

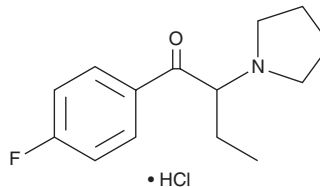
PRODUCT INFORMATION



4-fluoro- α -Pyrrolidinobutyphenone (hydrochloride)

Item No. 15516

CAS Registry No.:	2749302-50-7
Formal Name:	1-(4-fluorophenyl)-2-(pyrrolidin-1-yl)butan-1-one, monohydrochloride
Synonym:	4-fluoro PBP
MF:	C ₁₄ H ₁₈ FNO • HCl
FW:	271.8
Purity:	≥98%
UV/Vis.:	λ_{\max} : 253 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

Pyrrolidinophenones, including 4-methyl- α -PBP (4-MePBP; Item No. 9001190), are synthetic cathinones that have been identified in party pills.¹⁻³ Although not much is known regarding their pharmacological actions, pyrrolidinophenones have potential for abuse due to their close structural relation to pyrovalerone (Item No. 10817). 4-fluoro- α -PBP is an analog of 4-MePBP in which a fluorine is substituted for a methyl at the four position of the phenyl ring. Its physiological and toxicological properties have not been determined. This product is intended for forensic and research applications.

References

1. 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).
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. 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, 06/20/2022

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