

# Wallerstein Differential Biosart® 100 Nutrient Media



## Introduction

Wallerstein differential Biosart® 100 Nutrient Media is used for the isolation and enumeration of bacteria encountered in breweries and industrial fermentations.

The medium contains all the essential ingredients for the growth of lactic acid bacteria, yeasts, and molds, similar to the Wallerstein nutrient medium. Additionally, cycloheximide has been added to suppress the growth of *Saccharomyces cerevisiae*.

# Technical Specifications

<b>Order No.</b>	16400-02----WL-K (50 units)
<b>Media</b>	Wallerstein Differential
<b>Color</b>	Turquoise
<b>Storage</b>	Refrigerate (2 - 8 °C) after arrival*, dark and dry, use before expiry date on the label
<b>Shelf Life</b>	12 months
<b>For Use With</b>	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
Peptic digest of casein	5
Ferric chloride	0.0025
Yeast extract	4
Dextrose	50
Magnesium sulfate	0.125
Monopotassium phosphate	0.55
Potassium chloride	0.425
Calcium chloride	0.125
Bromocresol green	0.022
Manganese sulfate	0.0025
Cycloheximide	0.006
Water (AP-Quality)	Ad 1,000 mL

pH 5.5 ± 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

Yeast and mold: 3 - 5 days at 26 ± 2 °C, aerobic.

Lactic acid bacteria: 3 - 7 days at 26 ± 2 °C, anaerobic.

### Evaluation and Typical Results

*Lactobacillus brevis* forms small green colored colonies with sharp edge. *Lactobacillus plantarum* grow as small greenish or yellow colored and round colonies with sharp edge. *Escherichia coli* grow as white to cream-colored colonies with sharp edge.

The growth of a *Saccharomyces cerevisiae* is inhibited by cycloheximide. If cycloheximide resistant yeast grow, they are creamy, green-white.

## 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:

Yeast and mold: 3 - 5 days at 26 ± 2 °C, aerobic

Lactic acid bacteria: 3 - 7 days at 26 ± 2 °C, anaerobic

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

Microorganism	Test strain	Specification	Morphology
<i>Lactobacillus brevis</i>	WDCM 00099	$P_R \geq 0.85$	Green colonies
<i>Lactobacillus lindneri</i> ***	DSM 20690	$P_R \geq 0.85$	Very small, glassy green colonies
<i>Rhodotorula mucilaginosa</i> ***	DSM 70403	$P_R \geq 0.85$	Bright red to red brown colonies

\*\*\* Tested on a regular basis.

$P_R$  Productivity Ratio

### Selectivity: Qualitative

Inoculum: 10<sup>4</sup> CFU

Specification: No growth | total inhibition

Microorganism	Test strain	Specification
<i>Saccharomyces cerevisiae</i>	WDCM 00058	No growth   total inhibition

### Specificity: Qualitative

Inoculum: 10<sup>3</sup> CFU

TSA spread plate as control

Microorganism	Test strain	Specification   Morphology
<i>Penicillium commune</i> ***	DSM 2211	Grey blue, fluffy 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](http://www.sartorius.com)