

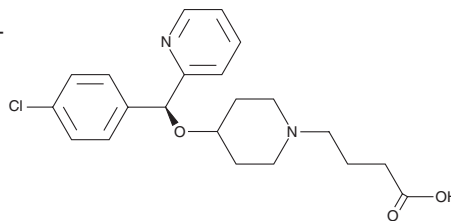
PRODUCT INFORMATION



Bepotastine

Item No. 39932

CAS Registry No.: 125602-71-3
Formal Name: 4-[(S)-(4-chlorophenyl)-2-pyridinylmethoxy]-1-piperidinebutanoic acid
MF: C₂₁H₂₅ClN₂O₃
FW: 388.9
Purity: ≥98%
Supplied as: A 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

Bepotastine is supplied as a solid. A stock solution may be made by dissolving the bepotastine in the solvent of choice, which should be purged with an inert gas. Bepotastine is soluble in organic solvents such as methanol and DMSO. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Bepotastine is an antagonist of the histamine H₁ receptor that is selective over H₃, α₁-, α₂-, and β-adrenergic, dopamine D_{2long}, serotonin 5-HT₂, muscarinic acetylcholine, and benzodiazepine receptors.¹ It reduces dye leakage from the nasal passages of rats acutely sensitized to an antigen (ED₅₀ = 0.03 mg/kg) and inhibits histamine-induced bronchoconstriction in the anesthetized dog (ED₅₀ = 3.2 μg/kg).^{2,3} Bepotastine prevents conjunctival vascular hyperpermeability in a guinea pig model of conjunctivitis in a dose-dependent manner.⁴ Formulations containing bepotastine have been used in the treatment of itching associated with allergic conjunctivitis.

References

1. Kato, M., Nishida, A., Aga, Y., *et al.* Pharmacokinetic and pharmacodynamic evaluation of central effect of the novel antiallergic agent bepotastine besilate. *Arzneimittelforschung*. **47(10)**, 1116-1124 (1997).
2. Matsubara, S., Ono, C., Yamazaki, N., *et al.* Inhibitory effects of bepotastine, a novel anti-allergic drug on histamine-induced bronchoconstriction in anesthetized dogs. *Yakuri to Chiryō* **25(4)**, 895-900 (1997).
3. Murata, T., Matsumoto, Y., Suzuki, T., *et al.* Effect of bepotastine besilate (TAU-284), a novel anti-allergic agent, on experimental allergic rhinitis. *Arerugi* **46(7)**, 576-584 (1997).
4. Kida, T., Fuji, A., Sakai, O., *et al.* Bepotastine besilate, a highly selective histamine H₁ receptor antagonist, suppresses vascular hyperpermeability and eosinophil recruitment in in vitro and in vivo experimental allergic conjunctivitis models. *Exp. Eye Res.* **91(1)**, 85-91 (2010).

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, 03/20/2024

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