

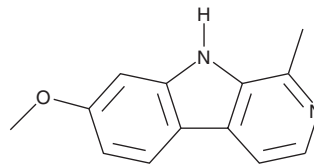
PRODUCT INFORMATION



Harmine

Item No. 35146

CAS Registry No.: 442-51-3
Formal Name: 7-methoxy-1-methyl-9H-pyrido[3,4-b]indole
MF: $C_{13}H_{12}N_2O$
FW: 212.2
Purity: $\geq 98\%$
UV/Vis.: λ_{\max} : 241, 301 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥ 5 years



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

Description

Harmine (Item No. 35146) is an analytical reference standard categorized as an alkaloid and monoamine oxidase inhibitor (MAOI).¹ Harmine is a component of ayahuasca, a botanical mixture that also contains DMT (Item Nos. 13959 | 33586 | 9003568) and has hallucinogenic properties. Harmine reduces certain symptoms of naloxone-induced opioid withdrawal in rats.² This product is intended for analytical forensic applications. This product is also available as a general research tool (Item No. 10010324).

References

1. Brierley, D.I. and Davidson, C. Developments in harmine pharmacology--implications for ayahuasca use and drug-dependence treatment. *Prog. Neuropsychopharmacol. Biol. Psychiatry* **39**(2), 263-272 (2012).
2. Aricioglu-Kartal, F., Kayir, H., and Uzbay, I.T. Effects of harman and harmine on naloxone-precipitated withdrawal syndrome in morphine-dependent rats. *Life Sci.* **73**(18), 2363-2371 (2003).

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, 07/06/2023

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