

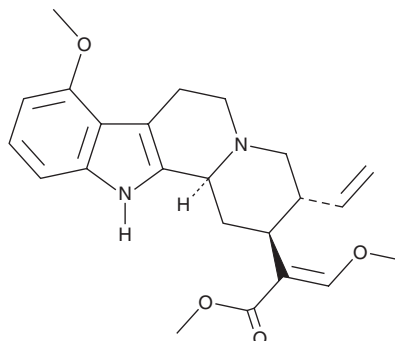
PRODUCT INFORMATION



Paynantheine

Item No. 21841

CAS Registry No.: 4697-66-9
Formal Name: (αE,2S,3R,12bS)-3-ethenyl-1,2,3,4,6,7,12,12b-octahydro-8-methoxy-α-(methoxymethylene)-indolo[2,3-a]quinolizine-2-acetic acid, methyl ester
Synonym: (+)-Paynantheine
MF: C₂₃H₂₈N₂O₄
FW: 396.5
Purity: ≥98%
UV/Vis.: λ_{max}: 226, 292 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

Paynantheine (Item No. 21841) is an analytical reference standard that is structurally similar to known opioids. Paynantheine is an alkaloid found in *M. speciosa* (Kratom in Thai).¹ It has antinociceptive activity and reduces alcohol intake in rats.^{2,3} This product is intended for research and forensic applications.

References

1. Philipp, A.A., Wissenbach, D.K., Weber, A.A., *et al.* Metabolism studies of the kratom alkaloids mitraciliatine and isopaynantheine, diastereomers of the main alkaloids mitragynine and paynantheine, in rat and human urine using liquid chromatography-linear ion trap-mass spectrometry. *J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.* **879**(15-16), 1049-1055 (2011).
2. León, F., Obeng, S., Mottinelli, M., *et al.* Activity of *Mitragyna speciosa* ("Kratom") Alkaloids at Serotonin Receptors. *J. Med. Chem.* **64**(18), 13510-13523 (2021).
3. Gutridge, A.M., Robins, M.T., Cassell, R.J., *et al.* G protein-biased kratom-alkaloids and synthetic carfentanilamide opioids as potential treatments for alcohol use disorder. *Br. J. Pharmacol.* **177**(7), 1497-1513 (2019).

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, 04/10/2025

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