
remel

**QC-Slide™ for Parasitology
Stain Controls**

INTENDED USE

Remel QC-Slide™ for Parasitology Stain Controls are specifically designed, unstained, ready-to-use microscope slides containing dried suspensions of known parasites. It is intended to be part of a quality control program to monitor microbiological stains and staining techniques used in parasite staining procedures.

PRINCIPLE

Accurate diagnosis of parasitic infections depends on proper collection, transport, detection, and identification of parasites in blood, fecal and other body fluid specimens. Permanently stained smears aid in detection and identification of parasites by providing detailed microscopic visualization of nuclear and cytoplasmic features.

Quality control programs are required by various clinical laboratory regulatory organizations and Good Laboratory Practice regulations. The performance of equipment, reagents, methods, and techniques must be monitored and documented.¹⁻⁵ QC-Slide™ Parasitology Stain Controls can be used as controls with commonly performed staining methods, eliminating the need to maintain samples for preparation of stain control slides.

PRECAUTIONS

This product is For Laboratory Use only and should be used by properly trained individuals. Precautions should be taken against the dangers of microbiological and sharps hazards by properly sterilizing and disposing of specimens, containers, and microscope slides after use. Directions should be read and followed carefully.

STORAGE

This product is ready for use and no further preparation is necessary. Store product in its original container at the temperature indicated on the package label.

PROCEDURE

Stain QC-Slide™ for Parasitology Stain Controls according to the procedures described by the manufacturer of the stains and reagents being used, or consult appropriate references.

INTERPRETATION

Staining reactions will be dictated by the method chosen. Refer to the instructions provided by the stain manufacturer or consult appropriate references.

QUALITY CONTROL

All lot numbers of QC-Slide™ for Parasitology Stain Controls have been analyzed to verify contents. Quality control testing should be performed in accordance with established laboratory procedures. If aberrant quality control results are noted, patient results should not be reported.

LIMITATIONS

Remel QC-Slide™ for Parasitology Stain Controls are intended as stain and staining technique controls, not as teaching slides.




BIBLIOGRAPHY

1. Garcia, L.S. 2001. Diagnostic Medical Parasitology. 4th ed. ASM, Washington, D.C.
2. Melvin, D. and M. Brooke. 1982. Laboratory Procedures for the Diagnosis of Intestinal Parasites. 3rd ed. U.S. Dept. of H.H.S. and CDC, Atlanta, GA.
3. 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.
4. Clinical and Laboratory Standards Institute (CLSI). 2005. Procedures for the Recovery and Identification of Parasites from the Intestinal Tract; Approved Guideline, 2nd ed. M28-A2. CLSI, Wayne, PA.
5. Clinical and Laboratory Standards Institute (CLSI). 2000. Laboratory Diagnosis of Blood-Borne Parasitic Diseases; Approved Guideline. M15-A. CLSI, Wayne, PA.

PACKAGING CONFIGURATIONS

QC-Slide™ for Parasitology Stain Controls are available in a variety of parasite/preservative combinations. Refer to individual product descriptions in the Remel catalog or visit www.remel.com for individual product descriptions.

Symbol Legend

REF	Catalog Number
IVD	In Vitro Diagnostic Medical Device
LAB	For Laboratory Use
	Consult Instructions for Use (IFU)
	Temperature Limitation (Storage Temp.)
LOT	Batch Code (Lot Number)
	Use By (Expiration Date)

QC-Slide™ is a trademark of Remel Inc.

IFU 2410010, Revised August 10, 2010

Printed in U.S.A.

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