

PRODUCT INFORMATION



7-hydroxy Mitragynine-d₃

Item No. 42292

Formal Name: (αE,2S,3S,7aS,12bS)-3-ethyl-1,2,3,4,6,7,7a,12b-octahydro-7a-hydroxy-8-(methoxy-d₃)-α-(methoxymethylene)-indolo[2,3-a]quinolizine-2-acetic acid, methyl ester

Synonyms: 9-methoxy Corynantheidine
Hydroxyindolenine-d₃, 7-HMG-d₃

MF: C₂₃H₂₇D₃N₂O₅

FW: 417.5

Chemical Purity: ≥98% (7-hydroxy mitragynine)

Deuterium

Incorporation: ≥99% deuterated forms (d₁-d₃); ≤1% d₀

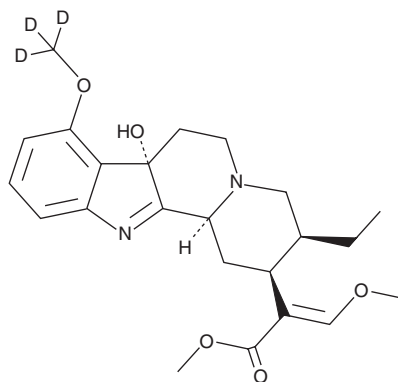
Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Special Conditions: Light sensitive, darkroom conditions, amber glassware

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



Description

7-hydroxy Mitragynine is intended for use as an internal standard for the quantification of 7-hydroxy mitragynine (Item Nos. 35468 | 13114) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

7-hydroxy Mitragynine is categorized as an opioid.¹ 7-hydroxy Mitragynine has been found in *M. speciosa* (Kratom in Thai) and is also an active metabolite of mitragynine (Item Nos. 18567 | 11151).^{2,3} This product is intended for research and forensic applications.

References

1. Takayama, H., Ishikawa, H., Kurihara, M., *et al.* Studies on the synthesis and opioid agonistic activities of mitragynine-related indole alkaloids: Discovery of opioid agonists structurally different from other opioid ligands. *J. Med. Chem.* **45**(9), 1949-1956 (2002).
2. Ponglux, D., Wongseripipatana, S., Takayama, H., *et al.* A new indole alkaloid, 7α-hydroxy-7H-mitragynine, from *Mitragyna speciosa* in thailand. *Planta Med.* **60**(6), 580-581 (1994).
3. Basiliere, S. and Kerrigan, S. Identification of metabolites and potential biomarkers of kratom in urine. *J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.* **1140**, 121971 (2020).

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, 04/06/2026

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