

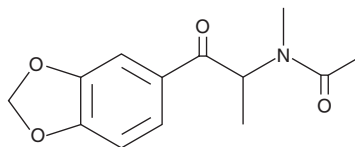
PRODUCT INFORMATION



N-acetyl-3,4-Methylenedioxymethcathinone

Item No. 9001492

CAS Registry No.: 1227293-15-3
Formal Name: N-[2-(1,3-benzodioxol-5-yl)-1-methyl-2-oxoethyl]-N-methyl-acetamide
Synonyms: N-acetyl-3,4-MDMC, N-acetyl Methylone
MF: C₁₃H₁₅NO₄
FW: 249.3
Purity: ≥95%
UV/Vis.: λ_{max}: 230, 276, 312 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

N-acetyl-3,4-Methylenedioxymethcathinone (N-acetyl-3,4-MDMC) is an acetylated analog of methylone (Item No. 10986), the MDMA-type designer drug that has been detected in products marketed as bath salts, plant food, and tablets.^{1,2} The biological actions of N-acetyl-3,4-MDMC are unknown. This product is intended for forensic and research purposes.

References

1. Kikura-Hanajiri, R., Uchiyama, N., and Goda, Y. Survey of current trends in the abuse of psychotropic substances and plants in Japan. *Leg. Med. (Tokyo)* **13(3)**, 109-115 (2011).
2. Bossong, M.G., Van Dijk, J.P., and Niesink, R.J. Methylone and mCPP, two new drugs of abuse? *Addict. Biol.* **10(4)**, 321-323 (2005).

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, 07/11/2023

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