

Biosart® 100 Nutrient Media

Microbiological Nutrient
Media for Filtration Units



Benefits

Some of the advantages you will benefit from when using Biosart® 100 Media:

Secure and Reliable

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

Economical

- Ready-to-use
- Long shelf life

Description

The membrane filtration method is the suitable technique for microbiological analysis of pharmaceuticals, water, cosmetics, foods and beverages. The use of ready-to-use disposable units is optimal for these applications. The membrane filtration method is worldwide accepted and complies with International Standards.

Biosart® 100 Monitors have been specifically designed for the detection and enumeration of microorganisms in pharmaceuticals, cosmetics, food, beverages, water 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.

Biosart® 100 Nutrient Media

Each box of Biosart® 100 Nutrient Media contains 50 ampoules with sterile media, each with 2.5 mL and a lot certificate. The media comply with international regulations and recommendations: International pharmacopoeias, DIN and ISO standards, the American Standards for Water and Foods, mineral water regulations, guidelines of the food and beverage industries.

Product Improvement

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.

Technical Data and Order Information

Biosart® 100 Nutrient Media, 2.5 mL, individually sterile filled in plastic ampoules, 50 pcs.

Detection Target	Media Type	Order No.	pH	Shelf Life (months)	Monitor Type ¹	Test Strains ²
Counting of total colony forming units						
Total count	Caso	16400-02----CA-K	7.3	12	2	01, 03, 05, 09, 18, 22, 25, 26
	R2A	16400-02----RA-K	7.2	12	2	01, 03, 05, 09, 18, 22, 26
	Total Count TGE	16400-02----TC-K	7.0	12	2	03, 07, 09, 18, 26
	Total Count TTC	16400-02----TZ-K	7.0	12	2	03, 07, 09, 18, 26
E. coli and coliforms, Enterobacteria						
E. coli and coliforms	m Endo	16400-02----EN-K	7.2	9	1	06, 07, 09, 21, 25, 28
	m FC	16400-02----MF-K	7.4	12	1	06, 07, 09, 11, 25
	Tergitol TTC	16400-02----TT-K	7.2	9	1	06, 07, 09, 11, 22
Other faecal bacteria						
Enterococci	Azide KF Strep	16400-02----KF-K	7.2	9	2	07, 08, 09, 22, 26
Non-faecal, pathogenic bacteria						
Pseudomonas aeruginosa	Cetrimide	16400-02----CE-K	7.2	12	1	04, 09, 21, 22, 26, 30
Yeasts and molds						
Yeasts and molds	m Green	16400-02----MG-K	4.6	12	1	03, 05, 20, 23, 24
	m Green Selective	16400-02----GS-K	4.6	12	3	03, 05, 09, 20, 23, 24
	Sabouraud	16400-02----SB-K	5.6	12	3	01, 05, 20, 23, 24
	Wallerstein WL Nutrient	16400-02----WN-K	5.5	12	1	05, 12, 19, 20, 23, 29
	Wort	16400-02----WZ-K	4.4	12	3	05, 20, 23, 24
Product-spoiling microorganisms						
Acid-tolerant microorganisms	Orange Serum	16400-02----OS-K	5.5	12	2	02, 05, 13, 14, 20, 23, 24
Bacteria in fermentation processes	Wallerstein Differential	16400-02----WL-K	5.5	12	1	12, 20, 23, 24, 29

¹The recommended Biosart® 100 Monitor Types are selected for optimum recovery of microorganisms on the corresponding nutrient media:

- 1) 16401: white membrane filter with black grid
- 2) 16402: green membrane filter with dark green grid
- 3) 16403: gray (after wetting black) membrane filter with white grid

²Test Strains [ATCC No.], [DSM No.]


01. *Aspergillus brasiliensis* 16404, 1988
02. *Bacillus cereus* 11778, 345
03. *Bacillus subtilis* subsp. *spizizenii* 6633, 347
04. *Brevundimonas diminuta* 19146, 1635
05. *Candida albicans* 10231, 1386
06. *Enterobacter aerogenes* 13048, 30053
07. *Enterococcus faecalis* 29212, 2570
08. *Enterococcus faecium* 19434, 20477
09. *Escherichia coli* 8739, 1576
10. *Geobacillus stearothermophilus* 7953, 5934
11. *Klebsiella pneumoniae* 13883, 30104
12. *Lactobacillus lindneri* DSM 20690
13. *Lactobacillus plantarum* subsp. *plantarum* 14917, 20174
14. *Leuconostoc mesenteroides* subsp. *mesenteroides* 8293, 20343
15. *Oenococcus oeni* 23279, 20252
16. *Lactobacillus acidophilus* 4356, 20079
17. Raw cane sugar solution (10%)
18. Tap water
19. *Pediococcus damnosus* 29358, 20331
20. *Penicillium commune* 10428, 2211
21. *Proteus mirabilis* 29906, 4479
22. *Pseudomonas aeruginosa* 9027, 1128
23. *Rhodotorula mucilaginosa* DSM 70403
24. *Saccharomyces cerevisiae* 9763, 1333
25. *Salmonella enterica* subsp. *enterica* serotype *typhimurium* 14028, 19587
26. *Staphylococcus aureus* subsp. *aureus* 6538, 799
27. *Staphylococcus epidermidis* 12228, 1798
28. *Escherichia coli* 25922, 1103
29. *Lactobacillus brevis* 14869, 20054
30. *Pseudomonas aeruginosa* 27853, 1117

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