

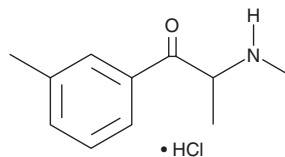
PRODUCT INFORMATION



3-Methylmethcathinone (hydrochloride)

Item No. 11224

CAS Registry No.: 1246816-62-5
Formal Name: 2-(methylamino)-1-(3-methylphenyl)-1-propanone, monohydrochloride
Synonyms: 3-methyl MC, 3-MMC
MF: C₁₁H₁₅NO • HCl
FW: 213.7
Purity: ≥97%
UV/Vis.: λ_{max}: 252, 292 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

4-Methylmethcathinone (4-methyl MC) is a psychoactive synthetic cathinone that has been identified in bath salts and powders.¹⁻³ 3-Methylmethcathinone (3-methyl MC) (hydrochloride) is a potential major impurity in preparations of 4-methyl MC. As an isomer of 4-methyl MC, it is likely to have psychoactive properties and may itself be marketed as a designer drug. The physiological and toxicological properties of this compound have not been determined. 3-methyl MC is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

References

1. Gibbons, S. and Zloh, M. An analysis of the 'legal high' mephedrone. *Bioorg. Med. Chem. Lett.* **20(14)**, 4135-4139 (2010).
2. Brandt, S.D., Freeman, S., Sumnall, H.R., et al. Analysis of NRG 'legal highs' in the UK: Identification and formation of novel cathinones. *Drug Test. Anal.* **3(9)**, 569-575 (2011).
3. 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).

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, 07/07/2023

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