

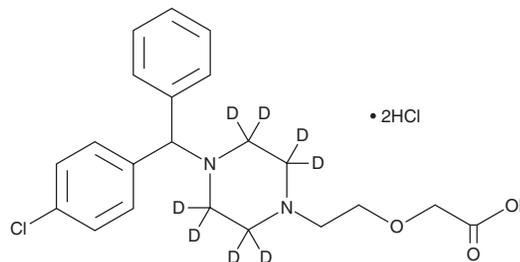
PRODUCT INFORMATION



Cetirizine-d₈ (hydrochloride)

Item No. 26445

CAS Registry No.: 2070015-04-0
Formal Name: 2-[2-[4-[(4-chlorophenyl)phenylmethyl]-1-piperazinyl-d₈]ethoxy]-acetic acid, dihydrochloride
MF: C₂₁H₁₇ClD₈N₂O₃ • 2HCl
FW: 469.9
Chemical Purity: ≥98% (Cetirizine)
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

Cetirizine-d₈ (hydrochloride) is intended for use as an internal standard for the quantification of cetirizine (Item No. 19686) 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).

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

Description

Cetirizine is a bioactive carboxylated metabolite of hydroxyzine (Item No. 24039) that acts as a selective histamine H₁ receptor antagonist (K_i = 10 nM).^{1,2} As a second generation antihistamine, it is non-sedating due to low lipophilicity, which prevents blood-brain barrier transit.³ Cetirizine is a racemic mixture composed of equal amounts of two enantiomers, (R)-cetirizine (Item No. 23992) and (S)-cetirizine, with pharmacological activity residing primarily in the (R) isomer.² Cetirizine inhibits eosinophil chemotaxis and leukotriene B₄ (LTB₄; Item No. 20110) release independent from H₁ antagonism.⁴ It inhibits aerosol histamine-induced bronchospasm in guinea pigs (ED₅₀ = 100 µg/kg, p.o.).⁵ Formulations containing cetirizine have been used in the treatment of allergic rhinitis and chronic urticaria.

References

1. Thurmond, R.L., Gelfand, E.W., and Dunford, P.J. The role of histamine H₁ and H₄ receptors in allergic inflammation: The search for new antihistamines. *Nat. Rev. Drug Discov.* **7**(1), 41-53 (2008).
2. Zhang, L., Cheng, L., and Hong, J. The clinical use of cetirizine in the treatment of allergic rhinitis. *Pharmacology* **92**(1-2), 14-25 (2013).
3. Tillement, J.P. The advantages for an H₁ antihistamine of a low volume of distribution. *Allergy* **55**(suppl 60), 17-21 (2000).
4. Köller, M., Hilger, R.A., Rihoux, J.P., et al. Cetirizine exerts anti-inflammatory effects on human neutrophils. *Int. Arch. Allergy. Immunol.* **110**(1), 52-56 (1996).
5. Llupia, J., Gras, J., and Llenas, J. Comparative anti-allergic effects of second-generation H₁-antihistamines ebastine, cetirizine and loratadine in preclinical models. *Arzneimittelforschung* **53**(2), 93-97 (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, 12/07/2022

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