

BRUCELLA AGAR (w/ 5% SHEEP BLOOD)

INTENDED USE

Remel Brucella Agar (w/ 5% Sheep Blood) is a solid medium recommended for use in qualitative procedures for primary isolation and cultivation of both fastidious aerobic and anaerobic microorganisms.

SUMMARY AND EXPLANATION

Brucella Agar was originally developed for isolation of *Brucella* spp. from potentially contaminated material such as dairy products.¹ It is prepared according to the formulation of the American Public Health Association for Albimi broth.^{2,3} Brucella Agar (w/ 5% Sheep Blood) is recommended for isolation of obligate and facultative anaerobes from clinical specimens.^{4,5}

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. Sheep blood provides growth factors required by certain anaerobic organisms and allows for demonstration of hemolysis by aerobic and anaerobic bacteria. 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	Sheep Blood.....	5 %
Yeast Extract.....	2.0 g	Agar.....	15.0 g
		Demineralized Water.....	1000.0 ml

pH 7.0 ± 0.2 @ 25°C

*Adjusted as required to meet performance standards.

PROCEDURE

1. Inoculate and streak the specimen as soon as possible after it is received in the laboratory.
2. If material is being cultured directly from a swab, roll the swab over a small area of the agar surface and streak for isolation.
3. Incubate aerobically or anaerobically at 33-37°C for up to 72 hours.
4. Confirm the growth of anaerobic organisms by subculture to aerobic blood agar plate.

QUALITY CONTROL

All lot numbers of Brucella Agar (w/ 5% Sheep Blood) have been tested using the following quality control organisms and have been found to be acceptable. This quality control testing meets or exceeds CLSI standards.⁶ Testing of control organisms should be performed in accordance with established laboratory quality control procedures. If aberrant quality control results are noted, patient results should not be reported.

CONTROL

- **Bacteroides fragilis* ATCC® 25285
- **Clostridium perfringens* ATCC® 13124
- **Escherichia coli* ATCC® 25922
- **Fusobacterium nucleatum* ATCC® 25586
- **Peptostreptococcus anaerobius* ATCC® 27337
- **Prevotella melaninogenica* ATCC® 25845
- **Staphylococcus aureus* ATCC® 25923
- **Streptococcus pneumoniae* ATCC® 6305
- **Streptococcus pyogenes* ATCC® 19615

®CLSI recommended organism

INCUBATION

- Anaerobic, up to 48 h @ 33-37°C
- Anaerobic, up to 48 h @ 33-37°C
- Aerobic, 18-24 h @ 33-37°C
- Anaerobic, up to 48 h @ 33-37°C
- Anaerobic, up to 48 h @ 33-37°C
- Anaerobic, up to 48 h @ 33-37°C
- Aerobic, 18-24 h @ 33-37°C
- Aerobic, 18-24 h @ 33-37°C
- Aerobic, 18-24 h @ 33-37°C

RESULTS

- Growth
- Growth, beta hemolysis
- Growth
- Growth
- Growth
- Growth
- Growth
- Growth, alpha hemolysis
- Growth, beta hemolysis

BIBLIOGRAPHY

1. Murray, P.R., E.J. Baron, J.H. Jorgensen, M.L. Landry, and M.A. Pfaller. 2007. Manual of Clinical Microbiology. 9th ed. ASM Press, Washington, D.C.
2. MacFaddin, J.F. 1985. Media for Isolation-Cultivation-Identification-Maintenance of Medical Bacteria. Vol. 1. Williams & Wilkins, Baltimore, MD.
3. Hausler, W.J. 1976. Standard Methods for the Examination of Dairy Products. 14th ed. American Public Health Association, Washington, D.C.
4. Summanen, P., E.J. Baron, D.M. Citron, C. Strong, H.M. Wexler, and S.M. Finegold. 1993. Wadsworth Anaerobic Bacteriology Manual. 5th ed. Star Publishing Co., Belmont, CA.
5. Engelkirk, P.G., J. Duben-Engelkirk, and V.R. Dowell. 1992. Principles and Practice of Clinical Anaerobic Bacteriology. Star Publishing Co., Belmont, CA.
6. Clinical and Laboratory Standards Institute (CLSI). 2004. Quality Control for Commercially Prepared Microbiological Culture Media; Approved Standard, 3rd ed. M22-A3. CLSI, Wayne, PA.

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

ATCC® is a registered trademark of American Type Culture Collection.
IFU 1252, Revised March 18, 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