

# LOWENSTEIN-JENSEN (L-J) BASE

## INTENDED USE

Remel Lowenstein-Jensen (L-J) Base is a solid medium recommended for use in qualitative procedures for the isolation and cultivation of *Mycobacterium* species.

## SUMMARY AND EXPLANATION

In 1931, Lowenstein formulated a medium for the cultivation of mycobacteria using congo-red and malachite-green dyes to partially inhibit a number of bacterial species.<sup>1</sup> In 1932, Jensen modified the medium by altering the citrate and phosphate contents, eliminating congo-red dye, and increasing the concentration of malachite-green dye.<sup>2</sup> In comparison studies of media for recovery of pathogenic *Mycobacterium* spp., L-J Medium was found to be superior to other media tested, including ATS (American Trudeau Society) Medium.<sup>3,4</sup> In 1960, Boisvert successfully used L-J Medium for niacin testing to identify *M. tuberculosis*.<sup>5</sup>

## PRINCIPLE

L-J Medium is a glycerated egg-potato medium. The glycerol and egg mixture provide fatty acids and protein essential for the metabolism of mycobacterium. Malachite-green dye is added to inhibit nonmycobacterial organisms and as a pH indicator.<sup>6</sup>

## REAGENTS (CLASSICAL FORMULA)\*

Potato Flour.....	30.0 g	Magnesium Citrate .....	0.6 g
Asparagine.....	3.6 g	Malachite Green .....	0.4 g
Monopotassium Phosphate.....	2.4 g	Magnesium Sulfate.....	0.24 g
		Deminerlized Water .....	600.0 ml

\*Adjusted as required to meet performance standards.

## PRECAUTIONS

This product is for Laboratory Use only. It is not intended for use in the diagnosis of disease.

## PREPARATION OF DEHYDRATED CULTURE MEDIUM

1. Suspend 37.2 g of medium in 600 ml of demineralized water, to which 12 ml of glycerol has been added.
2. Heat to boiling with agitation to completely dissolve.
3. Sterilize by autoclaving at 121°C for 15 minutes or following established laboratory procedures.
4. Cool to 50°C and add 1000 ml of homogenized whole egg and optional ingredients if desired. Mix until uniform.
5. Dispense into appropriate containers.
6. Inspissate or autoclave at 85°C for 45 minutes.

## PROCEDURE

1. Consult current editions of appropriate references for the recommended procedure for sample preparation, inoculation, incubation and testing.

## QUALITY CONTROL

Each lot number of L-J Medium has been manufactured, packaged, and processed in accordance with current Good Manufacturing Practice regulations. All lot numbers 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, sample results should not be reported.

### CONTROL

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

### INCUBATION

CO<sub>2</sub>, up to 21 days @ 35-37°C  
CO<sub>2</sub>, up to 21 days @ 35-37°C  
CO<sub>2</sub>, up to 21 days @ 35-37°C  
CO<sub>2</sub>, up to 21 days @ 35-37°C  
CO<sub>2</sub>, up to 21 days @ 35-37°C

### RESULTS

Growth  
Growth  
Growth  
Growth  
Growth

## LIMITATIONS

1. Media containing malachite green, such as L-J Medium, is photosensitive and should not be exposed to light during storage.<sup>6</sup>
2. Mycobacteria require moisture for growth. L-J tubes should be kept tightly sealed prior to inoculation to avoid dehydration of media.<sup>6</sup>
3. Lowenstein-Jensen Medium may display some variation in the light green color throughout the tube. This does not interfere with the growth of mycobacteria; however, color changes showing bright yellow or dark blue zones may indicate contamination.<sup>6</sup>
4. *Mycobacterium bovis* is inhibited by glycerol and will not grow on L-J Medium.<sup>6</sup>

## BIBLIOGRAPHY

1. Lowenstein, E. 1931. Zentralbl. Bakteriol. Parasitenkd. Infektionskr. Hyg. Abt. 1 Orig. 120:127-129.
2. Jensen, K.A. 1932. Zentralbl. Bakteriol. Parasitenkd. Infektionskr. Hyg. Abt. 1 Orig. 125:222-239.
3. Cummings, M.M. 1951. J. Clin. Pathol. 21:684-690.
4. Hardy, A.V., F.P. Dunbar, M.B. Jeffries, J.O. Bond, and A.G. Lewis. 1958. Am. J. Public Health. 48:754-757.
5. Boisvert, H. 1960. Ann. Inst. Pasteur. 99:600-607.
6. 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, sample collection, storage and transportation, materials required, quality control, and limitations.

ATCC® is a registered trademark of American Type Culture Collection.  
IFU 453751, Revised August 23, 2010

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