

CATARRHALIS SELECTIVE MEDIUM

INTENDED USE

Remel Catarrhalis Selective Medium is a solid medium recommended for use in the primary isolation of *Moraxella catarrhalis* from clinical specimens containing mixed flora.

SUMMARY AND EXPLANATION

Moraxella catarrhalis is now recognized as a significant pathogen associated with upper and lower respiratory tract infection.^{1,2} Timely isolation and identification of this organism is important because most strains produce beta-lactamase and are resistant to penicillin and ampicillin. Isolation of *M. catarrhalis* from respiratory specimens is complicated due to the high numbers of normal flora. Several selective and differential media formulations for *M. catarrhalis* have been proposed.³ These formulations allowed for growth of *Neisseria* spp. which made recognition and differentiation of *M. catarrhalis* difficult. In 1988, Vanechoutte et al. developed a highly selective medium which significantly increased the isolation of *M. catarrhalis* by decreasing the recovery of normal respiratory flora, especially *Neisseria* spp.⁴

PRINCIPLE

Casein and soy peptones are a source of nitrogen, amino acids, and peptides. Sheep blood supplies essential growth factors. Amphotericin B, trimethoprim, vancomycin, and sodium acetazolamide suppress the growth of normal respiratory flora. Amphotericin B is active against yeast and filamentous fungi. Vancomycin is active primarily against gram-positive bacteria, while trimethoprim inhibits gram-negative organisms. Sodium acetazolamide (a synthetic sulfonamide) is a potent inhibitor of carbonic anhydrase, a rare bacterial enzyme commonly produced by *Neisseria* spp.

REAGENTS (CLASSICAL FORMULA)*

Casein Peptone.....	15.0 g	Trimethoprim	5.0 mg
Soy Peptone.....	5.0 g	Amphotericin B	2.0 mg
Sodium Chloride.....	5.0 g	Sheep Blood.....	5 %
Sodium Acetazolamide	10.0 mg	Agar	15.0 g
Vancomycin.....	10.0 mg	Demineralized Water.....	1000.0 ml

pH 7.3 ± 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 a swab specimen is received, roll the swab over a small area of the agar surface and streak for isolation.
3. Incubate plate aerobically (not in CO₂) at 33-37°C for 24-48 hours.
4. Examine plate for typical colony morphology. *M. catarrhalis* forms a "hockey puck"-like colony which may be nudged across the plate intact with a bacteriological loop.²

QUALITY CONTROL

All lot numbers of Catarrhalis Selective Medium 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

Moraxella catarrhalis ATCC®25238
Candida albicans ATCC®10231
Neisseria sicca ATCC®9913
Proteus mirabilis ATCC®12453
Staphylococcus aureus ATCC®25923

INCUBATION

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

RESULTS

Good growth
Inhibition (marked to complete)
Inhibition (marked to complete)
Inhibition (marked to complete)
Inhibition (marked to complete)

PERFORMANCE CHARACTERISTICS

In a collaborative study, 792 sputum specimens were plated on routine primary isolation media and Catarrhalis Selective Medium. *M. catarrhalis* was isolated from 107 (14%) sputum specimens; of these, 34% were recovered only on Catarrhalis Selective Medium.⁵

LIMITATIONS

1. Other miscellaneous organisms, including pseudomonads, may grow on this medium. It is important to accurately identify colonies morphologically typical of *M. catarrhalis* with a Gram stain, oxidase, and traditional carbohydrate or enzyme tests.
2. This product is only part of the overall scheme for isolation and identification of *M. catarrhalis*. Further biochemical testing is required for identification of the test isolate. Consult appropriate references.^{2,6}
3. The inhibitory activity of sodium acetazolamide against carbonic anhydrase can be overcome by a carbon dioxide atmosphere.⁴

BIBLIOGRAPHY

1. Yuen, K.Y., W.H. Seto, and S.G. Ong. 1989. *J. Infect.* 19:251-256.
2. 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.
3. Soto-Hernandez, J.L., D. Nunley, S. Holtsclaw-Berk, and S.L. Berk. 1988. *J. Clin. Microbiol.* 26:405-408.
4. Vaneechoutte, M., G. Verschraegen, G. Claeys, and A. Van Den Abeele. 1988. *J. Clin. Microbiol.* 26:2544-2548.
5. Perry, J.L., S. Ainsworth, and G.R. Miller. 1990. 30th Interscience Conference on Antimicrobial Agents and Chemotherapy (ICAAC) Annual Meeting. Abstract 1066, p. 264.
6. Forbes, B.A., D.F. Sahm, and A.S. Weissfeld. 2007. *Bailey and Scott's Diagnostic Microbiology*. 12th ed. Mosby Elsevier, St. Louis, MO.

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 1575, Revised September 20, 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