

DERMATOPHYTE TEST MEDIUM (DTM)

INTENDED USE

Remel Dermatophyte Test Medium (DTM) is a solid medium recommended for use in qualitative procedures for selective isolation of pathogenic fungi (dermatophytes) from cutaneous sources.

SUMMARY AND EXPLANATION

The dermatophytes are fungi that possess keratinolytic properties that enable them to invade skin, nails, and hair.¹ The infections caused by these organisms are commonly referred to as ringworm and are classified by the Latin word *tinea* followed by the area of the body infected.² Dermatophyte Test Medium (DTM) was formulated by Taplin et al. for use in locations where specialized training and microscopic examination is not available.³ A pH indicator and three antimicrobial agents are incorporated into the agar to provide a differential and selective medium for isolation of dermatophytes belonging to the genera *Microsporum*, *Trichophyton*, or *Epidermophyton*.

PRINCIPLE

Soy peptone supplies the nitrogen and carbon compounds necessary for the growth of microorganisms. Dextrose is an energy source. Phenol red is a pH indicator which detects alkaline metabolites produced by dermatophytes, resulting in a red color development of the medium. Cycloheximide, chloramphenicol, and gentamicin are selective agents which inhibit most saprophytic fungi and many gram-positive and gram-negative bacteria, including some *Pseudomonas* spp.

REAGENTS (CLASSICAL FORMULA)*

Dextrose.....	10.0 g	Chloramphenicol	0.1 g
Soy Peptone.....	10.0 g	Gentamicin	0.1 g
Cycloheximide.....	0.5 g	Agar.....	20.0 g
Phenol Red	0.2 g	Deminerlized Water	1000.0 ml

pH 5.5 ± 0.2 @ 25°C

*Adjusted as required to meet performance standards.

PROCEDURE

1. Collect specimen following established procedures. Consult appropriate references for further instructions.^{1,2,4}
2. Allow DTM to equilibrate to room temperature prior to use. The agar surface should be dry before inoculation.
3. Place the specimen centrally on the surface of the medium and press into the agar to ensure firm contact.
4. Incubate in ambient air at 25-30°C for up to 14 days. Allow the cap on the tube to remain loose so that air may be exchanged during incubation.
5. Examine the medium at regular intervals for a red color development. On DTM, dermatophytes elaborate alkaline metabolites which elevate the pH of the medium and change the phenol red indicator from yellow to red.⁵ **Note:** Microscopic examination (e.g., wet-mount, KOH) is required for presumptive identification of an isolate as a dermatophyte.

QUALITY CONTROL

All lot numbers of Dermatophyte Test Medium (DTM) 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

Candida albicans ATCC® 10231
Trichophyton mentagrophytes ATCC® 9533
Aspergillus niger ATCC® 16404
Cryptococcus neoformans ATCC® 34877
Escherichia coli ATCC® 25922
Pseudomonas aeruginosa ATCC® 27853
Staphylococcus aureus ATCC® 25923

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

RESULTS

Good growth
Good growth, red zone
Inhibition (partial to complete)
Inhibition (partial to complete)
Inhibition (partial to complete)
Inhibition (partial to complete)
Inhibition (partial to complete)

LIMITATIONS

1. DTM is primarily used as a screening test for detection of dermatophytes. Additional testing may be required for definitive identification.²
2. Specimens from heavily soil-contaminated sources (e.g., feet and nails) may contain saprophytic fungi which redden the medium. Such cultures may require additional interpretation to distinguish contaminating fungi from dermatophytes.^{3,5}
3. Organisms other than dermatophytes (e.g., saprophytic fungi, yeast, and bacteria) may grow on DTM and produce alkaline metabolites. DTM should not be used as the only method for identification of dermatophytes.^{2,3}

BIBLIOGRAPHY

1. 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.
2. Forbes, B.A., D.F. Sahm, and A.S. Weissfeld. 2007. Bailey and Scott's Diagnostic Microbiology. 12th ed. Mosby Elsevier, St. Louis, MO.
3. Taplin, D., N. Zaias, G. Rebell, and H. Blank. 1969. Arch. Dermatol. 99:203-209.
4. Miller, J.M. 1999. A Guide to Specimen Management in Clinical Microbiology. 2nd ed. ASM Press, Washington, D.C.
5. 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 1365, Revised October 5, 2010

Printed in U.S.A.

remel

12076 Santa Fe Drive, Lenexa, KS 66215, USA

General Information: (800) 255-6730 Website: www.remel.com Email: remel@remel.com

Local/International Phone: (913) 888-0939 International Fax: (913) 895-4128