

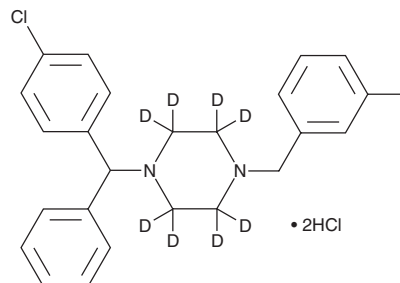
PRODUCT INFORMATION



Meclizine-d₈ (hydrochloride)

Item No. 26444

CAS Registry No.: 1432062-16-2
Formal Name: 1-[(4-chlorophenyl)phenylmethyl]-4-[[3-methylphenyl)methyl]-piperazine-2,2,3,3,5,5,6,6-d₈, dihydrochloride
Synonym: Meclozine-d₈
MF: C₂₅H₁₉ClD₈N₂ • 2HCl
FW: 471.9
Chemical Purity: ≥98% (Meclizine)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₈); ≤1% d₀
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

Meclizine-d₈ (hydrochloride) is intended for use as an internal standard for the quantification of meclizine (Item No. 14615) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated *versus* unlabeled).

Meclizine-d₈ (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the meclizine-d₈ (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Meclizine-d₈ (hydrochloride) is slightly soluble in chloroform and methanol.

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. U.S.A.* **75**(12), 6290-6294 (1978).
2. Pasquale, C.P., e Silva, P.M., 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, 12/22/2022

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