

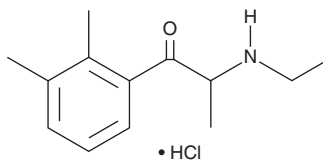
PRODUCT INFORMATION



2,3-Dimethylethcathinone (hydrochloride)

Item No. 11227

CAS Registry No.: 2705531-06-0
Formal Name: 1-(2,3-dimethylphenyl)-2-(ethylamino)propan-1-one, monohydrochloride
Synonym: 2,3-DMEC
MF: C₁₃H₁₉NO • HCl
FW: 241.8
Purity: ≥98%
UV/Vis.: λ_{max}: 252, 293 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

Substituted cathinones are psychoactive compounds commonly used as recreational drugs.^{1,2} Ethcathinone (Item No. 11241) is a substituted cathinone that produces amphetamine-like stimulus effects in rats trained to discriminate amphetamine from vehicle.³ It has been identified as a component of capsules or tablets.^{2,4} 2,3-Dimethylethcathinone is a disubstituted analog of ethcathinone. Its physiological and toxicological properties have not been investigated. This product is intended for forensic and research purposes.

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).
- Kikura-Hanajiri, R., Uchiyama, N., and Goda, Y. Survey of current trends in the abuse of psychotropic substances and plants in Japan. *Leg. Med. (Tokyo)* **13(3)**, 109-15 (2011).
- Dal Cason, T.A., Young, R., and Glennon, R.A. Cathinone: An investigation of several N-alkyl and methylenedioxy-substituted analogs. *Pharmacol. Biochem. Behav.* **58(4)**, 1109-1116 (1997).
- Camilleri, A., Johnston, M.R., Brennan, M., *et al.* Chemical analysis of four capsules containing the controlled substance analogues 4-methylmethcathinone, 2-fluoromethamphetamine, α-phthalimidopropiophenone and N-ethylcathinone. *Forensic Sci. Int.* **197(1-3)**, 59-66 (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, 07/13/2023

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