

Endo Biosart® 100 Nutrient Media



Introduction

Coliform bacteria serve as indicator organisms for fecal contamination.

Endo Biosart® 100 Nutrient Media is used for the detection and enumeration of *E. coli* and coliform bacteria. It is designed for cultivating these microorganisms from raw materials, water (general quality), natural water, wastewater, beverages, food and other products.

Sodium sulfite and fuchsin function as inhibitors, suppressing the growth of Gram-positive bacteria. Colonies of coliform bacteria fermenting lactose as a carbon source turn pink due to the reaction of acetaldehyde, a fermentation product, with sodium sulfite and fuchsin. In *E. coli* colonies, this reaction is very strong, resulting in a metallic sheen. Non-coliform bacteria grow colorless.

Technical Specifications

Order No.	16400-02----EN-K (50 units)
Media	Endo
Color	Pink
Storage	Refrigerate (2 - 8 °C) after arrival*, dark and dry, use before expiry date on the label
Shelf Life	9 months
For Use With	Biosart® 100 Monitor (16401) with white membrane filter and black grid, 0.45 µm

* Data have shown constant performance in microbiological tests after storage at 22 °C for 14 days.

Media Formulation**

Ingredients	g/L
Thiopeptone	5
Tryptose	10
Yeast extract	1.5
Sodium chloride	5
Casitone	5
Lactose	12.5
Monopotassium phosphate	1.375
Dipotassium phosphate	4.375
Sodium lauryl sulfate	0.05
Basic Fuchsin	1.05
Ethanol	20 mL
Water (AP-Quality)	Ad 1,000 mL

pH 7.3 ± 0.25 (at room temperature)

** Formula adjusted, standardized to suit performance parameters.

Instructions

The Biosart® 100 Monitor is a sterile, ready-to-use disposable unit featuring an integrated membrane filter and cellulose pad. After filtration, add the Biosart® 100 Nutrient Media from the ampoule and apply vacuum for 1 second. Remove the disposable unit from the manifold and seal the outlet. Finally, detach the funnel and close the lid to transform the monitor into a petri dish.

Incubation Conditions

18 - 24 hours at 36 ± 2 °C.

Evaluation and Typical Results

Escherichia coli forms red colonies with a metallic sheen and a red dot at the underside of the membrane. Other coliforms grow as dark to light red colonies without metallic sheen. Colorless colonies of lactose-negative bacteria are not counted.

Microbiological Quality Control

Sterility: Qualitative

Incubation conditions: 14 days at 30 - 35 °C

Specification: No growth or turbidity

Productivity: Quantitative

Inoculum: 50 - 150 CFU

Incubation conditions: 18 - 24 hours at 36 ± 2 °C

Specification: ≥ 85% membrane filtration on control agar as reference

Microorganism	Test strain	Specification	Morphology
<i>Escherichia coli</i>	WDCM 00012	$P_R \geq 0.85$	Colonies with metallic green sheen
<i>Escherichia coli</i> ***	WDCM 00013	$P_R \geq 0.85$	Colonies with metallic green sheen

*** Tested on a regular basis.

P_R Productivity Ratio

Selectivity: Qualitative

Inoculum: 10⁴ CFU

Specification: No growth to small glassy colonies

Microorganism	Test strain	Specification
<i>Enterococcus faecalis</i>	WDCM 00087	No growth to small glassy colonies

Specificity: Qualitative

Inoculum: 10³ CFU

TSA spread plate as control

Microorganism	Test strain	Specification Morphology
<i>Klebsiella aerogenes</i> ***	WDCM 00175	Red colonies (examination after 18 ± 1 hours), metallic sheen on control
<i>Proteus mirabilis</i> ***	WDCM 00023	Glassy, beige to bright pink colonies
<i>Salmonella enterica</i> ***	WDCM 00031	Glassy, beige to bright pink colonies

*** Tested on a regular basis.


The incubation conditions recommended by Sartorius can be adjusted based on the type of samples, in accordance with the reference standards or customer requirements. Descriptions of typical results illustrate the usual appearance of the specified microorganisms. However, in certain cases, the color and shape of the colonies may differ from the expected appearance. Additional tests may be required to confirm the results. Sartorius shall not be liable for any consequential or incidental damages incurred by customers from the use of its products.

Germany

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

USA

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

 For further contacts, visit
www.sartorius.com