
MANNITOL SALT AGAR

INTENDED USE

Remel Mannitol Salt Agar is a solid medium recommended for use in qualitative procedures for selective and differential isolation of staphylococci.

SUMMARY AND EXPLANATION

In 1942, Koch reported the use of 7.5% sodium chloride as a selective agent for the isolation of staphylococci.¹ Chapman confirmed the results of Koch and suggested the addition of 7.5% sodium chloride to phenol red mannitol agar.² Most strains of coagulase-positive staphylococci grow on Mannitol Salt Agar, producing yellow zones as a result of mannitol fermentation. Coagulase-negative strains of staphylococci produce small colonies with red-colored zones in the surrounding medium.

PRINCIPLE

Casein and meat peptones supply nitrogen, amino acids, and peptides necessary for bacterial growth. Sodium chloride in a concentration of 7.5% is a selective agent which inhibits many bacteria other than staphylococci. Phenol red is a pH indicator which causes a color change in the medium from red-orange to yellow when acid is produced. Staphylococci colonies that ferment mannitol will be surrounded by a yellow zone, while those that do not ferment mannitol will have a red zone.

REAGENTS (CLASSICAL FORMULA)*

Sodium Chloride.....	75.0 g	Beef Extract.....	1.0 g
D-Mannitol.....	10.0 g	Phenol Red.....	25.0 mg
Casein Peptone.....	5.0 g	Agar.....	15.0 g
Meat Peptone.....	5.0 g	Demineralized Water.....	1000.0 ml

pH 7.4 ± 0.2 @ 25°C

*Adjusted as required to meet performance standards.

PROCEDURE

1. Inoculate and streak the specimen as soon as possible after it is received in the laboratory.
2. If material is being cultured directly from a swab, roll the swab over a small area of the agar surface and streak for isolation.
3. Incubate the plate in ambient air at 33-37°C for 24-48 hours.
4. Examine plate for typical colony morphology and color change.

QUALITY CONTROL

All lot numbers of Mannitol Salt Agar have been tested using the following quality control organisms and have been found to be acceptable. This quality control testing meets or exceeds CLSI standards.³ 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

**Staphylococcus aureus* ATCC® 25923

**Proteus mirabilis* ATCC® 12453

**Staphylococcus epidermidis* ATCC® 12228

*CLSI recommended organism

INCUBATION

Ambient, 18-24 h @ 33-37°C

Ambient, 18-24 h @ 33-37°C

Ambient, 18-24 h @ 33-37°C

RESULTS

Growth, colonies with yellow zones

Inhibited swarming

Inhibition (partial), colonies with red zones

BIBLIOGRAPHY

1. Koch, F.D. 1942. Zentr. Bakt. Labt. Orig. 149:122.
2. Chapman, G.H. 1945. J. Bacteriol. 50:201.
3. Clinical and Laboratory Standards Institute (CLSI). 2004. Quality Control for Commercially Prepared Microbiological Culture Media; Approved Standard, 3rd ed. M22-A3. 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 1580, Revised March 10, 2011

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