

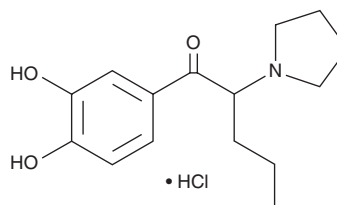
PRODUCT INFORMATION



3,4-Methylenedioxy Pyrovalerone metabolite 2 (hydrochloride)

Item No. 12004

CAS Registry No.: 2748289-38-3
Formal Name: 1-(3,4-dihydroxyphenyl)-2-(1-pyrrolidinyl)-1-pentanone, monohydrochloride
Synonyms: 3,4-Dihydroxypyrovalerone, 3,4-MDPV metabolite 2, 3,4-hydroxy- α -Pyrrolidinopentiophenone
MF: C₁₅H₂₁NO₃ • HCl
FW: 299.8
Purity: \geq 98%
UV/Vis.: λ_{max} : 209, 239, 289, 323 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Description

3,4-Methylenedioxy Pyrovalerone (3,4-MDPV, Item No. 10684) is an analog of pyrovalerone (Item No. 10817) that includes the 3,4-methylenedioxy moiety found on 3,4-methylenedioxymethamphetamine (Item No. 13971), a DEA Schedule I controlled substance. 3,4-MDPV is commonly abused as a recreational drug.¹⁻³ 3,4-MDPV metabolite 2 is produced during phase I metabolism of 3,4-MDPV.^{4,5} This compound can be further modified, e.g., by glucuronidation, before secretion in urine.⁴ This product is intended for forensic and research purposes.

References

1. Brandt, S.D., Sumnall, H.R., Measham, F., *et al.* Analyses of second-generation 'legal-highs' in the UK: Initial findings. *Drug Test. Anal.* **2(8)**, 377-382 (2010).
2. 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)**, 109-115 (2011).
3. 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).
4. Strano-Rossi, S., Cadwallader, A.B., de la Torre, X., *et al.* Toxicological determination and *in vitro* metabolism of the designer drug methylenedioxypyrovalerone (MPDV) by gas chromatography/mass spectrometry and liquid chromatography/quadrupole time-of-flight mass spectrometry. *Rapid Commun. Mass Spectrom.* **24(18)**, 2706-2714 (2010).
5. Meyer, M.R., Du, P., Schuster, F., *et al.* Studies on the metabolism of the α -pyrrolidinophenone designer drug methylenedioxy-pyrovalerone (MDPV) in rat and human urine and human liver microsomes using GC-MS and LC-high-resolution MS and its detectability in urine by GC-MS. *J. Mass Spectrom.* **45(12)**, 1426-1442 (2010).

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, 06/13/2023

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