

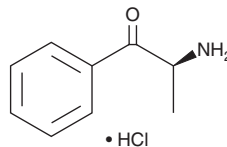
# PRODUCT INFORMATION



## (-)-(S)-Cathinone (hydrochloride)

Item No. 13869

**CAS Registry No.:** 72739-14-1  
**Formal Name:** 2-amino-1-phenyl-1-propanone, monohydrochloride  
**MF:** C<sub>9</sub>H<sub>11</sub>NO • HCl  
**FW:** 185.7  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 247 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

Cathinone, the main psychoactive alkaloid from the khat plant (*C. edulis*), is a potent central stimulant and a naturally occurring analog of amphetamine. (-)-(S)-Cathinone is several times more potent than its (+)-isomer, and its absolute stereochemistry is identical to that of (+)-(S)-amphetamine.<sup>1</sup> Like amphetamine, (-)-(S)-cathinone is a potent releaser of norepinephrine and dopamine from their intracellular stores.<sup>2,3</sup> With an ED<sub>50</sub> value of 0.42 μmol/kg, (-)-(S)-cathinone produces amphetamine-like stimulus effects in rats trained to discriminate 1 mg/kg of (+)-(S)-amphetamine from vehicle.<sup>1</sup> 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. Be.* **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).

#### 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