

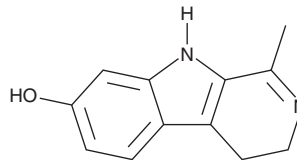
PRODUCT INFORMATION



Harmalol

Item No. 35145

CAS Registry No.: 525-57-5
Formal Name: 4,9-dihydro-1-methyl-3H-pyrido[3,4-b]indol-7-ol
Synonyms: 11-hydroxy Harmalan, Harmidol, NSC 72293
MF: C₁₂H₁₂N₂O
FW: 200.2
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥5 years



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

Description

Harmalol (Item No. 35145) is an analytical reference standard categorized as a β -carboline alkaloid.³ Harmalol is an active metabolite of harmaline (Item No. 10995).¹ It is a minor component of ayahuasca, a botanical mixture that has hallucinogenic properties.² This product is intended for research and forensic applications.

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. Navickiene, S., Santos, L.F.S., Santos, M.C., *et al.* Use of coconut charcoal and menthone-thiosemicarbazone polymer as solid phase materials for the determination of N,N-dimethyltryptamine, harmine, harmaline, harmalol, and tetrahydroharmine in Ayahuasca beverage by liquid chromatography-tandem mass spectrometry. *J. Braz. Chem. Soc.* **30(1)**, 180-187 (2019).
3. Tarpley, M., Oladapo, H.O., Strepay, D., *et al.* Identification of harmine and β -carboline analogs from a high-throughput screen of an approved drug collection; profiling as differential inhibitors of DYRK1A and monoamine oxidase A and for *in vitro* and *in vivo* anti-cancer studies. *Eur. J. Pharm. Sci.* **162**, 105821 (2021).

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