

PRODUCT INFORMATION



4-acetoxy DMT (hydrochloride)

Item No. 14056

Formal Name: 3-[2-(dimethylamino)ethyl]-1H-indol-4-yl-4-acetate, monohydrochloride

Synonyms: Acetylpsilocin, 4-acetoxy-N,N-Dimethyltryptamine, 4-AcO DMT

MF: C₁₄H₁₈N₂O₂ • HCl

FW: 282.8

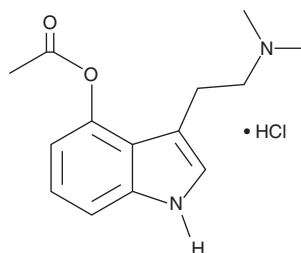
Purity: ≥90%

UV/Vis.: λ_{max}: 219, 277 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥2 years



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

Laboratory Procedures

4-acetoxy DMT (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the 4-acetoxy DMT (hydrochloride) in the solvent of choice. 4-acetoxy DMT (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of 4-acetoxy DMT (hydrochloride) in these solvents is approximately 20, 10, and 30 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 4-AcO DMT (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 4-AcO DMT (hydrochloride) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

4-AcO DMT is an O-acetylated psilocin that is believed to be a synthetic prodrug of psilocin, the psychedelic hallucinogen derived from various mushrooms.¹ Although its pharmacological and toxicological properties are not well known, its activity has been proposed to be somewhat similar to that of psilocybin and psilocin.¹ This product is intended for forensic and research applications.

Reference

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).

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, 05/21/2018

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