

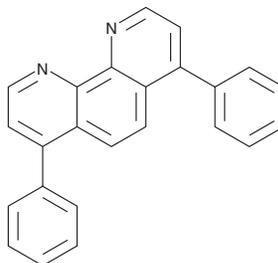
PRODUCT INFORMATION



Bathophenanthroline

Item No. 34811

CAS Registry No.: 1662-01-7
Formal Name: 4,7-diphenyl-1,10-phenanthroline
Synonyms: BPhen, NSC 637659
MF: C₂₄H₁₆N₂
FW: 332.4
Purity: ≥98%
UV/Vis.: λ_{max}: 220, 274 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

Bathophenanthroline is supplied as a solid. A stock solution may be made by dissolving the bathophenanthroline in the solvent of choice, which should be purged with an inert gas. Bathophenanthroline is soluble in DMSO.

Description

Bathophenanthroline is a colorimetric reagent for the detection of iron.^{1,2} It forms a complex with ferrous iron that can be quantified by colorimetric detection at 535 nm as a measure of iron concentration.

References

1. Ghosh, M.M., O'Connor, J.T., and Engelbrecht, R.S. Bathophenanthroline method for the determination of ferrous iron. *J. Am. Water Work. Assoc.* **59(7)**, 897-905 (1967).
2. Patel, M. and Ramavataram, D.V.S.S. A simple, rapid and improved colorimetric assay for non transferrin thalassemia patients bound iron estimation in thalassemia patients. *Int. J. Pharm. Life Sci.* **4(1)**, 2294-2305 (2013).

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, 11/22/2022

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