

PRODUCT INFORMATION



dihydro-7-hydroxy Mitragynine

Item No. 43910

CAS Registry No.: 1158901-38-2
Formal Name: (αE,2S,3S,7aS,12aR,12bS)-3-ethyl-1,2,3,4,6,7,7a,12,12a,12b-decahydro-7a-hydroxy-8-methoxy-α-(methoxymethylene)-indolo[2,3-a]quinolizine-2-acetic acid, methyl ester

Synonym: MGM-15

MF: C₂₃H₃₂N₂O₅

FW: 416.5

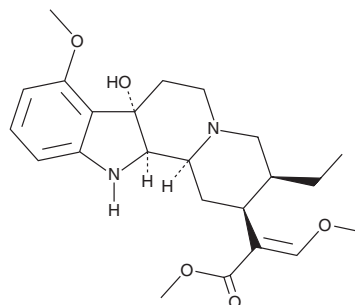
Purity: ≥98%

Supplied as: A crystalline solid

Storage: -80°C

Stability: ≥4 years

Special Conditions: Light sensitive



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

Description

dihydro-7-hydroxy Mitragynine (Item No. 43910) is an analytical reference standard categorized as an opioid.^{1,2} This product is intended for research and forensic applications.

References

1. Matsumoto, K., Narita, M., Muramatsu, N., *et al.* Orally active opioid μ/δ dual agonist MGM-16, a derivative of the indole alkaloid mitragynine, exhibits potent antiallodynic effect on neuropathic pain in mice. *J. Pharmacol. Exp. Ther.* **348(3)**, 383-392 (2014).
2. Gour, A., Mukhopadhyay, S., Henderson, A., *et al.* From kratom to semi-synthetic opioids: The rise and risks of MGM- 15. *Drug Test Anal.* (2025).

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, 11/21/2025

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