

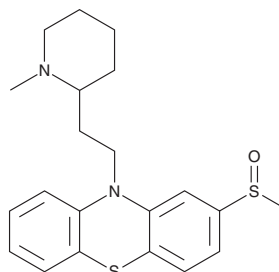
PRODUCT INFORMATION



Mesoridazine

Item No. 31370

CAS Registry No.: 5588-33-0
Formal Name: 10-[2-(1-methyl-2-piperidyl)ethyl]-2-(methylsulfinyl)-10H-phenothiazine
Synonym: NSC 186066
MF: C₂₁H₂₆N₂OS₂
FW: 386.6
Purity: ≥95%
Supplied as: A solution in acetonitrile
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Mesoridazine is supplied as a solution in acetonitrile. To change the solvent, simply evaporate the acetonitrile under a gentle stream of nitrogen and immediately add the solvent of choice. Mesoridazine is slightly soluble in DMSO (warmed) and methanol.

Description

Mesoridazine is an active metabolite of the typical antipsychotic thioridazine (Item No. 14400).^{1,2} It is formed via sulfoxidation of thioridazine by the cytochrome P450 (CYP) isoform CYP2D6 in human liver microsomes.³ It binds to histamine H₁, dopamine D₂, muscarinic acetylcholine, and α₁- and α₂-adrenergic receptors (K_ds = 1.8, 19, 69, 2, and 1,600 nM, respectively).² Mesoridazine (10, 30, and 45 mg/kg, i.p.) prevents mescaline-induced hyperactivity in mice.⁴

References

1. von Bahr, C., Movin, G., Nordin, C., *et al.* Plasma levels of thioridazine and metabolites are influenced by the debrisoquin hydroxylation phenotype. *Clin. Pharmacol. Ther.* **49(3)**, 234-240 (1991).
2. Richelson, E. and Nelson, A. Antagonism by neuroleptics of neurotransmitter receptors of normal human brain in vitro. *Eur. J. Pharmacol.* **103(3-4)**, 197-204 (1984).
3. Wójcikowski, J., Maurel, P., and Daniel, W.A. Characterization of human cytochrome P450 enzymes involved in the metabolism of the piperidine-type phenothiazine neuroleptic thioridazine. *Drug Metab. Dispos.* **34(3)**, 471-476 (2006).
4. Shah, N.S. Influence of psychotropic drugs and β-diethylaminoethyl-diphenylpropylacetate (SKF 525-A) on mescaline-induced behavior and on tissue levels of mescaline in mice. *Biochem. Pharmacol.* **25(5)**, 591-597 (1976).

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, 11/08/2022

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