

Version 2a Last updated 31 January 2024

# ab245888 Wright-Giemsa Stain Kit

For the differential staining of blood smears, bone marrow and blood parasites.

View Alcian Blue Stain Kit (pH 0.5) datasheet:

<https://www.abcam.com/ab245888>

[use [www.abcam.cn/ab245888](http://www.abcam.cn/ab245888) for China, or [www.abcam.co.jp/ab245888](http://www.abcam.co.jp/ab245888) 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. General guidelines, precautions, and troubleshooting	3
4. Assay Procedure	4
5. Notes	5

# 1. Overview

The Wright-Giemsa Stain Kit (ab245888) is intended to be used for differential staining of blood smears, bone marrow and blood parasites.

## **Staining Interpretation:**

Erythrocytes: Pink-Tan

Leukocytes: Blue-Purple

Neutrophils: Light Purple or lavender granules in cytoplasm

Eosinophils: Bright Red or Red-Orange granules in cytoplasm

Basophils: Deep Purple or Violet-Black granules in cytoplasm

Platelets: Violet-Purple granules in light blue cytoplasm

## **Control Tissue:**

Blood smear on clean 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)	Storage temperature (after prep)
Wright-Giesma Solution	500 mL	RT	RT
Phosphate Buffer Solution (pH 6.8)	2 x 500 mL	RT	RT

## 3. Materials required but not supplied

- Methanol, 500 ml
- Distilled water, 1 Gallon

## 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](http://www.abcam.com/assaykitguidelines)

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

## 5. Assay Procedure

- Equilibrate all materials and prepared reagents to room temperature just prior to use and gently agitate.
  - Prepare working Wright-Giemsa Solution by mixing 25 mL of Wright-Giemsa Solution with 25 mL of phosphate Buffer Solution, pH 6.8.
- 1.1 Smear a small drop of blood on a clean microscope slide and allow to air dry.
  - 1.2 Fix by placing in absolute Methanol for 5 minutes.
  - 1.3 Place slide in staining tray and flood with Working Wright-Giemsa Solution for 5 minutes. Note: Agitate slide occasionally to insure proper staining.
  - 1.4 Rinse slide in deionized/distilled water.
  - 1.5 Flood slide with Phosphate Buffer Solution, pH 6.8 until no stain runs off.
  - 1.6 Allow slide to remain in Phosphate Buffer Solution, pH 6.8 for an additional 1 minute.
  - 1.7 Dip slide in distilled water and air dry at room temperature.
  - 1.8 Dip slide several times in Xylene or Xylene Substitute.
  - 1.9 Mount in synthetic resin.

## 6. Notes

# Technical Support

Copyright © 2024 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

**For all technical or commercial enquiries please go to:**

[www.abcam.com/contactus](http://www.abcam.com/contactus)

[www.abcam.cn/contactus](http://www.abcam.cn/contactus) (China)

[www.abcam.co.jp/contactus](http://www.abcam.co.jp/contactus) (Japan)