

A-8 AGAR SELECTIVE and NONSELECTIVE

INTENDED USE

Remel A-8 Agar, Selective and Nonselective, are solid media recommended for use in qualitative procedures for the cultivation, identification, and differentiation of *Ureaplasma urealyticum* and *Mycoplasma hominis*.

SUMMARY AND EXPLANATION

The discovery of a urease enzyme system in *U. urealyticum* made methods for identification of ureaplasmas possible.¹ Urease enzymes hydrolyze urea with the production of ammonia and resultant alkalinity. A differential medium for *Ureaplasma* (A-7 Agar) was developed by Shepard and Lunceford which contained urea with manganous sulfate as an indicator. The ammonia produced during urea hydrolysis results in the precipitation of a metallic oxide reaction product within and on the surface of *Ureaplasma* colonies.² In 1979, Shepard and Combs added putrescine dihydrochloride to A-7 differential medium to enhance colony size. This resulted in a significant increase in the mean colony size of *Ureaplasma* isolates and intensified visualization of ammonia diffusion in the agar surrounding the colonies.^{3,4}

PRINCIPLE

Casein and soy peptones supply nitrogenous substances and amino acids necessary for bacterial growth. Sodium chloride is a source of essential electrolytes and maintains osmotic equilibrium. Dipotassium phosphate provides a buffering action. Dextrose is added as an energy source. L-cysteine hydrochloride is a reducing agent which minimizes the formation of inhibitory substances. Normal horse serum is added as a source of sterols. Yeast extract supplies B-complex vitamins and serves as a growth enhancer. Putrescine is a polyamine, which enhances growth and intensifies ammonia diffusion in the agar. Urea is required for the growth of ureaplasmas. GHL triptide is a liver cell growth factor and calcium chloride is a precipitating agent. Cefoperazone is a broad-spectrum antibiotic which inhibits many gram-negative and gram-positive organisms; amphotericin B inhibits yeast and molds.

REAGENTS (CLASSICAL FORMULA)*

Casein Peptone.....	17.0 g	GHL Tripeptide	0.02 mg
Sodium Chloride.....	5.0 g	Normal Horse Serum.....	200.0 ml
Soy Peptone.....	3.0 g	Urea 10%.....	10.0 ml
Dextrose.....	2.5 g	Yeast Extract	10.0 ml
Dipotassium Phosphate	2.5 g	L-Cysteine Hydrochloride	1.0 ml
Putrescine Dihydrochloride	1.7 g	Agar.....	11.8 g
Calcium Chloride.....	0.5 g	Deminerlized Water	770.0 ml

pH 6.0 +/- 0.2 @ 25°C

The following ingredients are added per liter of A-8 Selective Agar:

Amphotericin B.....	2.5 mg	Cefoperazone	20.0 mg
---------------------	--------	--------------------	---------

The following ingredient is added per liter of A-8 Nonselective Agar:

Penicillin G	1,000,000 U
--------------------	-------------

*Adjusted as required to meet performance standards.

PROCEDURE

1. Prepare a 1:10 dilution of the specimen in a suitable transport medium, such as 10B Broth (REF 20302) or TSB w/ 0.5 Bovine Albumin (REF 065006), following recommended guidelines.
2. Using a sterile pipette, transfer an aliquot from the broth to the A-8 Agar, Selective or Nonselective.
3. Streak the plate for isolation and seal closed to restrict dehydration.
4. Incubate the plate in 5% CO₂ or in an anaerobic atmosphere at 35-37°C.
5. Examine microscopically for growth and typical colonial morphology for up to 7 days. *U. urealyticum* produces golden brown colonies on A-8 Agar. Colonies of *M. hominis* may exhibit a typical "fried egg" appearance.^{4,5}

NOTE: Serial dilutions to 10⁻³ of any specimen will optimize recovery and is recommended to overcome potential inhibitory substances that may be present in the medium or in the specimen.^{3,4}

QUALITY CONTROL

All lot numbers A-8 Agar, Selective and Nonselective, 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

A-8 Nonselective:

Ureaplasma urealyticum ATCC® 27618
Mycoplasma hominis ATCC® 23114
Staphylococcus aureus ATCC® 25923

INCUBATION

CO₂, 72 h @ 33-37°C
CO₂, 72 h @ 33-37°C
Ambient, 18-24 h @ 33-37°C

RESULTS

Good growth, golden brown colonies
Good growth
Inhibition (partial to complete)

A-8 Selective:

Ureaplasma urealyticum ATCC® 27618
Mycoplasma hominis ATCC® 23114
Candida albicans ATCC® 10231
Escherichia coli ATCC® 25922
Staphylococcus aureus ATCC® 25923

CO₂, 72 h @ 33-37°C
CO₂, 72 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

Good growth, golden brown colonies
Good growth
Inhibition (partial to complete)
Inhibition (partial to complete)
Inhibition (partial to complete)

LIMITATIONS

1. Ureaplasmas are extremely susceptible to a rapid, steep death phase in culture due to urea depletion and elevated pH and must be monitored and subcultured to appropriate media on a regular basis to obtain maximum viability.⁵

BIBLIOGRAPHY

1. Shepard, M.C. and C.D. Lunceford. 1976. J. Clin. Microbiol. 3:613-625.
2. Shepard, M.C. and R.C. Combs. 1979. J. Clin. Microbiol. 10:931-933.
3. Versalovic, J., K.C. Carroll, G. Funke, J.H. Jorgensen, M.L. Landry, and D.W. Warnock. 2011. Manual of Clinical Microbiology. 10th ed. ASM Press, Washington, D.C.
4. Winn, W.C., S.D. Allen, W.M. Janda, E.W. Koneman, G.W. Procop, P.C. Schreckenberger, and G.L. Woods. 2006. Koneman's Color Atlas and Textbook of Diagnostic Microbiology. 6th ed. Lippincott Williams & Wilkins, Philadelphia, PA.
5. 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 20204, Revised July 10, 2012

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