

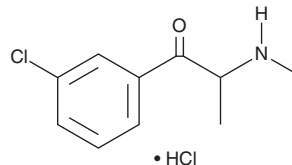
PRODUCT INFORMATION



3-Chloromethcathinone (hydrochloride)

Item No. 44462

CAS Registry No.: 1607439-32-6
Formal Name: 1-(3-chlorophenyl)-2-(methylamino)-1-propanone, monohydrochloride
Synonym: 3-CMC
MF: C₁₀H₁₂ClNO • HCl
FW: 234.1
Purity: ≥98%
Supplied as: A 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-Chloromethcathinone (hydrochloride) (Item No. 44462) is an analytical reference standard categorized as a cathinone. 3-Chloromethcathinone increases spontaneous locomotor activity in mice.¹ It has been associated with fatality and has been found in seized illicit drug samples.^{2,3} This product is intended for research and forensic applications.

References

1. Wojcieszak, J., Kuczyńska, K., and Zawilska, J.B. Four synthetic cathinones: 3-Chloromethcathinone, 4-chloromethcathinone, 4-fluoro- α -pyrrolidinopentiophenone, and 4-methoxy- α -pyrrolidinopentiophenone produce changes in the spontaneous locomotor activity and motor performance in mice with varied profiles. *Neurotox. Res.* **38**(2), 536-551 (2020).
2. Romańczuk, A., Rojek, S., Synowiec, K., *et al.* The stability of synthetic cathinones and the study of potential intake biomarkers in the biological material from a case of 3-CMC poisoning. *J. Anal. Toxicol.* **47**(5), 470-480 (2023).
3. Piorunska-Sedlak, K. and Stypulkowska, K. Strategy for identification of new psychoactive substances in illicit samples using attenuated total reflectance infrared spectroscopy. *Forensic Sci. Int.* **312**, 110262 (2020).

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, 09/10/2025

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