

PRODUCT INFORMATION



Speciogynine

Item No. 25794

CAS Registry No.: 4697-67-0
Formal Name: (αE,2S,3R,12bS)-3-ethyl-1,2,3,4,6,7,12,12b-octahydro-8-methoxy-α-(methoxymethylene)-indolo[2,3-a]quinolizine-2-acetic acid, methyl ester

Synonym: (+)-Speciogynine

MF: C₂₃H₃₀N₂O₄

FW: 398.5

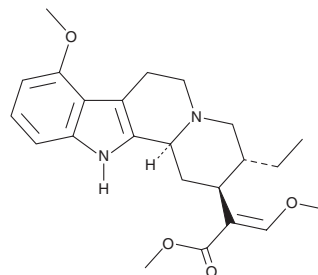
Purity: ≥98%

UV/Vis.: λ_{max}: 226, 293 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥2 years



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

Description

Speciogynine (Item No. 25794) is an analytical reference standard that is structurally similar to known opioids.^{1,2} Speciogynine is an alkaloid found in *M. speciosa* (Kratom in Thai). It has antinociceptive activity and reduces alcohol intake in rats.^{3,4} This product is intended for research and forensic applications.

References

1. Philipp, A.A., Wissenbach, D.K., Weber, A.A., *et al.* Phase I and II metabolites of speciogynine, a diastereomer of the main *Kratom* alkaloid mitragynine, identified in rat and human urine by liquid chromatography coupled to low- and high-resolution linear ion trap mass spectrometry. *J. Mass. Spectrom.* **45(11)**, 1344-1357 (2010).
2. Kruegel, A.C., Gassaway, M.M., Kapoor, A., *et al.* Synthetic and receptor signaling explorations of the *Mitragyna* alkaloids: Mitragynine as an atypical molecular framework for opioid receptor modulators. *J. Am. Chem. Soc.* **138(21)**, 6754-6764 (2016).
3. 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).
4. Guttridge, 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, 06/15/2023

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