

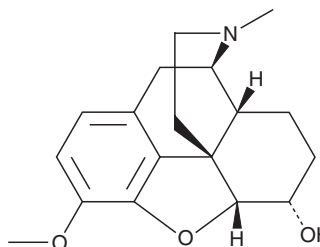
# PRODUCT INFORMATION



## Dihydrocodeine

Item No. 15460

**CAS Registry No.:** 125-28-0  
**Formal Name:** 4,5 $\alpha$ -epoxy-3-methoxy-17-methyl-morphinan-6 $\alpha$ -ol  
**Synonyms:** DH-Codeine, Dihydrin, Drocode, 6 $\alpha$ -Hydrocodol, NSC 231319, Paracodine  
**MF:** C<sub>18</sub>H<sub>23</sub>NO<sub>3</sub>  
**FW:** 301.4  
**Purity:**  $\geq$ 98%  
**Supplied as:** A neat solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 9 years



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

### Description

Dihydrocodeine (Item No. 15460) is an analytical reference material categorized as an opioid.<sup>1</sup> Dihydrocodeine is an active metabolite of hydrocodone (Item Nos. ISO60144 | 15461 | 22166).<sup>2</sup> Formulations containing dihydrocodeine have been used in the treatment of severe pain and have been associated with misuse and polypharmacy-related overdose.<sup>3</sup> Dihydrocodeine 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 for reference materials.

### 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  $\mu$ - and  $\delta$ -opioid receptors. *J. Pharmacol. Exp. Ther.* **308**(2), 547-554 (2004).
2. Cone, E.J., Darwin, W.D., Gorodetzky, C.W., *et al.* Comparative metabolism of hydrocodone in man, rat, guinea pig, rabbit, and dog. *Drug Metab. Dispos.* **6**(4), 488-493 (1978).
3. Rock, K.L., Reynolds, L.M., Rees, P., *et al.* Highlighting the hidden dangers of a 'weak' opioid: Deaths following use of dihydrocodeine in England (2001-2020). *Drug Alcohol Depend.* **233**, 109376 (2022).

#### 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, 11/08/2022

#### 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