

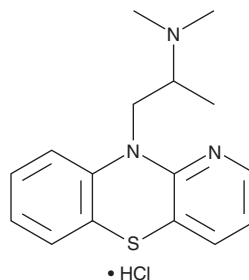
PRODUCT INFORMATION



Isothipendyl (hydrochloride)

Item No. 33582

CAS Registry No.: 1225-60-1
Formal Name: N,N,α-trimethyl-10H-pyrido[3,2-b]
[1,4]benzothiazine-10-ethanamine,
monohydrochloride
Synonym: NSC 169186
MF: C₁₆H₁₉N₃S • HCl
FW: 321.9
Purity: ≥98%
UV/Vis.: λ_{max}: 250 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

Isothipendyl is supplied as a crystalline solid. A stock solution may be made by dissolving the isothipendyl in the solvent of choice, which should be purged with an inert gas. Isothipendyl (hydrochloride) is soluble in the organic solvent DMSO.

Description

Isothipendyl (hydrochloride) is an antihistamine.¹ It protects against seizures induced by supramaximal electroshock (MES) in rats (ED₅₀ = 57.9 mg/kg). Formulations containing isothipendyl (hydrochloride) have been used in the treatment of itching associated with insect bites.

Reference

1. Saxena, V.C., Bapat, S.K., and Dhawan, B.N. An experimental evaluation of the anticonvulsant activity of some antihistaminic drugs. *Jpn. J. Pharmacol.* **19(4)**, 477-484 (1969).

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, 12/05/2022

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