

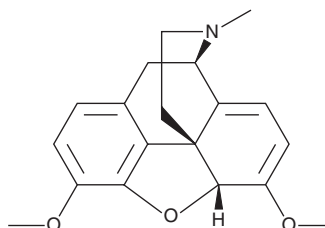
PRODUCT INFORMATION



Thebaine

Item No. 36540

CAS Registry No.: 115-37-7
Formal Name: 6,7,8,14-tetrahydro-4,5 α -epoxy-3,6-dimethoxy-17-methyl-morphinan
Synonyms: Paramorphine, (-)-Thebaine
MF: C₁₉H₂₁NO₃
FW: 311.4
Purity: $\geq$ 98%
Supplied as: A crystalline 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

Thebaine (Item No. 36540) is an analytical reference standard categorized as a precursor in the synthesis of morphine (Item Nos. ISO60147 | 15464) and codeine (Item Nos. ISO60140 | 15459).¹ Thebaine is a potential impurity in the illicit synthesis of heroin (Item Nos. ISO60187 | 9001543 | 9003076).² Thebaine is regulated as a Schedule II compound in the United States. This product is intended for research and forensic applications.

References

1. Dekeyser, S., van Tomme, M., Boone, B., *et al.* Occupational allergic contact dermatitis caused by thebaine and oripavine. *Contact Dermatitis* **76**(2), 120-122 (2017).
2. Schwemer, T., Rössler, T., Ahrens, B., *et al.* Characterization of a heroin manufacturing process based on acidic extracts by combining complementary information from two-dimensional gas chromatography and high resolution mass spectrometry. *Forensic Chem.* **4**, 9-18 (2017).

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, 07/12/2023

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