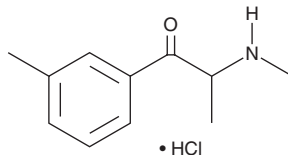


PRODUCT INFORMATION

3-Methylmethcathinone (hydrochloride) (exempt preparation)

Item No. 38141

CAS Registry No.: 1246816-62-5
Formal Name: 2-(methylamino)-1-(3-methylphenyl)-1-propanone, monohydrochloride
Synonyms: 3-methyl MC, 3-MMC
MF: C₁₁H₁₅NO • HCl
FW: 213.7
Purity: ≥97%
UV/Vis.: λ_{max}: 252, 292 nm
Supplied as: A 1 mg/ml solution in methanol
Storage: -20°C
Stability: ≥3 years



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

Description

3-Methylmethcathinone (hydrochloride) (exempt preparation) (Item No. 38141) is an analytical reference standard categorized as a cathinone. 3-Methylmethcathinone is a potential impurity in the synthesis of 4-methylmethcathinone. It has been associated with hospitalizations for mono-intoxication and overdose fatality in combination with other compounds.¹ 3-Methylmethcathinone is regulated as a Schedule I compound in the United States. 3-Methylmethcathinone (hydrochloride) (exempt preparation) (Item No. 38141) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

Reference

1. Bäckberg, M., Lindeman, E., Beck, O., *et al.* Characteristics of analytically confirmed 3-MMC-related intoxications from the Swedish STRIDA project. *Clin. Toxicol. (Phila)* **53**(1), 46-53 (2015).

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, 01/17/2025

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