

# PSEUDOMONAS ISOLATION AGAR

## INTENDED USE

Remel Pseudomonas Isolation Agar is a solid medium recommended for use in qualitative procedures for the isolation and identification of *Pseudomonas* species from clinical and nonclinical specimens.

## SUMMARY AND EXPLANATION

Pseudomonas Isolation Agar is a modification of the Medium A formulation of King, Ward, and Raney.<sup>1</sup> Irgasan<sup>®</sup> is added as a potent broad-spectrum antimicrobial. The formulation of this medium was developed to enhance the production of pyocyanin which aids in the presumptive identification of *Pseudomonas aeruginosa*. Pseudomonas Isolation Agar is especially useful in isolating *Pseudomonas* from industrial materials such as cosmetics and lotion.

## PRINCIPLE

*P. aeruginosa* is the only species of bacteria known to produce pyocyanin, a blue-green pigment which diffuses into the medium surrounding the growth. Gelatin and meat peptones provide the nutrients necessary for bacterial growth. The medium has a low phosphorous concentration which facilitates pyocyanin production. Magnesium chloride and potassium sulfate incorporated in the medium also enhance pyocyanin production. Glycerol added to the prepared medium also enhances pyocyanin production and is an energy source. Irgasan<sup>®</sup> is the selective agent which inhibits many gram-positive and gram-negative bacteria other than *Pseudomonas* spp.

## REAGENTS (CLASSICAL FORMULA)\*

Gelatin Peptone .....	10.0 g	Magnesium Chloride .....	1.4 g
Meat Peptone .....	10.0 g	Irgasan <sup>®</sup> .....	25.0 mg
Potassium Sulfate .....	10.0 g	Agar .....	13.6 g
		Demineralized Water .....	1000.0 ml

pH 7.0 ± 0.2 @ 25°C

\*Adjusted as required to meet performance standards.

## PRECAUTIONS

This product is For Laboratory Use only. It is not intended for use in the diagnosis of disease or other conditions.

## PREPARATION OF DEHYDRATED CULTURE MEDIUM

1. Suspend 45 g of medium in 980 ml of demineralized water.
2. Add 20 ml of glycerol.
3. Heat to boiling with agitation to completely dissolve.
4. Sterilize by autoclaving at 121°C for 15 minutes or following established laboratory procedures.
5. Dispense into appropriate containers.

## PROCEDURE

1. Consult current editions of appropriate references for the recommended procedure for sample preparation, inoculation, testing, and interpretation.

## QUALITY CONTROL

Each lot number of Pseudomonas Isolation 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. Testing of control organisms should be performed in accordance with established laboratory quality control procedures. If aberrant quality control results are noted, sample results should not be reported.

### CONTROL

*Pseudomonas aeruginosa* ATCC<sup>®</sup> 27853  
*Pseudomonas fluorescens* ATCC<sup>®</sup> 13525  
*Escherichia coli* ATCC<sup>®</sup> 25922  
*Staphylococcus aureus* ATCC<sup>®</sup> 25923

### 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

Growth, blue-green pigment  
Growth, no pigment  
Inhibition (partial to complete)  
Inhibition (partial to complete)

## LIMITATIONS

1. Occasional strains of *Pseudomonas aeruginosa* fail to produce pyocyanin.<sup>3</sup>

## BIBLIOGRAPHY

1. King, E.O., M.K. Ward, and D.E. Raney. 1954. J. Lab. Clin. Med. 44:301-306.
2. Brown, V.I. and E.J.L. Lowbury. 1965. J. Clin. Pathol. 18:752-756.
3. Gaby, W.L. and E. Free. 1953. J. Bacteriol. 65:746.

Refer to the front of Remel *Technical Manual of Microbiological Media* for **General Information** regarding precautions, product storage and deterioration, sample collection, storage and transportation, materials required, quality control, and limitations.

Irgasan<sup>®</sup> is a registered trademark of Ciba-Geigy for 2,4,4'-Trichloro-2-Hydroxydiphenol-ether.  
ATCC<sup>®</sup> is a registered trademark of American Type Culture Collection.

IFU 454391, Revised October 7, 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) [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