

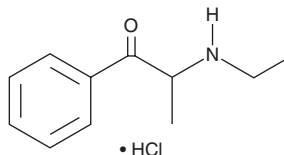
# PRODUCT INFORMATION



## Ethcathinone (hydrochloride)

Item No. 11241

**CAS Registry No.:** 51553-17-4  
**Formal Name:** 2-(ethylamino)-1-phenyl-1-propanone, monohydrochloride  
**Synonyms:** N-Ethylcathinone, RMI 8201A  
**MF:** C<sub>11</sub>H<sub>15</sub>NO • HCl  
**FW:** 213.7  
**Purity:** ≥97%  
**UV/Vis.:** λ<sub>max</sub>: 204, 248 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.<sup>1,2</sup> Ethcathinone (hydrochloride) is a substituted cathinone that produces amphetamine-like stimulus effects in rats trained to discriminate amphetamine from vehicle.<sup>3</sup> It has been identified as a component of capsules or tablets.<sup>2,4</sup> This product is intended for research and forensic applications.

### References

- Schifano, F., Albanese, A., Fergus, S., *et al.* Mephedrone (4-methylmethcathinone; 'meow meow'): Chemical, pharmacological and clinical issues. *Psychopharmacology (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, 07/07/2023

#### 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](http://WWW.CAYMANCHEM.COM)