

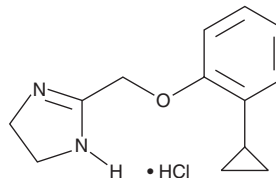
PRODUCT INFORMATION



Cirazoline (hydrochloride)

Item No. 21791

CAS Registry No.: 40600-13-3
Formal Name: 2-[(2-cyclopropylphenoxy)methyl]-4,5-dihydro-1H-imidazole, monohydrochloride
MF: C₁₃H₁₆N₂O • HCl
FW: 252.7
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Cirazoline (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the cirazoline (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Cirazoline (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of cirazoline (hydrochloride) in these solvents is approximately 3, 33, and 16 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of cirazoline (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of cirazoline (hydrochloride) in PBS, pH 7.2, is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Cirazoline is an α_1 -adrenergic receptor (α_1 -AR) agonist (K_{iS} = 120, 960, and 660 nM for recombinant α_{1A} -, α_{1B} -, and α_{1D} -ARs, respectively, in CHO cell membranes).¹ It acts as a full agonist at α_{1A} - and a partial agonist at α_{1B} - and α_{1D} -ARs *in vitro* (EC_{50S} = 70.7, 79.4, and 239.8 nM, respectively). It also acts as an antagonist at α_2 -ARs with a pA_2 value of 7.56 to inhibit the norepinephrine-induced twitch response in isolated pig ileum.² Cirazoline (0.01-1 μ g/kg) decreases blood pressure when administered *via* microinjection to the nucleus reticularis lateralis (NRL) of anesthetized normotensive cats.³ It enhances spatial memory and reduces depressive- and anxiety-like behavior in mice when administered at a concentration of 10 mg/L in drinking water for 2-9 months.⁴ It also enhances performance in the variable delayed response task in aged rhesus monkeys at high doses of 1-10 μ g/kg but impairs performance at lower doses of 0.01-1 μ g/kg.⁵

References

1. Horie, K., Obika, K., Foglar, R., *et al.* *Br. J. Pharmacol.* **116**(1), 1611-1618 (1995).
2. Ruffolo, R.R., Jr. and Waddell, J.E. *Br. J. Pharmacol.* **77**(1), 169-176 (1982).
3. Bousquet, P., Feldman, J., and Schwartz, J. *J. Pharmacol. Exp. Ther.* **230**(1), 232-236 (1984).
4. Doze, V.A., Papay, R.S., Goldenstein, B.L., *et al.* *Mol. Pharmacol.* **80**(4), 747-758 (2011).
5. Arnsten, A.F.T. and Jentsch, J.D. *Pharmacol. Biochem. Behav.* **58**(1), 55-59 (1997).

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, 12/16/2022

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