

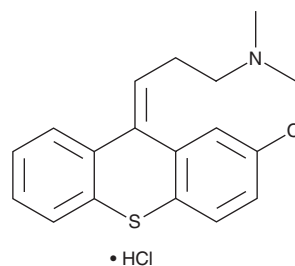
PRODUCT INFORMATION



Chlorprothixene (hydrochloride)

Item No. 20772

CAS Registry No.: 6469-93-8
Formal Name: (3Z)-3-(2-chloro-9H-thioxanthen-9-ylidene)-N,N-dimethyl-1-propanamine, monohydrochloride
Synonyms: cis-Chlorprothixene, NSC 169899
MF: C₁₈H₁₈ClNS • HCl
FW: 352.3
Purity: ≥98%
UV/Vis.: λ_{max}: 230, 270, 329 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

Chlorprothixene (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the chlorprothixene (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Chlorprothixene (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of chlorprothixene (hydrochloride) in these solvents is approximately 30 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of chlorprothixene (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of chlorprothixene (hydrochloride) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Chlorprothixene is a thioxanthine antipsychotic that functions by antagonizing dopamine D₂ receptors.¹ It can block a subset of GABA_A receptors in rat cortex that is also blocked by clozapine (Item No. 12059).² It has also been shown to be effective against *P. falciparum* growth with an EC₅₀ value of 1.7 μM.³

References

1. Froimowitz, M., and Cody, V. Biologically active conformers of phenothiazines and thioxanthenes. Further evidence for a ligand model of dopamine D₂ receptor antagonists. *J. Med. Chem.* **36(15)**, 2219-2227 (1993).
2. Squires, R.F., and Saederup, E. Clozapine and several other antipsychotic/antidepressant drugs preferentially block the same 'core' fraction of GABA_A receptors. *Neurochem. Res.* **23(10)**, 1283-1290 (1998).
3. Weisman, J.L., Liou, A.P., Shelat, A.A., et al. Searching for new antimalarial therapeutics amongst known drugs. *Chem. Biol. Drug Des.* **67(6)**, 409-416 (2006).

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, 12/01/2022

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