

PRODUCT INFORMATION



3-Acetylmorphine

Item No. 9001954

CAS Registry No.: 5140-28-3

Formal Name: (4R,7S,12bS)-7-hydroxy-3-methyl-2,3,4,4aR,7,7aR-hexahydro-1H-4,12-methanobenzofuro[3,2-e]isoquinolin-9-yl acetate

Synonyms: 3-MAM, O³-Monoacetylmorphine

MF: C₁₉H₂₁NO₄

FW: 327.4

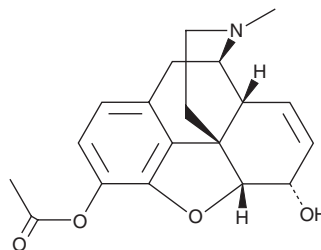
Purity: ≥98%

UV/Vis.: λ_{max}: 278 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥4 years



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

Description

3-Acetylmorphine (3-MAM) is a less active metabolite of heroin (Item No. 9001543) compared to the more active 6-MAM (Item Nos. 19418 | 15880) and morphine (Item No. ISO60147).^{1,2} 3-MAM has relatively weak affinity for μ-opioid receptors due to the placement of the acetyl group in the 3 position. This product is intended only for forensic and research purposes.

References

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).
2. Cone, E.J., Jufer, R., Darwin, W.D., *et al.* Forensic drug testing for opiates. VII. Urinary excretion profile of intranasal (snorted) heroin. *J. Anal. Toxicol.* **20(6)**, 379-392 (1996).

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, 07/25/2023

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