

---

# LAURYL TRYPTOSE BROTH (LST BROTH)

---

## INTENDED USE

Remel Lauryl Tryptose Broth (LST Broth) (Lauryl Sulfate Tryptose) is a liquid medium recommended for use in qualitative procedures for the detection of coliforms in water, wastewater, and foods.

## SUMMARY AND EXPLANATION

Lauryl Tryptose Broth was developed by Mallman and Darby for the detection of coliform organisms in water, wastewater, and foods.<sup>1,2</sup> They evaluated a variety of wetting agents and determined sodium lauryl sulfate added to selective media produced the best results for the detection of the coliform group. The production of gas in Lauryl Tryptose Broth serves as a presumptive test for the presence of coliforms in water and foods. Lauryl Tryptose Broth is recommended by the American Public Health Association, AOAC International (AOAC), and the U.S. Food and Drug Administration for detection of coliforms in water, wastewater, and foods.<sup>3-6</sup>

## PRINCIPLE

Casein peptone supplies nitrogen, carbon, and sulfur required for bacterial growth. Lactose is the carbohydrate source fermented by coliforms. Dipotassium phosphate and monopotassium phosphate serve as buffering agents. Sodium chloride supplies essential electrolytes and maintains osmotic equilibrium. Sodium lauryl sulfate is a selective agent that inhibits many organisms other than coliforms. A positive fermentation reaction is indicated by turbidity in the broth and gas bubbles in the inverted Durham tube.

## REAGENTS (CLASSICAL FORMULA)\*

Casein Peptone.....	20.0 g	Dipotassium Phosphate .....	2.75 g
Lactose.....	5.0 g	Monopotassium Phosphate.....	2.75 g
Sodium Chloride.....	5.0 g	Sodium Lauryl Sulfate .....	0.1 g
		Demineralized Water .....	1000.0 ml

pH 6.8 ± 0.2 @ 25°C

\*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 or other conditions.

## PREPARATION OF DEHYDRATED CULTURE MEDIUM

1. Suspend 35.6 g of medium in 1000 ml of demineralized water.
2. Warm slightly to dissolve completely.
3. Dispense as required into test tubes or bottles containing inverted Durham tubes.
4. Sterilize at 121°C for 15 minutes or following established laboratory procedures.

## PROCEDURE

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

## INTERPRETATION OF RESULTS

Positive Test - Turbidity with gas production\* within 48 hours

Negative Test - Turbidity with no gas production within 48 hours

\*Gas production is indicated by the appearance of bubbles in the Durham tube.

## QUALITY CONTROL

Each lot number of Lauryl Tryptose Broth 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.

### CONTROL

*Escherichia coli* ATCC® 25922

*Salmonella enterica* serovar Typhimurium ATCC® 14028

*Staphylococcus aureus* ATCC® 25923

### INCUBATION

Ambient, 18-24 h @ 33-37°C

Ambient, 18-24 h @ 33-37°C

Ambient, 18-24 h @ 33-37°C

### RESULTS

Growth w/ gas production

Growth w/o gas production

Marked to complete inhibition

## LIMITATIONS

1. Media may be cloudy or form a precipitate when refrigerated, but clears at room temperature.<sup>7</sup>
2. A bubble in the Durham tube with no turbidity in the broth does not indicate a positive test.

## BIBLIOGRAPHY

1. Darby, C.W. and W.L. Mallman. 1939. J. Am. Water Works Assoc. 31:689-706.
2. Mallman, W.L. and C.W. Darby. 1941. Am. J. Public Health. 31:127-134.
3. Clesceri, L.S., A.E. Greenberg, and A.D. Eaton. 1998. Standard Methods for the Examination of Water and Wastewater. 20<sup>th</sup> ed. APHA, Washington, D.C.
4. Downes, F.P. and K. Ito. 2001. Compendium of Methods for the Microbiological Examination of Foods. 4<sup>th</sup> ed. APHA, Washington, D.C.
5. Food and Drug Administration. 2001. Bacteriological Analytical Manual Online. Chapter 4, Revised September 2002. Authors: P. Feng, S.D. Weagant, and M.A. Grant. AOAC International, Gaithersburg, MD.
6. Horwitz, W. 2010. Official Methods of Analysis Online. 18<sup>th</sup> ed. AOAC International, Gaithersburg, MD.
7. MacFaddin, J.F. 1985. Media for Isolation-Cultivation-Identification-Maintenance of Medical Bacteria. Vol. 1. Williams and 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 9446, Revised April 12, 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