

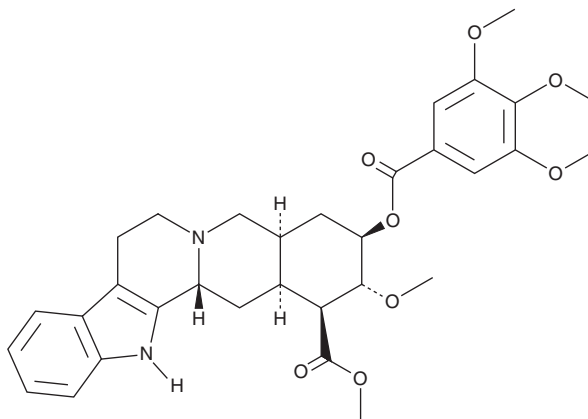
PRODUCT INFORMATION



Deserpidine

Item No. 44222

CAS Registry No.: 131-01-1
Formal Name: (3 β ,16 β ,17 α ,18 β ,20 α)-17-methoxy-18-[(3,4,5-trimethoxybenzoyl)oxy]-yohimban-16-carboxylic acid, methyl ester
Synonyms: Canescine, 11-Desmethoxyreserpine, NSC 72138, Raunormine, Recanescine
MF: C₃₂H₃₈N₂O₈
FW: 578.7
Purity: $\geq 98\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Synthetic



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

Laboratory Procedures

Deserpidine is supplied as a solid. A stock solution may be made by dissolving the deserpidine in the solvent of choice, which should be purged with an inert gas. Deserpidine is soluble (≥ 10 mg/ml) in DMSO.

Description

Deserpidine is an alkaloid that has been found in *R. canescens* and a derivative of reserpine (Item No. 16474) that has diverse biological activities.^{1,2} Deserpidine (5 mg/kg) induces cardiac norepinephrine depletion and sedation in mice.¹ It decreases blood pressure in dogs when administered at a dose of 0.5 mg/kg.² Deserpidine also decreases heart rate and body temperature and induces miosis and diarrhea in rabbits and induces emesis in pigeons.

References

1. Creveling, C.R., Daly, J.W., Parfitt, R.T., *et al.* The chemorelease of norepinephrine from mouse hearts. IV. Structure-activity relationships. Reserpines and yohimbines. *J. Med. Chem.* **11**(3), 596-598 (1968).
2. Schneider, J.A., Plummer, A.J., Earl, A.E., *et al.* Pharmacological studies with deserpidine, a new alkaloid from *Rauwolfia canescens*. *J. Pharmacol. Exp. Ther.* **114**(1), 10-13 (1955).

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, 10/21/2025

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