

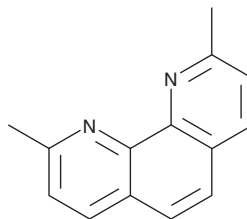
PRODUCT INFORMATION



Neocuproine

Item No. 20874

CAS Registry No.: 484-11-7
Formal Name: 2,9-dimethyl-1,10-phenanthroline
Synonyms: NSC 4280, VUF 7738
MF: $C_{14}H_{12}N_2$
FW: 208.3
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 231, 269 nm
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

Neocuproine is supplied as a crystalline solid. A stock solution may be made by dissolving the neocuproine in the solvent of choice, which should be purged with an inert gas. Neocuproine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of neocuproine in ethanol is approximately 10 mg/ml and approximately 25 mg/ml in DMSO and DMF.

Neocuproine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, neocuproine should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Neocuproine has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Neocuproine is a phenanthroline compound that is used as a chelating agent in a wide array of transition metal-catalyzed reactions.¹ It can also be used in the synthesis of aryl ketones.

Reference

1. Gouda, A.A. and Amin, A.S. Copper(II)-neocuproine reagent for spectrophotometric determination of captopril in pure form and pharmaceutical formulations. *Arabian J. Chem.* **3**, 159-165 (2010).

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, 12/01/2022

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