

PRODUCT INFORMATION



Codeine (phosphate hydrate)

Item No. 26514

CAS Registry No.: 41444-62-6
Formal Name: (5 α ,6 α)-7,8-didehydro-4,5-epoxy-3-methoxy-17-methyl-morphinan-6-ol, phosphate, hydrate (2:2:1)
Synonyms: Codicept, Coducept, Methyilmorphine, Morphine 3-methyl ether

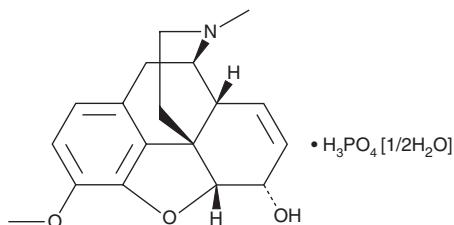
MF: C₁₈H₂₁NO₃ • H₃PO₄ [1/2H₂O]
FW: 406.4

Purity: ≥98%
UV/Vis.: λ_{max}: 212 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥5 years



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

Description

Codeine (phosphate hydrate) (Item No. 26514) is an analytical reference standard categorized as an opioid.¹ Codeine is an analgesic and formulations containing it have commonly been abused.² Codeine is regulated as a Schedule II compound in the United States. This product is intended for research and forensic applications.

References

1. Codd, E.E., Shank, R.P., Schupsky, J.J., *et al.* Serotonin and norepinephrine uptake inhibiting activity of centrally acting analgesics: Structural determinants and role in antinociception. *J. Pharmacol. Exp. Ther.* **274**(3), 1263-1270 (1995).
2. Cone, E.J., Darwin, W.D., and Buchwald, W.F. Assay for codeine, morphine and ten potential urinary metabolites by gas chromatography--mass fragmentography. *J. Chromatogr.* **275**(2), 307-318 (1983).

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, 07/07/2023

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