

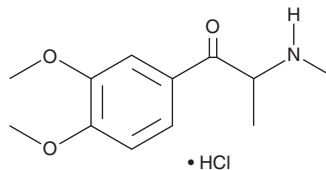
PRODUCT INFORMATION



3,4-Dimethoxymethcathinone (hydrochloride)

Item No. 17447

CAS Registry No.: 22930-82-1
Formal Name: 1-(3,4-dimethoxyphenyl)-2-(methylamino)-1-propanone, monohydrochloride
Synonyms: 3-methoxy Methedrone, 3,4-DMMC
MF: C₁₂H₁₇NO₃ • HCl
FW: 259.7
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

Mephedrone (Item No. 10801) is a designer drug of the phenethylamine class that shares substantial structural similarities with methcathinone (Item No. 11709) and methamphetamine (Item No. 13997). Compounds with these structural characteristics have psychoactive effects and have been identified in products sold as bath salts and plant food.¹⁻³ 3,4-Dimethoxymethcathinone is an analog of mephedrone with potential for abuse. Its biological properties have not been elucidated. This product is intended for research and forensic applications.

References

1. 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).
2. Gibbons, S. and Zloh, M. An analysis of the 'legal high' mephedrone. *Bioorg. Med. Chem. Lett.* **20(14)**, 4135-4139 (2010).
3. Drugs and chemicals of concern. *Drug Enforcement Administration* (2010).

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, 05/09/2023

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