

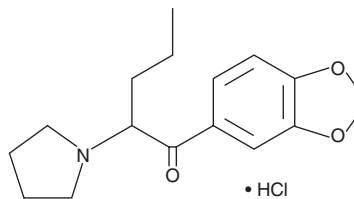
PRODUCT INFORMATION



3,4-Methylenedioxy Pyrovalerone (hydrochloride)

Item No. 10684

CAS Registry No.: 24622-62-6
Formal Name: 1-(1,3-benzodioxol-5-yl)-2-(1-pyrrolidinyl)-1-pentanone, monohydrochloride
Synonym: 3,4-MDPV
MF: C₁₆H₂₁NO₃ • HCl
FW: 311.8
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥7 years



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

Description

3,4-Methylenedioxy pyrovalerone (hydrochloride) (3,4-MDPV) (Item No. 10684) is an analytical reference material categorized as a cathinone. Pyrovalerone (Item No. 10817) and its analogs are inhibitors of the transporters for certain monoamine neurotransmitters, including dopamine and norepinephrine, preventing their uptake.^{1,2} 3,4-MDPV is an analog of pyrovalerone which includes the 3,4-methylenedioxy moiety found on 3,4-methylenedioxymethamphetamine, a DEA Schedule I controlled substance. While its physiological, neurological, and toxicological actions have not been characterized, 3,4-MDPV has been reported by the DEA to be abused as a central nervous system stimulant.³ Its effective dose and chemical interactions are unknown, but it has been used alone and in combination with other stimulating compounds. Products containing 3,4-MDPV have been marketed in Europe and Australia; they have also been seized by law enforcement in several states.³ 3,4-MDPV and some of its metabolites have recently been characterized by spectroscopic analysis.⁴⁻⁷ 3,4-MDPV is to be used in the forensic analysis of samples that may contain this compound. This product, the hydrochloride salt of 3,4-MDPV, has superior solubility in aqueous solvents, compared to the free base (Item No. 10624). This product is intended for research and forensic applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards for reference materials.

References

1. Heron, C., Costentin, J., and Bonnet, J.J. *Eur. J. Pharmacol.* **264**(3), 391-398 (1994).
2. Meltzer, P.C., Butler, D., Deschamps, J.R., et al. *J. Med. Chem.* **49**(4), 1420-1432 (2006).
3. U.S. Drug Enforcement Administration DEA, Diversion Control Division. In: Drug and Chemical Evaluation Section [Internet]. Available from: https://www.deadiversion.usdoj.gov/drugs_concern/mdpv.pdf.
4. Westphal, F., Junge, T., Rosner, P., et al. *Forensic Sci. Int.* **190**(1-3), 1-8 (2009).
5. Meyer, M.R., Du, P., Schuster, F., et al. *J. Mass Spectrom.* **45**(12), 1426-1442 (2010).
6. Yohannan, J.C. and Bozenko, J.S. *Microgram J.* **7**(1), 12-15 (2010).
7. Strano-Rossi, S., Cadwallader, A.B., de la Torre, X., et al. *Rapid Commun. Mass Spectrom.* **24**(18), 2706-2714 (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, 02/22/2021

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