

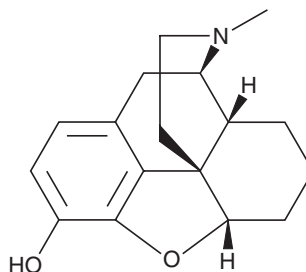
# PRODUCT INFORMATION



## Desomorphine (exempt preparation)

Item No. 15714

**CAS Registry No.:** 427-00-9  
**Formal Name:** 4,5 $\alpha$ -epoxy-17-methyl-morphinan-3-ol  
**Synonyms:** Dihydrodesoxymorphine, Krokodil, Permonid  
**MF:** C<sub>17</sub>H<sub>21</sub>NO<sub>2</sub>  
**FW:** 271.4  
**Purity:**  $\geq 98\%$   
**Supplied as:** A 1 mg/ml solution in methanol  
**Storage:** -20°C  
**Stability:**  $\geq 3$  years



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

### Description

Desomorphine is a morphine derivative in which the 6-hydroxyl group and the 7,8 double bond have been reduced. Compared to morphine, it demonstrates similar opioid-like effects with rapid onset, 10 times greater potency, and short duration of action.<sup>1</sup> This product is intended for forensic and research applications.

### Reference

1. Su, C.-J., Srimurugan, S., Chen, C., et al. Sol-gel titania-coated needles for solid phase dynamic extraction-GC/MS analysis of desomorphine and desocodeine. *Anal. Sci.* **27(11)**, 1107-1113 (2011).

#### 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, 05/19/2025

#### 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](http://WWW.CAYMANCHEM.COM)