
POTATO DEXTROSE AGAR

INTENDED USE

Remel Potato Dextrose Agar is a solid medium recommended for use in qualitative procedures for the identification, cultivation, and enumeration of yeast and molds.

SUMMARY AND EXPLANATION

Potato Dextrose Agar is recommended by the American Public Health Association (APHA) and AOAC International (AOAC) for plate counts of yeast and molds in the examination of dairy products and foods.¹⁻³ This medium is commonly used in slide-culture preparations of fungi to stimulate sporulation.^{4,5} It is also used to maintain stock cultures of dermatophytes. Rebell and Taplin used Potato Dextrose Agar for differentiation of atypical dermatophytes based on pigment production.⁶

PRINCIPLE

Infusion from potatoes provides a nutritious base for luxuriant development of most fungi. Dextrose is incorporated in the medium as a growth stimulant. Sterile tartaric acid may be added to this medium to lower the pH to 3.5 to inhibit bacterial growth.

REAGENTS (CLASSICAL FORMULAE)*

Dextrose.....	20.0 g	Agar.....	15.0 g
Potato Infusion Solids.....	4.0 g	Demineralized Water.....	1000.0 ml

pH 5.6 ± 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 39 g of medium in 1000 ml of demineralized water.
2. Heat to boiling with agitation to completely dissolve.
3. Sterilize by autoclaving at 121°C for 15 minutes or following established laboratory procedures.
4. For a medium with a pH of 3.5, add 1.8 ml of sterile 10% tartaric acid to 100 ml of cooled, sterilized base.
5. Mix thoroughly and 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 Potato Dextrose 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

Candida albicans ATCC® 10231
Cryptococcus neoformans ATCC® 34877
Trichophyton mentagrophytes ATCC® 9533

INCUBATION

Ambient, up to 72 h @ 25-30°C
Ambient, up to 72 h @ 25-30°C
Ambient, up to 72 h @ 25-30°C

RESULTS

Good growth
Good growth
Good growth

BIBLIOGRAPHY

1. Wehr, H.M. and J.F. Frank. 2004. Standard Methods for the Examination of Dairy Products. 17th ed. APHA, Washington, D.C.
2. Downes, F.P. and K. Ito. 2001. Compendium of Methods for the Microbiological Examination of Foods. 4th ed. APHA, Washington, D.C.
3. Food and Drug Administration. 2001. Bacteriological Analytical Manual Online. Chapter 18, Revised January 2001. Authors: Tournas, V., M.E. Stack, P.B. Mislivec, H.A. Koch, and R. Bandler. AOAC International, Gaithersburg, MD.
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.
5. Forbes, B.A., D.F. Sahm, and A.S. Weissfeld. 2007. Bailey and Scott's Diagnostic Microbiology. 12th ed. Mosby Elsevier, St. Louis, MO.
6. Rebell, G. and D. Taplin. 1970. Dermatophytes, Their Recognition and Identification. University of Miami Press, Coral Gables, FL.

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.

ATCC® is a registered trademark of American Type Culture Collection.
IFU 454311, Revised October 21, 2010

Printed in U.S.A.

remel

12076 Santa Fe Drive, Lenexa, KS 66215, USA

General Information: (800) 255-6730 Local/International Phone: (913) 888-0939 International Fax: (913) 895-4128
Website: www.remel.com www.remel.com Email: remel@remel.com