

---

# YM AGAR BASE

---

## INTENDED USE

Remel YM Agar Base is a medium recommended for use in qualitative procedures for the isolation and cultivation of yeasts and molds.

## SUMMARY AND EXPLANATION

YM Agar is prepared according to the formulation of Wickerham.<sup>1</sup> It is a nonselective growth medium for yeasts and molds. It can also serve as the basal medium for selective agar with the addition of antibiotics or by acidification.

## PRINCIPLE

Yeast and malt extracts and peptone provide necessary nutrients for growth. Dextrose provides a source of energy. Antibiotics may be incorporated or the pH may be reduced by the addition of acid to increase selectivity.

## REAGENTS (CLASSICAL FORMULA)\*

Dextrose.....	10.0 g	Malt Extract.....	3.0 g
Meat Peptone.....	5.0 g	Yeast Extract.....	3.0 g
		Agar.....	20.0 g

pH 6.2 ± 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 41 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. 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 YM Agar Base 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

*Aspergillus brasiliensis* ATCC® 16404  
*Candida albicans* ATCC® 10231  
*Candida glabrata* ATCC® 2001  
*Cryptococcus neoformans* ATCC® 34877  
*Escherichia coli* ATCC® 25922

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

### RESULTS

Growth  
Growth  
Growth  
Growth  
Growth

## BIBLIOGRAPHY

1. Wickerham, L.J. 1951. U.S. Dept. Agric. Tech. Bull. No. 1029.

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 456032, Revised August 11, 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