

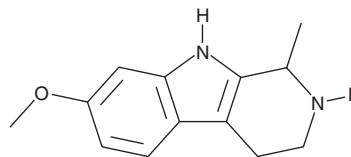
PRODUCT INFORMATION



Tetrahydroharmine

Item No. 35037

CAS Registry No.: 17019-01-1
Formal Name: 2,3,4,9-tetrahydro-7-methoxy-1-methyl-1H-pyrido[3,4-b]indole
Synonyms: Leptaflorine, THH
MF: C₁₃H₁₆N₂O
FW: 216.3
Purity: ≥98%
UV/Vis.: λ_{max}: 224 nm
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

Tetrahydroharmine (Item No. 35037) is an analytical reference standard categorized as an alkaloid and monoamine oxidase inhibitor (MAOI).¹ Tetrahydroharmine is a component of ayahuasca, a botanical mixture that also contains DMT (Item Nos. 13959 | 33586 | 9003568) and has hallucinogenic properties. This product is intended for analytical forensic applications. This product is also available as a general research tool (Item No. 14449).

Reference

1. 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