

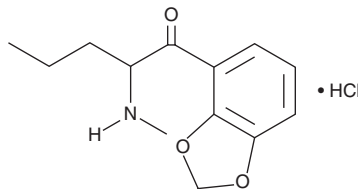
# PRODUCT INFORMATION



## 2,3-Pentylone isomer (hydrochloride)

Item No. 11463

**CAS Registry No.:** 2749428-57-5  
**Formal Name:** 1-(1,3-benzodioxol-4-yl)-2-(methylamino)-1-pentanone, monohydrochloride  
**MF:** C<sub>13</sub>H<sub>17</sub>NO<sub>3</sub> • HCl  
**FW:** 271.7  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 231, 260, 345 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

Pentylone is a synthetic cathinone that has been detected in party pills and powders.<sup>1,2</sup> Structurally, pentylone features a 3,4-methylenedioxy group attached to the phenone moiety which characterizes cathinones. 2,3-Pentylone isomer (hydrochloride) is a structural isomer of pentylone, having the methylenedioxy group attached at carbons 2 and 3 of the terminal phenyl group. The physiological and toxicological properties of this compound have not been characterized. This product is intended for forensic and research applications.

### References

1. Analysis of NRG 'legal highs' in the UK: Identification and formation of novel cathinones. Brandt, S.D., et al. Retrieved August 18, 2010, from [www.drugtestinganalysis.com](http://www.drugtestinganalysis.com).
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).

#### 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](http://WWW.CAYMANCHEM.COM)