

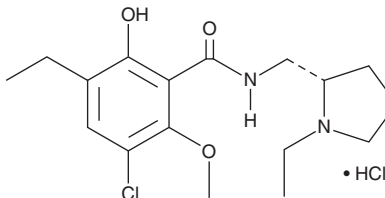
PRODUCT INFORMATION



Eticlopride (hydrochloride)

Item No. 29112

CAS Registry No.: 97612-24-3
Formal Name: 3-chloro-5-ethyl-N-[[[(2S)-1-ethyl-2-pyrrolidinyl]methyl]-6-hydroxy-2-methoxybenzamide, monohydrochloride
Synonyms: S-(-)-Eticlopride, (-)-Eticlopride
MF: C₁₇H₂₅ClN₂O₃ • HCl
FW: 377.3
Purity: ≥98%
UV/Vis.: λ_{max}: 213 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Eticlopride (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the eticlopride (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Eticlopride (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of eticlopride (hydrochloride) in ethanol is approximately 30 mg/ml and approximately 25 mg/ml in DMSO and DMF.

Description

Eticlopride is a dopamine D₂ and D₃ receptor antagonist (K_is = 0.029 and 0.46 nM in MN9D mouse neuronal cells expressing rat D₂ and human D₃ receptors, respectively).¹ It is selective for dopamine D₂ and D₃ receptors (IC₅₀s = 1 and 113 nM, respectively) over dopamine D₁, α₂-adrenergic, and β-adrenergic, histamine H₁, and muscarinic receptors (IC₅₀s = 700->100,000 nM), as well as the serotonin (5-HT) receptor subtypes 5-HT₁ and 5-HT₂ (IC₅₀s = 6,200 and 830, respectively), but does bind α₁-adrenergic receptors (α₁-ARs; IC₅₀ = 110 nM) in radioligand binding assays.² Eticlopride (10 μg/kg) inhibits stereotyped behavior in rats induced by 7-hydroxy-N,N-di-n-propyl-2-aminotetralin (7-OH-DPAT).³ It also inhibits ketamine- and cocaine-induced hypermotility in rats when administered at doses of 20 and 50 μg/kg, respectively.

References

1. Tang, L., Todd, R.D., Heller, A., *et al.* Pharmacological and functional characterization of D₂, D₃ and D₄ dopamine receptors in fibroblast and dopaminergic cell lines. *J. Pharmacol. Exp. Ther.* **268**(1), 495-502 (1994).
2. Hall, H., Köhler, C., and Gawell, L. Some in vitro receptor binding properties of [³H]eticlopride, a novel substituted benzamide, selective for dopamine-D₂ receptors in the rat brain. *Eur. J. Pharmacol.* **111**(2), 191-199 (1985).
3. Giuliani, D., and Ferrari, F. Involvement of dopamine receptors in the antipsychotic profile of (-) eticlopride. *Physiol. Behav.* **61**(4), 563-567 (1997).

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, 12/03/2019

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