

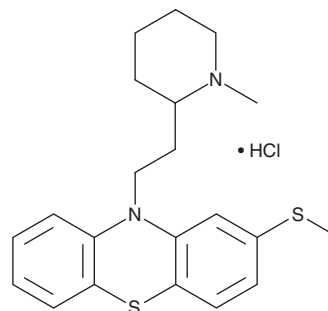
PRODUCT INFORMATION



Thioridazine (hydrochloride)

Item No. 14400

CAS Registry No.: 130-61-0
Formal Name: 10-[2-(1-methyl-2-piperidinyl)ethyl]-2-(methylthio)-10H-phenothiazine, monohydrochloride
Synonyms: Aldazine, NSC 186060
MF: C₂₁H₂₆N₂S₂ • HCl
FW: 407.0
Purity: ≥98%
UV/Vis.: λ_{max}: 230, 263, 316 nm
Supplied as: A crystalline 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

Thioridazine (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the thioridazine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Thioridazine (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of thioridazine (hydrochloride) in ethanol is approximately 10 mg/ml and approximately 25 mg/ml in DMSO and DMF.

Description

Thioridazine is a typical antipsychotic.¹ It binds to dopamine D₂, histamine H₁, M₃ muscarinic, and α₁- and α₂-adrenergic receptors (K_is = 5-341.3 nM), as well as the serotonin (5-HT) receptor subtypes 5-HT_{1A}, 5-HT_{2A}, 5-HT_{2C}, 5-HT₆, and 5-HT₇ (K_is = 10-180.7 nM). Thioridazine (5 mg/kg) reduces amphetamine-induced repetitive head bobbing and oral behavior in rats.² It reduces conditioned fear stress-induced freezing behavior in rats when administered at doses ranging from 3 to 100 mg/kg.³ Thioridazine is also active against multidrug-resistant tuberculosis *in vitro* and *in vivo*.⁴

References

1. Kroeze, W.K., Hufeisen, S.J., Popadak, B.A., *et al.* H1-histamine receptor affinity predicts short-term weight gain for typical and atypical antipsychotic drugs. *Neuropsychopharmacology* **28**(3), 519-526 (2003).
2. Tschanz, J.T. and Rebec, G.V. Atypical antipsychotic drugs block selective components of amphetamine-induced stereotypy. *Pharmacol. Biochem. Behav.* **31**(3), 519-522 (1988).
3. Ishida-Tokuda, K., Ohno, Y., Sakamoto, H., *et al.* Evaluation of perospirone (SM-9018), a novel serotonin-2 and dopamine-2 receptor antagonist, and other antipsychotics in the conditioned fear stress-induced freezing behavior model in rats. *Jpn. J. Pharmacol.* **72**(2), 119-126 (1996).
4. Amaral, L. Thioridazine: An old neuroleptic effective against totally drug resistant tuberculosis. *Acta Med. Port.* **25**(2), 118-121 (2012).

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, 11/15/2022

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