

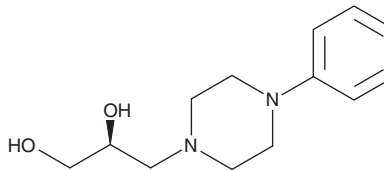
PRODUCT INFORMATION



Levodropropizine

Item No. 32954

CAS Registry No.: 99291-25-5
Formal Name: (2S)-3-(4-phenyl-1-piperazinyl)-1,2-propanediol
Synonym: (-)-(S)-Dropropizine
MF: C₁₃H₂₀N₂O₂
FW: 236.3
Purity: ≥95%
UV/Vis.: λ_{max}: 248 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

Levodropropizine is supplied as a crystalline solid. A stock solution may be made by dissolving the levodropropizine in the solvent of choice, which should be purged with an inert gas. Levodropropizine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of levodropropizine in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

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 levodropropizine can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of levodropropizine in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Levodropropizine is the levo isomer of the antitussive agent dropropizine (Item No. 31110) and a peripherally acting antitussive agent.¹ It reduces the frequency of citric acid-induced cough in guinea pigs when administered at doses of 10 and 20 mg/kg.²

References

1. Lavezzo, A., Melillo, G., Clavenna, G., *et al.* Peripheral site of action of levodropropizine in experimentally-induced cough: Role of sensory neuropeptides. *Pulm. Pharmacol.* **5(2)**, 143-147 (1992).
2. Clay, E.H., Patacchini, R., Trevisani, M., *et al.* Ozone-induced hypertussive responses in rabbits and guinea pigs. *J. Pharmacol. Exp. Ther.* **357(1)**, 73-83 (2016).

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, 11/10/2022

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