

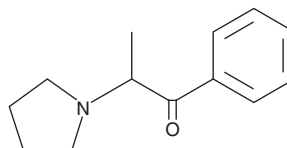
PRODUCT INFORMATION



α -Pyrrolidinopropiophenone (hydrochloride)

Item No. 10445

CAS Registry No.: 92040-10-3
Formal Name: 1-phenyl-2-(1-pyrrolidinyl)-1-propanone, monohydrochloride
Synonym: α -PPP
MF: C₁₃H₁₇NO • HCl
FW: 239.7
Purity: $\geq 97\%$
UV/Vis.: λ_{\max} : 250 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥ 5 years



• HCl

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

Description

α -Pyrrolidinopropiophenone (hydrochloride) (Item No. 10445) is an analog of the appetite suppressant diethylcathinone and is related to the designer drug 4'-methyl- α -pyrrolidinopropiophenone (MPPP). In male Wistar rats, it is metabolized to a variety of products, including cathinone, the main psychoactive alkaloid of *Catha edulis*.^{1,2} α -Pyrrolidinopropiophenone (hydrochloride) is intended to be used for forensic applications.

References

1. Springer, D., Fritschi, G., and Maurer, H. Metabolism of the new designer drug α -pyrrolidinopropiophenone (PPP) and the toxicological detection of PPP and 4'-methyl- α -pyrrolidinopropiophenone (MPPP) studied in rat urine using gas chromatography-mass spectrometry. *J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.* **796(2)**, 253-266 (2003).
2. Staack, R.F. and Maurer, H.H. Metabolism of designer drugs of abuse. *Curr. Drug Metab.* **6(3)**, 259-274 (2005).

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, 07/18/2023

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