

PIKA Weihenstephan™ FastOrange™ B Agar

Catalog Number 4488196

Pub. no. MAN0009371 Rev. 1.0

Description	Amount	Storage
Agar for detection of beer-spoiling bacteria	12 X 170 mL	Store at room temperature. Protect from light. Refer to the expiration date on the label.

WARNING! Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from www.lifetechnologies.com/support.

IMPORTANT! Follow all applicable local, state/provincial, and/or national regulations for handling of biological samples.

Product description

PIKA Weihenstephan™ FastOrange™ B Agar is a nutrient medium developed for the detection of beer-spoiling bacteria in brewery samples. Positive results due to growth of acid-producing microorganisms, such as *Lactobacilli* or *Pediococci*, are detected by a violet-to-yellow color change of the medium.

The nutrient medium is not suitable for consumption.

Guidelines for use

Prepare agar

- Heat bottle in a 90°C water bath to melt the agar.
Alternatively, heat uncapped bottle in a microwave on a low setting until agar has melted.

IMPORTANT! Remove bottle caps before microwaving the bottle to avoid a possible explosion. Monitor the heating process carefully to make sure the agar does not boil over in the microwave.

- Prepare plates:

Sample type	Agar plates
Clear liquids and filtered samples	Streak plates: pour liquid agar into sterile, vented Petri dishes and let cool to solidify.
Yeast-containing and turbid, non-filterable liquids	Pour plates: store liquid agar at 55°C until ready to plate samples.

Avoid long holding times for liquid agar, and avoid repeatedly melting the agar.

IMPORTANT! Do not attempt to resterilize the agar by autoclaving or with other heat treatments, because the agar will lose its growth-supporting characteristics.

Follow these recommended guidelines for your sample type.

Clear samples (e.g., beer, water, or filtered sample)

- Add sample to FastOrange™ B Agar plates:
 - Membrane filters: place directly on agar plate.
 - Liquids: streak directly on agar plate.
- Incubate plates at the appropriate temperature for 2–7 days.

Yeast-containing and turbid, non-filterable samples

- Pour liquid sample into a sterile Petri dish.
The volume should not exceed 20–30% of the total plate volume.
- Add at least the same volume of FastOrange™ B Agar that is at 55°C.
- Mix by careful swirling and let cool to solidify.
- Incubate plates at the appropriate temperature for 2–7 days.

Incubation conditions

The choice of the oxygen content during incubation will lead to different specificity:

Oxygen content	Microorganisms
Anaerobic	Exclusive support of growth of beer-spoiling lactic acid bacteria, such as <i>Lactobacilli</i> and <i>Pediococci</i>
Aerobic	<ul style="list-style-type: none"> Lactic acid bacteria Acetic acid bacteria Indicator bacteria

Results

Sample type	Visual observations
Clear liquids and filtered samples	<ul style="list-style-type: none"> Growth of bacterial colonies on the medium. For acid-producing microorganisms, violet-to-yellow color change of the medium.
Yeast-containing and turbid, non-filterable liquids	<ul style="list-style-type: none"> Growth of bacterial colonies on or in the medium. For acid-producing microorganisms, violet-to-yellow color change of the medium.

Rarely, growth of yeast (for example, *Brettanomyces*) may occur, because of a resistance to Cycloheximide in the medium. Yeast and bacteria can be clearly differentiated by microscopic examination of colonies.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

For testing of Food and Environmental samples only.

The information in this guide is subject to change without notice.

DISCLAIMER: LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

Limited Use Label License No. 492: Environmental Testing, Quality Control/Quality Assurance Testing, Food and Agricultural Testing

Notice to Purchaser: The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product (a) to perform internal research for the sole benefit of the purchaser; and (b) for environmental testing, quality control/quality assurance testing, food and agricultural testing, including reporting results of purchaser's activities in environmental testing, quality control/quality assurance testing, food and agricultural testing for a fee or other commercial consideration. No other right is hereby granted expressly, by implication, or by estoppel. This product is for environmental testing, quality control/ quality assurance testing, food and agricultural testing and research purposes only.

The purchase of this product does not grant the purchaser any additional rights, including (without limitation) the right to transfer or resell the product in any form or the right to use the product as a therapeutic agent or diagnostics test component. For information on obtaining additional rights, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

© 2013 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation and/or its affiliates or their respective owners. FastOrange is a trademark of Gudrun Vogeser. PIKA Weihenstephan is a trademark of PIKA Weihenstephan GmbH.

Headquarters

5791 Van Allen Way | Carlsbad, CA 92008 USA | Phone +1 760 603 7200 | Toll Free in USA 800 955 6288

For support visit lifetechnologies.com/support or email techsupport@lifetech.com

lifetechnologies.com

10 September 2013

