

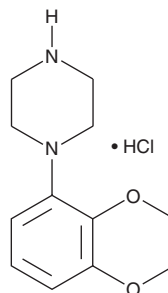
PRODUCT INFORMATION



Etoprazine (hydrochloride)

Item No. 18428

CAS Registry No.: 98206-09-8
Formal Name: 1-(2,3-dihydro-1,4-benzodioxin-5-yl)-piperazine, monohydrochloride
Synonyms: DU-28,853, DU-28,893
MF: C₁₂H₁₆N₂O₂ • HCl
FW: 256.7
Purity: ≥98%
UV/Vis.: λ_{max}: 216 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

Etoprazine (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the etoprazine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Etoprazine (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of etoprazine (hydrochloride) in these solvents is approximately 25, 20, 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 etoprazine (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of etoprazine (hydrochloride) in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Etoprazine is a phenylpiperazine compound that acts as a partial agonist at serotonin 5-HT_{1A}, 5-HT_{1B}, and 5-HT_{2B} receptors (K_is = 40, 52, and 81 nM, respectively).¹ It exerts a dose-dependent decrease in aggressive behavior in resident-intruder tests with rats (ID₅₀ = 0.24 mg/kg).²

References

- Schipper, J., Tulp, M.T., and Sijbesma, H. Neurochemical profile of etoprazine. *Drug Metabol. Drug Interact.* **8(102)**, 85-114 (1990).
- de Boer, S.F., Lesourd, M., Mocaer, E., et al. Selective antiaggressive effects of alnespirone in resident-intruder test are mediated via 5-hydroxytryptamine1A receptors: A comparative pharmacological study with 8-hydroxy-2-dipropylaminotetralin, ipsapirone, buspirone, etoprazine, and WAY-100635. *J. Pharmacol. Exp. Ther.* **288(3)**, 1125-1133 (1999).

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, 12/14/2022

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