

Version 3 Last updated 19 December 2018

ab150670

Giemsa Stain Kit

For the histological visualization of Cells present in Hematopoietic Tissues and Certain Microorganisms.

[View kit datasheet: www.abcam.com/ab150670](http://www.abcam.com/ab150670)
(use www.abcam.cn/ab150670 for China, or www.abcam.co.jp/ab150670 for Japan)

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. Staining Protocol	5
6. FAQs / Troubleshooting	7
7. Typical Data	8
8. Notes	9

1. Overview

Giemsa Stain Kit (ab150670) is intended for use in the visualization of cells present in hematopoietic tissues and certain microorganisms. This kit may be used on formalin-fixed, paraffin-embedded or frozen sections.

Staining Interpretation:

Nuclei	Blue/Violet
Cytoplasm	Light Blue
Collagen	Pale Pink
Muscle Fibers	Pale Pink
Erythrocytes	Gray, Yellow or Pink
Rickettsia	Reddish-Purple
<i>Helicobacter pylori</i>	Blue
Mast Cells	Dark Blue with Red Granules

Control Tissue: Blood film. Bone Marrow. Spleen. Any well fixed tissue.

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.

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

Item	Quantity	Storage temperature (before prep)
May-Grunwald Stock Solution	500 mL	RT
Giemsa Stock Solution	500 mL	RT
Phosphate Buffer Solution (pH 6.8)	500 mL	RT

3. Materials Required, Not Supplied

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

- Xylene or xylene substitute.
- Acetic Acid Solution (0.25%).
- Distilled water.

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.

5.1 Preparation of Solutions:

1. Prepare Working May-Grunwald Solution. Mix.

Component	Volume (mL)
May-Grunwald Stock Solution	25
Phosphate Buffer Solution (pH6.8)	25

2. Prepare Working Giemsa Solution. Mix.

Component	Volume (mL)
Giemsa Stock Solution	2.5
Phosphate Buffer Solution (pH6.8)	50

5.2 Standard Procedure:

1. Deparaffinize sections if necessary and hydrate in distilled water.
2. Place slide in staining tray and flood with Working May-Grunwald solution for 5-7 minutes. Agitate slide occasionally to ensure proper staining.
3. Carefully flood slide with Phosphate Buffer Solution (pH6.8) until no stain runs off.
4. Flood with Working Giemsa Solution for 10-15 minutes. Agitate slide occasionally to ensure proper staining.
5. Carefully flood slide with Phosphate Buffer Solution (pH6.8) until no stain runs off.
6. Allow slide to remain in Phosphate Buffer Solution, pH 6.8 for an additional 3 minutes.
7. Dip slide quickly in distilled water and air dry at room temperature.
8. Dip slide in Xylene or Xylene substitute.
9. Mount in synthetic resin.

5.3 Mast Cell Procedure:

1. Deparaffinize sections if necessary and hydrate in distilled water.
2. Place slide in staining tray and flood with Working May-Grunwald solution for 5-7 minutes. Agitate slide occasionally to ensure proper staining.
3. Carefully flood slide with Phosphate Buffer solution (pH6.8) until no stain runs off.
4. Flood with Working Giemsa Solution for 10-15 minutes. Agitate slide occasionally to ensure proper staining.
5. Carefully flood slide with Phosphate Buffer Solution (pH6.8) until no stain runs off.
6. Differentiate by dipping slide in Acetic Acid Solution (0.25%) until background is desired intensity.
7. Dip slide for 10 seconds in Phosphate Buffer Solution, pH 6.8 while agitating gently.
8. Dip slide quickly in distilled water and air dry at room temperature.
9. Dip slide in Xylene or Xylene substitute.
10. Mount in synthetic resin

6. FAQs / Troubleshooting

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

7. Notes

Technical Support

Copyright © 2012-2018 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