

---

# McCARTHY AGAR

---

## INTENDED USE

Remel McCarthy Agar is a solid medium recommended for use in qualitative procedures for the selective isolation and presumptive differentiation of *Gardnerella vaginalis*.

## SUMMARY AND EXPLANATION

Gardner and Duker first isolated *G. vaginalis* from women with nonspecific vaginitis in 1955.<sup>1</sup> Dunkelberg et al. developed peptone starch dextrose (PSD) agar which, unlike previously described media, did not contain blood.<sup>2</sup> Smith modified PSD by adding a pH indicator to the medium, allowing detection of *G. vaginalis* by production of acid from starch.<sup>3</sup> In 1977, Mickelsen et al. modified Smith's formulation by addition of 1% cornstarch, which makes the medium opaque.<sup>4</sup> Colonies of *G. vaginalis* are characterized by zones of clearing that appear around the colonies, presumably because of an ability to solubilize large aggregates of cornstarch.

## PRINCIPLE

Casein and meat peptones supply nutrients necessary to support bacterial growth. Phosphate buffer controls pH changes that may occur as result of amine production. Sodium chloride maintains osmotic equilibrium. GCHI Enrichment is a defined supplement which provides vitamins, amino acids, coenzymes, and other factors that stimulate growth. Colistin and nalidixic acid are selective agents, inhibitory to most *Enterobacteriaceae*. Cornstarch neutralizes fatty acids that may be toxic to *G. vaginalis* and is the fermentable substrate in this medium. *G. vaginalis* is recognized by zones of clearing in the opaque medium around the colonies.

## REAGENTS (CLASSICAL FORMULAE)\*

Cornstarch.....	11.0 g	Monopotassium Phosphate.....	1.0 g
Casein Peptone.....	7.5 g	Colistin.....	10.0 mg
Meat Peptone.....	7.5 g	Nalidixic Acid.....	10.0 mg
Sodium Chloride.....	5.0 g	●GCHI Enrichment.....	5.0 ml
Dipotassium Phosphate.....	4.0 g	Agar.....	12.0 g
		Deminerlized Water.....	1000.0 ml

pH 7.2 ± 0.2 @ 25°C

### ●GCHI Enrichment:

Glucose.....	100.0 g	Coccarboxylase.....	0.1 g
Cysteine Hydrochloride.....	25.9 g	Guanine Hydrochloride.....	0.03 g
L-Glutamine.....	10.0 g	Ferric Nitrate.....	0.02 g
L-Cystine.....	1.1 g	p-Aminobenzoic Acid.....	0.013 g
Adenine.....	1.0 g	Vitamin B12.....	0.01 g
NAD.....	0.25 g	Thiamine Hydrochloride.....	0.003 g
		Deminerlized Water.....	1000.0 ml

\*Adjusted as required to meet performance standards.

## PROCEDURE

1. Inoculate and streak the specimen as soon as possible after it is received in the laboratory.
2. If the material is being cultured directly from a swab, roll the swab over a small area of the agar surface and streak for isolation.
3. Incubate plates in 5-10% CO<sub>2</sub> at 33-37°C for up to 48 hours.
4. Examine plate daily for typical colony morphology. On McCarthy Agar, colonies of *G. vaginalis* are surrounded by zones of clearing due to the hydrolysis of cornstarch.

## QUALITY CONTROL

All lot numbers of McCarthy Agar 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

*Gardnerella vaginalis* ATCC® 14018  
*Escherichia coli* ATCC® 25922  
*Proteus mirabilis* ATCC® 12453

### INCUBATION

CO<sub>2</sub>, up to 48 h @ 33-37°C  
Aerobic, 18-24 h @ 33-37°C  
Aerobic, 18-24 h @ 33-37°C

### RESULTS

Growth, zone of clearing around colonies  
Inhibition (partial to complete)  
Inhibition (partial to complete)

## LIMITATIONS

1. Some *Streptococcus* spp. may hydrolyze cornstarch on McCarthy Agar. Additional tests are required for presumptive identification of *G. vaginalis* including, a Gram-stain demonstrating characteristic morphology (i.e., gram-positive, gram-negative, or gram-variable coccobacilli), beta-hemolysis on human blood agar, and negative catalase and oxidase.<sup>7</sup>
2. Obligately anaerobic strains of *G. vaginalis* have been reported; however, studies have demonstrated that anaerobic isolates occur with such infrequency that routine anaerobic incubation of genital cultures is unwarranted.<sup>8</sup>

## BIBLIOGRAPHY

1. Gardner, H.L. and C.D. Dukes. 1955. Am. J. Obstet. Gynecol. 69:962-976.
2. Dunkelberg, W.E., Jr., R. Skaggs, and D.S. Kellogg, Jr. 1970. Appl. Microbiol. 19:47-52.
3. Smith, R.F. 1975. Health Lab Science. 12:219-224.
4. Mickelson, P.A., L.R. McCarthy, and M.E. Mangum. 1977. J. Clin. Microbiol. 5:488-489.
5. MacFaddin, J.F. 1985. Media for Isolation-Cultivation-Identification-Maintenance of Medical Bacteria. Vol. 1. Williams & Wilkins, Baltimore, MD.
6. Bailey, R.K., J.L. Voss, and R.F. Smith. 1979. J. Clin. Microbiol. 9:65-71.
7. Winn, W.C., Jr., S.D. Allen, W.M. Janda, E.W. Koneman, G.W. Procop, P.C. Schreckenberger, and G.L. Woods. 2006. Koneman's Color Atlas and Textbook of Diagnostic Microbiology. 6<sup>th</sup> ed. Lippincott Williams and Wilkins, Philadelphia, PA.

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 1590, Revised February 11, 2008

Printed in U.S.A.

# remel

12076 Santa Fe Drive, Lenexa, KS 66215, USA

General Information: (800) 255-6730    Technical Service: (800) 447-3641    Order Entry: (800) 447-3635

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

Website: [www.remel.com](http://www.remel.com)    Email: [remel@remel.com](mailto:remel@remel.com)