

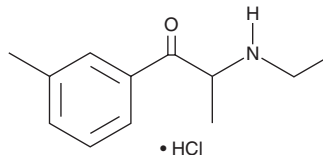
PRODUCT INFORMATION



3-Methylethcathinone (hydrochloride)

Item No. 11222

CAS Registry No.: 2493976-59-1
Formal Name: 2-(ethylamino)-1-(3-methylphenyl)-1-propanone, monohydrochloride
Synonym: 3-MEC
MF: C₁₂H₁₇NO • HCl
FW: 227.7
Purity: ≥98%
UV/Vis.: λ_{max}: 206, 252, 291 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Description

Substituted cathinones, like 4-methylethcathinone (Item No. 9001069) and 3,4-dimethylmethcathinone (Item No. 9001098) are psychoactive compounds that may be used as designer drugs.¹ 3-Methylethcathinone (hydrochloride) is a substituted cathinone with potential for abuse. Its biological properties have not been elucidated. 3-Methylethcathinone is regulated as a Schedule I compound in the United States. This product is intended for forensic and research applications.

Reference

1. Jankovics, P., Váradi, A., Tölgyesi, L., *et al.* Identification and characterization of the new designer drug 4'-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, 005/14/2021

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