

PRODUCT INFORMATION



Basic Blue 7 Item No. 44934

CAS Registry No.: 2390-60-5

Formal Name: N-[4-[[4-(diethylamino)phenyl]
[4-(ethylamino)-1-naphthalenyl]methylene]-
2,5-cyclohexadien-1-ylidene]-N-ethyl-
ethanaminium, monochloride

Synonyms: C.I. 42595, Solvent Blue 5, Victoria Pure Blue

MF: $C_{33}H_{40}N_3 \cdot Cl$

FW: 514.2

UV/Vis.: λ_{max} : 334, 590 nm

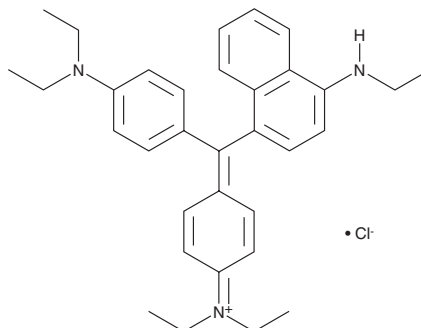
Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Special Conditions: Light sensitive; Store in amber vials

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



Laboratory Procedures

Basic blue 7 is supplied as a solid. A stock solution may be made by dissolving the basic blue 7 in the solvent of choice, which should be purged with an inert gas. Basic blue 7 is soluble (≥ 10 mg/ml) in ethanol and DMSO.

Description

Basic blue 7 is a cationic textile dye.¹ It displays an absorption maximum of 619 nm.² Basic blue 7 has been found as a contaminant in textile industry wastewater effluents.¹

References

1. Tunali Akar, S., Koc, E., Sayin, F., *et al.* Design and modeling of the decolorization characteristics of a regenerable and eco-friendly geopolymer: Batch and dynamic flow mode treatment aspects. *J. Environ. Manage.* **298**, 113548 (2021).
2. Sekine, Y., Nankawa, T., Yunoki, S., *et al.* Eco-friendly carboxymethyl cellulose nanofiber hydrogels prepared via freeze cross-linking and their applications. *App. Polymer Mater.* **2(12)**, 5482-5491 (2020).

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, 02/16/2026

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