

Version 3c Last updated 14 June 2021

ab150665 Combined Eosinophil- Mast Cell (Hematologic Stain)

For the simultaneous visualization of eosinophils and mast cells.

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

Table of Contents

1. Overview	1
2. Materials Supplied and Storage	2
3. Materials Required, Not Supplied	3
4. General guidelines, precautions, and troubleshooting	4
5. Assay Procedure	5
6. FAQs / Troubleshooting	6
7. Typical Data	7
8. Notes	8

1. Overview

The Combined Eosinophil-Mast Cell (Hematologic Stain) is intended for the simultaneous visualization of eosinophils and mast cells.

Staining Interpretation:

Mast cells	Bright Blue
Eosinophils	Bright Red
Nuclei	Blue

Control Tissue:

Any well-fixed section (3-5 microns) or any air-dried smear on a clear slide.

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)
Astra Blue Solution	125 ml	RT
Vital New Red Solution	125 ml	RT
Hematoxylin (Modified Mayer's 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:

- 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 Procedure

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

- 5.1** Deparaffinize sections if necessary and hydrate in distilled water.
- 5.2** Stain tissue section with Astra Blue Solution for 1-3 minutes.
- 5.3** Rinse quickly in 2 changes of distilled water.
- 5.4** Stain tissue section with Vital New Red Solution for 30 minutes with frequent agitation.
- 5.5** Rinse quickly in 2 changes of distilled water.
Optional:
 - a. Counterstain in Hematoxylin (Modified Mayer's solution) for 15-30 seconds.
 - b. Rinse slide in running tap water for 2 minutes followed by one dip in distilled water.
- 5.6** Blot dry.
- 5.7** Dehydrate quickly through 3 changes of absolute alcohol (about 20 seconds each).
- 5.8** Shake off excess and blot dry.
- 5.9** Air dry for about 1 minute.
- 5.10** Clean and mount in synthetic resin.

6. FAQs / Troubleshooting

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

7. Notes

Technical Support

Copyright © 2021 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