

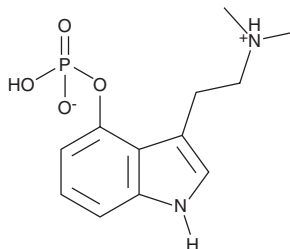
PRODUCT INFORMATION



Psilocybin Analytical Reference Standard

Item No. 39698

CAS Registry No.: 520-52-5
Formal Name: 3-[2-(dimethylamino)ethyl]-1H-indol-4-ol, 4-(dihydrogen phosphate)
MF: C₁₂H₁₇N₂O₄P
FW: 284.3
Purity: ≥99%
Supplied as: A crystalline solid
Storage: 2-8°C
Stability: ≥30 months



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

Description

Cayman Chemical's psilocybin analytical reference standard (Item No. 39698) is a highly purified and characterized crystalline solid ideally suited as a reference material for analytical testing of GMP psilocybin (Item No. 36463). Limit atmospheric exposure and maintain the material in a tightly sealed container. Psilocybin is regulated as a Schedule I compound in the United States.

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, 01/22/2025

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