

BISMUTH SULFITE AGAR BASE

INTENDED USE

Remel Bismuth Sulfite Agar Base is a medium recommended for use in qualitative procedures for the isolation and differentiation of *Salmonella* species from clinical specimens and other materials.

SUMMARY AND EXPLANATION

Bismuth Sulfite Agar was introduced by Wilson and Blair in 1927 for selective isolation of typhoid and paratyphoid groups of bacteria from stool specimens.¹ The formulation was modified by Hajna in 1951, and by Hajna and Damon in 1956, by substituting bismuth ammonium citrate and ferric ammonium citrate for bismuth citrate and ferrous sulfate, respectively.^{2,3} Bismuth Sulfite Agar has been recommended by the *United States Pharmacopeia* (USP) for use in microbial limit testing.⁴ It is recommended by the Food and Drug Administration (FDA) and the American Public Health Association (APHA) for isolation of *Salmonella* and other pathogens from foods and dairy products.⁵⁻⁸

PRINCIPLE

Beef extract and meat peptone supply amino acids, peptides, carbohydrates, vitamins, and nitrogenous compounds required for the growth of microorganisms. Dextrose provides a ready source of energy. Ferrous sulfate is an indicator of hydrogen sulfide production. Bismuth is a heavy metal having inhibitory properties for certain microorganisms. Brilliant green dye is a selective agent. The characteristic black or green coloration of *Salmonella* species is a metallic precipitate that forms when hydrogen sulfide, produced by *Salmonella* species from sulfur compounds in the medium, reacts with ferrous sulfate.

REAGENTS (CLASSICAL FORMULA)*

Meat Peptone.....	10.0 g	Disodium Phosphate.....	4.0 g
Sodium Sulfite.....	6.0 g	Bismuth Ammonium Citrate.....	2.0 g
Beef Extract.....	5.0 g	Ferrous Sulfate.....	0.3 g
Dextrose.....	5.0 g	Brilliant Green.....	25.0 mg
		Agar.....	20.0 g

pH 7.6 ± 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 52 g of medium in 1000 ml of demineralized water.
2. Heat to boiling with agitation to completely dissolve. **Do not autoclave.**
3. Cool medium to 50-55°C and mix to evenly disperse the precipitate.
4. Dispense into plates. (**Note:** Medium should be used the day it is prepared or within user's validated parameters for use.)

PROCEDURE

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

QUALITY CONTROL

Each lot number of Bismuth Sulfite Agar Base 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

Salmonella enterica serovar Arizonae ATCC® 13314
Salmonella enterica serovar Choleraesuis ATCC® 13312
Salmonella enterica serovar Typhimurium ATCC® 14028
Salmonella enterica serovar Typhimurium ATCC® 19430
Escherichia coli ATCC® 25922
Pseudomonas aeruginosa ATCC® 27853
Enterococcus faecalis ATCC® 29212
Shigella flexneri ATCC® 12022

INCUBATION

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

RESULTS

Growth, black colonies w/ metallic sheen
Growth, black colonies w/ metallic sheen
Growth, black colonies w/ metallic sheen
Growth, black colonies w/ metallic sheen
Marked inhibition
Marked inhibition
No growth
No growth

BIBLIOGRAPHY

1. Wilson, W.J. and E.M. McV. Blair. 1927. J. Hyg. 26:374-391.
2. Hajna, A.A. 1951. Public Health Lab. 9:23-28.
3. Hajna, A.A. and S.R. Damon. 1956. Appl. Microbiol. 4:341-345.
4. The United States Pharmacopeia. 2009. 32nd ed. United States Pharmacopeial Convention, Rockville, MD.
5. Wehr, H.M. and J.F. Frank. 2004. Standard Methods for the Examination of Dairy Products. 17th ed. APHA, Washington, D.C.
6. Eaton, A.D., L.S. Clesceri, E.W. Rice, and A.E. Greenberg. 2005. Standard Methods for the Examination of Water and Wastewater. 21st ed. APHA, Washington, D.C.
7. Downes, F.P. and K. Ito. 2001. Compendium of Methods for the Microbiological Examination of Foods. 4th ed. APHA, Washington, D.C.
8. Food and Drug Administration. 2000. Bacteriological Analytical Manual Online. AOAC International, Gaithersburg, MD. <http://www.fda.gov/Food/ScienceResearch/LaboratoryMethods/BacteriologicalAnalyticalManualBAM/ucm055778.htm>.

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 452402, Revised August 23, 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