

# MIDDLEBROOK 7H9 BROTH w/ GLYCEROL

## INTENDED USE

Remel Middlebrook 7H9 Broth w/ Glycerol is a liquid medium recommended for use in qualitative procedures for the isolation and cultivation of *Mycobacterium* species.

## SUMMARY AND EXPLANATION

Middlebrook 7H9 Broth was first described by Middlebrook and Cohn in 1958 for the cultivation of mycobacteria.<sup>1</sup> The supplemental nutrients, glycerol, oleic acid, albumin, and dextrose were found to enhance the growth of most mycobacteria.<sup>2,3</sup> This medium is used as a basal medium to prepare inocula for biochemical identification tests and antibiotic susceptibility testing procedures.<sup>4,5</sup>

## PRINCIPLE

Middlebrook 7H9 Broth w/ Glycerol contains inorganic compounds to supply potassium, sulfur, magnesium, and phosphorus which are necessary for the growth of mycobacteria. Inorganic copper, iron, zinc, and calcium are growth stimulators. Sodium citrate, when converted to citric acid, holds certain inorganic cations in solution. Ammonium sulfate is a nitrogen source and dextrose and glycerol are carbon sources. Albumin binds free fatty acids which may be toxic to *Mycobacterium* species. Catalase destroys toxic peroxides which may be present in the medium and catalyzes the reaction of iron with molecular oxygen. Biotin plays an important role in carboxylation and decarboxylation reactions, while pyridoxine is a precursor of enzymatic activity.

## REAGENTS (CLASSICAL FORMULAE)\*

Disodium Phosphate.....	2.5 g	Copper Sulfate.....	1.0 mg
Monopotassium Phosphate.....	1.0 g	Pyridoxine.....	1.0 mg
Ammonium Sulfate.....	0.5 g	Zinc Sulfate.....	1.0 mg
Monosodium Glutamate.....	0.5 g	Biotin.....	0.5 mg
Sodium Citrate.....	0.1 g	Calcium Chloride.....	0.5 mg
Magnesium Sulfate.....	0.05 g	Glycerol.....	2.0 ml
Ferric Ammonium Citrate.....	0.04 g	•ADC Enrichment.....	100.0 ml
		Deminerlized Water.....	900.0 ml

pH 6.6 ± 0.2 @ 25°C

### •ADC Enrichment:

Bovine Albumin (Fraction V).....	50.0 g	Sodium Chloride.....	8.5 g
Dextrose.....	20.0 g	Catalase.....	30.0 mg
		Deminerlized Water.....	1000.0 ml

\*Adjusted as required to meet performance standards.

## PROCEDURE

1. Follow established laboratory procedures for sample preparation, media inoculation, and incubation. Consult appropriate references for further instructions.<sup>3-5</sup>

## QUALITY CONTROL

All lot numbers of Middlebrook 7H9 Broth w/ Glycerol 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

*Mycobacterium fortuitum* ATCC® 6841  
*Mycobacterium kansasii* ATCC® 12478  
*Mycobacterium intracellulare* ATCC® 13950  
*Mycobacterium scrofulaceum* ATCC® 19981  
*Mycobacterium tuberculosis* ATCC® 25177

### INCUBATION

CO<sub>2</sub>, up to 14 days @ 33-37°C  
CO<sub>2</sub>, up to 14 days @ 33-37°C  
CO<sub>2</sub>, up to 14 days @ 33-37°C  
CO<sub>2</sub>, up to 14 days @ 33-37°C  
CO<sub>2</sub>, up to 14 days @ 33-37°C

### RESULTS

Growth  
Growth  
Growth  
Growth  
Growth

## BIBLIOGRAPHY

1. Middlebrook, G. and M.L. Cohn. 1958. Am. J. Public Health. 48:844.
2. Vestal, A.L. 1981. Procedures for the Isolation and Identification of Mycobacteria. U.S. Dept. of H.H.S. and CDC, Atlanta, GA.
3. Kent, P.T. and G.P. Kubica. 1985. Public Health Mycobacteriology, A Guide for the Level III Laboratory. U.S. Dept. of H.H.S. and CDC, Atlanta, GA.
4. Howard, B.J., J.F. Keiser, A.S. Weissfeld, T.F. Smith, R.C. Tilton, and J. Comerford. 1994. Clinical and Pathogenic Microbiology. 2<sup>nd</sup> ed. Mosby, St. Louis, MO.
5. Clinical and Laboratory Standards Institute (CLSI). 2003. Susceptibility Testing of Mycobacteria, Nocardiae, and Other Aerobic Actinomycetes; Approved Standard, M24-A. 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.

IFU 61344, Revised February 3, 2011

Printed in U.S.A.

**remel**

12076 Santa Fe Drive, Lenexa, KS 66215, USA  
General Information: (800) 255-6730 Website: [www.remel.com](http://www.remel.com) Email: [remel@remel.com](mailto:remel@remel.com)  
Local/International Phone: (913) 888-0939 International Fax: (913) 895-4128