

PRODUCT INFORMATION



Meclizine N-oxide

Item No. 35054

CAS Registry No.: 114624-69-0

Formal Name: 1-[(4-chlorophenyl)phenylmethyl]-4-[(3-methylphenyl)methyl]-piperazine 4-oxide

Synonym: Meclozine N-oxide

MF: $C_{25}H_{27}ClN_2O$

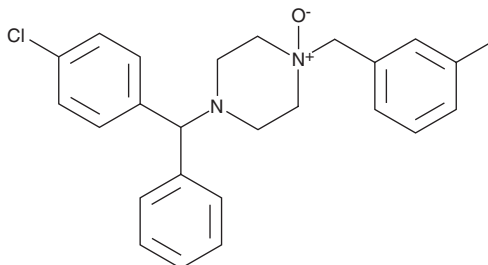
FW: 407.0

Purity: $\geq 90\%$

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

Meclizine N-oxide is supplied as a solid. A stock solution may be made by dissolving the meclizine N-oxide in the solvent of choice, which should be purged with an inert gas. Meclizine N-oxide is soluble in the organic solvent DMSO.

Description

Meclizine N-oxide is a metabolite of the histamine H_1 receptor antagonist meclizine (Item No. 14615).¹ It is also a degradation product and potential impurity in commercial preparations of meclizine.²

References

1. Goenechea, V.S., Rücker, G., Brzezinka, H., *et al.* Biotransformation of meclizine in the human body. *J. Clin. Chem. Clin. Biochem.* **26**(2), 105-115 (1988).
2. Peraman, R., Manikala, M., Kondreddy, V.K., *et al.* A stability-indicating RP-HPLC method for the quantitative analysis of meclizine hydrochloride in tablet dosage form. *J. Chromatogr. Sci.* **53**(5), 793-799 (2015).

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, 11/19/2024

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