

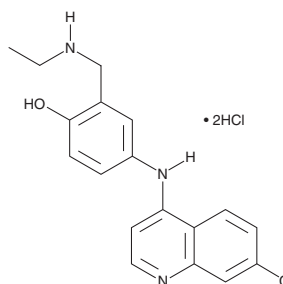
PRODUCT INFORMATION



N-desethyl Amodiaquine (hydrochloride)

Item No. 20822

CAS Registry No.: 79049-30-2
Formal Name: 4-[(7-chloro-4-quinolinyl)amino]-2-[(ethylamino)methyl]-phenol, dihydrochloride
MF: C₁₈H₁₈ClN₃O • 2HCl
FW: 400.7
Purity: ≥95%
UV/Vis.: λ_{max}: 222, 254, 337 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

N-desethyl Amodiaquine (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the N-desethyl amodiaquine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. N-desethyl Amodiaquine (hydrochloride) is soluble in methanol. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

N-desethyl Amodiaquine is a primary metabolite of the antimalarial compound amodiaquine (Item No. 15954), produced by the action of cytochrome P450 isoform 2C8.¹ N-desethyl Amodiaquine is highly active against *P. falciparum* and can synergize with amodiaquine.²

References

1. Li, X.-Q., Björkman, A., Andersson, T.B., *et al.* Amodiaquine clearance and its metabolism to N-desethylamodiaquine is mediated by CYP2C8: A new high affinity and turnover enzyme-specific probe substrate. *J. Pharmacol. Exp. Ther.* **300(2)**, 399-407 (2002).
2. Mariga, S.T., Gil, J.P., Sisowath, C., *et al.* Synergism between amodiaquine and its major metabolite, desethylamodiaquine, against *P. falciparum* *in vitro*. *Antimicrob. Agents Chemother.* **48(11)**, 4089-4096 (2004).

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