

PRODUCT INFORMATION



Reserpiline

Item No. 31332

CAS Registry No.: 131-02-2
Formal Name: (4S,4aS,13bR,14aS)-4a,5,7,8,13,13b,14,14a-octahydro-10,11-dimethoxy-4-methyl-4H-indolo[2,3-a]pyrano[3,4-g]quinolizine-1-carboxylic acid, methyl ester

MF: C₂₃H₂₈N₂O₅

FW: 412.5

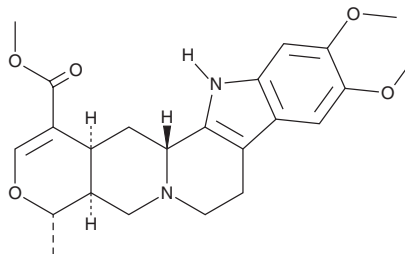
Purity: ≥95%

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

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

Description

Reserpiline is an indole alkaloid that has been found in *R. tetraphylla*.¹ Oral administration of reserpiline (12.5 and 25 mg/kg) decreases amphetamine-induced hyperactivity, indicating antipsychotic potential, in mice.

Reference

1. 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, 11/04/2022

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