

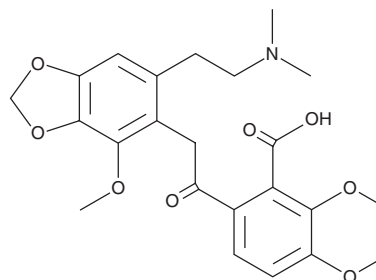
# PRODUCT INFORMATION



## Narceine

Item No. 45067

**CAS Registry No.:** 131-28-2  
**Formal Name:** 6-[2-[6-[2-(dimethylamino)ethyl]-4-methoxy-1,3-benzodioxol-5-yl]acetyl]-2,3-dimethoxy-benzoic acid  
**MF:** C<sub>23</sub>H<sub>27</sub>NO<sub>8</sub>  
**FW:** 445.5  
**Purity:** ≥98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Description

Narceine (Item No. 45067) is an analytical reference standard categorized as an alkaloid. Narceine has been found in *P. somniferum* and as an impurity in seized heroin samples.<sup>1,2</sup> This product is intended for research and forensic applications.

### References

1. C.C., Presley, and Lindsley, C.W. DARK Classics in chemical neuroscience: Opium, a historical perspective. *ACS Chem. Neurosci.* **9**(10), 2503-2518 (2018).
2. Cui, X., Lian, R., Chen, J., *et al.* Source identification of heroin by rapid detection of organic impurities using direct analysis in real time with high-resolution mass spectrometry and multivariate statistical analysis. *Microchem. J.* **147**, 121-126 (2019).

#### 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, 03/04/2026

#### 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