

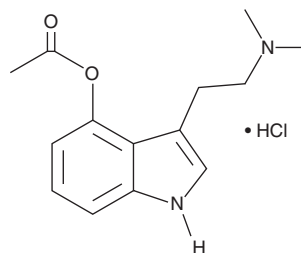
PRODUCT INFORMATION



4-acetoxy DMT (hydrochloride)

Item No. 14056

CAS Registry No.: 2748484-99-1
Formal Name: 3-[2-(dimethylamino)ethyl]-1H-indol-4-ol-4-acetate, monohydrochloride
Synonyms: 4-acetoxy-N,N-Dimethyltryptamine, O-Acetylpsilocin, Psilacetin
MF: C₁₄H₁₈N₂O₂ • HCl
FW: 282.8
Purity: ≥90%
UV/Vis.: λ_{max}: 219, 277 nm
Supplied as: A solution in acetonitrile
Storage: -20°C
Stability: ≥3 years



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

Description

4-acetoxy DMT (hydrochloride) (Item No. 14056) is an analytical reference standard categorized as a tryptamine.^{1,2} 4-acetoxy DMT is a prodrug form of psilocin (Item Nos. 9003135 | 11864 | 36971), a psychedelic hallucinogen derived from various mushrooms. 4-acetoxy DMT induces the head-twitch response (HTR) in mice, indicating hallucinogenic potential.² This product is intended for research and forensic applications.

References

1. Nichols, D.E. and Frescas, S. Improvements to the synthesis of psilocybin and a facile method for preparing the O-acetyl prodrug of psilocin. *Synthesis* **6**, 935-938 (1999).
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).

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, 07/25/2023

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