

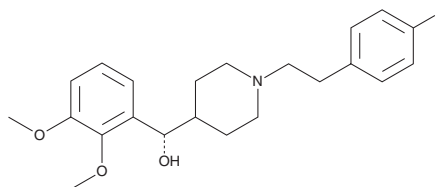
PRODUCT INFORMATION



(-)-Volinanserin

Item No. 42955

CAS Registry No.: 175673-57-1
Formal Name: α S-(2,3-dimethoxyphenyl)-1-[2-(4-fluorophenyl)ethyl]-4-piperidinemethanol
Synonyms: (-)-MDL 100907, (-)-M100907, (S)-MDL 100907
MF: C₂₂H₂₈FNO₃
FW: 373.5
Purity: ≥98%
Supplied as: A 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

(-)-Volinanserin is supplied as a solid. A stock solution may be made by dissolving the (-)-volinanserin in the solvent of choice, which should be purged with an inert gas. (-)-Volinanserin is soluble (≥10 mg/ml) in ethanol.

Description

(-)-Volinanserin is an antagonist of the serotonin (5-HT) receptor subtype 5-HT_{2A} (IC₅₀ = 15.6 nM).¹ It is selective for 5-HT_{2A} over the 5-HT_{2C}, dopamine D₂, and α_1 -adrenergic receptors (IC₅₀s = 750, >10,000, and 1,000 nM, respectively)

Reference

1. Heinrich, T., Böttcher, H., Prücher, H., *et al.* 1-(1-Phenethylpiperidin-4-yl)-1-phenylethanols as potent and highly selective 5-HT_{2A} antagonists. *ChemMedChem* **1**(2), 245-255 (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, 04/01/2025

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