

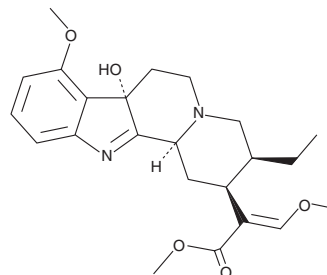
PRODUCT INFORMATION



7-hydroxy Mitragynine

Item No. 13114

CAS Registry No.: 174418-82-7
Formal Name: (αE,2S,3S,7aS,12bS)-3-ethyl-1,2,3,4,6,7,7a,12b-octahydro-7a-hydroxy-8-methoxy-α-(methoxymethylene)-indolo[2,3-a]quinolizine-2-acetic acid, methyl ester
Synonyms: 7-Hydroxymitragynine, 9-methoxy Corynantheidine Hydroxyindolenine
MF: C₂₃H₃₀N₂O₅
FW: 414.5
Purity: ≥98%
UV/Vis.: λ_{max}: 221 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Description

7-hydroxy Mitragynine (Item No. 13114) is an analytical reference standard 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, 06/13/2023

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