

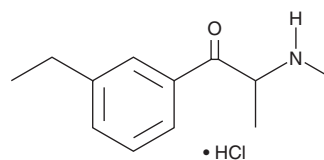
PRODUCT INFORMATION



3-Ethylmethcathinone (hydrochloride)

Item No. 9001082

CAS Registry No.: 2446466-60-8
Formal Name: 1-(3-ethylphenyl)-2-(methylamino)-1-propanone, monohydrochloride
Synonym: 3-EMC
MF: C₁₂H₁₇NO • HCl
FW: 227.7
Purity: ≥98%
UV/Vis.: λ_{max}: 208, 253, 292 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

3-Ethylmethcathinone (hydrochloride) is a potential stimulant/entactogen designer drug that is of the amphetamine, phenethylamine, and cathinone chemical classes. It is structurally related to 4-methylmethcathinone (Item No. 10801), a designer drug that has been detected in products marketed as bath salts, plant food, and tablets.¹⁻³ The physiological and toxicological properties of this compound are not known. This product is intended for forensic purposes.

References

1. Brandt, S.D., Sumnall, H.R., Measham, F., *et al.* Analyses of second-generation 'legal-highs' in the UK: Initial findings. *Drug Test. Anal.* **2(8)**, 377-382 (2010).
2. Gibbons, S. and Zloh, M. An analysis of the 'legal high' mephedrone. *Bioorg. Med. Chem. Lett.* **20(14)**, 4135-4139 (2010).
3. James, D., Adams, R.D., Spears, R., *et al.* Clinical characteristics of mephedrone toxicity reported to the UK National Poisons Information Service. *Emerg. Med. J.* **28(8)**, 686-689 (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, 06/23/2023

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