

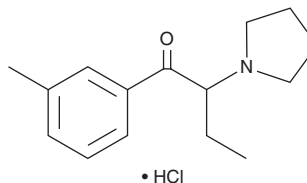
PRODUCT INFORMATION



3-methyl- α -Pyrrolidinobutiophenone (hydrochloride)

Item No. 9001189

Formal Name: 2-(pyrrolidin-1-yl)-1-(m-tolyl)butan-1-one, monohydrochloride
Synonym: 3-methyl PBP
MF: C₁₅H₂₁NO • HCl
FW: 267.8
Purity: ≥98%
UV/Vis.: λ_{max} : 255, 295 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

4-Methyl- α -pyrrolidinobutiophenone (4-methyl PBP) (Item No. 9001190) is a synthetic cathinone with psychoactive properties which has recently been identified in party pills and powders.^{1,2} The metabolism of methyl PBP has also been characterized.^{3,4} 3-methyl PBP (hydrochloride) is an isomer of 4-methyl PBP having the methyl group attached at the 3 position of the phenyl ring, instead of the 4 position. The biological and toxicological properties of this compound are unknown. This product is intended for forensic and research applications.

References

1. Analysis of NRG 'legal highs' in the UK: Identification and formation of novel cathinones. Brandt, S.D., *et al.* Retrieved August 18, 2010, from www.drugtestinganalysis.com.
2. Westphal, F., Junge, T., Rösner, P., *et al.* Mass spectral and NMR spectral data of two new designer drugs with an α -aminophenone structure: 4'-Methyl- α -pyrrolidinohexanophenone and 4'-methyl- α -pyrrolidinobutyrophenone. *Forensic Sci. Int.* **169(1)**, 32-42 (2007).
3. Meyer, M.R. and Maurer, H.H. Metabolism of designer drugs of abuse: An updated review. *Curr. Drug Metab.* **11(5)**, 468-482 (2010).
4. Peters, F.T., Meyer, M.R., Theobald, D.S., *et al.* Identification of cytochrome P450 enzymes involved in the metabolism of the new designer drug 4'-methyl- α -pyrrolidinobutyrophenone. *Drug Metab. Dispos.* **36(1)**, 163-168 (2008).

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, 07/06/2023

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