PEPTONE WATER BASE

INTENDED USE

Remel Peptone Water Base is a medium recommended for use in qualitative procedures for the cultivation of nonfastidious organisms and for carbohydrate fermentation studies.

SUMMARY AND EXPLANATION

This medium is based on the formulation of Shread, Donovan, and Lee.¹ It is a minimal medium for all-purpose growth and nonselective enrichment. It may be used for isolation of *Enterobacteriaceae* from foods.²

PRINCIPLE

Peptone provides a source of necessary nutrients. The peptone in this medium does not contain any fermentable carbohydrates, making this medium suitable for carbohydrate fermentation tests. Sodium chloride is a source of essential electrolytes and maintains osmotic equilibrium.

REAGENTS (CLASSICAL FORMULA)*

Meat Peptone......10.0 g

Sodium Chloride......5.0 g

pH 7.2 ± 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 15 g of medium in 1000ml of demineralized water.
- 2. Mix well and dispense into appropriate containers.
- 3. Sterilize at 121°C for 15 minutes or following established laboratory procedures.
- 4. For fermentation studies, dispense into tubes containing inverted fermentation vials (Durham tubes). When adding carbohydrate solutions, the volume of water used to rehydrate the medium should be reduced by an equivalent amount.

PROCEDURE

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

QUALITY CONTROL

Each lot number of Peptone Water Base 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.

INCUBATION

Ambient, 18-24 h @ 33-37°C

CONTROL

Escherichia coli ATCC[®] 25922 Salmonella enterica serovar Enteritidis ATCC[®] 13076 Salmonella enterica serovar Typhi ATCC[®] 19430 Salmonella enterica serovar Typhimurium ATCC[®] 14028

BIBLIOGRAPHY

- 1. Shread, P., T.J. Donovan, and J.V. Lee. 1981. Soc. Gen. Microbiol. Q. 8:184.
- 2. 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.

 $ATCC^{\otimes}$ is a registered trademark of American Type Culture Collection. IFU 454242, Revised August 23, 2010

Printed in U.S.A.

RESULTS

Growth

Growth

Growth

Growth



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