

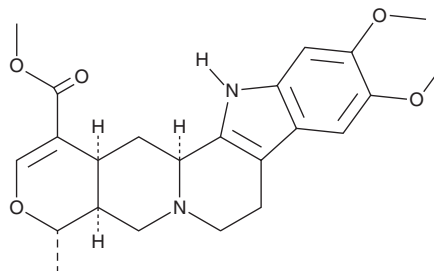
PRODUCT INFORMATION



Isoreserpiline

Item No. 31641

CAS Registry No.: 572-67-8
Formal Name: (19 α ,20 α)-16,17-didehydro-10,11-dimethoxy-19-methyl-oxayohimban-16-carboxylic acid, methyl ester
Synonyms: 3-Isoreserpilinic Acid, methyl ester, Neoreserpiline
MF: C₂₃H₂₈N₂O₅
FW: 412.5
Purity: $\geq 95\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Plant/*Neisosperma poweri*



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

Laboratory Procedures

Isoreserpiline is supplied as a solid. A stock solution may be made by dissolving the isoreserpiline in the solvent of choice, which should be purged with an inert gas. Isoreserpiline is soluble in DMSO.

Description

Isoreserpiline is an indole alkaloid that has been found in *R. tetraphylla* and has diverse biological activities.^{1,2} It has a synergistic effect on the antibacterial activity induced by nalidixic acid (Item No. 19807) in nalidixic acid-sensitive and -resistant strains of *E. coli*.¹ It also reduces amphetamine-induced hyperactivity in mice when administered at a dose of 25 mg/kg.²

References

1. Dwivedi, G.R., Gupta, S., Maurya, A., *et al.* Synergy potential of indole alkaloids and its derivative against drug-resistant *Escherichia coli*. *Chem. Biol. Drug Des.* **86**(6), 1471-1481 (2015).
2. Gupta, S., Khanna, V.K., Maurya, A., *et al.* Bioactivity guided isolation of antipsychotic constituents from the leaves of *Rauwolfia tetraphylla* L. *Fitoterapia* **83**(6), 1092-1099 (2012).

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, 10/25/2022

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