

PRODUCT INFORMATION

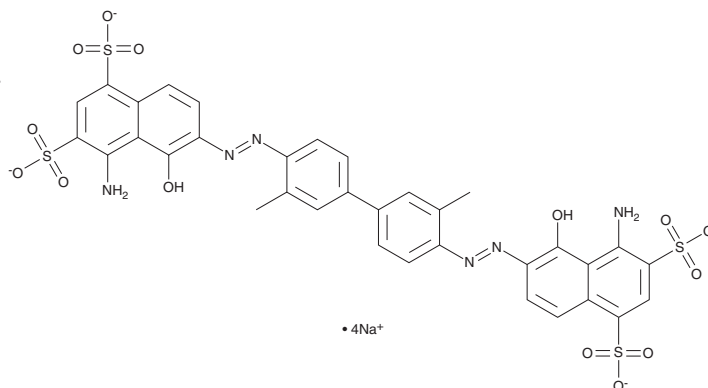


Evans Blue

Item No. 44513

CAS Registry No.: 314-13-6
Formal Name: 6,6'-[(3,3'-dimethyl[1,1'-biphenyl]-4,4'-diyl)bis(2,1-diazenediyl)]bis[4-amino-5-hydroxy-1,3-naphthalenedisulfonic acid, tetrasodium salt

Synonyms: Direct Blue 53, T-1824
MF: C₃₄H₂₄N₆O₁₄S₄ • 4Na
FW: 960.8
UV/Vis.: λ_{max}: 236, 316, 606 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

Evans blue is supplied as a solid. A stock solution may be made by dissolving the Evans blue in the solvent of choice, which should be purged with an inert gas. Evans blue is soluble (≥10 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 Evans blue can be prepared by directly dissolving the solid in aqueous buffers. Evans blue is soluble (≥10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Evans blue is a synthetic bis-azo dye.¹ It displays an absorbance maximum at 620 nm. Evans blue has been used for the estimation of blood volume and determination of blood-brain barrier integrity, among other applications.^{1,2}

References

1. Yao, L., Xue, X., Yu, P., *et al.* Evans blue dye: A revisit of its applications in biomedicine. *Contrast Media & Mol. Imaging* **2018**, 7628037 (2018).
2. Goldim, M.P.S., Della Giustina, A., and Petronilho, F. Using evans blue dye to determine blood-brain barrier integrity in rodents. *Curr. Protec. Immunol.* **126(1)**, e83 (2019).

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, 12/09/2025

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