacteria        | Orange Serum, Wallerstein Differential  |
|  | Enterobacteria, E. coli and coliforms                     | Endo  |
|  | Total colony count  | TGE   Tryptone Glucose Extract, Total Count TTC   |
|  | Yeasts and molds  | m Green yeast and mold   Schaufus Pottinger, m Green yeast and mold selective, Wallerstein Nutrient, Wort |
| Sugar, sugar products  | E. coli and coliforms                                     | Endo  |
|  | Total colony count  | Total Count TTC   |
|  | Yeasts and molds  | m Green yeast and mold   Schaufus Pottinger, m Green yeast and mold selective, Wort                       |
| Water (general quality), mineral water, natural water, waste water | Acid-tolerant microorganisms, Lactic-acid bacteria        | Orange Serum  |
|  | Enterobacteria, E. coli and coliforms                     | Endo, m FC, Teepol   Lauryl Sulphate, Tergitol TTC  |
|  | Enterococci, Enterococcus faecalis                        | Azide   KF Strep  |
|  | Pseudomonas aeruginosa                                    | Cetrimide   |
|  | Total colony count  | Caso, R2A, TGE   Tryptone Glucose Extract   |
|  | Yeasts and molds, Candida albicans                        | Sabouraud   |
| Wine   | Acetobacter   | Orange Serum (by adding 5 - 8% ethanol)   |
|  | Acid-tolerant microorganisms, Lactic-acid bacteria        | Orange Serum, Wallerstein Differential  |
|  | Yeasts and molds  | m Green yeast and mold   Schaufus Pottinger, Wallerstein Nutrient, Wort                                   |

\* These Biosart® 100 Media types are suitable for the determination of the mentioned microorganisms, although the media are not explicit declared in references.

### Easy work flow - reliable results

- 1 | Pour the sample**
- 2 | Apply vacuum and filter the sample**
- 3 | Add the Biosart® 100 Nutrient Media**
- 4 | Close the outlet**
- 5 | Remove the funnel**
- 6 | Incubate the petri dish**

**Remarks**  
The pictures show typical appearance of the mentioned microorganisms. In particular cases, color and shape of the colonies could vary from the expected habitus. Further tests may be necessary to validate the result.

Sartorius Stedim Biotech shall not be liable for consequential and/or incidental damage sustained by any customer from the use of its products.


Biosart® 100 Nutrient Media are subject to continuous product improvement as part of our product development program to align our products with changing application requirements. For current specifications and lot release criteria please visit our homepage under: [www.sartorius-stedim.com/BiosartMediaSearch](http://www.sartorius-stedim.com/BiosartMediaSearch).

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