

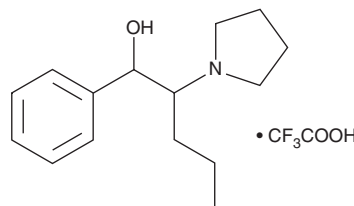
PRODUCT INFORMATION



α -Pyrrolidinopentiophenone metabolite 1 (trifluoroacetate salt)

Item No. 14093

CAS Registry No.: 1797986-63-0
Formal Name: α -phenyl- β -propyl-1-pyrrolidineethanol, 2,2,2-trifluoroacetate salt
Synonyms: α -PVP metabolite 1, α -Pyrrolidinovalerophenone, 2-(1-pyrrolidinyl)-Valerophenone metabolite 1
MF: $C_{15}H_{23}NO \cdot CF_3COOH$
FW: 347.4
Purity: $\geq 98\%$
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 5 years



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

Description

α -Pyrrolidinopentiophenone (α -PVP) is an analog of the psychoactive stimulant pyrovalerone (Item No. 10836). α -PVP metabolite 1 is an analog of α -PVP in which the keto group is reduced. The biological properties of this compound have not been evaluated. This product is intended for forensic and research purposes.

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