

PRODUCT INFORMATION



10-fluoro dihydro-7-hydroxy Mitragynine

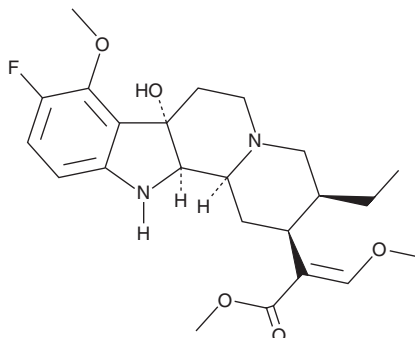
Item No. 43911

CAS Registry No.: 1158901-40-6
Formal Name: 3-ethyl-9-fluoro-1,2S,3S,4,6,7,7aS,12,12aR,12bS-decahydro-7a-hydroxy-8-methoxy- α E-(methoxymethylene)-indolo[2,3-a]quinolizine-2-acetic acid, methyl ester

Synonym: MGM-16
MF: $C_{23}H_{31}FN_2O_5$
FW: 434.5
Purity: $\geq 95\%$
Supplied as: A solid
Storage: -80°C
Stability: ≥ 2 years

Special Conditions: Light sensitive; avoid protic solvents

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



Description

10-fluoro dihydro-7-hydroxy Mitragynine (Item No. 43911) is an analytical reference standard categorized as an opioid.¹ 10-fluoro dihydro-7-hydroxy Mitragynine has antinociceptive and antiallodynic effects in mice. This product is intended for research and forensic applications.

Reference

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).

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, 06/30/2026

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