

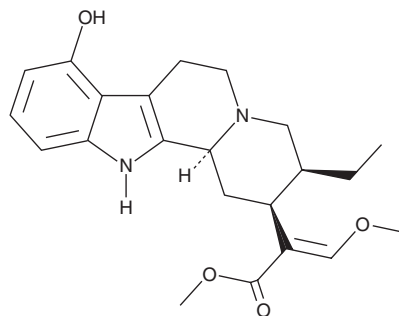
PRODUCT INFORMATION



9-O-desmethyl Mitragynine

Item No. 27540

CAS Registry No.: 425623-55-8
Formal Name: (αE,2S,3S,12bS)-3-ethyl-1,2,3,4,6,7,12,12b-octahydro-8-hydroxy-α-(methoxymethylene)-indolo[2,3-a]quinolizine-2-acetic acid, methyl ester
Synonym: 9-hydroxy Corynantheidine
MF: C₂₂H₂₈N₂O₄
FW: 384.5
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Description

9-O-desmethyl Mitragynine (Item No. 27540) is an analytical reference standard categorized as an opioid.^{1,2} 9-O-Mitragynine is an active metabolite of mitragynine (Item Nos. 18567 | 11151). This product is intended for research and forensic applications.

References

1. 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).
2. 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).

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, 05/05/2026

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