

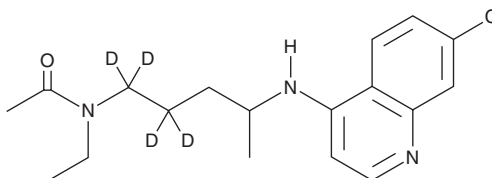
PRODUCT INFORMATION



N-acetyl Desethylchloroquine-d₄

Item No. 30908

CAS Registry No.: 1216619-20-3
Formal Name: N-[4-[(7-chloro-4-quinolinyl)amino]pentyl-1,1,2,2-d₄]-N-ethyl-acetamide
Synonym: N-acetyl-(mono) Desethylchloroquine
MF: C₁₈H₂₀ClD₄N₃O
FW: 337.9
Chemical Purity: ≥95% (N-acetyl Desethylchloroquine)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₄); ≤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

N-acetyl Desethylchloroquine-d₄ is intended for use as an internal standard for the quantification of N-acetyl desethylchloroquine 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).

N-acetyl Desethylchloroquine-d₄ is supplied as a solid. A stock solution may be made by dissolving the N-acetyl desethylchloroquine-d₄ in the solvent of choice, which should be purged with an inert gas. N-acetyl Desethylchloroquine-d₄ is slightly soluble in DMSO, chloroform, and methanol.

Description

N-acetyl Desethylchloroquine is a minor metabolite of chloroquine (Item No. 14194).^{1,2}

References

1. Brown, N.D., Stermer-Cox, M.G., Poon, B.T., *et al.* Separation and identification of a plasma and urinary mono-acetylated conjugate of chloroquine in man by ion-pair high-performance liquid chromatography. *J. Chromatogr.* **309**(2), 426-430 (1984).
2. Kotecka, B.M., Edstein, M.D., and Rieckmann, K.H. Chloroquine bioassay of plasma specimens obtained from soldiers on chloroquine plus doxycycline for malaria prophylaxis. *Int. J. Parasitol.* **26**(11), 1325-1329 (1996).

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, 11/04/2022

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