

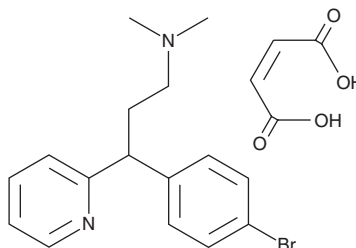
PRODUCT INFORMATION



Brompheniramine (maleate)

Item No. 20996

CAS Registry No.: 980-71-2
Formal Name: γ -(4-bromophenyl)-N,N-dimethyl-2-pyridinepropanamine, 2Z-butenedioate
Synonym: (\pm)-Brompheniramine
MF: C₁₆H₁₉BrN₂ • C₄H₄O₄
FW: 435.3
Purity: \geq 95%
UV/Vis.: λ_{max} : 261 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Brompheniramine (maleate) is supplied as a crystalline solid. A stock solution may be made by dissolving the brompheniramine (maleate) in the solvent of choice. Brompheniramine (maleate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of brompheniramine (maleate) in these solvents is approximately 5, 10, and 15 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of brompheniramine (maleate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of brompheniramine (maleate) in PBS, pH 7.2, is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Brompheniramine is a first generation antihistamine of the propylamine class. It antagonizes histamine H₁ receptors (K_i = 6.06 nM) and also inhibits the reuptake of serotonin and norepinephrine.^{1,2}

References

1. Cusack, B., Nelson, A., and Richelson, E. Binding of antidepressants to human brain receptors: Focus on newer generation compounds. *Psychopharmacology (Berl)* **114(4)**, 559-565 (1994).
2. Högberg, T., Ulf, B., Renyi, A.L., et al. Synthesis of pyridylallylamines related to zimelidine and their inhibition of neuronal monoamine uptake. *J. Med. Chem.* **24(12)**, 1499-1507 (1981).

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, 12/19/2022

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