

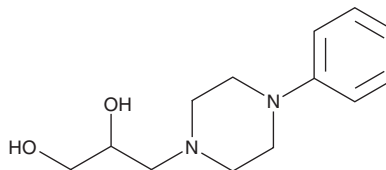
PRODUCT INFORMATION



Dropropizine

Item No. 31110

CAS Registry No.: 17692-31-8
Formal Name: 3-(4-phenyl-1-piperazinyl)-1,2-propanediol
Synonym: (±)-Dropropizine
MF: $C_{13}H_{20}N_2O_2$
FW: 236.3
Purity: $\geq 95\%$
UV/Vis.: λ_{\max} : 247 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

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

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 dropropizine can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of dropropizine in PBS, pH 7.2, is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Dropropizine is an antitussive agent.^{1,2} It reduces the number and intensity of mechanical irritation-induced coughs in conscious cats when administered at doses of 100 and 30 mg/kg, respectively.

References

1. Nosál'ová, G., Mokry, J., and Hassan, K.M.T. Antitussive activity of the fruit extract of *Emblca officinalis* Gaertn. (Euphorbiaceae). *Phytomedicine* **10(6-7)**, 583-589 (2003).
2. Franova, S. and Nosalova, G. Antitussive effect of diltiazem in experimental conditions. *Bratisl. Lek. Listy* **105(5-6)**, 203-206 (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

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, 12/09/2022

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