

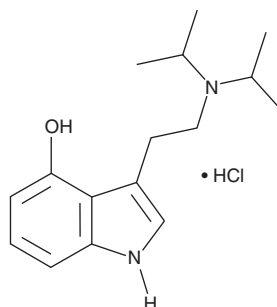
PRODUCT INFORMATION



4-hydroxy DiPT (hydrochloride)

Item No. 11312

CAS Registry No.: 63065-90-7
Formal Name: 3-[2-[bis(1-methylethyl)amino]ethyl]-1H-indol-4-ol, monohydrochloride
Synonym: 4-hydroxy-N,N-Diisopropyltryptamine
MF: C₁₆H₂₄N₂O • HCl
FW: 296.8
Purity: ≥95%
UV/Vis.: λ_{max}: 222, 267, 293 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥3 years



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

Description

4-hydroxy DiPT (hydrochloride) (Item No. 11312) is an analytical reference standard categorized as a tryptamine.¹ 4-hydroxy DiPT induces the head-twitch response (HTR) in mice, indicating hallucinogenic potential, and substitutes for the amphetamine DOM in a two-lever drug discrimination test in rats.^{2,3} This product is intended for research and forensic applications.

References

1. Rickli, A., Moning, O.D., Hoener, M.C., *et al.* Receptor interaction profiles of novel psychoactive tryptamines compared with classic hallucinogens. *Eur. Neuropsychopharmacol.* **26(8)**, 1327-1337 (2016).
2. Klein, A.K., Chatha, M., Laskowski, L.J., *et al.* Investigation of the structure-activity relationships of psilocybin analogues. *ACS Pharmacol. Transl. Sci.* **4(2)**, 533-542 (2020).
3. Gatch, M.B., Hoch, A., and Carbonaro, T.M. Discriminative stimulus effects of substituted tryptamines in rats. *ACS Pharmacol. Transl. Sci.* **4(2)**, 467-471 (2020).

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, 05/21/2024

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