

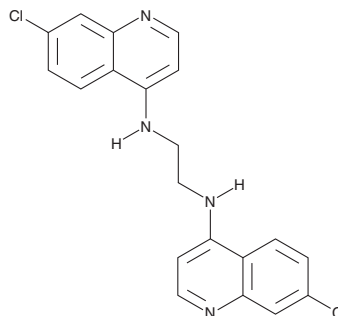
# PRODUCT INFORMATION



**NSC 5844**

Item No. 23443

**CAS Registry No.:** 140926-75-6  
**Formal Name:** N<sup>1</sup>,N<sup>2</sup>-bis(7-chloro-4-quinolinyl)-  
1,2-ethanediamine  
**MF:** C<sub>20</sub>H<sub>16</sub>Cl<sub>2</sub>N<sub>4</sub>  
**FW:** 383.3  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 219, 254, 336 nm  
**Supplied as:** A crystalline 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

NSC 5844 is supplied as a crystalline solid. A stock solution may be made by dissolving the NSC 5844 in the solvent of choice, which should be purged with an inert gas. NSC 5844 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of NSC 5844 in these solvents is approximately 0.25 and 0.33 mg/ml, respectively. NSC 5844 is also slightly soluble in ethanol.

## Description

NSC 5844 is a bis-quinoline with diverse biological activities.<sup>1-3</sup> It inhibits the growth of *P. falciparum* strains that are sensitive (D-6) and resistant (W-2) to chloroquine (Item No. 14194) *in vitro* (IC<sub>50</sub>s = 17 and 27 nM, respectively) but lacks activity against *P. berghei* *in vivo*.<sup>1</sup> NSC 5844 inhibits the growth of MDA-MB-468 and MCF-7 breast cancer cells with GI<sub>50</sub> values of 7.35 and 14.80 μM, respectively.<sup>2</sup> It also inhibits 24% of botulinum neurotoxin serotype A light chain (BoNT/A LC) metalloprotease activity at a concentration of 20 μM.<sup>3</sup>

## References

1. Vennerstrom, J.L., Ellis, W.Y., Ager, A.L., Jr., *et al.* Bisquinolines. 1. N,N-bis(7-chloroquinolin-4-yl)alkane-diamines with potential against chloroquine-resistant malaria. *J. Med. Chem.* **35**(11), 2129-2134 (1992).
2. Zhang, H., Solomon, V.R., Hu, C., *et al.* Synthesis and *in vitro* cytotoxicity evaluation of 4-aminoquinoline derivatives. *Biomed. Pharmacother.* **62**(2), 65-69 (2008).
3. Burnett, J.C., Schmidt, J.J., Stafford, R.G., *et al.* Novel small molecule inhibitors of botulinum neurotoxin A metalloprotease activity. *Biochem. Biophys. Res. Commun.* **310**(1), 84-93 (2003).

### 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, 11/08/2022

## 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](http://WWW.CAYMANCHEM.COM)