

PRODUCT INFORMATION



Basic Blue 26

Item No. 44826

CAS Registry No.: 2580-56-5

Formal Name: N-[4-[[4-(dimethylamino)phenyl]
[4-(phenylamino)-1-naphthalenyl]methylene]-
2,5-cyclohexadien-1-ylidene]-N-methyl-
methanaminium, monochloride

Synonyms: C.I. 44045, Calcozine Blue B, Victoria Blue B

MF: $C_{33}H_{32}N_3 \cdot Cl$

FW: 506.1

Abs. Max: 605 nm

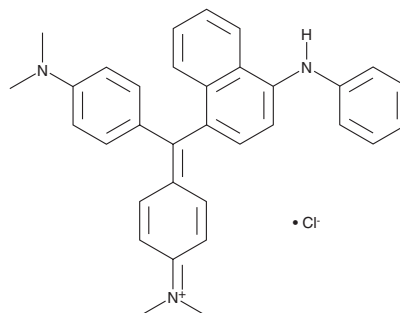
UV/Vis.: λ_{max} : 210, 252, 600 nm

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Special Conditions: Light sensitive



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

Laboratory Procedures

Basic blue 26 is supplied as a solid. A stock solution may be made by dissolving the basic blue 26 in the solvent of choice, which should be purged with an inert gas. Basic blue 26 is soluble (≥ 10 mg/ml) in ethanol and DMSO.

Description

Basic blue 26 is a cationic textile dye.¹ It displays an absorption maximum of 605 nm. Basic blue 26 has been found as a contaminant in textile industry wastewater effluents.²

References

1. Eryomin, V.A. and Poznyakov, S.P. Quantitative determination of phospholipids using the dyes Victoria blue R and B. *Anal. Biochem.* **180(1)**, 186-191 (1989).
2. Belaid, K.D. and Seddiki, H. Influence of anions and cations on the oxidation of basic blue 26 and basic violet 3 dyes by the Fenton process. *SEES* **5(1)**, 2423-2440 (2024).

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, 12/17/2025

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