

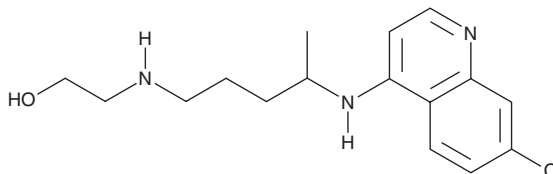
PRODUCT INFORMATION



Desethyl Hydroxychloroquine

Item No. 31152

CAS Registry No.: 4298-15-1
Formal Name: 2-[[4-[(7-chloro-4-quinolinyl)amino]pentyl]amino]-ethanol
Synonyms: DHCQ, Cletoquine, (±)-Desethylhydroxychloroquine
MF: $C_{16}H_{22}ClN_3O$
FW: 307.8
Purity: ≥95%
Supplied as: A 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

Desethyl hydroxychloroquine is supplied as a solid. A stock solution may be made by dissolving the desethyl hydroxychloroquine in the solvent of choice, which should be purged with an inert gas. Desethyl hydroxychloroquine is soluble in methanol, DMSO, and dimethyl formamide.

Description

Desethyl hydroxychloroquine is an active metabolite of hydroxychloroquine (Item No. 17911).¹ It is formed via the desethylation of hydroxychloroquine in the liver, a process mediated by the cytochrome P450 (CYP) isoforms CYP2D6, CYP3A4, CYP3A5, and CYP2C8.

Reference

- Charlier, B., Pigeon, M., Dal Piaz, F., *et al.* Development of a novel ion-pairing HPLC-FL method for the separation and quantification of hydroxychloroquine and its metabolites in whole blood. *Biomed. Chromatogr.* **32**(8), e4258 (2018).

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, 10/24/2022

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