

PRODUCT INFORMATION

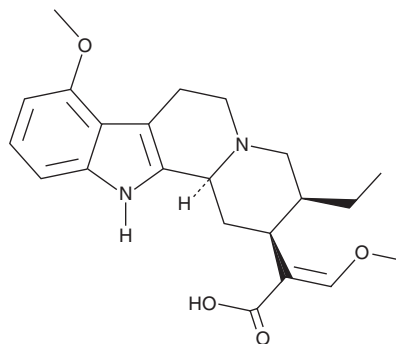


16-carboxy Mitragynine

Item No. 27538

CAS Registry No.: 1224429-90-6
Formal Name: (αE,2S,3S,12bS)-3-ethyl-1,2,3,4,6,7,12,12b-octahydro-8-methoxy-α-(methoxymethylene)-indolo[2,3-a]quinolizine-2-acetic acid

MF: C₂₂H₂₈N₂O₄
FW: 384.5
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥3 years



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

Description

16-carboxy Mitragynine (Item No. 27538) is an analytical reference standard categorized as an opioid.¹ 16-carboxy Mitragynine is a metabolite of mitragynine (Item Nos. 18567 | 11151). This product is intended for research and forensic applications.

Reference

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).

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, 02/26/2024

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