

BACTEROIDES FRAGILIS ISOLATION AGAR (Bacteroides Bile Esculin) (BBE)

INTENDED USE

Remel Bacteroides Fragilis Isolation Agar is a solid medium recommended for use in qualitative procedures for the isolation and presumptive identification of *Bacteroides fragilis* group.

SUMMARY AND EXPLANATION

Members of the *B. fragilis* group are frequently isolated from human clinical infections.¹ Because clinical specimens for anaerobic culture frequently contain mixed flora, the use of selective media is recommended.² Livingston et al. formulated a selective and differential medium, Bacteroides Esculin Agar, which provided recovery and presumptive identification of the *Bacteroides fragilis* group.³

PRINCIPLE

The selectivity of this medium is provided by addition of gentamicin and bile. Gentamicin inhibits most facultative anaerobes and bile inhibits anaerobic gram-negative bacilli with the exception of *B. fragilis* group and other bile resistant *Bacteroides* and *Fusobacteria*. Differentiation is facilitated by the hydrolysis of esculin to esculetin by the *B. fragilis* group. Ferric ammonium citrate is added to the medium to react with esculetin and produce a brown-black complex surrounding the colony. Vitamin K is a growth factor. Hemin also is a growth factor and allows testing for catalase production.⁴

REAGENTS (CLASSICAL FORMULA)*

Bile (Oxgall)	20.0 g	Ferric Ammonium Citrate	0.5 g
Casein Peptone	15.0 g	Gentamicin	0.1 g
Meat Peptone	5.0 g	Vitamin K	10.0 mg
Sodium Chloride	5.0 g	Hemin	5.0 mg
Esculin	1.0 g	Agar	15.0 g
		Deminerlized Water	1000.0 ml

pH 7.0 ± 0.2 @ 25°C

*Adjusted as required to meet performance standards.

PROCEDURE

1. Prior to use, reduce the plates for a minimum of 24 hours by placing them in an anaerobic environment at room temperature.
2. Inoculate specimens for anaerobic culture on both selective and non-selective media as soon as possible after receipt in the laboratory; streak plates for isolation.
3. Incubate anaerobically at 33-37°C for 48-72 hours.
4. Confirm anaerobic growth by subculture to an aerobic blood agar plate.

QUALITY CONTROL

All lot numbers of Bacteroides Fragilis Isolation Agar 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, patient results should not be reported.

CONTROL

Bacteroides fragilis ATCC® 25285
Bacteroides thetaiotaomicron ATCC® 29148
Clostridium perfringens ATCC® 13124
Escherichia coli ATCC® 25922
Pseudomonas aeruginosa ATCC® 27853
Staphylococcus aureus ATCC® 25923

INCUBATION

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, blackening of media
Growth, blackening of media
Inhibition (partial to complete)
Inhibition (partial to complete)
Inhibition (partial to complete)
Inhibition (partial to complete)

LIMITATIONS

1. Fusobacteria are resistant to 20% bile and may grow on Bacteroides Fragilis Isolation Agar; however, *Fusobacterium* spp. are catalase-negative.⁴
2. Some enterococci and enteric gram-negative bacilli may grow, but the colonial morphology produced is not characteristic of *B. fragilis* group.⁴
3. Some penicillin-resistant strains of *Bacteroides* are inhibited by 20% bile and may not grow on this medium.⁴
4. Yeast is not inhibited on this medium; however, if growth occurs the colonies will not cause blackening of the agar.⁴

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. Dowell, V.R., Jr., G.L. Lombard., F.S. Thompson, and A.Y. Armfield. 1977. CDC Laboratory Manual. CDC, Atlanta, GA.
3. Livingston, S.J., S.D. Kominos, and R.B. Yee. 1978. J. Clin. Microbiol. 7:448-453.
4. MacFaddin, J.F. 1985. Media for Isolation-Cultivation-Identification-Maintenance of Medical Bacteria. Vol. 1. Williams & Wilkins, Baltimore, MD.

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 1104, Revised September 18, 2007

Printed in U.S.A.

remel

12076 Santa Fe Drive, Lenexa, KS 66215, USA

General Information: (800) 255-6730 Technical Service: (800) 447-3641 Order Entry: (800) 447-3635

Local/International Phone: (913) 888-0939 International Fax: (913) 895-4128

Website: www.remel.com Email: remel@remel.com