

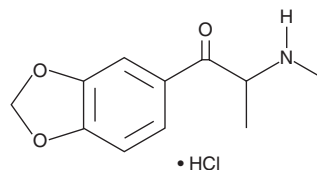
PRODUCT INFORMATION



Methylone (hydrochloride)

Item No. 10986

CAS Registry No.: 186028-80-8
Formal Name: 1-(1,3-benzodioxol-5-yl)-2-(methylamino)-1-propanone, monohydrochloride
Synonyms: M₁, bk-MDMA, β-keto Methyleneedioxy Methamphetamine, 3,4-Methyleneedioxy-N-methylcathinone
MF: C₁₁H₁₃NO₃ • HCl
FW: 243.7
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥3 years



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

Description

Methylone (hydrochloride) (Item No. 10986) is an analytical reference material that is categorized as a cathinone. Methylone blocks reuptake of dopamine, 5-HT, and norepinephrine in rat synaptosomes (IC₅₀s = 2.9, 2.3, and 0.74 μM, respectively).¹ It also induces release of the same neurotransmitters from synaptosomes (EC₅₀s = 0.38, 0.22, and 0.093 μM, respectively). It has been detected in products marketed as bath salts, plant food, and tablets.^{2,3} Methylone is regulated as a Schedule I 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.

References

1. Nagai, F., Nonaka, R., and Satoh Hisashi Kamimura, K. The effects of non-medically used psychoactive drugs on monoamine neurotransmission in rat brain. *Eur. J. Pharmacol.* **559(2-3)**, 132-137 (2007).
2. 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)**, 109-15 (2011).
3. 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, 10/12/2021

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