

Product datasheet

Fontana-Masson Stain Kit (Melanin Stain) ab150669

★★★★★ 2 Abreviews 14 References 1 Image

Overview

Product name Fontana-Masson Stain Kit (Melanin Stain)

Product overview The Fontana-Masson Stain is intended for use in the histological visualization of Argentaffin cells and Melanin in paraffin or frozen sections. In addition, the Fontana-Masson stain has been reported to be useful in identifying Capsule-Deficient Cryptococcus Neoformans and typical Cryptococcus Neoformans.

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

Argentaffin Cells	Black
Melanin	Black
Nuclei	Red
Cytoplasm	Light Pink
Cryptococci Cell Wall	Black

Control Tissue: Any paraffin-embedded tissue that contains Hair Follicles or Skin for Melanin. Small intestine for Argentaffin.

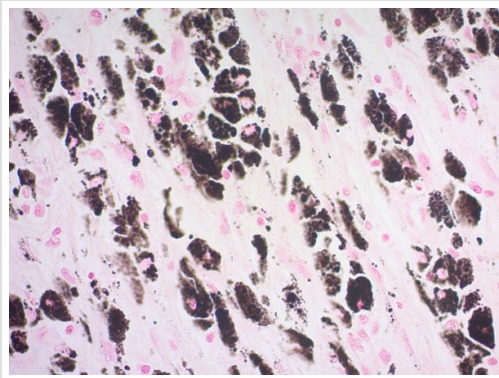
Properties

Storage instructions Store at +4°C. Please refer to protocols.

Components	100 tests
Gold Chloride Solution (0.2%)	1 x 125ml

Components	100 tests
Nuclear Fast Red Solution	1 x 125ml
Silver Nitrate Solution (10%)	5 x 9ml
Sodium Thiosulfate Solution (5%)	1 x 125ml

Images



ab150669 Fontana-Masson Stain Kit (Melanin Stain) staining formalin-fixed-paraffin embedded human skin melanoma.

ab150669 - Fontana-Masson Stain Kit (Melanin Stain)

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