

MACCONKEY AGAR w/ SORBITOL

INTENDED USE

Remel MacConkey Agar w/ Sorbitol is a solid medium recommended for use in qualitative procedures as a differential and selective medium for the detection of sorbitol-negative *Escherichia coli*.

SUMMARY AND EXPLANATION

Escherichia coli serotype O157:H7 (VTEC) produces a verotoxin which causes hemorrhagic colitis outbreaks, with symptoms ranging from mild diarrhea to idiopathic hemolytic uremic syndrome.^{1,2} On routine isolation media, colonies of VTEC are indistinguishable from those of non-VTEC *E. coli*. In 1986, March and Ratnam introduced MacConkey Agar w/ Sorbitol for isolation of verotoxin-producing *E. coli* O157:H7 from clinical specimens. Unlike most serotypes of *E. coli*, VTEC does not ferment sorbitol. Haldane et al. implemented the use of MacConkey Agar w/ Sorbitol as a primary isolation medium to distinguish verotoxin-producing strains of *E. coli* which are sorbitol-negative from non-VTEC *E. coli* which is sorbitol-positive.³

PRINCIPLE

Gelatin, casein, and meat peptones provide nitrogen, amino acids, and peptides essential for bacterial growth. Sodium chloride is a source of essential electrolytes and maintains osmotic equilibrium. Bile salts and crystal violet are selective agents which inhibit gram-positive organisms. Lactose is omitted from this formulation and is substituted with D-sorbitol. Verotoxin-producing strains of *E. coli* ferment lactose but not D-sorbitol and form colorless colonies on MacConkey Agar w/ Sorbitol.

REAGENTS (CLASSICAL FORMULA)*

Gelatin Peptone	17.0 g	Meat Peptone	1.5 g
D-Sorbitol	10.0 g	Neutral Red	30.0 mg
Sodium Chloride.....	5.0 g	Crystal Violet	1.0 mg
Bile Salts	1.5 g	Agar	13.5 g
Casein Peptone	1.5 g	Demineralized Water	1000.0 ml

pH 7.1 ± 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. Selective and nonselective media should be inoculated to increase the chance of recovering enteric pathogens when the number of gram-negative organisms is low and to provide isolation of other organisms present in the specimen.
4. Incubate plates in ambient air at 33-37°C for 18-24 hours.
5. Observe plates for typical colonial morphology.

Pour Tube: Melt the pour tube in a boiling water bath and cool to 45-50°C. Mix and dispense into a sterile petri dish and proceed with the instructions above.

QUALITY CONTROL

All lot numbers of MacConkey Agar w/ Sorbitol 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, patient results should not be reported.

CONTROL

Escherichia coli ATCC® 43888
Escherichia coli ATCC® 25922
Proteus mirabilis ATCC® 12453
Salmonella enterica serovar Typhimurium ATCC® 14028
Enterococcus faecalis ATCC® 29212
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
Ambient, 18-24 h @ 33-37°C
Ambient, 18-24 h @ 33-37°C
Ambient, 18-24 h @ 33-37°C

RESULTS

Growth, colorless colonies
Growth, pink colonies
Growth, colorless colonies, swarming inhibited
Growth, pink colonies
Inhibition (partial to complete)
Inhibition (partial to complete)

LIMITATIONS

1. Sorbitol-negative isolates should be confirmed as *E. coli* by current accepted methodologies. Confirmed *E. coli* isolates should be serotyped using an *E. coli* O157 and H7 antiserum. These strains are reportable to State Public Health Departments.⁴

BIBLIOGRAPHY

1. March, S.B. and S. Ratnam. 1986. J. Clin. Microbiol. 23:869-872.
2. Krishnan, C., A. Fitzgerald, S.A. Dakin, and R.J. Behme. 1987. J. Clin. Microbiol. 25:1043-1047.
3. Haldane, D.J. M., M.A.S. Dam, and J.D. Anderson. 1986. J. Clin. Microbiol. Oct: 652-653.
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.

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 1556, Revised March 23, 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