

---

# CETRIMIDE SELECTIVE AGAR

---

## INTENDED USE

Remel Cetrimide Selective Agar is a solid medium recommended for use in qualitative procedures for selective isolation and presumptive identification of *Pseudomonas aeruginosa*.

## SUMMARY AND EXPLANATION

King et al. devised a medium called Tech Agar which was used to enhance pyocyanin production by *Pseudomonas*.<sup>1</sup> In 1951 and 1955, Lowbury described the use of cetrimide in a selective medium for *P. aeruginosa*.<sup>2,3</sup> The formula of Cetrimide Selective Agar is identical to that of Tech Agar with one exception, it is modified by addition of cetrimide. Cetrimide Selective Agar is formulated in conformance with harmonized *United States Pharmacopeia* (USP)/*European Pharmacopeia* (EP) guidelines.<sup>4</sup>

## PRINCIPLE

*P. aeruginosa* is characterized by production of pyocyanin, a blue-green, water-soluble, nonfluorescent phenazine pigment. Potassium sulfate and magnesium chloride stimulate the production of pyocyanin and fluorescein. An ultraviolet light is used to visualize fluorescein production. Cetrimide inhibits bacteria other than *P. aeruginosa*, by causing nitrogen and phosphorus to be released from bacterial cells.

## REAGENTS (CLASSICAL FORMULA)\*

Gelatin Peptone .....	20.0 g	Cetrimide .....	0.3 g
Potassium Sulfate .....	10.0 g	Glycerol .....	10.0 ml
Magnesium Chloride .....	1.4 g	Agar .....	13.6 g
		Demineralized Water .....	1000.0 ml

pH 7.2 ± 0.2 @ 25°C

\*Adjusted as required to meet performance standards.

## PREPARATION OF DEHYDRATED CULTURE MEDIUM

1. Suspend 45.3 g of medium in 1000 ml of demineralized water.
2. Add 10 ml of glycerol.
3. Heat to boiling with agitation to completely dissolve.
4. Sterilize by autoclaving at 121°C for 15 minutes or according to established laboratory procedures.
5. Dispense into appropriate containers.

## PROCEDURE

1. Consult current editions of appropriate references for the recommended procedure for sample preparation, inoculation, and testing.<sup>4,7</sup>

## INTERPRETATION OF THE TEST

### Pyocyanin Production:

Positive Test - Blue-green pigmentation surrounding growth

Negative Test - No color development

### Fluorescein Production (requires use of an ultraviolet light source):

Positive Test - Yellow-green fluorescence

Negative Test - No fluorescence

## QUALITY CONTROL

Each lot number of Cetrimide Selective Agar has been manufactured, packaged, and processed in accordance with current Good Manufacturing Practice regulations. All lot numbers have been tested using the following quality control organisms and have been found to be acceptable. This quality control testing meets or exceeds harmonized USP/EP guidelines for microbial limits testing. Testing of control organisms should be performed in accordance with established laboratory quality control procedures.

## CONTROL

*Pseudomonas aeruginosa* ATCC® 9027  
*Pseudomonas aeruginosa* ATCC® 27853  
*Escherichia coli* ATCC® 8739  
*Escherichia coli* ATCC® 25922

## INCUBATION

Aerobic, 18-24 h @ 30-35°C  
Aerobic, 18-24 h @ 30-35°C  
Aerobic, 18-24 h @ 30-35°C  
Aerobic, 18-24 h @ 30-35°C

## RESULTS

Growth, blue-green pigmentation, yellow-green fluorescence  
Growth, blue-green pigmentation, yellow-green fluorescence  
Inhibition (partial to complete)  
Inhibition (partial to complete)

## LIMITATIONS

1. Some enteric gram-negative bacilli may exhibit growth on Cetrimide Selective Agar and produce slight yellowing of the medium. This yellow color is easily distinguished from fluorescein, by its lack of fluorescence.<sup>5</sup>
2. Of the pseudomonads, only *P. aeruginosa* is known to excrete pyocyanin as well as produce pyorubin simultaneously with pyocyanin and/or fluorescein. Pyorubin is a pink to red or dark maroon pigment.<sup>5</sup>
3. Occasional strains of *P. aeruginosa* may fail to produce pyocyanin.<sup>5</sup>
4. *P. aeruginosa* may lose its fluorescence under ultraviolet light if the cultures are left at room temperature for a short time. Fluorescence reappears when plates are reincubated.<sup>5</sup>
5. If Cetrimide Agar slants are inoculated, tubes should be incubated with caps loosened.<sup>5</sup>
6. Cetrimide Selective Agar is both selective and differential. As such, low levels of the target organism may not be recoverable when inoculated directly onto this medium. Consult appropriate references to determine methods for recovery optimization.

## BIBLIOGRAPHY

1. King, E.O., M.K. Ward, and D.E. Raney. 1954. J. Lab. Clin. Med. 44:301-307.
2. Lowbury, E.J. 1951. J. Clin. Pathol. 4:66-72.
3. Lowbury, E.J. and A.G. Collins. 1955. J. Clin. Pathol. 8:47-48.
4. The United States Pharmacopeia. United States Pharmacopeial Convention, Rockville, MD.
5. MacFaddin, J.F. 1985. Media for Isolation-Cultivation-Identification-Maintenance of Medical Bacteria. Vol. 1. Williams & Wilkins, Baltimore, MD.
6. Gaby, W.L. and E. Free. 1953. J. Bacteriol. 65:746.
7. Food and Drug Administration. 2001. Bacteriological Analytical Manual Online. AOAC International, Gaithersburg, 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 1292-MLT, Revised August 25, 2009

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