

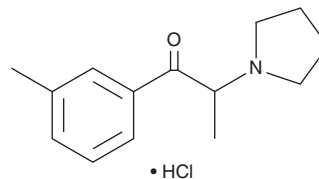
PRODUCT INFORMATION



3-methyl- α -Pyrrolidinopropiophenone (hydrochloride)

Item No. 11485

CAS Registry No.: 2749302-38-1
Formal Name: 1-(3-methylphenyl)-2-(1-pyrrolidinyl)-1-propanone, monohydrochloride
Synonyms: 3-methyl- α -PPP, 3-methyl PPP
MF: C₁₄H₁₉NO • HCl
FW: 253.8
Purity: \geq 98%
UV/Vis.: λ_{\max} : 254, 293 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 2 years



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

Description

The α -pyrrolidinophenones (α -PPPs) are a class of designer drugs, typified by α -PPP (Item No. 10445) and 4-methyl- α -PPP (Item No. 10446).¹⁻³ Although structurally related to amphetamines and cathinones, the pharmacology and toxicology of members of this class have not been elucidated.⁴ 3-methyl- α -PPP is a structural isomer of 4-methyl- α -PPP, having the methyl group at the 3 position, rather than the 4 position, of the phenyl group. This product is intended for forensic and research 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. 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. 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).
4. 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, 05/23/2022

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