

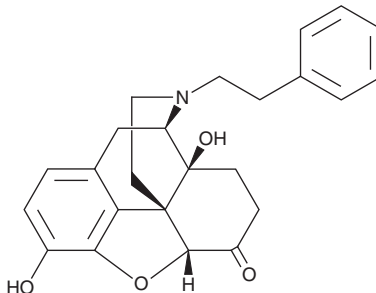
PRODUCT INFORMATION



N-Phenethylnoroxymorphone

Item No. 35677

CAS Registry No.: 4778-94-3
Formal Name: 4,5 α -epoxy-3,14-dihydroxy-17-(2-phenylethyl)-morphinan-6-one
Synonym: (+)-Phenethylnoroxymorphone
MF: C₂₄H₂₅NO₄
FW: 391.5
Purity: $\geq$ 98%
Supplied as: A neat solid
Storage: -20°C
Stability: $\geq$ 4 years



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

Description

N-Phenethylnoroxymorphone (Item No. 35677) is an analytical reference standard categorized as an opioid.¹ N-Phenethylnoroxymorphone potentiates morphine-induced analgesia in rats.² This product is intended for research and forensic applications.

References

1. Ben Haddou, T., Béni, S., Hosztafi, S., *et al.* Pharmacological investigations of N-substituent variation in morphine and oxymorphone: Opioid receptor binding, signaling and antinociceptive activity. *PLoS One* **9(6):e99231**, (2014).
2. Selfridge, B.R., Wang, X., Zhang, Y., *et al.* Structure-activity relationships of (+)-naltrexone-inspired toll-like receptor 4 (TLR4) antagonists. *J. Med. Chem.* **58(12)**, 5038-5052 (2015).

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, 02/14/2024

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