

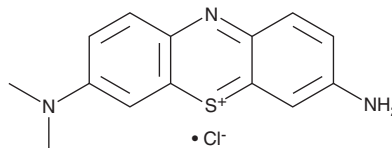
PRODUCT INFORMATION



Azure A

Item No. 44525

CAS Registry No.: 531-53-3
Formal Name: 3-amino-7-(dimethylamino)-
phenothiazin-5-ium, monochloride
MF: $C_{14}H_{14}N_3S \cdot Cl$
FW: 291.8
Absorbance: 620-634 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

Azure A is supplied as a solid. A stock solution may be made by dissolving the azure A in the solvent of choice, which should be purged with an inert gas. Azure A is slightly soluble (0.1-1 mg/ml) in DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of azure A can be prepared by directly dissolving the solid in aqueous buffers. Azure A is slightly soluble (0.1-1 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Azure A is a phenothiazine dye.¹ It displays an absorbance maximum between 620-634 nm. Azure A has been used as a starting material in the synthesis of DNA damage-sensing electrochemical probes and NADH-detecting electrodes.^{2,3}

References

1. Sabnis, R.W. Handbook of biological dyes and stains: Synthesis and industrial applications. John Wiley & Sons, Inc. 1st. (2010).
3. Porfireva, A., Platinina, K., Evtugyn, V., et al. Electrochemical DNA sensor based on poly(Azure A) obtained from the buffer saturated with chloroform. *Sensors (Basel)* **21**(9), 2949 (2021).
4. Revenga-Parra, M., Gómez-Anquela, C., García-Mendiola, T., et al. Grafted Azure A modified electrodes as disposable β -nicotinamide adenine dinucleotide sensors. *Anal. Chim. Acta* **747**, 84-91 (2012).

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, 11/04/2025

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