

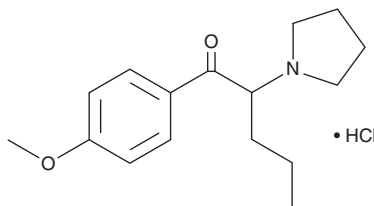
PRODUCT INFORMATION



4-methoxy- α -Pyrrolidinopentiophenone (hydrochloride)

Item No. 14097

CAS Registry No.: 5537-19-9
Formal Name: 1-(4-methoxyphenyl)-2-(1-pyrrolidinyl)-1-pentanone, monohydrochloride
Synonyms: 4-MeO- α -PVP, 4-methoxy- α -Pyrrolidinovalerophenone, 4'-methoxy-2-(1-pyrrolidinyl)-Valerophenone
MF: C₁₆H₂₃NO₂ • HCl
FW: 297.8
Purity: ≥98%
UV/Vis.: λ_{max} : 225, 291 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 are a family of compounds that feature a pyrrolidinyl group affixed to the narcotic cathinone. Many have been identified as components of designer drugs. 4-methoxy- α -Pyrrolidinopentiophenone (4-MeO- α -PVP) (hydrochloride) shares structural features with the stimulant α -PVP. The mechanism of action is unknown for α -PVP, but is believed to be similar to that of the designer drug MDPV (Item No. 10624), which acts as a norepinephrine-dopamine reuptake inhibitor.^{1,2} This product is intended to be used for forensic and research applications.

References

1. Heron, C., Costentin, J., and Bonnet, J.J. Evidence that pure uptake inhibitors including cocaine interact slowly with the dopamine neuronal carrier. *Eur. J. Pharmacol.* **264**(3), 391-398 (1994).
2. Meltzer, P.C., Butler, D., Deschamps, J.R., *et al.* 1-(4-methylphenyl)-2-pyrrolidin-1-yl-pentan-1-one (pyrovalerone) analogs. A promising class of monoamine uptake inhibitors. *J. Med. Chem.* **49**(4), 1420-1432 (2006).

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, 07/14/2023

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