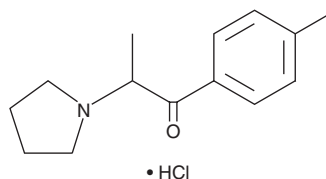


PRODUCT INFORMATION

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%
UV/Vis.: λ_{max} : 202, 261 nm
Supplied as: A solution in methanol
Storage: -20°C
Stability: ≥5 years



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

Laboratory Procedures

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.

Description

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), 109-15 (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.* (2010).
3. 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. Springer, D., Fritschi, G., and Maurer, H.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, 10/01/2025

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