

BIBLIOGRAPHY

1. Cary SG, Blair EB. New Transport Medium for Shipment of Clinical Specimens, I. Fecal specimens. J Bacteriol 1964;88:96-98.
2. Christian DL, Ederer GM. Evaluation of Bacteriological Transport Media. Am J Technol 1973;39:12-16.
3. Edwards RR, Ewing WH. Identification of Enterobacteriaceae, 3rd ed. Minneapolis: Burgess Publishing Co., 1972;28, 337-338.
4. Eisenberg HD, Schoenkecht FD, Von Gravenitz A. Collection and Processing of Bacterial Specimens. Cumitech 9. Washington, D.C.: Amer. Soc. Microbio., 1979.
5. Ewing WH. Transport Methods for Enterobacteriaceae and Allied Bacteria. Public Health Lab 1971;29:8-23.
6. Ewing WH, McWhorter AC, Montague TS. Transport Media in the Detection of Salmonella typhi in Carriers. Public Health Lab. 24:63-65.
7. Gaines S, Haque SU, Paniom W, Duongmani C, Cary SG, Blair EB. A Field Trial of a New Transport Medium for Collection of Feces for Bacteriologic Examination. Am J Trop Med Hyg 1965;14:136-140.
8. Kelly MT, Brenner DJ, Farmer III JJ. "Enterobacteriaceae" Manual of Clinical Microbiology, 4th ed. Washington, D.C.: Am. Soc. Microbiol, 1985:263-277.
9. Lior H, Krol P. Abstr. Annu. Meet. Am. Soc. Microbiol. 1979;C:174-339.
10. Leuchtefeld NW, Wang WL, Blaser MJ, Reller LB. Evaluation of Transport and Storage Techniques for Isolation of Campylobacter fetus spp. jejuni from Turkey Fecal Specimen. J clin Microbiol 1981;13:438-443.
11. Monsur KA. Bacteriological Diagnosis of Cholera under Field Conditions. Bul WHO 1963;28:387-389.
12. Morris GK, Patton CM. "Campylobacter" Manual of Clinical Microbiology 4th ed. Washington, D.C.: Am Soc Microbiol. 1985:302-308.
13. Garcia MS, Bruckner D. Diagnostic Medical Parasitology, 3rd ed. Washington D.C.: ASM Press, 1997: 594-822.

Protocol Cary Blair Vials

Protocol Cary Blair Vials, 20/pk

Other Protocol Parasitology Products

Protocol LV-PVA Vials, 20/pk
Protocol SAF Vials, 20/pk
Protocol Clean Vials, 20/pk
Protocol 10% Buff. Formalin Vials, 20/pk
Protocol Modified (Cu) PVA Vials, 20/pk
Protocol C & S Vials, 20/pk
Protocol MIF Vials, 20/pk
Protocol Zn-PVA Vials, 20/pk
Protocol LV-PVA/ Formalin Kit, 2x10 vl/pk
Protocol SAF / Clean Kit, 2x10 vl/pk
Protocol LV-PVA / Clean Kit, 2x10 vl/pk
Protocol SAF / C&S Kit, 2x10 vl/pk
Protocol Modified (Cu) PVA/Formalin, 2x10 vl/pk
Protocol Zn-PVA/10%Formalin Kits, 2x10 vl/pk
Protocol PVA/Formalin/Clean, 3x10 vl/pk
Protocol PVA/Formalin/C&S, 3x10 vl/pk
Protocol SAF/Clean/C&S, 3x10 vl/pk
Protocol Zn-PVN Formalin/ Clean, 3x10 vl/pk
Protocol Mod. (Cu) PVA/Formalin/Clean, 3x10 vl/pk
Protocol Zn-PVA Formalin/ C&S, 3x10 vl/pk
Protocol MIF Kit
Protocol MIF Bulk Kit
Protocol Trichrome, 500 mL
Protocol LV-PVA Bulk, 500 mL
Protocol Zn-PVA Bulk, 500 mL
Protocol Bulk Tubes and caps, 15 mL
Protocol Iodine, 50 mL
Protocol 50 mL Concentration System, 120 ea.
Protocol 15 mL Concentration System, 50 ea.
Protocol 15 mL Concentration System, 4x50 ea.
Protocol Ethyl Acetate

Item

23-005-47

Item

23-005-29
23-005-41
23-005-31
23-005-46
23-005-33
23-005-43
23-005-35
23-005-37
23-005-23
23-005-44
23-005-25
23-005-39
23-005-27
23-005-45
23-005-22
23-005-24
23-005-26
23-005-28
23-005-30
23-005-32
23-005-34
23-005-36
23-005-38
23-005-49
23-005-40
23-005-42
23-005-48
23-005-50
23-005-51
23-005-52
23-005-68

Protocol™

Application for Cary Blair Medium

PURPOSE:

Protocol Parasitology Systems provide standardized procedures for the routine collection, transportation, preservation and examination of stool specimens for enteric pathogens. Protocol Systems are for use by patient or healthcare worker and help minimize the need for immediate transport and processing of specimens.

EXPLANATION

Diagnosis of enteric bacterial disease is confirmed by isolation and identification of pathogenic organisms in stool specimens. Procedures such as freezing, incubation and refrigeration do not insure recovery and identification of enteric bacterial pathogens. Cary Blair transport medium has been formulated to facilitate collection and transportation while maintaining the bacterial population for optimum recovery up to 96 hours after, passage. Proper use of the Protocol Cary Blair System assures the parasitologist that enteric pathogens such as Yersinia, Salmonella, Shigella, Campylobacter, Cholera Vibrio, if present, will be preserved.

CONTENTS:

Each vial contains 15 mL of Cary Blair Medium. Directions for use are also included.

SPECIMEN COLLECTION

1. Caution: Patient should not use antacids, barium, bismuth, antibiotics, anti malarial agents, antidiarrheal medication or oily laxatives prior to specimen collection. After administration of any of these compounds, specimen collection should be delayed for 5 to 10 days or at least two weeks after barium or antibiotics.
2. Several specimens, collected intermittently over several days, should be examined to insure recovery of organisms.
3. Specimens must be collected properly to avoid contamination with urine or water (see collection instructions). Specimens are best collected in., a bedpan. A clean dry container such as a milk container may be used by removing the top and washing thoroughly. Another option is to place plastic wrap over the toilet seat opening.
4. A suitable area (i.e. bloody, slimy, watery) from the sides, ends and middle of the stool should be selected using the collection spoon provided. Fill with sufficient stool to bring the liquid level up to the "Fill" line. This will result in approximately 5 mL of sample.
5. Stir each specimen with the spoon provided, tighten the cap and shake firmly until the specimen is adequately mixed. When mixing is complete the specimen should appear uniform.
6. Complete the label on each vial and replace the vials in the plastic bag. Transport the specimen to the laboratory. Specimen may be refrigerated or kept at room temperature.

SPECIMEN PROCESSING

Protocol Cary Blair Medium contains a buffering system that prevents development of an acid pH shift which may adversely affect recovery of Shigella and Salmonella species. If the buffering capacity of this system is exceeded, it will be indicated by the presence of a yellow color. If this yellow color is present, the medium has reached an adversely acidic pH and should not be used. To insure optimum recovery of organisms, the medium should be red in color. Store Protocol Cary Blair System vials at 20-25°C. The Protocol System allows for a variety of procedures to be utilized and is only part of a complete examination.

Before Processing:

1. Add sufficient sample to bring the level to the Fill line.
2. Maintain the vial at room temperature, avoid extreme temperatures.
3. Mix sample thoroughly using the spoon provided on the cap. The mixture should appear homogeneous.
4. Important: The pH indicator must be red in color. If the medium has turned yellow, this indicates an acid pH and is unsuitable for use. A new specimen must be collected.

Processing:

1. Mix the contents of the vial thoroughly.
2. Remove the vial cap and lay on a paper towel saturated with disinfectant.
3. Inoculate medium from the Cary Blair vial as indicated by the laboratory standard operating procedure.
4. Reseal the vial and dispose of according to your laboratory infectious waste policy. Autoclaving/sterilization prior to disposal is recommended.

PRECAUTIONS

Occupational Safety and Health Act regulations (including Universal Precautions) should be used for handling all specimens.

For assistance please call our Technical Service Department toll-free at 1-800-528-0494 between the hours of 8 A.M. and 5 P.M. Eastern Standard Time.

CAS NUMBERS:

Cary Blair Medium

Agar	9002-18-0
DI Water	7732-18-5
KH ₂ PO ₄	7778-77-0
Na ₂ HPO ₄	7558-79-4
Phenol Red	143-74-8
Sodium Chloride	7647-14-5
Sodium Thioglycolate	367-51-1

Fisher Diagnostics
a division of Fisher Scientific Company, LLC
a subsidiary of Thermo Fisher Scientific Inc.
Middletown, VA 22645-1905 USA
Phone: 1-800-528-0494
1-540-869-3200
Fax: 1-540-869-8132

Thermo
SCIENTIFIC