

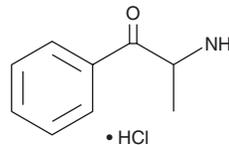
PRODUCT INFORMATION



DL-Cathinone (hydrochloride)

Item No. 14138

CAS Registry No.: 16735-19-6
Formal Name: 2-amino-1-phenyl-1-propanone, monohydrochloride
Synonyms: α -Aminopropiophenone, NSC 166465
MF: C₉H₁₁NO • HCl
FW: 185.7
Purity: \geq 98%
UV/Vis.: λ_{max} : 247 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 5 years



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

Description

Cathinone, the main psychoactive alkaloid from the khat plant (*C. edulis*), is a potent central stimulant and a naturally occurring analog of amphetamine. DL-Cathinone is a racemic mixture of (-)-(S)-cathinone (Item No. 13869) and its less potent (+)-isomer.¹ Like amphetamine, DL-cathinone and (-)-(S)-cathinone are potent releasers of norepinephrine and dopamine.²⁻⁴ With an ED₅₀ value of 0.24 mg/kg, DL-cathinone produces amphetamine-like effects on the operant behavior of rats trained to discriminate 0.2-0.8 mg/kg of (D)-amphetamine.⁴ This product is intended for research and forensic applications.

References

1. 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).
2. Kalix, P. and Glennon, R.A. Further evidence for an amphetamine-like mechanism of action of the alkaloid cathinone. *Biochem. Pharmacol.* **35(18)**, 3015-3019 (1986).
3. Glennon, R.A., and Liebowitz, S.M. Serotonin receptor affinity of cathinone and related analogues. *J. Med. Chem.* **25(4)**, 393-397 (1982).
4. Schechter, M.D., Rosecrans, J.A., and Glennon, R.A. Comparison of behavioral effects of cathinone, amphetamine and apomorphine. *Pharmacol. Biochem. Behav.* **20(2)**, 181-184 (1984).

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, 07/25/2023

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