

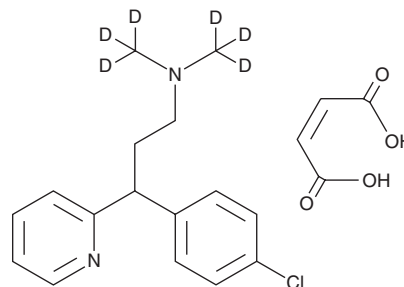
# PRODUCT INFORMATION



## Chlorpheniramine-d<sub>6</sub> (maleate)

Item No. 35767

**CAS Registry No.:** 1219806-45-7  
**Formal Name:** 3-(4-chlorophenyl)-N,N-bis(methyl-d<sub>3</sub>)-3-(pyridin-2-yl)propan-1-amine maleate  
**Synonyms:** (±)-Chlorpheniramine-d<sub>6</sub>, Chlorprophenpyridamine-d<sub>6</sub>  
**MF:** C<sub>16</sub>H<sub>13</sub>ClD<sub>6</sub>N<sub>2</sub> • C<sub>4</sub>H<sub>4</sub>O<sub>4</sub>  
**FW:** 396.9  
**Chemical Purity:** ≥98% (Chlorpheniramine)  
**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>6</sub>); ≤1% d<sub>0</sub>  
**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

Chlorpheniramine-d<sub>6</sub> (maleate) is intended for use as an internal standard for the quantification of chlorpheniramine (Item No. 21253) 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).

Chlorpheniramine-d<sub>6</sub> (maleate) is supplied as a solid. A stock solution may be made by dissolving the chlorpheniramine-d<sub>6</sub> (maleate) in the solvent of choice, which should be purged with an inert gas. Chlorpheniramine-d<sub>6</sub> (maleate) is slightly soluble in methanol.

### Description

Chlorpheniramine is a histamine H<sub>1</sub> receptor antagonist with an IC<sub>75</sub> value of 0.0016 µg/ml for reversal of histamine-induced spasms in isolated guinea pig ileum.<sup>1</sup> It protects against intravenous histamine-induced death (PD<sub>50</sub> = 0.15 mg/kg) and delays induction of aerosolized histamine-induced coughing (ED<sub>100sec</sub> = 0.44 mg/kg) in guinea pigs. Chlorpheniramine (20 mg/kg, i.p.) prevents histamine-induced passive cutaneous anaphylaxis (PCA) in rabbits.<sup>2</sup> It also reduces respiratory resistance and hypersecretion of tracheobronchial fluid in a dog model of histamine-induced asthma.<sup>3</sup> Formulations containing chlorpheniramine have been used in the treatment of seasonal allergies.

### References

1. Lish, P.M., Robbins, S.I., and Peters, E.L. Specificity of antihistamine drugs and involvement of the adrenergic system in histamine deaths in the guinea pig. *J. Pharmacol. Exp. Ther.* **153**(3), 538-543 (1966).
2. Henson, P.M. and Cochrane, C.G. Immunological induction of increased vascular permeability. I. A rabbit passive cutaneous anaphylactic reaction requiring complement, platelets, and neutrophils. *J. Exp. Med.* **129**(1), 153-165 (1969).
3. Yamatake, Y., Sasagawa, S., Yanaura, S., et al. Involvement of histamine H<sub>1</sub>- and H<sub>2</sub>-receptors in induced asthmas in dogs. *Jpn. J. Pharmacol.* **27**(6), 791-797 (1977).

#### 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, 12/13/2022

#### 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