

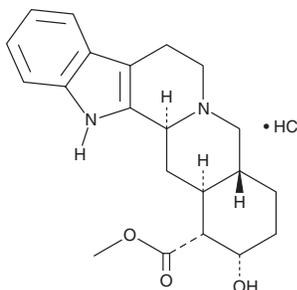
PRODUCT INFORMATION



Yohimbine (hydrochloride)

Item No. 19869

CAS Registry No.: 65-19-0
Formal Name: 17 α -hydroxy-yohimban-16 α -carboxylic acid, methyl ester, monohydrochloride
Synonyms: NIH 9689, NSC 19509, Yohimbe
MF: C₂₁H₂₆N₂O₃ • HCl
FW: 390.9
Purity: \geq 98%
UV/Vis.: λ_{max} : 220, 271, 289 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Description

Yohimbine is a natural alkaloid that has highest affinity for α_2 -adrenergic receptors (pK_i s = 8.2, 8.7, and 9.6 for human α_{2A} , α_{2B} , and α_{2C} , respectively), with antagonist activity at these receptors.^{1,2} It less potently antagonizes serotonin 5-HT₁ receptors (pK_i s = 7.3, 6.8, and 7.6 for human HT_{1A}, HT_{1B}, and HT_{1D}, respectively) and α_1 -adrenergic receptors (pK_i s = 6.7, 6.8, and 6.8 for human α_{1A} , α_{1B} , and α_{1D} , respectively).¹ Yohimbine is used in veterinary medicine to reverse the sedative effects of α_2 -agonists.

References

1. Millan, M.J., Newman-Tancredi, A., Audinot, V., *et al.* Agonist and antagonist actions of yohimbine as compared to fluparoxan at α_2 -adrenergic receptors (AR)s, serotonin (5-HT)_{1A}, 5-HT_{1B}, 5-HT_{1D} and dopamine D₂ and D₃ receptors. Significance for the modulation of frontocortical monoaminergic transmission and depressive states. *Synapse* **35**(2), 79-95 (2000).
2. Schwartz, D.D., and Clark, T.P. Selectivity of atipamezole, yohimbine and tolazoline for alpha-2 adrenergic receptor subtypes: Implications for clinical reversal of alpha-2 adrenergic receptor mediated sedation in sheep. *J. Vet. Pharmacol. Ther.* **21**(5), 342-347 (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, 11/29/2022

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