

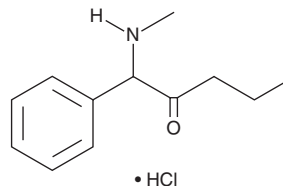
PRODUCT INFORMATION



Isopentedrone (hydrochloride)

Item No. 11563

CAS Registry No.: 1429402-13-0
Formal Name: 1-(methylamino)-1-phenyl-2-pentanone, monohydrochloride
MF: C₁₂H₁₇NO • HCl
FW: 227.7
Purity: ≥98%
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

Pentedrone (Item No. 11011) is a substituted cathinone and designer drug that is similar to methcathinone, a controlled psychoactive stimulant. Isopentedrone is a variant of pentedrone in which the α-propyl and β-keto groups have switched positions. This by-product of the synthesis of pentedrone has been detected in seized designer drug powders.¹ The physiological and toxicological effects of this compound have not been characterized. This product is intended for forensic and research purposes.

Reference

1. Westphal, F., Junge, T., Girreser, U., *et al.* Mass, NMR and IR spectroscopic characterization of pentedrone and pentylone and identification of their isocathinone by-products. *Forensic Sci. Int.* **217(1-3)**, 157-167 (2012).

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, 07/18/2023

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