PRODUCT INFORMATION



Opipramol

Item No. 29471

CAS Registry No.: Formal Name:	315-72-0 4-[3-(5H-dibenz[b,f]azepin-5-yl) propyl]-1-piperazineethanol	
Synonyms:	G-33040, GR-33040, NSC 169867	
MF:	C ₂₃ H ₂₉ N ₃ O	
FW:	363.5	
Purity:	≥98%	
UV/Vis.:	λ _{max} : 215, 257 nm	Ň
Supplied as:	A crystalline solid	
Storage:	-20°C	ОН
Stability:	≥4 years	

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

Laboratory Procedures

Opipramol is supplied as a crystalline solid. A stock solution may be made by dissolving the opipramol in the solvent of choice, which should be purged with an inert gas. Opipramol is slightly soluble in acetonitrile, chloroform, and methanol.

Description

Opipramol is a sigma-1 (σ_1) and σ_2 receptor ligand.¹ It binds to σ_1 and σ_2 receptors in guinea pig brain membrane preparations (IC₅₀s = 7 and 56 nM, respectively). It is selective for σ_1 and σ_2 over histamine H₂, dopamine D₁, α_1 - and α_2 -adrenergic, and muscarinic M₁ receptors (IC₅₀s = 4,300, 900, 200, 6,100, and 3,300 nM, respectively) and has no effect on serotonin (5-HT) or norepinephrine uptake ($IC_{50}s = >10,000$ nM for both), but does bind to histamine H_1 and dopamine D_2 receptors ($IC_{50}s = 12$ and 120 nM, respectively), as well as the 5-HT receptor subtype 5-HT₂ ($IC_{50} = 120$ nM). Opipramol (0.01 mg/kg) increases social interaction time in a social exploration test in rats, indicating anxiolytic-like activity. It also decreases immobility time in the forced swim test in rats, indicating antidepressant-like activity, when administered at a dose of 10 mg/kg.

Reference

1. Müller, W.E., Siebert, B., Holoubek, G., et al. Neuropharmacology of the anxiolytic drug opipramol, a sigma site ligand. Pharmacopsychiatry 37(Suppl 3), S189-S197 (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

uyer 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, 12/09/2022

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