

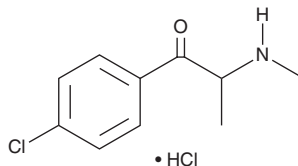
PRODUCT INFORMATION



4-Chloromethcathinone (hydrochloride)

Item No. 16436

CAS Registry No.: 2319878-22-1
Formal Name: 1-(4-chlorophenyl)-2-(methylamino)-1-propanone, monohydrochloride
Synonyms: Clephedrone, 4-CMC
MF: C₁₀H₁₂ClNO • HCl
FW: 234.1
Purity: ≥98%
UV/Vis.: λ_{max}: 260 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-Chloromethcathinone is an analog of methcathinone that has a chlorine atom inserted at the 4 position of the phenyl ring. While the physiological and toxicological properties of this compound are not known, a similar compound, 4-bromomethcathinone (Item No. 12089) inhibits norepinephrine reuptake, like methcathinone, but fails to stimulate spontaneous locomotor activity in rats.^{1,2} This product is intended for forensic and research applications.

References

1. Foley, K.F. and Cozzi, N.V. Inhibition of transport function and desipramine binding at the human noradrenaline transporter by N-ethylmaleimide and protection by substrate analogs. *Naunyn Schmiedebergs Arch. Pharmacol.* **365**(6), 457-461 (2002).
2. Foley, K.F. and Cozzi, N.V. Novel aminopropiophenones as potential antidepressants. *Drug Develop. Res.* **60**(4), 252-260 (2003).

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, 06/30/2023

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