

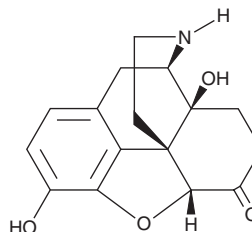
PRODUCT INFORMATION



Noroxymorphone

Item No. 15903

CAS Registry No.: 33522-95-1
Formal Name: 4,5 α -epoxy-3,14-dihydroxy-morphinan-6-one
Synonyms: EN 3169, (-)-Noroxymorphone
MF: C₁₆H₁₇NO₄
FW: 287.3
Purity: $\geq$ 98%
UV/Vis.: λ_{max} : 283 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: $\geq$ 5 years



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

Description

Noroxymorphone (Item No. 15903) is an analytical reference standard categorized as an opioid.¹ Noroxymorphone is an active metabolite of oxycodone (Item Nos. ISO60148 | 15465), noroxycodone (Item Nos. 35521 | 27217), and oxymorphone (Item Nos. ISO60171 | 15911 | 29435).¹⁻³ It is a precursor in the synthesis of naltrexone (Item Nos. ISO60192 | 15520) and naloxone (Item Nos. ISO60191 | 15594).⁴ Noroxymorphone is regulated as a Schedule II compound in the United States. This product is intended for research and forensic applications.

References

1. Thompson, C.M., Wojno, H., Greiner, E., *et al.* Activation of G-proteins by morphine and codeine congeners: Insights to the relevance of O- and N-demethylated metabolites at μ - and δ -opioid receptors. *J. Pharmacol. Exp. Ther.* **308**(2), 547-554 (2004).
2. Kokki, M., Heikkinen, M., Valitalo, P., *et al.* Maturation of oxycodone pharmacokinetics in neonates and infants: Oxycodone and its metabolites in plasma and urine. *Br. J. Clin. Pharmacol.* (2016).
3. Korjamo, T., Tolonen, A., Ranta, V.-P., *et al.* Metabolism of oxycodone in human hepatocytes from different age groups and prediction of hepatic plasma clearance. *Front. Pharmacol.* **82**(87), 1-7 (2012).
4. Lalovic, B., Phillips, B., Risler, L.L., *et al.* Quantitative contribution of CYP2D6 and CYP3A to oxycodone metabolism in human liver and intestinal microsomes. *Drug Metab. Dispos.* **32**(4), 447-454 (2004).
5. Gutmann, B., Weigl, U., Cox, D.P., *et al.* Batch- and continuous-flow aerobic oxidation of 14-hydroxy opioids to 1,3-oxazolidines-A concise synthesis of noroxymorphone. *Chemistry* **22**(30), 10393-10398 (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, 03/20/2025

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