

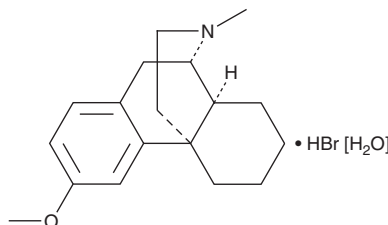
# PRODUCT INFORMATION



## Dextromethorphan (hydrobromide hydrate)

Item No. 13950

**CAS Registry No.:** 6700-34-1  
**Formal Name:** (9 $\alpha$ ,13 $\alpha$ ,14 $\alpha$ )-3-methoxy-17-methylmorphinan, hydrobromide monohydrate  
**Synonyms:** DXM, D-Methorphan  
**MF:** C<sub>18</sub>H<sub>25</sub>NO • HBr [H<sub>2</sub>O]  
**FW:** 370.3  
**Purity:**  $\geq$ 98%  
**UV/Vis.:**  $\lambda_{\max}$ : 227, 280 nm  
**Supplied as:** A crystalline 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

Dextromethorphan (hydrobromide hydrate) (Item No. 13950) is an analytical reference standard categorized as an antitussive.<sup>1</sup> Dextromethorphan has analgesic and anticonvulsant activities in rodents.<sup>2,3</sup> It has been found as an adulterant in 3,4-MDMA and has potential for abuse.<sup>4,5</sup> 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<sub>B</sub> 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 pentylentetrazole- 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