

CAMPY CEFEX AGAR

INTENDED USE

Remel Campy Cefex Agar is a solid medium recommended for use in qualitative procedures for selective and differential isolation of *Campylobacter* species from food and poultry.

SUMMARY AND EXPLANATION

Campylobacter is recognized as a major cause of acute diarrheal disease in humans. Consumption of contaminated food and water accounts for 70% of *Campylobacter*-related illnesses each year, with poultry identified as a primary reservoir.¹ In 1992, Stern developed Campy Cefex Agar for the isolation of *Campylobacter* from poultry.² Oyarzabal et al. reported Campy Cefex Agar demonstrated the best results of the six agars tested for recovery of *Campylobacter* spp. from poultry carcass rinses.³ In 2007, Campy Cefex Agar was recommended by the National Advisory Committee on Microbiological Criteria for Foods (NACMCF) for isolation of *Campylobacter* spp. from poultry.⁴ Campy Cefex Agar is also recommended by the American Public Health Association (APHA) and United States Department of Agriculture (USDA) for the isolation of *Campylobacter* spp. from foods and poultry.^{5,6}

PRINCIPLE

Casein and meat peptones supply nitrogenous substances, carbon, and sulfur required for the growth of *Campylobacter* spp. Sodium chloride is a source of essential electrolytes and maintains osmotic equilibrium. Yeast extract provides B-complex vitamins. Dextrose is a ready source of energy. Ferrous sulfate, sodium pyruvate, and sodium bisulfite enhance the aerotolerance of *Campylobacter* and facilitate its growth. Cefoperazone is a selective agent which inhibits commensal microbial fecal flora and cycloheximide inhibits the growth of fungi. Laked horse blood provides hemin and other essential growth factors. Agar is a solidifying agent.

REAGENTS (CLASSICAL FORMULA)*

Casein Peptone.....	10.0 g	Sodium Pyruvate	0.5 g
Meat Peptone.....	10.0 g	Sodium Bisulfite.....	0.3 g
Sodium Chloride.....	5.0 g	Cycloheximide	0.2 g
Yeast Extract.....	2.0 g	Cefoperazone	33.0 mg
Dextrose.....	1.0 g	Laked Horse Blood.....	5 %
Ferrous Sulfate.....	0.5 g	Agar.....	15.0 g
		Deminerlized 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.

PROCEDURE

Consult current editions of appropriate references for the recommended procedure for sample preparation, inoculation, testing, and interpretation.^{4,5,6}

QUALITY CONTROL

Each lot number of Campy Cefex 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

Campylobacter jejuni ATCC® 33291
Cryptococcus neoformans ATCC® 34877
Pseudomonas aeruginosa ATCC® 27853

INCUBATION

Microaerophilic, up to 72 h @ 40-42°C
Ambient, up to 72 h @ 25-30°C
Ambient, 18-24 h @ 33-37°C

RESULTS

Growth
No growth
No growth

LIMITATIONS

1. Cephalothin-sensitive *Campylobacter* spp. may not be recovered on Campy Cefex Agar because it contains cefoperazone.
2. Colonies isolated on Campy Cefex Agar require additional biochemical testing for definitive identification as *Campylobacter* spp.

BIBLIOGRAPHY

1. Food and Drug Administration. 2000. Bacteriological Analytical Manual Online. Retrieved, January 17, 2011 from: <http://www.fda.gov/Food/ScienceResearch/LaboratoryMethods/BacteriologicalAnalyticalManualBAM/UCM072616>.
2. Stern, N.J., B. Wojton, and K. Kwiatek. 1992. J. Food Prot. 55:514-517.
3. Oyarzabal, O.A., K.S. Macklin, J.M. Barbaree, and R.S. Miller. 2005. Appl. Environ. Microbiol. 71:3351-3354.
4. NACMCF. 2007. Analytical Utility of *Campylobacter* Methodologies. J. Food Prot. 70:241-250.
5. United States Department of Agriculture. 2010. Compliance Guideline for Controlling Salmonella and *Campylobacter* in Poultry. 3rd ed. Retrieved, January 12, 2011 from: http://www.fsis.usda.gov/PDF/Compliance_Guide_Controling_Salmonella_Campylobacter_Poultry_0510.pdf.
6. Downes, F.P. and K. Ito. 2001. Compendium of Methods for the Microbiological Examination of Foods. 4th ed. APHA, Washington, D.C.

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 110138, Revised March 22, 2011

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