

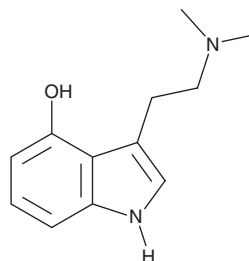
PRODUCT INFORMATION



Psilocin

Item No. 11864

CAS Registry No.: 520-53-6
Formal Name: 3-[2-(dimethylamino)ethyl]-1H-indol-4-ol
Synonyms: 4-hydroxy-N,N-Dimethyltryptamine,
4-hydroxy DMT
MF: C₁₂H₁₆N₂O
FW: 204.3
Purity: ≥95%
UV/Vis.: λ_{max}: 223 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

Psilocin (Item No. 11864) is an analytical reference standard categorized as a tryptamine.¹ Psilocin is the active metabolite of the psychedelic compound psilocybin (Item Nos. 9003134 | 14041) and the synthetic tryptamine 4-acetoxy DMT (psilacetin; Item Nos. 14056 | 35188).^{2,3} It induces the head-twitch response (HTR) in mice, indicating hallucinogenic potential.⁴ Psilocin is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

References

1. Blair, J.B., Kurrasch-Orbaugh, D., Marona-Lewicka, D., *et al.* Effect of ring fluorination on the pharmacology of hallucinogenic tryptamines. *J. Med. Chem.* **43(24)**, 4701-4710 (2000).
2. Halberstadt, A.L., Koedood, L., Powell, S.B., *et al.* Differential contributions of serotonin receptors to the behavioral effects of indoleamine hallucinogens in mice. *J. Psychopharmacol.* **25(11)**, 1548-1561 (2011).
3. Geiger, H.A., Wurst, M.G., and Daniels, R.N. DARK Classics in Chemical Neuroscience: Psilocybin. *ACS Chem. Neurosci.* **9(10)**, 2438-2447 (2018).
4. Glatfelter, G.C., Pottier, E., Partilla, J.S., *et al.* Structure-activity relationships for psilocybin, baeocystin, aeruginascin, and related analogues to produce pharmacological effects in mice. *ACS Pharmacol. Transl. Sci.* **5(11)**, 1181-1196 (2022).

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, 07/24/203

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