

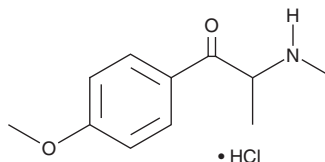
PRODUCT INFORMATION



Methedrone (hydrochloride)

Item No. 10529

CAS Registry No.: 879665-92-6
Formal Name: 1-(4-methoxyphenyl)-2-(methylamino)-1-propanone, monohydrochloride
Synonyms: *para*-Methoxymethcathinone, PMMC, Methoxyphedrine, 4-Methedrone
MF: C₁₁H₁₅NO₂ • HCl
FW: 229.7
Purity: ≥98%
UV/Vis.: λ_{max}: 223, 286 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

Methedrone is a phenethylamine, amphetamine, and cathinone derivative that acts as a triple reuptake/release/reversible monoamine oxidase inhibitor. Its use as a recreational drug and its toxicity has recently been reported.^{1,2} Methedrone is closely related to mephedrone, a psychoactive compound which is regulated in many countries, including the United States. The hydrochloride formulation of this compound is intended for use as a standard for the forensic analysis of samples that may contain methedrone.

References

1. 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).
2. Prosser, J.M. and Nelson, L.S. The toxicology of bath salts: A review of synthetic cathinones. *J. Med. Toxicol.* **8(1)**, 33-42 (2012).

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, 07/18/2023

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