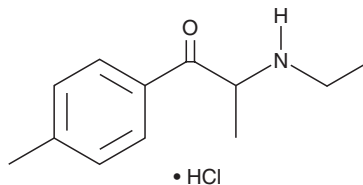


PRODUCT INFORMATION

4-Methylethcathinone (hydrochloride)

Item No. 9001069

CAS Registry No.: 1266688-86-1
Formal Name: 2-(ethylamino)-1-(4-methylphenyl)-1-propanone, monohydrochloride
MF: C₁₂H₁₇NO • HCl
FW: 227.7
Purity: ≥98%
UV/Vis.: λ_{max}: 260 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

4-Methylethcathinone (4-MEC) is a cathinone derivative identified in several designer, recreational drugs that are sold as “legal high” replacements for controlled stimulants such as methamphetamine and 3,4-methylenedioxymethamphetamine.¹⁻³ The hydrochloride formulation of this compound is intended for use as a standard for the forensic analysis of samples that may contain 4-MEC.

References

- Schifano, F., Albanese, A., Fergus, S., *et al.* Mephedrone (4-methylmethcathinone; ‘meow meow’): Chemical, pharmacological and clinical issues. *Psychopharmacol. (Berl)*. **214(3)**, 593-602 (2011).
- Santali, E.Y., Cadogan, A.-K., Nic Daeid, N., *et al.* Synthesis, full chemical characterisation and development of validated methods for the quantification of (±)-4'-methylmethcathinone (mephedrone): A new “legal high”. *J. Pharm. Biomed. Anal.* **56(2)**, 246-255 (2011).
- Gibbons, S. and Zloh, M. An analysis of the ‘legal high’ mephedrone. *Bioorg. Med. Chem. Lett.* **20(14)**, 4135-39 (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, 06/22/2023

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