

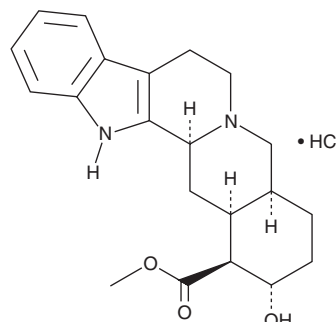
# PRODUCT INFORMATION



## Rauwolscine (hydrochloride)

Item No. 15596

**CAS Registry No.:** 6211-32-1  
**Formal Name:** (16 $\beta$ ,17 $\alpha$ ,20 $\alpha$ )-17-hydroxy-yohimban-16-carboxylic acid, methyl ester, monohydrochloride  
**Synonyms:** Isoyohimbine, NSC 407307,  $\alpha$ -Yohimbine  
**MF:** C<sub>21</sub>H<sub>26</sub>N<sub>2</sub>O<sub>3</sub> • HCl  
**FW:** 390.9  
**Purity:**  $\geq$ 98%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 223, 281 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.

### Laboratory Procedures

Rauwolscine (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the rauwolscine (hydrochloride) in the solvent of choice. Rauwolscine (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of rauwolscine (hydrochloride) in these solvents is approximately 2 and 0.1 mg/ml, respectively.

Rauwolscine (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, rauwolscine (hydrochloride) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Rauwolscine (hydrochloride) has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Rauwolscine is a natural alkaloid that acts as a selective and reversible  $\alpha_2$ -adrenergic receptor antagonist ( $K_i = 12$  nM).<sup>1</sup> It is a stereoisomer of yohimbine, which potently antagonizes both  $\alpha_1$ - and  $\alpha_2$ -adrenergic receptors.<sup>1</sup> Rauwolscine also acts as a receptor antagonist at the serotonin 5-HT<sub>2B</sub> receptor ( $K_i = 14.3$  nM) and as a weak partial agonist at 5-HT<sub>1A</sub> ( $IC_{50} = 1.3$   $\mu$ M).<sup>2-4</sup> The  $\alpha_2$ -adrenergic receptor has diverse physiological functions and antagonists like rauwolscine have numerous applications, including the modulation of mood and behavior.<sup>5,6</sup>

### References

1. Perry, B.D. and U'Prichard, D.C. *Eur. J. Pharmacol.* **76(4)**, 461-464 (1981).
2. Kaumann, A.J. *Naunyn Schmiedebergs Arch. Pharmacol.* **323(2)**, 149-154 (1983).
3. Arthur, J.M., Casañas, S.J., and Raymond, J.R. *Biochem. Pharmacol.* **45(11)**, 2337-2341 (1993).
4. Wainscott, D.B., Sasso, D.A., Kursar, J.D., et al. *Naunyn Schmiedebergs Arch. Pharmacol.* **357(1)**, 17-24 (1998).
5. Brede, M., Philipp, M., Knaus, A., et al. *Biol. Cell* **96(5)**, 343-348 (2004).
6. Cottingham, C. and Wang, Q. *Neurosci. Biobehav. Rev.* **36(10)**, 2214-2225 (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