

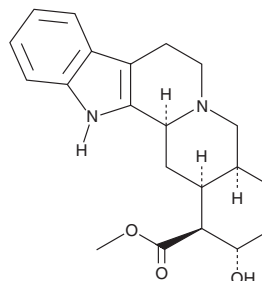
PRODUCT INFORMATION



Rauwolscline

Item No. 46130

CAS Registry No.: 131-03-3
Formal Name: (16 β ,17 α ,20 α)-17-hydroxy-yohimban-16-carboxylic acid, methyl ester
Synonyms: Isoyohimbine, Corynanthidine, α -Yohimbine
MF: C₂₁H₂₆N₂O₃
FW: 354.4
Purity: $\geq 95\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Plant/*Rauwolfia vomitoria*



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

Laboratory Procedures

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

Description

Rauwolscline is a monoterpene indole alkaloid that has been found in *R. tetraphylla* and has cardiovascular activities.¹⁻⁵ It is an antagonist of the α_2 -adrenergic receptor and serotonin (5-HT) receptor subtype 5-HT_{2B} (K_i s = 4 and 14.3 nM, respectively) and a partial agonist of the 5-HT_{1A} receptor (IC_{50} = 1.5 μ M).³⁻⁵ Rauwolscline (1 and 3 μ g/kg) restores electrically induced tachycardia after inhibition by clonidine (Item No. 15949) and inhibits hypertension induced by phenylephrine in anesthetized dogs.²

References

1. Stander, E.A., Lehka, B., Carqueijeiro, I., *et al.* The Rauwolfia tetraphylla genome suggests multiple distinct biosynthetic routes for yohimbane monoterpene indole alkaloids. *Commun. Biol.* **6**(1), 1197 (2023).
2. Shepperson, N.B., Duval, N., Massingham, R., *et al.* Pre- and postsynaptic alpha adrenoceptor selectivity studies with yohimbine and its two diastereoisomers rauwolscline and corynanthine in the anesthetized dog. *J. Pharmacol. Exp. Ther.* **219**(2), 540-546 (1981).
3. Perry, B.D. and U'Prichard, D.C. [³H]rauwolscline (α -yohimbine): A specific antagonist radioligand for brain α_2 -adrenergic receptors. *Eur. J. Pharmacol.* **76**(4), 461-464 (1981).
4. Kaumann, A.J. Yohimbine and rauwolscline inhibit 5-hydroxytryptamine-induced contraction of large coronary arteries of calf through blockade of 5 HT₂ receptors. *Naunyn Schmiedeberg's Arch. Pharmacol.* **323**(2), 149-154 (1983).
5. Wainscott, D.B., Sasso, D.A., Kursar, J.D., *et al.* [³H]Rauwolscline: An antagonist radioligand for the cloned human 5-hydroxytryptamine_{2B} (5-HT_{2B}) receptor. *Naunyn Schmiedeberg's Arch. Pharmacol.* **357**(1), 17-24 (1998).

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, 06/17/2026

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