

BLOOD AGAR (TSA w/ 5% Sheep Blood)

INTENDED USE

Remel Blood Agar (TSA w/ 5% Sheep Blood) is a solid medium recommended for use in qualitative procedures for the cultivation of a wide variety of microorganisms and visualization of hemolytic reactions.

SUMMARY AND EXPLANATION

Tryptic Soy Agar is based on the soybean casein digest agar formula described in the U.S. Pharmacopeia.¹ Blood is incorporated in the medium to facilitate growth of more fastidious bacteria and to observe hemolytic reactions. The absence of carbohydrates and reducing sugars permits the demonstration of hemolysis which is an important differentiating characteristic for bacteria, especially *Streptococcus* species.

PRINCIPLE

Blood Agar (TSA w/ 5% Sheep Blood) contains casein and soy peptones which provide nitrogen, amino acids, and peptides necessary for bacterial growth. Sodium chloride supplies essential electrolytes and maintains osmotic equilibrium. Sheep blood enriches the medium by providing essential growth factors and allows hemolytic reactions to be demonstrated. Group B streptococci produce a protein-like compound called the CAMP factor which acts synergistically with beta toxin, produced by some strains of *Staphylococcus aureus*.² This reaction occurs when a streak of β -lysin-producing *S. aureus* is inoculated perpendicular to a streak of group B *Streptococcus* resulting in an area of complete lysis in the shape of an arrowhead or crescent. Blood Agar (TSA w/ 5% Sheep Blood) is suitable for performing the CAMP test.

REAGENTS (CLASSICAL FORMULA)*

Casein Peptone	15.0 g	Sheep Blood	5 %
Sodium Chloride	5.0 g	Agar	15.0 g
Soy Peptone	5.0 g	Demineralized Water	1000.0 ml

pH 7.3 $\pm$ 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 plate aerobically, or in 5-10% CO₂, for 18-24 hours at 33-37°C.
4. Examine plate for typical colony morphology and hemolytic reactions.

QUALITY CONTROL

All lot numbers of Blood Agar (TSA w/ 5% Sheep Blood) 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

Enterococcus faecalis ATCC® 29212
**Escherichia coli* ATCC® 25922
Listeria monocytogenes ATCC® 7646
Moraxella catarrhalis ATCC® 25238
**Staphylococcus aureus* ATCC® 25923
Streptococcus agalactiae ATCC® 12386
**Streptococcus pneumoniae* ATCC® 6305
**Streptococcus pyogenes* ATCC® 19615

INCUBATION

Ambient, 18-24 h @ 33-37°C
Ambient, 18-24 h @ 33-37°C
Ambient, 18-24 h @ 33-37°C
CO₂, 18-24 h @ 33-37°C
Ambient, 18-24 h @ 33-37°C
Ambient, 18-24 h @ 33-37°C
CO₂, 18-24 h @ 33-37°C
Ambient, 18-24 h @ 33-37°C

RESULTS

Growth, gamma hemolysis
Growth
Growth, beta hemolysis
Growth
Growth
Growth, beta hemolysis
Growth, alpha hemolysis
Growth, beta hemolysis

CAMP Test**

**Streptococcus agalactiae* ATCC® 12386
**Streptococcus pyogenes* ATCC® 19615

Ambient, 18-24 h @ 33-37°C
Ambient, 18-24 h @ 33-37°C

Arrowhead-shaped zone of increased beta hemolysis
No arrowhead formation

*CLSI recommended organism

**CAMP Test performed with *S. aureus* ATCC® 33862

BIBLIOGRAPHY

1. The United States Pharmacopeia. 2006. 24th ed., Sup. 2., 29th rev. United States Pharmacopeial Convention, Rockville, MD.
2. Christie, R., N.E. Atkins, and E. Munch-Peterson. 1944. Aust. J. Exp. Biol. 22:197-200.
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 1200, Revised September 18.2014

Printed in U.S.A.

remel

12076 Santa Fe Trail 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