
SALINE (0.85%)

INTENDED USE

Remel Saline (0.85%) is a liquid medium recommended for use in qualitative procedures that require the use of an isotonic diluent.

SUMMARY AND EXPLANATION

Normal Saline (0.85 %) has been recommended by various manufacturers and authors for preparing test suspensions of organisms to help maintain bacterial cell integrity and viability.¹ Normal saline is recommended by Clinical and Laboratory Standards Institute (CLSI) for preparation of bacterial suspensions used for susceptibility testing.^{2,3}

PRINCIPLE

This medium lacks properties that may interfere with biochemical reactions and/or antibiotic susceptibility tests. The sodium chloride concentration in Saline (0.85%) provides osmotic protection for microbial cells.

REAGENTS (CLASSICAL FORMULA)*

Sodium Chloride..... 8.5 g Demineralized Water 1000.0 ml

*Adjusted as required to meet performance standards.

PROCEDURE

Consult appropriate references for the recommended procedures or test methods requiring 0.85% Saline.²⁻⁵

QUALITY CONTROL

All lot numbers of Saline (0.85%) have been tested for microbial content and aesthetic value only and have been found to be acceptable. To ensure efficacy of the product, functional testing should be performed by the user in accordance with established laboratory quality control procedures. If aberrant quality control results are noted, patient results should not be reported.

BIBLIOGRAPHY

1. Bauer, A.W., W.M. Kirby, J.C. Sherris, and M. Turck. 1966. Am. J. Clin. Pathol. 45:493-496.
2. Clinical and Laboratory Standards Institute (CLSI). 2009. Performance Standards for Antimicrobial Disk Susceptibility Tests; Approved Standard, 10th ed. M2-A10. CLSI, Wayne, PA.
3. Clinical and Laboratory Standards Institute (CLSI). 2006. Methods for Dilution Antimicrobial Susceptibility Tests for Bacteria that Grow Aerobically; Approved Standard, 7th ed. M7-A7. CLSI, Wayne, PA.
4. Murray, P.R., E.J. Baron, J.H. Jorgensen, M.L. Landry, and M.A. Pfaller. 2007. Manual of Clinical Microbiology. 9th ed. ASM Press, Washington, D.C.
5. Forbes, B.A., D.F. Sahm, and A.S. Weissfeld. 2007. Bailey and Scott's Diagnostic Microbiology. 12th ed. Mosby Elsevier, St. Louis, MO.

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 64430, Revised February 1, 2010

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