

PRODUCT INFORMATION



Giemsa Stain

Item No. 38336

CAS Registry No.: 51811-82-6

Synonym: Azure Stain

Supplied as: A mixture of methylene blue, azure B, and eosin Y

Storage: -20°C

Stability: ≥4 years

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Giemsa stain is supplied as a solid. A stock solution may be made by dissolving the giemsa stain in the solvent of choice, which should be purged with an inert gas. Giemsa stain is soluble in methanol and acetonitrile.

Description

Giemsa stain is a nucleic acid stain that has commonly been used in the detection of malaria and other parasites, as well as in cytogenetic applications.^{1,2}

References

1. Barcia, J.J. The Giemsa stain: Its history and applications. *Int. J. Surg. Pathol.* **15**(3), 292-296 (2007).
2. Dolan, M. The role of the Giemsa stain in cytogenetics. *Biotech. Histochem.* **86**(2), 94-97 (2011).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/13/2025

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM