

Version 3 Last updated 11 November 2019

ab150669 Fontana-Masson (Melanin Stain)

For the histological visualization of Argentaffin cells and melanin.

[View kit datasheet: www.abcam.com/\[\]](http://www.abcam.com/)
(use [www.abcam.cn/\[\]](http://www.abcam.cn/) for China, or [www.abcam.co.jp/\[\]](http://www.abcam.co.jp/) for Japan)

This product is for research use only and is not intended for diagnostic use.

Table of Contents

1. Overview	3
2. Materials Supplied and Storage	4
3. Materials Required, Not Supplied	4
4. General guidelines, precautions, and troubleshooting	5
5. Staining Protocol	6
6. FAQs / Troubleshooting	7

1. Overview

The Fontana-Masson (Melanin Stain) (ab150669) 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.

Staining Interpretation:

Argentaffin cells	Black
Melanin	Black
Cryptococci Cell Wall	Black
Nuclei	Red
Cytoplasm	Light pink

Control Tissue:

Any paraffin-embedded tissue that contains hair follicles or skin for melanin.

Small intestine for Argentaffin.

2. Materials Supplied and Storage

Store kit at room temperature immediately on receipt and check below for storage for individual components. Kit can be stored for 1 year from receipt, if components have not been reconstituted.

Keep away from open flame and refer to the safety datasheet.

Item	Quantity	Storage temperature (before prep)
Gold Chloride Solution (0.2%)	125 mL	4°C
Silver Nitrate Solution (10%)	5 x 9 mL	4°C
Sodium Thiosulfate Solution (5%)	125 mL	RT
Nuclear Fast Red Solution	125 mL	RT

3. Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully perform this assay:

- Concentrate ammonium hydroxide
- Absolute alcohol

4. General guidelines, precautions, and troubleshooting

Please observe safe laboratory practice and consult the safety datasheet.

For general guidelines, precautions, limitations on the use of our assay kits and general assay troubleshooting tips, particularly for first time users, please consult our guide:

www.abcam.com/assaykitguidelines

For typical data produced using the assay, please see the assay kit datasheet on our website.

5. Staining Protocol

- Equilibrate all materials and prepared reagents to room temperature just prior to use and gently agitate.

Prepare Ammoniacal Silver Solution immediately prior to use:

In new or chemically cleaned glassware, mix 27 ml distilled/deionized water with one vial of Silver Nitrate Solution and blend completely. Carefully add concentrated Ammonium Hydroxide one drop at a time, swirling gently after each drop.

Initially the mixture will turn dark brown and then gradually become transparent with a fine layer of sediment. The solution is ready for immediate use when all sediment dissolves.

- 5.1** Deparaffinize sections if necessary and hydrate in distilled water.
- 5.2** Place freshly mixed Ammoniacal Silver solution in a 58-60°C water bath and allow adequate time for the temperature to equilibrate.
- 5.3** Incubate slide in warmed Ammoniacal Silver solution for 30-60 minutes or until tissue section becomes yellowish/brown in color.

ΔNote: Melanin typically stains in 30 minutes while Argentaffin stains in 50-60 minutes.

- 5.4** Rinse slide in several changes of distilled water.
- 5.5** Incubate slide in Gold Chloride Solution (0.2%) at room temperature for 30 seconds.
- 5.6** Rinse slide in several changes of distilled water.
- 5.7** Incubate slide in Sodium Thiosulfate Solution (5%) at room temperature for 1-2 minutes.
- 5.8** Rinse for 2 minutes in running tap water for followed by 2 changes of distilled water.
- 5.9** Incubate slide in Nuclear Fast Red Solution for 5 minutes.
- 5.10** Rinse for 2 minutes in running tap water for followed by 2 changes of distilled water.
- 5.11** Dehydrate very quickly in 3 changes of fresh Absolute Alcohol.
- 5.12** Clear and mount in synthetic resin (or organic/ permanent mounting medium).

6. FAQs / Troubleshooting

General troubleshooting points are found at www.abcam.com/assaykitguidelines.

Technical Support

Copyright © 2012-2019 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

Austria

wissenschaftlicherdienst@abcam.com | 019-288-259

France

supportscientifique@abcam.com | 01.46.94.62.96

Germany

wissenschaftlicherdienst@abcam.com | 030-896-779-154

Spain

soportecientifico@abcam.com | 91-114-65-60

Switzerland

technical@abcam.com

Deutsch: 043-501-64-24 | Français: 061-500-05-30

UK, EU and ROW

technical@abcam.com | +44(0)1223-696000

Canada

ca.technical@abcam.com | 877-749-8807

US and Latin America

us.technical@abcam.com | 888-772-2226

Asia Pacific

hk.technical@abcam.com | (852) 2603-6823

China

cn.technical@abcam.com | 400 921 0189 / +86 21 2070 0500

Japan

technical@abcam.co.jp | +81-(0)3-6231-0940

Singapore

sg.technical@abcam.com | 800 188-5244

Australia

au.technical@abcam.com | +61-(0)3-8652-1450

New Zealand

nz.technical@abc.com | +64-(0)9-909-7829