

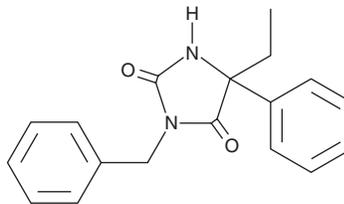
PRODUCT INFORMATION



(±)-N-3-Benzylirvanol

Item No. 27298

CAS Registry No.: 93879-40-4
Formal Name: 5-ethyl-5-phenyl-3-(phenylmethyl)-
2,4-imidazolidinedione
MF: C₁₈H₁₈N₂O₂
FW: 294.4
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

(±)-N-3-Benzylirvanol is supplied as a solid. A stock solution may be made by dissolving the (±)-N-3-benzylirvanol in the solvent of choice, which should be purged with an inert gas. (±)-N-3-Benzylirvanol is slightly soluble in DMSO and methanol.

Description

(±)-N-3-Benzylirvanol is a racemic mixture of (+)-N-3-benzylirvanol and (-)-N-3-benzylirvanol.^{1,2} (+)-N-3-benzylirvanol and (-)-N-3-benzylirvanol are inhibitors of the cytochrome P450 (CYP) isoform CYP2C19 (K_is = 0.25 and 5.3 μM, respectively).¹ (±)-N-3-Benzylirvanol (20 mg/kg, i.p.) lowers plasma levels of cholesterol in mice.²

References

1. Suzuki, H., Kneller, M.B., Haining, R.L., *et al.* (+)-N-3-Benzyl-nirvanol and (-)-N-3-benzyl-phenobarbital: New potent and selective in vitro inhibitors of CYP2C19. *Drug Metab. Dispos.* **30(3)**, 235-239 (2002).
2. Maguire, J.H., Murthy, A.R., and Hall, I.H. Hypolipidemic activity of antiepileptic 5-phenylhydantoin in mice. *Eur. J. Pharmacol.* **117(1)**, 135-138 (1985).

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, 11/16/2022

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