

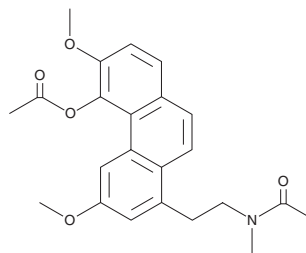
PRODUCT INFORMATION



ATM4 4-acetoxy analog

Item No. 15498

CAS Registry No.: 91295-74-8
Formal Name: N-[2-[5-(acetyloxy)-3,6-dimethoxy-1-phenanthrenyl]ethyl]-N-methyl-acetamide
MF: C₂₃H₂₅NO₅
FW: 395.5
Purity: ≥98%
UV/Vis.: λ_{max}: 259, 284, 303, 315 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

ATM4 4-acetoxy analog (Item No. 15498) is an analytical reference standard categorized as a potential impurity in the synthesis of heroin (Item Nos. ISO60187 | 9001543).¹ ATM4 4-acetoxy analog has been found in seized heroin samples. This product is intended for research and forensic applications.

Reference

1. 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, 06/20/2023

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