

Product datasheet

Fite Stain Kit (Microorganism Stain) ab150668

1 Image

Overview

Product name Fite Stain Kit (Microorganism Stain)

Product overview Fite Stain Kit ab150668 is intended for use in the histological visualization of Mycobacterium *Lepra bacillus* and Nocardia. This kit may be used on formalin-fixed, paraffin-embedded or frozen sections.

Other products for staining tissue sections

Find more kits and reagents in the [special stains guide](#), or products for antigen retrieval, blocking, signal amplification, visualization, counterstaining, and mounting in the [IHC kits and reagents guide](#).

Notes

Staining Interpretation:

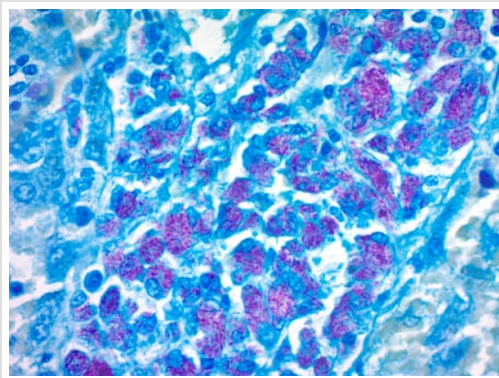
<i>Lepra bacillus</i>	Red
Nocardia	Red
Background	Blue

Control Tissue: Any well fixed paraffin-embedded tissue.

Properties

Storage instructions Store at room temperature. Please refer to protocols.

Components	100 tests
Acid Alcohol Solution (1%)	1 x 500ml
Carbol Fuchsin Solution	1 x 125ml
Methylene Blue solution	1 x 125ml
Xylene-Peanut Oil Solution	1 x 125ml



ab150668 - Fite's Stain Kit (Microorganism Stain)

ab150668 Fite's Stain Kit (Microorganism Stain) staining formalin-fixed-paraffin embedded acid fast bacteria.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours

- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors