

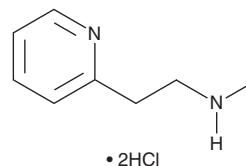
PRODUCT INFORMATION



Betahistine (hydrochloride)

Item No. 27602

CAS Registry No.: 5579-84-0
Formal Name: N-methyl-2-pyridineethanamine, dihydrochloride
MF: C₈H₁₂N₂ • 2HCl
FW: 209.1
Purity: ≥98%
UV/Vis.: λ_{max}: 261, 268 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

Betahistine (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the betahistine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Betahistine (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of betahistine (hydrochloride) in ethanol is approximately 2.5 mg/ml and approximately 2 mg/ml in DMSO and DMF.

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 betahistine (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of betahistine (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

Betahistine is a histamine H₃ receptor antagonist and histamine H₁ receptor agonist.¹ *In vivo*, betahistine dilates pre-capillary arterioles and increases blood flow to the stria vascularis in guinea pigs.^{2,3} Betahistine (8 mg/kg) reduces weight gain and increased feeding behavior induced by olanzapine (Item No. 11937) in female rats.⁴ Formulations containing betahistine have been used in the treatment of balance disorders and vertigo symptoms associated with Meniere's disease.

References

1. Timmerman, H. Histamine agonists and antagonists. *Acta Otolaryngol. Suppl.* **479**, 5-11 (1991).
2. Bertlich, M., Ihler, F., Sharaf, K., *et al.* Betahistine metabolites, aminoethylpyridine, and hydroxyethylpyridine increase cochlear blood flow in guinea pigs *in vivo*. *Int. J. Audiol.* **53(10)**, 753-759 (2014).
3. Bertlich, M., Ihler, F., Weiss, B.G., *et al.* Role of capillary pericytes and precapillary arterioles in the vascular mechanism of betahistine in a guinea pig inner ear model. *Life Sci.* **187**, 17-21 (2017).
4. Lian, J., Huang, X.-F., Pai, N., *et al.* Ameliorating antipsychotic-induced weight gain by betahistine: Mechanisms and clinical implications. *Pharmacol. Res.* **106**, 51-63 (2016).

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, 11/18/2022

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