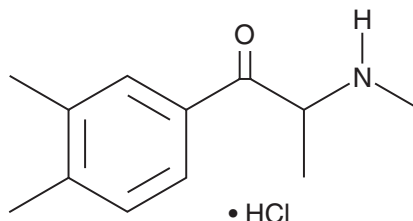


PRODUCT INFORMATION

3,4-Dimethylmethcathinone (hydrochloride)

Item No. 9001098

CAS Registry No.: 1081772-06-6
Formal Name: 1-(3,4-dimethylphenyl)-2-(methylamino)-1-propanone, monohydrochloride
Synonym: 3,4-DMMC
MF: C₁₂H₁₇NO • HCl
FW: 227.7
Purity: ≥98%
UV/Vis.: λ_{max}: 205, 262 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

3,4-DMMC (hydrochloride) is a potential designer drug with combined features of amphetamines, cathinones, and phenethylamines. It is structurally related to 4-methylmethcathinone, a psychoactive compound that has been identified in products sold as bath salts and plant food.¹⁻³ The physiological and toxicological properties of this compound have not been elucidated. 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), 4135-39 (2010).
2. James, D., Adams, R.D., Spears, R., *et al.* Clinical characteristics of mephedrone toxicity reported to the UK National Poisons Information Service. *Emerg. Med. J.* **28**(8), 686-689 (2011).
3. Jankovics, P., Váradi, A., Tölgyesi, L., *et al.* Identification and characterization of the new designer drug 40-methylethcathinone (4-MEC) and elaboration of a novel liquid chromatography-tandem mass spectrometry (LC-MS/MS) screening method for seven different methcathinone analogs. *Forensic Sci. Int.* **210**(1-3), 213-220 (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, 06/27/2023

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