

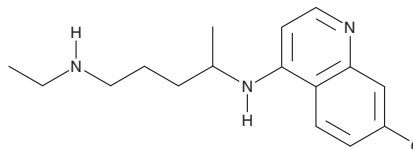
PRODUCT INFORMATION



Desethylchloroquine

Item No. 30113

CAS Registry No.: 1476-52-4
Formal Name: N⁴-(7-chloro-4-quinolinyl)-N¹-ethyl-1,4-pentanediamine
Synonyms: N-Desethylchloroquine, NSC 13254
MF: C₁₆H₂₂ClN₃
FW: 291.8
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Desethylchloroquine is supplied as a solid. A stock solution may be made by dissolving the desethylchloroquine in the solvent of choice, which should be purged with an inert gas. Desethylchloroquine is slightly soluble in chloroform and methanol.

Description

Desethylchloroquine is a major active metabolite of chloroquine (Item No. 14194).¹ Desethylchloroquine is formed when chloroquine undergoes dealkylation, primarily by the cytochrome P450 (CYP) isoforms CYP2C8 and CYP3A4 and to a lesser extent by CYP2D6. It inhibits the growth of the *P. falciparum* strain LA136 *in vitro* (IC₅₀ = 9.9 ng/ml).²

References

1. Projean, D., Baune, B., Farinotti, R., *et al.* In vitro metabolism of chloroquine: identification of CYP2C8, CYP3A4, and CYP2D6 as the main isoforms catalyzing N-desethylchloroquine formation. *Drug Metab. Dispos.* **31(6)**, 748-754 (2003).
2. Aderounmu, F.A. In vitro assessment of the antimalarial activity of chloroquine and its major metabolites. *Ann. Trop. Med. Parasitol.* **78(6)**, 581-585 (1984).

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, 04/01/2020

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