

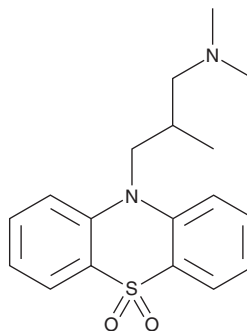
PRODUCT INFORMATION



(±)-Oxomemazine

Item No. 39511

CAS Registry No.: 3689-50-7
Formal Name: N,N,β-trimethyl-10H-phenothiazine-10-propanamine, 5,5-dioxide
Synonym: RP 6847
MF: C₁₈H₂₂N₂O₂S
FW: 330.4
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

(±)-Oxomemazine is supplied as a solid. A stock solution may be made by dissolving the (±)-oxomemazine in the solvent of choice, which should be purged with an inert gas. (±)-Oxomemazine is soluble in organic solvents such as chloroform. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

(±)-Oxomemazine is a phenothiazine antihistamine.¹ It is protective against anaphylactic microshock in guinea pigs when administered at doses ranging from 5 to 10 µg/kg.

Reference

1. Herxheimer, H. and Stresemann, E. The effect of some new antihistamines on the anaphylactic microshock of the guinea-pig. *Br. J. Pharmacol. Chemother.* **21(3)**, 414-418 (1963).

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, 01/26/2024

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