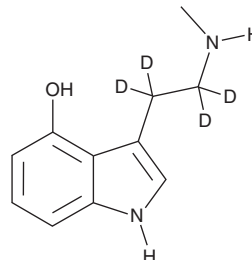


PRODUCT INFORMATION



Norpsilocin-d₄ Item No. 37127

CAS Registry No.: 2742678-87-9
Formal Name: 3-[2-(methylamino)ethyl-1,1,2,2-d₄]-1H-indol-4-ol
Synonyms: ω-N-methyl-4-Hydroxytryptamine-d₄,
4-hydroxy-N-Methyltryptamine-d₄,
4-hydroxy NMT-d₄
MF: C₁₁H₁₀D₄N₂O
FW: 194.3
Chemical Purity: ≥95% (Norpsilocin)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₄); ≤1% d₀
UV/Vis.: λ_{max}: 223, 267, 294 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥3 years



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

Description

Norpsilocin-d₄ (Item No. 37127) is intended for use as an internal standard for the quantification of norpsilocin (Item Nos. 37943 | 23586) 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).

Norpsilocin is categorized as a tryptamine.¹ It has been found in *P. cubensis*, a mushroom with hallucinogenic properties. Norpsilocin is a potential metabolite of baeocystin (Item Nos. 37763 | 31583) based on the published metabolism of psilocybin (Item Nos. 9003134 | 14041).² At the time norpsilocin-d₄ (Item No. 37127) was made available for purchase, specific metabolism data had not been published. Contact us if updated information on this molecule is now available. This product is intended for research and forensic applications.

References

1. Lenz, C., Wick, J., and Hoffmeister, D. Identification of ω-N-methyl-4-hydroxytryptamine (Norpsilocin) as a *Psilocybe* natural product. *J. Nat. Prod.* **80**(10), 2835-2838 (2017).
2. Dinis-Oliveira, R.J. Metabolism of psilocybin and psilocin: Clinical and forensic toxicological relevance. *Drug Metab. Rev.* **49**(1), 84-91 (2017).

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, 07/11/2023

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