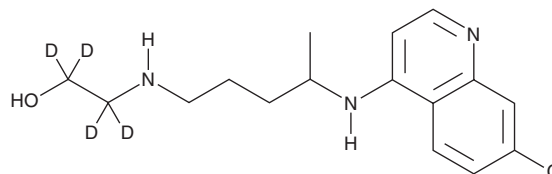


# PRODUCT INFORMATION



## Desethyl Hydroxychloroquine-d<sub>4</sub> Item No. 31162

**CAS Registry No.:** 1854126-47-8  
**Formal Name:** 2-[[4-[(7-chloro-4-quinoliny)amino]pentyl]amino]-ethan-1,1,2,2-d<sub>4</sub>-ol  
**Synonyms:** Cletoquine-d<sub>4</sub>, (±)-Desethylhydroxychloroquine-d<sub>4</sub>, DHCQ-d<sub>4</sub>  
**MF:** C<sub>16</sub>H<sub>18</sub>ClD<sub>4</sub>N<sub>3</sub>O  
**FW:** 311.8  
**Chemical Purity:** ≥95% (Desethyl Hydroxychloroquine)  
**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>4</sub>); ≤1% d<sub>0</sub>  
**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-d<sub>4</sub> is intended for use as an internal standard for the quantification of desethyl hydroxychloroquine (Item No. 31152) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Desethyl hydroxychloroquine-d<sub>4</sub> is supplied as a solid. A stock solution may be made by dissolving the desethyl hydroxychloroquine-d<sub>4</sub> in the solvent of choice, which should be purged with an inert gas. Desethyl hydroxychloroquine-d<sub>4</sub> is soluble in methanol, DMSO, and dimethyl formamide.

### Description

Desethyl hydroxychloroquine is an active metabolite of hydroxychloroquine (Item No. 17911).<sup>1</sup> 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, 10/25/2022

### 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