

THAYER MARTIN AGAR (IMPROVED)

INTENDED USE

Remel Thayer Martin Agar (Improved) is a solid medium recommended for use in qualitative procedures for isolation of *Neisseria gonorrhoeae* and *Neisseria meningitidis*.

SUMMARY AND EXPLANATION

In 1945, Johnston introduced a medium to grow *N. gonorrhoeae* in 24 hours.¹ This was the advent of GC Agar Base. In 1964, Thayer and Martin formulated a selective medium for isolation of *N. gonorrhoeae* and *N. meningitidis* that contained antibiotics capable of suppressing the commensal microbial flora present in the anal canal, vagina, pharynx, and other sites.² The antibiotics used initially, ristocetin and polymyxin B, were replaced in 1966 by vancomycin, colistimethate, and nystatin.³ Since 1970, the addition of trimethoprim to Thayer-Martin media to suppress swarming of *Proteus* has gained wide-spread acceptance.⁴ In subsequent years, multiple studies have demonstrated improved recovery rates for pathogenic *Neisseria* spp. using selective media.⁵⁻⁷ Martin and Lewis recommended the use of anisomycin, considered a more stable antifungal agent, in place of nystatin to increase the shelf life of the media.⁸

PRINCIPLE

Thayer Martin Agar (Improved) contains hemoglobin, which provides the X factor (hemin), and GCHI Enrichment, which provides the V factor, vitamins, amino acids, coenzymes, and dextrose. The vancomycin is increased over earlier Thayer Martin formulas to provide improved inhibition of gram-positive cocci. Anisomycin is added to suppress the growth of *Candida albicans* and trimethoprim lactate effectively suppresses the swarming of *Proteus* spp. The medium also contains colistin to inhibit most gram-negative organisms, including *Pseudomonas* spp. and an increased concentration of dextrose which allows for better growth of gonococci. The tablet provided in the JEMBEC™ and Pili-Pocket systems liberates carbon dioxide into the enclosed pack when the tablet comes in contact with the water vapor present in the medium.

REAGENTS (CLASSICAL FORMULAE)*

Casein Peptone.....	7.5 g	Anisomycin	20.0 mg
Meat Peptone.....	7.5 g	Colistin.....	7.5 mg
Sodium Chloride.....	5.0 g	Trimethoprim Lactate.....	6.25 mg
Dipotassium Phosphate	4.0 g	Vancomycin	4.0 mg
Dextrose.....	1.5 g	Hemoglobin Solution	350.0 ml
Corn Starch.....	1.0 g	●GCHI Enrichment	10.0 ml
Monopotassium Phosphate.....	1.0 g	Agar.....	10.0 g
		Demineralized Water.....	650.0 ml

pH 7.2 ± 0.2 @ 25°C

●GCHI Enrichment:

Glucose.....	100.0 g	Coccarboxylase.....	0.1 g
Cysteine Hydrochloride.....	25.9 g	Guanine Hydrochloride.....	0.03 g
L-Glutamine.....	10.0 g	Ferric Nitrate.....	0.02 g
L-Cystine.....	1.1 g	p-Aminobenzoic Acid.....	0.013 g
Adenine.....	1.0 g	Vitamin B12.....	0.01 g
NAD.....	0.25 g	Thiamine Hydrochloride.....	0.003 g
		Demineralized Water.....	1000.0 ml

*Adjusted as required to meet performance standards.

PROCEDURE

Note: Inoculate the specimen as soon as possible after receipt in the laboratory. Inoculation of selective and nonselective media from specimens likely to contain commensal microbial flora has been shown to increase the recovery of pathogenic *Neisseria* spp.^{9,10}

Transgrow Procedure:

1. Hold Transgrow medium in an upright position during inoculation to minimize the loss of carbon dioxide. Remove the bottle cap for only the shortest period necessary. Inoculate the medium by placing the swab at the bottom of the bottle and crossing from side to side over the surface of the medium. Remove and discard the swab.
2. Tighten cap securely after inoculation.
3. If the bottle is to be transported after inoculation, viability of the organism may be extended by incubating the culture for 16-20 hours at 33-37°C prior to shipment.

Plate Procedure:

1. Inoculate the plate by rolling the swab over the agar surface in a large "Z" pattern to sufficiently transfer the specimen.
2. Cross-streak the plate using a sterile inoculating loop to achieve isolated colonies.
3. Incubate plate in 3-7% CO₂ at 33-37°C and examine after 24-48 hours. Continue incubation of negative plates for 72-96 hours before reporting as negative.

JEMBEC™ / Pili-Pocket Plate Procedure for Carbon Dioxide Generation:

1. Using forceps, remove the CO₂ tablet from the foil pouch and place in the well of the plate.
2. Inoculate the plate by rolling the swab over the agar surface in a large "Z" pattern to sufficiently transfer the specimen.
3. Cross-streak the plate using a sterile inoculating loop to achieve isolated colonies.
4. Secure the top of the plate tightly and label with the patient information.
5. Place the plate in an environmental maintenance pouch, seal securely, and incubate in an inverted position.

EXPECTED RESULTS

Colonies of *N. gonorrhoeae* are translucent, raised, gray, and may be mucoid. *N. meningitidis* colonies are larger than *N. gonorrhoeae*, bluish-gray in color, and may be mucoid. Test isolates should be examined by Gram stain to verify they are gram-negative diplococci. Additional biochemical and/or serological testing is required for definitive identification, following established laboratory guidelines. Consult appropriate references for further instructions.^{9,10}

QUALITY CONTROL

All lot numbers of Thayer Martin Agar (Improved) 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.¹¹ 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 ORGANISM

Neisseria gonorrhoeae ATCC® 43070
**Neisseria gonorrhoeae* ATCC® 43069
**Neisseria meningitidis* ATCC® 13090
**Candida albicans* ATCC® 60193
**Escherichia coli* ATCC® 25922
**Neisseria sicca* ATCC® 9913
**Proteus mirabilis* ATCC® 43071
**Staphylococcus epidermidis* ATCC® 12228

INCUBATION

CO₂, 24-48 h @ 33-37°C
CO₂, 24-48 h @ 33-37°C
CO₂, 24-48 h @ 33-37°C
CO₂, 24-48 h @ 33-37°C
CO₂, 24-48 h @ 33-37°C
CO₂, 24-48 h @ 33-37°C
CO₂, 24-48 h @ 33-37°C
CO₂, 24-48 h @ 33-37°C

RESULTS

Growth
Growth
Growth
Inhibition (partial)
Inhibition (partial)
Inhibition (complete)
Inhibition (partial)
Inhibition (partial)

*CLSI recommended organism

LIMITATIONS

1. Cultures for pathogenic *Neisseria* should be incubated in 3-7% CO₂. Higher concentrations of CO₂ may be inhibitory to some strains.⁹
2. Organisms other than *N. gonorrhoeae* and *N. meningitidis* may grow on Thayer Martin Agar (Improved). Further testing is required for identification confirmation following established laboratory procedures. Consult appropriate references for further instructions.^{9,10}
3. Media containing antibiotics may be too selective for some strains of pathogenic *Neisseria*. Even though the clinical symptoms or microscopic examination may suggest infection with *Neisseria* spp., recovery of the pathogen may fail due to overgrowth with commensal microbial flora.^{9,10}

BIBLIOGRAPHY

1. Johnston, J. 1945. J. Vener. Dis. Inform. 26:239.
2. Thayer, J.D. and J.E. Martin, Jr. 1964. Public Health Rep. 79:49.
3. Thayer, J.D. and J.E. Martin, Jr. 1966. Public Health Rep. 81:559-562.
4. Riddell, R.H. and A.C. Buck. 1970. J. Clin. Pathol. 23:481-483.
5. Cross, R.C., M.B. Hoger, R. Neibaur, B. Pasternack, and F.J. Brady. 1971. Public Health Rep. 86:990-992.
6. Chen, N.C., S.S. Hipp, W.D. Lawton, and H.A. Gaafar. 1974. Health Lab. Sci. 11:173-177.
7. Faur, Y. C., M.H. Weisburd, and M.E. Wilson. 1978. Health Lab. Sci. 15:22-27.
8. Mirrett, S., L.B. Reller, and J.S. Knapp. 1981. J. Clin. Microbiol. 14:94-99.
9. 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.
10. Forbes, B.A., D.F. Sahm, and A.S. Weissfeld. 2007. Bailey and Scott's Diagnostic Microbiology. 12th ed. Mosby Elsevier, St. Louis, MO.
11. 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.
JEMBEC™ is a trademark of Miles Scientific.

IFU 1884, Revised March 31, 2008

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