

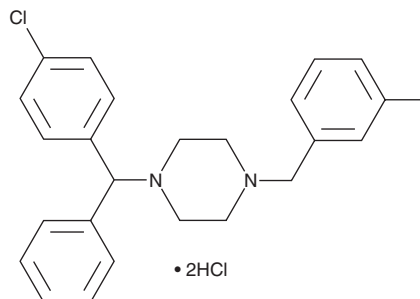
PRODUCT INFORMATION



Meclizine (hydrochloride)

Item No. 14615

CAS Registry No.: 1104-22-9
Formal Name: 1-[(4-chlorophenyl)phenylmethyl]-4-[(3-methylphenyl)methyl]piperazine, dihydrochloride
Synonyms: Meclozine, NSC 28728
MF: C₂₅H₂₇ClN₂ • 2HCl
FW: 463.9
Purity: ≥98%
UV/Vis.: λ_{max}: 229 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

Meclizine (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the Meclizine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Meclizine (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of Meclizine (hydrochloride) in these solvents is approximately 0.11, 0.17, and 5 mg/ml, respectively.

Description

Meclizine is a histamine H₁ receptor antagonist (K_i = 250 nM).¹ Pretreatment with meclizine (2.5-30 mg/kg, i.p.) reduces ovalbumin-induced mast cell degranulation, edema, and leukocyte influx in rat lung.² Formulations containing meclizine have been used in the treatment of motion sickness and vertigo.

References

1. Tran, V.T., Chang, R.S.L., and Snyder, S.H. Histamine H₁ receptors identified in mammalian brain membranes with [³H]mepyramine. *Proc. Natl. Acad. Sci. USA* **75(12)**, 6290-6294 (1978).
2. Pasquale, C.P., Silva, P.M.R.e., Lima, M.C., et al. Suppression by cetirizine of pleurisy triggered by antigen in actively sensitized rats. *Eur. J. Pharmacol.* **223(1)**, 9-14 (1992).

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, 12/12/2022

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