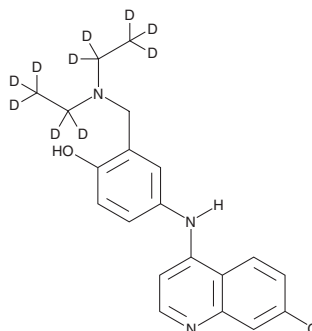


PRODUCT INFORMATION



Amodiaquine-d₁₀ Item No. 28525

CAS Registry No.: 1189449-70-4
Formal Name: 2-((bis(ethyl-d₅)amino)methyl)-4-((7-chloroquinolin-4-yl)amino)phenol
Synonyms: Camoquine-d₁₀, Flavoquine-d₁₀
MF: C₂₀H₁₂ClD₁₀N₃O
FW: 365.9
Chemical Purity: ≥95% (Amodiaquine)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₁₀); ≤1% d₀
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

Amodiaquine-d₁₀ is intended for use as an internal standard for the quantification of amodiaquine (Item No. 15954) 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).

Amodiaquine-d₁₀ is supplied as a solid. A stock solution may be made by dissolving the amodiaquine-d₁₀ in the solvent of choice, which should be purged with an inert gas. Amodiaquine-d₁₀ is soluble in organic solvents such as methanol and chloroform.

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, 12/20/2022

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