

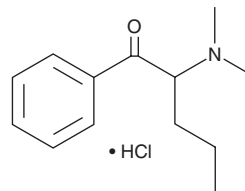
PRODUCT INFORMATION



α -Dimethylaminopentiophenone (hydrochloride)

Item No. 9001579

Formal Name:	2-(dimethylamino)-1-phenyl-1-pentanone, monohydrochloride
Synonym:	Dimethylpentedrone, N,N-dimethyl Pentedrone
MF:	C ₁₃ H ₁₉ NO • HCl
FW:	241.8
Purity:	≥98%
UV/Vis.:	λ_{\max} : 252 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

α -Pyrrolidinopentiophenone (α -PVP; Item No. 16058) is an analog of pyrovalerone (Item No. 10817), a psychoactive compound that is regulated as Schedule V in the United States. Pyrrolidinophenone-derived designer drugs, like α -PVP, are emerging drugs of interest.^{1,2} α -Dimethylaminopentiophenone is a novel α -PVP analog in which a dimethylamino group takes the place of the pyrrolidine. The biochemical, physiological, and toxicological properties of α -dimethylaminopentiophenone have not been determined. This product is intended for forensic and research applications.

References

1. Staack, R.F. and Maurer, H.H. Metabolism of designer drugs of abuse. *Curr. Drug Metab.* **6(3)**, 259-274 (2005).
2. Meyer, M.R. and Maurer, H.H. Metabolism of designer drugs of abuse: An updated review. *Curr. Drug Metab.* **11(5)**, 468-482 (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, 07/25/2023

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