

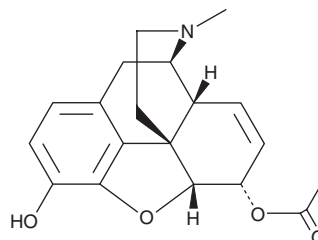
PRODUCT INFORMATION



6-Acetylmorphine

Item No. 15880

CAS Registry No.: 2784-73-8
Formal Name: (5 α ,6 α)-7,8-didehydro-4,5-epoxy-17-methyl-morphinan-3,6-diol 6-acetate
Synonyms: 6-AM, 6-MAM, 6-Monoacetylmorphine
MF: C₁₉H₂₁NO₄
FW: 327.4
Purity: \geq 98%
Supplied as: A neat solid
Storage: -20°C
Stability: \geq 2 years



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

Description

6-Acetylmorphine (Item No. 15880) is an analytical reference material categorized as an opioid metabolite.¹ It is a major metabolite of heroin (Item Nos. 9001543 | 9003076 | ISO60187). 6-Acetylmorphine is regulated as a Schedule II 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 for reference materials.

Reference

1. Zuccaro, P., Ricciarello, R., Pichini, S., *et al.* Simultaneous determination of heroin 6-monoacetylmorphine, morphine, and its glucuronides by liquid chromatography-atmospheric pressure ionspray-mass spectrometry. *J. Anal. Toxicol.* **21(4)**, 268-277 (1997).

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, 04/27/2022

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