

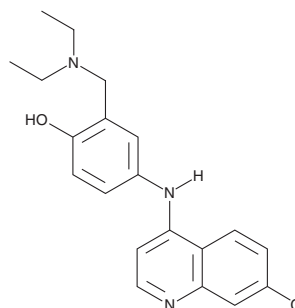
PRODUCT INFORMATION



Amodiaquine

Item No. 15954

CAS Registry No.: 86-42-0
Formal Name: 4-[(7-chloro-4-quinolinyl)amino]-2-[(diethylamino)methyl]-phenol
Synonyms: Camoquine, Flavoquine, NSC 13453, SN 10751
MF: C₂₀H₂₂ClN₃O
FW: 355.9
Purity: ≥95%
UV/Vis.: λ_{max}: 224, 343 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

Amodiaquine is supplied as a crystalline solid. A stock solution may be made by dissolving the amodiaquine in the solvent of choice, which should be purged with an inert gas. Amodiaquine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of amodiaquine in these solvents is approximately 2, 5, and 2.5 mg/ml, respectively.

Description

Amodiaquine is a prodrug form of the antimalarial compound N-desethyl amodiaquine (Item No. 20822).^{1,2} It is active against several strains of *P. falciparum* *in vitro* (EC₅₀s = 0.23-0.52 nM) and exhibits a synergistic effect when used in combination with N-desethyl amodiaquine.¹ Amodiaquine dose-dependently inhibits development of parasitemia in a mouse model of *P. berghei* infection.³

References

1. Mariga, S.T., Gil, J.P., Sisowath, C., *et al.* Synergism between amodiaquine and its major metabolite, desethylamodiaquine, against *Plasmodium falciparum* *in vitro*. *Antimicrob. Agents Chemother.* **48(11)**, 4089-4096 (2004).
2. Sá, J.M., Chong, J.L., and Wellems, T.E. Malaria drug resistance: New observations and developments. *Essays Biochem.* **51**, 137-160 (2014).
3. Jacobs, R.L., Alling, D.W., and Cantrell, W.F. An evaluation of antimalarial combinations against *Plasmodium berghei* in the mouse. *J. Parasitol.* **49(6)**, 920-925 (1963).

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, 12/15/2022

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