

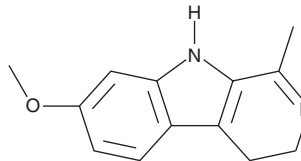
PRODUCT INFORMATION



Harmaline

Item No. 10995

CAS Registry No.: 304-21-2
Formal Name: 4,9-dihydro-7-methoxy-1-methyl-3H-pyrido[3,4-b]indole
MF: C₁₃H₁₄N₂O
FW: 214.3
Purity: ≥98%
UV/Vis.: λ_{max}: 205, 220, 335 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

Harmaline (Item No. 10995) is an analytical reference standard categorized as a β-carboline alkaloid and monoamine oxidase (MAO) inhibitor.^{1,2} Harmaline has been found as an additive in herbal mixtures containing synthetic cannabinoids and described as contributing to poisoning when combined with tryptamines.²⁻⁴ It is a minor component of ayahuasca, a botanical mixture that has hallucinogenic properties.¹ This product is intended for research and forensic applications.

References

1. 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).
2. Liu, C.-H., Chu, W.-L., Liao, S.-C., *et al.* Syrian rue seeds interacted with acacia tree bark in an herbal stew resulted in *N,N*-dimethyltryptamine poisoning. *Clin. Toxicol. (Phila)* **57(10)**, 867-869 (2019).
3. Dresen, S., Ferreirós, N., Pütz, M., *et al.* Monitoring of herbal mixtures potentially containing synthetic cannabinoids as psychoactive compounds. *J. Mass Spectrom.* **45(10)**, 1186-94 (2010).
4. Brush, D.E., Bird, S.B., and Boyer, E.W. Monoamine oxidase inhibitor poisoning resulting from internet misinformation on illicit substances. *J. Toxicol. Clin. Toxicol.* **42(2)**, 191-195 (2004).

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, 06/15/2023

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