

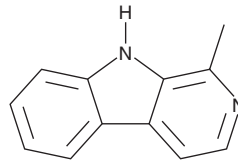
PRODUCT INFORMATION



Harmane

Item No. 33843

CAS Registry No.: 486-84-0
Formal Name: 1-methyl-9H-pyrido[3,4-b]indole
Synonyms: 1-Methyl- β -Carboline, NSC 54439
MF: C₁₂H₁₀N₂
FW: 182.2
Purity: \geq 98%
UV/Vis.: λ_{max} : 214, 235, 289 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 5 years



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

Description

Harmane (Item No. 33843) is an analytical reference standard categorized as a β -carboline alkaloid.¹ Harmane has been found in *Psilocybe* mushrooms. This product is intended for analytical forensic applications. This product is also available as a general research tool (Item No. 29613).

Reference

1. Blei, F., Dörner, S., Fricke, J., *et al.* Simultaneous production of psilocybin and a cocktail of β -carboline monoamine oxidase inhibitors in "magic" mushrooms. *Chemistry* **26(3)**, 729-734 (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, 07/05/2023

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