

# PRODUCT INFORMATION

## 2,3-Methylenedioxy Pyrovalerone (hydrochloride)

Item No. 9001051

**Formal Name:** 1-(1,3-benzodioxol-4-yl)-2-(1-pyrrolidinyl)-1-pentanone, monohydrochloride

**MF:** C<sub>16</sub>H<sub>21</sub>NO<sub>3</sub> • HCl

**FW:** 311.8

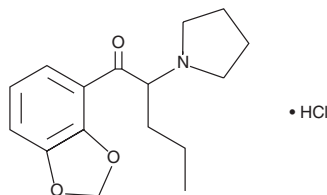
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 231, 261, 348 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

2,3-Methylenedioxy Pyrovalerone (2,3-MDPV) is a potential psychoactive designer drug. It is a structural isomer of 3,4-MDPV, a designer drug that has been detected in products marketed as bath salts, plant food, and tablets.<sup>1-3</sup> This product is intended for forensic 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), 109-15 (2011).
3. Yohannan, J.C. and Bozenko, J.S. The characterization of 3,4-methylenedioxypyrovalerone (MDPV). *Microgram Journal* **7**(1), 12-15 (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, 07/26/2023

#### 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