

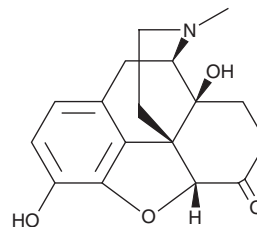
PRODUCT INFORMATION



Oxymorphone

Item No. 15911

CAS Registry No.: 76-41-5
Formal Name: (5 α)-4,5-epoxy-3,14-dihydroxy-17-methyl-morphinan-6-one
Synonyms: 14-Hydroxydihydromorphinone, NIH 10323, NSC 19045
MF: C₁₇H₁₉NO₄
FW: 301.3
Purity: \geq 98%
Supplied as: A neat 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

Oxymorphone (Item No. 15911) is an analytical reference material categorized as an opioid.^{1,2} Oxymorphone is a metabolite of oxycodone (Item Nos. ISO60148 | 15465 | 26513) that has been abused recreationally.³ Oxymorphone is regulated as a Schedule II compound in the United States. This product is intended for research and forensic applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards.

References

1. Fang, W.B., Lofwall, M.R., Walsh, S.L., *et al.* Determination of oxycodone, noroxycodone and oxymorphone by high-performance liquid chromatography-electrospray ionization-tandem mass spectrometry in human matrices: *In vivo* and *in vitro* applications. *J. Anal. Toxicol.* **37(6)**, 337-344 (2013).
2. 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)**, (2014).
3. Babalonis, S., Lofwall, M.R., Nuzzo, P.A., *et al.* Pharmacodynamic effects of oral oxymorphone: Abuse liability, analgesic profile and direct physiological effects in human. *Addict. Biol.* **21(1)**, 146-158 (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, 01/23/2020

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