

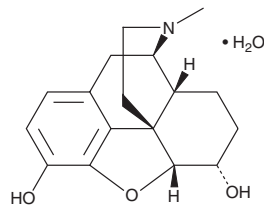
PRODUCT INFORMATION



Dihydromorphine (hydrate)

Item No. 21536

CAS Registry No.: 65943-40-0
Formal Name: (5 α ,6 α)-4,5-epoxy-17-methyl-morphinan-3,6-diol, monohydrate
Synonym: 6 α -Hydromorphol
MF: C₁₇H₂₁NO₃ • H₂O
FW: 305.4
Purity: $\geq$ 98%
Supplied as: A neat 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

Dihydromorphine (hydrate) (Item No. 21536) is an analytical reference material categorized as an opioid metabolite.¹ It is an active metabolite of hydromorphone (Item No. ISO60146).² Dihydromorphine (hydrate) is regulated as a Schedule I 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.

References

1. Pfeiffer, A. and Herz, A. Different types of opiate agonists interact distinguishably with mu, delta and kappa opiate binding sites. *Life Sci.* **31**(12-13), 1355-1358 (1982).
2. Bogusz, M.J., Maier, R.D., Krüger, K.D., *et al.* Determination of common drugs of abuse in body fluids using one isolation procedure and liquid chromatography--atmospheric-pressure chemical-ionization mass spectrometry. *J. Anal. Toxicol.* **22**(7), 549-558 (1998).

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, 07/16/2024

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