

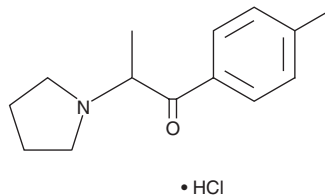
PRODUCT INFORMATION



4'-methyl- α -Pyrrolidinopropiophenone (hydrochloride)

Item No. 10446

CAS Registry No.: 1313393-58-6
Formal Name: 1-(4-methylphenyl)-2-(1-pyrrolidinyl)-1-propanone, monohydrochloride
Synonyms: 4'-methyl- α -PPP, 4'-methyl PPP
MF: C₁₄H₁₉NO • HCl
FW: 253.8
Purity: \geq 97%
UV/Vis.: λ_{max} : 202, 261 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

Pyrrolidinophenones (PPPs) are a family of compounds that feature a pyrrolidinyl group affixed to the narcotic cathinone. Many have been identified as components of designer drugs.¹ 4'-methyl- α -PPP (hydrochloride) shares structural features with the stimulant α -PPP and has been detected in bath salts and other formulations.^{1,2} The metabolism of 4'-methyl- α -PPP has been characterized.^{3,4} This product is intended to be used for forensic applications.

References

1. 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)**, 109-115 (2011).
2. Brandt, S.D., Freeman, S., Sumnall, H.R., et al. Analysis of NRG 'legal highs' in the UK: Identification and formation of novel cathinones. *Drug Test. Anal.* **3(9)**, 569-575 (2011).
3. Springer, D., Staack, R.F., Paul, L.D., et al. Identification of cytochrome P450 enzymes involved in the metabolism of 4'-methyl-alpha-pyrrolidinopropiophenone, a novel scheduled designer drug, in human liver microsomes. *Xenobiotica* **31(8)**, 979-982 (2003).
4. 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).

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