

# VOGEL-JOHNSON AGAR

## INTENDED USE

Remel Vogel-Johnson Agar is a solid medium recommended for use in qualitative procedures for the early detection of coagulase-positive, mannitol-fermenting strains of *Staphylococcus aureus* from contaminated foods and clinical specimens.

## SUMMARY AND EXPLANATION

In 1955, Zebovitz et al. developed Tellurite-Glycine Agar as a selective medium for the detection of coagulase-positive staphylococci.<sup>1</sup> Vogel and Johnson modified Tellurite-Glycine Agar by increasing the mannitol content and adding phenol red as a pH indicator.<sup>2</sup> Vogel-Johnson Agar is recommended by the Association of Official Analytical Chemists (AOAC) for food bacteriology.<sup>3,4</sup>

## PRINCIPLE

Casein peptone supplies nitrogen, amino acids, and peptides necessary for bacterial growth. Yeast extract provides B-complex vitamins. The fermentation of mannitol by coagulase-positive staphylococci is indicated by development of yellow zones surrounding the colonies. This is caused by a pH change that results in the phenol red indicator turning from red to yellow. *S. aureus* reduces potassium tellurite to metallic tellurium resulting in black colonies. Potassium tellurite, lithium chloride, and glycine are selective agents which inhibit non-staphylococcal organisms, including many gram-negative and gram-positive bacteria.

## REAGENTS (CLASSICAL FORMULA)\*

Casein Peptone.....	10.0 g	Yeast Extract .....	5.0 g
Mannitol.....	10.0 g	Phenol Red.....	25.0 mg
Glycine .....	10.0 g	Potassium Tellurite 1%.....	20.0 ml
Dipotassium Phosphate .....	5.0 g	Agar.....	15.0 g
Lithium Chloride .....	5.0 g	Deminerlized Water.....	1000.0 ml

pH 7.2 ± 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. The selective qualities of this medium are such that the inoculum may be applied heavily.
3. Incubate plates in ambient air at 33-37°C or at 30-35°C for 24-48 hours.
4. After 24 hours incubation, examine the plate for black colonies which may or may not be surrounded by a yellow zone. After 48 hours incubation, other organisms may exhibit slight growth.
5. Perform a Gram stain and coagulase test on all black colonies. Additional testing may be required for definitive identification of *S. aureus*. Consult appropriate references for further instructions.<sup>3</sup>

## QUALITY CONTROL

All lot numbers of Vogel-Johnson Agar 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

*Staphylococcus aureus* ATCC® 6538  
*Staphylococcus aureus* ATCC® 25923  
*Escherichia coli* ATCC® 25922  
*Staphylococcus epidermidis* ATCC® 12228

## INCUBATION

Ambient, up to 48 h @ 33-37°C  
Ambient, up to 48 h @ 33-37°C  
Ambient, 18-24 h @ 33-37°C  
Ambient, 18-24 h @ 33-37°C

## RESULTS

Black colonies with yellow zones  
Black colonies with yellow zones  
Inhibition (partial to complete)  
Inhibition (partial to complete)

## LIMITATIONS

1. If tellurite is reduced but mannitol is not fermented, the medium surrounding the colonies may be a deeper red color due to the utilization of proteins with resultant alkalinity.<sup>4</sup>
2. Direct inoculation of Vogel-Johnson Agar with a light inoculum is not recommended. Microscopic examination of plates may be useful to assist in early visualization of characteristic halo and color of colony.<sup>4</sup>

## BIBLIOGRAPHY

1. Zebovitz, E., J.B. Evans, and, C.F. Niven. 1955. J. Bacteriol. 70:686-690.
2. Vogel, R.A. and M.J. Johnson. 1960. Public Health Lab. 18:131.
3. Food and Drug Administration. 2000. Bacteriological Analytical Manual Online. Media Index, Updated 07/17/2009. AOAC International, Gaithersburg, MD.
4. 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, specimen collection, storage and transportation, materials required, quality control, and limitations.

ATCC® is a registered trademark of American Type Culture Collection.

IFU 1970, Revised October 6, 2010

Printed in U.S.A.

**remel**

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