

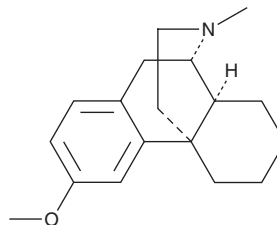
PRODUCT INFORMATION



Dextromethorphan

Item No. 11157

CAS Registry No.: 125-71-3
Formal Name: (9 α ,13 α ,14 α)-3-methoxy-17-methyl-morphinan
Synonyms: DXM, D-Methorphan
MF: C₁₈H₂₅NO
FW: 271.4
Purity: \geq 98%
UV/Vis.: λ_{max} : 280 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

Dextromethorphan (Item No. 11157) is an analytical reference standard categorized as an antitussive.¹ Dextromethorphan has analgesic and anticonvulsant activities in rodents.^{2,3} It has been found as an adulterant in 3,4-MDMA and has potential for abuse.^{4,5} This product is intended for research and forensic applications.

References

1. Bolser, D.C., Aziz, S.M., DeGennaro, F.C., *et al.* Antitussive effects of GABA_B agonists in the cat and guinea-pig. *Br. J. Pharmacol.* **110(1)**, 491-495 (1993).
2. Chen, Y.W., Chu, K.S., Lin, C.N., *et al.* Dextromethorphan or dextrorphan have a local anesthetic effect on infiltrative cutaneous analgesia in rats. *Anesth. Analg.* **104(5)**, 1251-1255 (2007).
3. Jamali, H. and Heydari, A. Effect of dextromethorphan/quinidine on pentylenetetrazole- induced clonic and tonic seizure thresholds in mice. *Neurosci. Lett.* **729**, 134988 (2020).
4. Cole, C., Jones, L., McVeigh, J., *et al.* CUT: A guide to adulterants, bulking agents and other contaminants found in illicit drugs. Liverpool John Moores University (2010).
5. Chomchai, C. and Manaboriboon, B. Stimulant methamphetamine and dextromethorphan use among Thai adolescents: Implications for health of women and children. *J. Med. Toxicol.* **8(3)**, 291-294 (2012).

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, 03/01/2024

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