

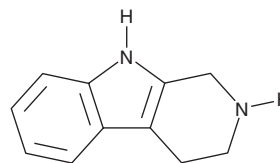
PRODUCT INFORMATION



1,2,3,4-Tetrahydro- β -carboline

Item No. 35038

CAS Registry No.: 16502-01-5
Formal Name: 2,3,4,9-tetrahydro-1H-pyrido[3,4-b]indole
Synonyms: Tetrahydronorharman, Tetrahydro- β -carboline, THBC, Tryptoline
MF: C₁₁H₁₂N₂
FW: 172.2
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 223 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

1,2,3,4-Tetrahydro- β -carboline (Item No. 35038) is an analytical reference standard categorized as a β -carboline alkaloid and a tryptamine metabolite.¹ 1,2,3,4-Tetrahydro- β -carboline is a metabolite of N,N-DMT (Item Nos. 13959 | 9003568 | 33586) and tryptamine (Item No. 33682).^{1,2} It is a minor component of ayahuasca, a botanical mixture that has hallucinogenic properties.¹ This product is intended for analytical forensic applications. This product is also available as a general research tool (Item No. 31220).

References

1. Hamill, J., Hallak, J., Dursun, S.M., et al. Ayahuasca: Psychological and physiologic effects, pharmacology and potential uses in addiction and mental illness. *Curr. Neuropsychopharmacol.* **17**(2), 108-128 (2019).
2. Rommelspacher, H., Bade, P., Coper, H., et al. Inhibition of the reuptake of serotonin by tryptoline. *Naunyn Schmiedeberg's Arch. Pharmacol.* **292**(1), 93-95 (1976).

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