

Caso Biosart® 100 Nutrient Media



Introduction

Soybean-Casein Digest Biosart® 100 Nutrient Media is used for isolating microorganisms and for determining the total colony forming unit (CFU) count. The culture medium is designed for cultivating microorganisms in pharmaceuticals, cosmetics, raw materials, water (general quality), waste water, foods and other products.

Technical Specifications

Order No.	16400-02----CA-K (50 units)
Media	Caso
Color	Beige
Storage	Refrigerate (2 - 8 °C) after arrival*, dark and dry, use before expiry date on the label
Shelf Life	12 months
For Use With	Biosart® 100 Monitor (16402) with green membrane filter and dark green 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
Pancreatic digest of casein	17
Papaic digest of soybean meal	3
D(+)-Glucose monohydrate	2.5
Sodium chloride	5
Dipotassium phosphate	2.5
Water (AP-Quality)	Ad 1,000 mL

pH 7.4 ± 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 and for yeasts, molds and tap water up to 5 days at 30 - 35 °C.

Evaluation and Typical Results

Predominantly bacteria of different sizes, shapes and colors.

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 and for yeasts, molds and tap water up to 5 days at 30 - 35 °C

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

Microorganism	Test strain	Specification	Morphology
<i>Pseudomonas aeruginosa</i>	WDCM 00026	$P_R \geq 0.85$	Beige colonies
<i>Bacillus subtilis</i>	WDCM 00003	$P_R \geq 0.85$	Beige colonies, matt, wrinkled
<i>Candida albicans</i>	WDCM 00054	$P_R \geq 0.85$	Cream-colored colonies
<i>Staphylococcus aureus</i> ***	WDCM 00032	$P_R \geq 0.85$	Yellow colonies
<i>Salmonella enterica</i> ***	WDCM 00031	$P_R \geq 0.85$	Beige, glossy colonies
<i>Escherichia coli</i> ***	WDCM 00012	$P_R \geq 0.85$	Beige colonies
Tap water***	N/A	$P_R \geq 0.85$	Beige, yellow, bright red colonies

*** Tested on a regular basis.

P_R Productivity Ratio

Specificity: Qualitative

Inoculum: 10³ CFU

TSA spread plate as control

Microorganism	Test strain	Specification Morphology
<i>Aspergillus brasiliensis</i>	DSM 1988	Strong growth, black-haired colonies


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