

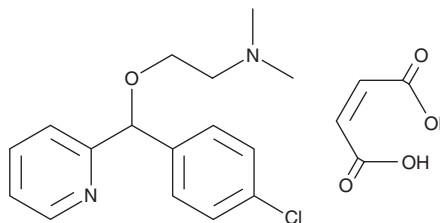
PRODUCT INFORMATION



Carbinoxamine (maleate)

Item No. 23937

CAS Registry No.: 3505-38-2
Formal Name: 2-[(4-chlorophenyl)-2-pyridinylmethoxy]-N,N-dimethylethanamine, (2Z)-2-butenedioate
Synonym: NSC 62362
MF: C₁₆H₁₉ClN₂O • C₄H₄O₄
FW: 406.9
Purity: ≥98%
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

Carbinoxamine (maleate) is supplied as a solid. A stock solution may be made by dissolving the carbinoxamine (maleate) in the solvent of choice, which should be purged with an inert gas. Carbinoxamine (maleate) is soluble in organic solvents such as chloroform and methanol.

Description

Carbinoxamine is a competitive histamine H₁ receptor antagonist (K_i = 2.3 nM) and first generation antihistamine.¹ It also competes with [³H]diltiazem, an L-type calcium channel blocker, for binding to the benzothiazepine site on rat cardiomyocytes (K_i = 1.08 nM).² Carbinoxamine decreases negative inotropic activity in isolated guinea pig left atria by 76% (EC₅₀ = 250 nM) and negative chronotropic activity in guinea pig spontaneously beating isolated right atria (EC₅₀s = 250 and 480 nM, respectively) by 48% compared with control. Formulations containing carbinoxamine have been used in the treatment of allergic rhinitis.

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. Carosati, E., Budriesi, R., Ioan, P., et al. Discovery of novel and cardioselective diltiazem-like calcium channel blockers via virtual screening. *J. Med. Chem.* **51**(18), 5552-5565 (2008).

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, 11/15/2022

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