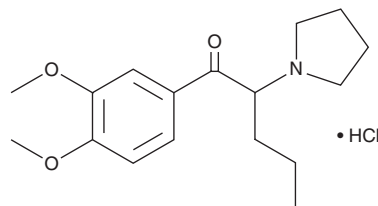


PRODUCT INFORMATION

3,4-dimethoxy- α -Pyrrolidinopentiophenone (hydrochloride)

Item No. 14106

CAS Registry No.: 850351-99-4
Formal Name: 1-(3,4-dimethoxyphenyl)-2-(1-pyrrolidinyl)-1-pentanone, monohydrochloride
Synonyms: 3,4-dimethoxy- α -PVP, 3,4-dimethoxy- α -Pyrrolidinovalerophenone, 3,4-dimethoxy-2-(1-pyrrolidinyl)-Valerophenone
MF: C₁₇H₂₅NO₃ • HCl
FW: 327.9
Purity: ≥98%
UV/Vis.: λ_{max} : 235, 285, 317 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

Methylenedioxy Pyrovalerone (MDPV) (Item No. 10624) is a cathinone commonly found in products sold as bath salts or plant food.^{1,2} It is regulated as a Schedule 1 compound in the United States.³ 3,4-dimethoxy- α -Pyrrolidinopentiophenone is an analog of MDPV, differing only by an opening of the ring structure on the phenyl group. The physiological and toxicological properties of this compound have not been investigated. This product is intended for research and forensic 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. Prosser, J.M. and Nelson, L.S. The toxicology of bath salts: A review of synthetic cathinones. *J. Med. Toxicol.* **8**(1), 33-42 (2012).
3. Methylenedioxy pyrovalerone (MDPV) (1-(1,3-Benzodioxol-5-yl)-2-(1-pyrrolidinyl)-1-pentanone]. In: Drug Enforcement Administration. Diversion Control Division (2010). Available from: http://www.deadiversion.usdoj.gov/drug_chem_info/mdpv.pdf

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