
BRUCELLA AGAR

INTENDED USE

Remel Brucella Agar is a solid medium recommended for use in qualitative procedures for isolation and cultivation of *Brucella* spp. and other fastidious microorganisms.

SUMMARY AND EXPLANATION

Brucella Agar was originally developed for isolation of *Brucella* spp. from potentially contaminated material such as dairy products. It is prepared according to the formulation for Albimi broth.¹ It is recommended by the American Public Health Association (APHA) and the United States Department of Agriculture (USDA) for recovery of fastidious microorganisms from foods and dairy products.^{2,3} With the addition of blood, Brucella Agar may be used to determine the hemolytic reactions of aerobic and anaerobic microorganisms.

PRINCIPLE

Casein and meat peptones supply nitrogen, amino acids, and peptides necessary for bacterial growth. Dextrose is a ready source of energy. Yeast extract provides B-complex vitamins and enhances growth. Sodium chloride is a source of essential electrolytes and maintains osmotic equilibrium. Sodium bisulfite is a reducing agent. Agar is a solidifying agent.

REAGENTS (CLASSICAL FORMULA)*

Casein Peptone.....	15.0 g	Dextrose.....	1.0 g
Meat Peptone.....	5.0 g	Sodium Bisulfite.....	0.1 g
Sodium Chloride.....	5.0 g	Agar.....	15.0 g
Yeast Extract.....	2.0 g	Demineralized Water.....	1000.0 ml

pH 7.0 ± 0.2 @ 25°C

*Adjusted as required to meet performance standards.

PRECAUTIONS

This product is For Laboratory Use only. It is not intended for use in the diagnosis of disease or other conditions.

PREPARATION OF DEHYDRATED CULTURE MEDIUM

1. Suspend 43 g of medium in 1000 ml of demineralized water.
2. Heat to boiling with agitation to completely dissolve.
3. Sterilize by autoclaving at 121°C for 15 minutes or following established laboratory guidelines.
4. If blood is to be added, cool medium to 45-50°C and aseptically add defibrinated blood to a final concentration of 5%.
5. Dispense into appropriate containers.

PROCEDURE

CAUTION: *Brucella* is considered a Biosafety Level 2 (BSL-2) pathogen. Use appropriate personal protective equipment and devices, including a biological safety cabinet for processing samples and performance of aerosol-generating procedures. Refer to appropriate references for specific instructions regarding handling of samples suspected to contain *Brucella* spp.⁴

1. Consult current editions of appropriate references for the recommended procedure for sample preparation, inoculation, testing, and interpretation.

QUALITY CONTROL

Each lot number of Brucella Agar has been manufactured, packaged, and processed in accordance with current Good Manufacturing Practice regulations. All lot numbers have been tested using the following quality control organisms and have been found to be acceptable. Testing of control organisms should be performed in accordance with established laboratory quality control procedures. If aberrant quality control results are noted, sample results should not be reported.

CONTROL

Escherichia coli ATCC® 25922
Pseudomonas aeruginosa ATCC® 27853
Staphylococcus aureus ATCC® 25923

INCUBATION

Aerobic, 18-24 h @ 33-37°C
Aerobic, 18-24 h @ 33-37°C
Aerobic, 18-24 h @ 33-37°C

RESULTS

Growth
Growth
Growth

BIBLIOGRAPHY

1. MacFaddin, J.F. 1985. Media for Isolation-Cultivation-Identification-Maintenance of Medical Bacteria. Vol. 1. Williams & Wilkins, Baltimore, MD.
2. Hausler, W.J. 1976. Standard Methods for the Examination of Dairy Products. 14th ed. American Public Health Association, Washington, D.C.
3. United States Department of Agriculture. 2008. Laboratory Guidebook. Retrieved 03/03/2010 from: http://origin-www.fsis.usda.gov/Science/Microbiological_Lab_Guidebook/index.asp.
4. Chosewood, L.C. and D.E. Wilson. 2007. Biosafety in Microbiological and Biomedical Laboratories. 5th ed. U.S. Dept. of H.H.S., CDC, and National Institutes of Health, Washington, D.C.

Refer to the front of Remel *Technical Manual of Microbiological Media* for **General Information** regarding precautions, product storage and deterioration, sample collection, storage and transportation, materials required, quality control, and limitations.

ATCC® is a registered trademark of American Type Culture Collection.
IFU 1250, Revised March 24, 2010

Printed in U.S.A.

remel

12076 Santa Fe Drive, Lenexa, KS 66215, USA
General Information: (800) 255-6730 Website: www.remel.com Email: remel@remel.com
Local/International Phone: (913) 888-0939 International Fax: (913) 895-4128