

Wort Biosart® 100 Nutrient Media



Introduction

Wort Biosart® 100 Nutrient Media is used for the detection and enumeration of yeasts and molds. It is designed for cultivating these microorganisms in raw materials, beverages, beer, wine, soft drinks, concentrates, foods, and other products.

The medium contains carbon sources for an optimal nutrient supply and the low pH value support the growth of yeasts and molds while inhibiting bacterial contaminants.

Technical Specifications

Order No.	16400-02----WZ-K (50 units)
Media	Wort
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 (16403) with black membrane filter and white 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
Malt extract	15
D-Maltose	12.5
Dextrin	2.5
Casein peptone	1
Potassium hydrogen phosphate	1
Ammonium chloride	1
Glycerin	2.5 mL
Water (AP-Quality)	Ad 1,000 mL

pH 4.4 ± 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 to transform the monitor into a petri dish.

Incubation Conditions

3 - 5 days at 20 - 25 °C.

Evaluation and Typical Results

Yeasts usually develop smooth white or colored colonies. Molds generally form velvety or fluffy cottonlike colonies that look white in the early growth phase and may take various colors after conidiospore production.

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: 3 - 5 days at 20 - 25 °C

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

Microorganism	Test strain	Specification	Morphology
<i>Candida albicans</i>	WDCM 00054	$P_R \geq 0.85$	Creamy-white colonies
<i>Saccharomyces cerevisiae</i> ***	WDCM 00058	$P_R \geq 0.85$	Creamy-white colonies
<i>Rhodotorula mucilaginosa</i> ***	DSM 70403	$P_R \geq 0.85$	Orange-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>Penicillium commune</i>	DSM 2211	White-grey, fluffy 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