

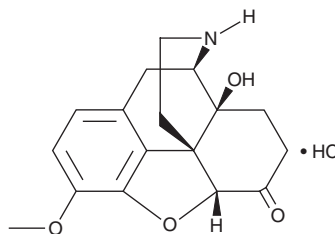
PRODUCT INFORMATION



Noroxycodone (hydrochloride)

Item No. 27217

CAS Registry No.: 52446-25-0
Formal Name: 4,5 α -epoxy-14-hydroxy-3-methoxy-morphinan-6-one, monohydrochloride
MF: C₁₇H₁₉NO₄ • HCl
FW: 337.8
Purity: \geq 98%
Supplied as: A neat solid
Storage: -20°C
Stability: \geq 6 years



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

Description

Noroxycodone (hydrochloride) (Item No. 27217) is an analytical reference material categorized as an opioid.^{1,2} Noroxycodone is a metabolite of oxycodone (Item Nos. ISO60148 | 15465).^{3,4} 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 for reference materials.

References

1. Leow, K.P. and Smith, M.T. The antinociceptive potencies of oxycodone, noroxycodone and morphine after intracerebroventricular administration to rats. *Life Sci.* **54(17)**, 1229-1236 (1994).
2. Lalovic, B., Kharasch, E., Hoffer, C., *et al.* Pharmacokinetics and pharmacodynamics of oral oxycodone in healthy human subjects: Role of circulating active metabolites. *Clin. Pharmacol. Ther.* **79(5)**, 461-479 (2006).
3. Romand, S., Spaggiari, D., Marsousi, N., *et al.* Characterization of oxycodone in vitro metabolism by human cytochromes P450 and UDP-glucuronosyltransferases. *J. Pharm. Biomed. Anal.* **144**, 129-137 (2017).
4. 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).

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, 04/29/2019

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