

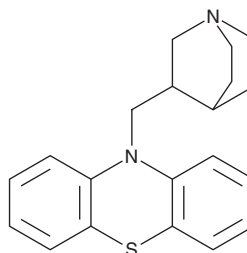
PRODUCT INFORMATION



Mequitazine

Item No. 45325

CAS Registry No.: 29216-28-2
Formal Name: 10-(1-azabicyclo[2.2.2]oct-3-ylmethyl)-10H-phenothiazine
Synonym: NSC 303612
MF: C₂₀H₂₂N₂S
FW: 322.5
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

Mequitazine is supplied as a solid. A stock solution may be made by dissolving the mequitazine in the solvent of choice, which should be purged with an inert gas. Mequitazine is sparingly soluble (1-10 mg/ml) in ethanol and slightly soluble (0.1-1 mg/ml) in DMSO.

Description

Mequitazine is a dual antagonist of histamine H₁ and muscarinic acetylcholine receptors (K_s = 4.4 and 5 nM, respectively).¹ *In vivo*, oral or subcutaneous administration of mequitazine increases survival in a rat model of mast cell degranulation induced by compound 48/80 (Item No. 22173; ED₅₀ = 1.02 and 0.26 mg/kg, respectively).² Formulations containing mequitazine have been used in the treatment of urticaria and allergic rhinitis.

References

1. Kubo, N., Shirakawa, S., Kuno, T., *et al.* Antimuscarinic effects of antihistamines: Quantitative evaluation by receptor-binding assay. *Jpn. J. Pharmacol.* **43(3)**, 277-282 (1987).
2. Neimegeers, C.J.E., Awouters, F.H.L., and Janssen, P.A.J. The *in vivo* pharmacological profile of histamine (H₁) antagonists in the rat. *Drug Dev. Res.* **2**, 559-566 (1982).

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, 03/23/2026

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