Product Information

4'-methyl- α -Pyrrolidinopropiophenone (hydrochloride) (exempt preparation)

Item No. 16051

CAS Registry No.: 1313393-58-6

Formal Name: 2-(pyrrolidin-1-yl)-1-(p-tolyl)propan-1-

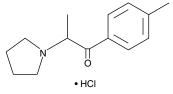
one, monohydrochloride

Synonyms: 4'-methyl-α-PPP, 4'-methyl PPP

MF: C₁₄H₁₉NO • HCl

FW: 253.8 **Purity:** ≥97%

Stability: ≥2 years at -20°C Supplied as: A solution in methanol λ_{max} : 202, 261 nm UV/Vis.:



Laboratory Procedures

For long term storage, we suggest that 4'-methyl-\alpha-pyrrolidinopropiophenone (4'-methyl-\alpha-PPP) (hydrochloride) (exempt preparation) be stored as supplied at -20°C. It should be stable for at least two years.

4'-methyl-α-PPP (hydrochloride) (exempt preparation) is supplied as a solution in methanol. To change the solvent, simply evaporate the methanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of 4'-methyl-α-PPP (hydrochloride) (exempt preparation) in these solvents is approximately 20, 10, and 3 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. If an organic solvent-free solution of 4'-methyl-α-PPP (hydrochloride) (exempt preparation) is needed, it can be prepared by evaporating the methanol and directly dissolving the neat oil in aqueous buffers. The solubility of 4'-methyl-α-PPP (hydrochloride) (exempt preparation) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

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. 14'-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), 109-15 (2011).
- 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. (2010).
- Springer, D., Staack, R.F., Paul, L.D., et al. Identification of cytochrome P450 enzymes involved in the metabolism of 4'-methyl-α-pyrrolidinopropiophenone, a novel scheduled designer drug, in human liver microsomes. Xenobiotica **31(8)**, 979-982 (2003).
- 4. Dietmar, S., Fritschi, G., and Maurer, H. Metabolism of the new designer drug α-pyrrolidinoproppiophenone (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).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10446

WARNING: This product is for laboratory research only: not for administration to humans. Not for human or veterinary DIAGNOSTIC OR THERAPEUTIC USE.

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will <u>meet our specifications</u>

at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitity does not apply in the case of intentional acts or negligence of Cayman, is directors or its employees.

Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that

Buyers exclusive remedy and Caymans sole hability increments man be immed to a teams of the purchase of the material at its destination. Failure of Buyer to give said notice within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 09/25/2014

Cayman Chemical

Mailing address

1180 E. Ellsworth Road Ann Arbor, MI 48108 USA

Phone

(800) 364-9897 (734) 971-3335

(734) 971-3640

custserv@caymanchem.com

www.caymanchem.com