

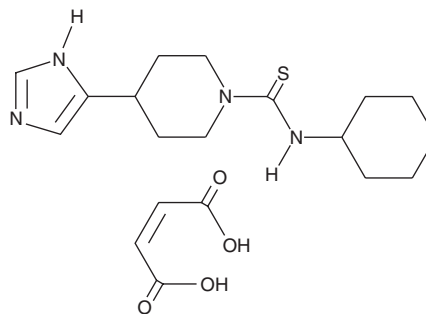
PRODUCT INFORMATION



Thioperamide (maleate)

Item No. 10011127

CAS Registry No.: 148440-81-7
Formal Name: N-cyclohexyl-4-(1H-imidazol-5-yl)-1-piperidinecarbothioamide, 2Z-butenedioate
MF: C₁₅H₂₄N₄S • C₄H₄O₄
FW: 408.5
Purity: ≥98%
UV/Vis.: λ_{max}: 249 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

Thioperamide (maleate) is supplied as a crystalline solid. A stock solution may be made by dissolving the thioperamide (maleate) in an organic solvent purged with an inert gas. Thioperamide (maleate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of thioperamide (maleate) in these solvents is approximately 10, 25, and 30 mg/ml, respectively.

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 thioperamide (maleate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of thioperamide (maleate) in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Thioperamide (maleate) is a selective histamine H₃ receptor antagonist that crosses the blood-brain barrier. This compound binds to rat cerebral cortical cells *in vitro* with a pK_i value of 8.4 and inhibits histamine binding *in vivo* with an ED₅₀ value of 1 mg/kg.¹ At a dose of 5 mg/kg thioperamide (maleate) inhibits kindled seizures in rats by decreasing histamine and gamma-aminobutyric acid.²

References

1. Meier, G., Apelt, J., Reichert, U., *et al.* Influence of imidazole replacement in different structural classes of histamine H₃-receptor antagonists. *Eur. J. Pharmac. Sci.* **13(3)**, 249-259 (2001).
2. Harada, C., Fujii, Y., Hirai, T., *et al.* Inhibitory effect of iodophenprofit, a selective histamine H₃ antagonist, on amygdaloid kindled seizures. *Brain Res. Bull.* **63(2)**, 143-146 (2004).

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, 01/31/2024

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