

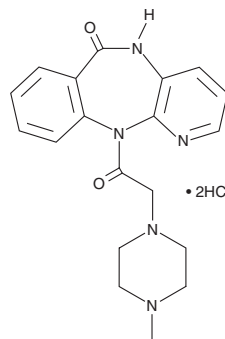
PRODUCT INFORMATION



Pirenzepine (hydrochloride)

Item No. 29527

CAS Registry No.: 29868-97-1
Formal Name: 5,11-dihydro-11-[2-(4-methyl-1-piperazinyl)acetyl]-6H-pyrido[2,3-b][1,4]benzodiazepin-6-one, dihydrochloride
Synonym: Gastrozepin
MF: C₁₉H₂₁N₅O₂ • 2HCl
FW: 424.3
Purity: ≥98%
UV/Vis.: λ_{max}: 280 nm
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

Pirenzepine (hydrochloride) is supplied as a crystalline solid. Aqueous solutions of pirenzepine (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of pirenzepine (hydrochloride) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Pirenzepine is an antagonist of M₁ muscarinic acetylcholine receptors (K_i = 11.48 nM).¹ It is selective for M₁ over M₂, M₃, and M₄ receptors (K_is = 602.56, 151.36, and 199.53 nM, respectively). Pirenzepine inhibits ascending reflex contraction of the circular smooth muscle in isolated guinea pig ileal segments induced by intraluminal balloon inflation (IC₅₀ = 501.19 nM). It inhibits methacholine-induced increases in ileal pressure in guinea pigs (ID₅₀ = 724.44 nmo/kg). Pirenzepine inhibits oxotremorine-induced gastric ulcer, gastric acid secretion, and salivation in rats (ED₅₀s = 13, 37.5, and 620 µg/kg i.v., respectively).² It prevents form-deprivation myopia (FDM) in a chick model of experimental myopia.³

References

1. Doods, H.N., Entzeroth, M., Ziegler, H., *et al.* Pharmacological profile of selective muscarinic receptor antagonists on guinea-pig ileal smooth muscle. *Eur. J. Pharmacol.* **253(3)**, 275-281 (1994).
2. Del Soldato, P. and Pagani, F. Pharmacodynamic evaluation of selective antimuscarinic properties of pirenzepine in the rat. *Pharmacol. Res. Commun.* **14(3)**, 279-287 (1982).
3. Luft, W.A., Ming, Y., and Stell, W.K. Variable effects of previously untested muscarinic receptor antagonists on experimental myopia. *Invest. Ophthalmol. Vis. Sci.* **44(3)**, 1330-1338 (2003).

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