

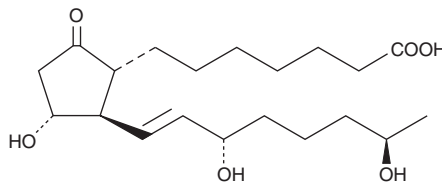
PRODUCT INFORMATION



PV9 (hydrochloride)

Item No. 15062

CAS Registry No.: 2749897-19-4
Formal Name: 1-phenyl-2-(pyrrolidin-1-yl)octan-1-one, monohydrochloride
Synonym: α -POP
MF: C₁₈H₂₇NO • HCl
FW: 309.9
Purity: \geq 98%
UV/Vis.: λ_{max} : 252 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 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. 9001083) is an analog of pyrovalerone (Item No. 10817), a known monoamine transport inhibitor that prevents the reuptake of dopamine and norepinephrine.^{1,2} PV9 is a novel α -PVP analog in which the pentyl side chain is elongated by three carbons. The biochemical, physiological, and toxicological properties of PV9 have not been determined. This product is intended for forensic and research applications.

References

1. Heron, C., Costentin, J., and Bonnet, J.J. Evidence that pure uptake inhibitors including cocaine interact slowly with the dopamine neuronal carrier. *Eur. J. Pharmacol.* **264**(3), 391-398 (1994).
2. Meltzer, P.C., Butler, D., Deschamps, J.R., *et al.* 1-(4-methylphenyl)-2-pyrrolidin-1-yl-pentan-1-one (pyrovalerone) analogs. A promising class of monoamine uptake inhibitors. *J. Med. Chem.* **49**(4), 1420-1432 (2006).

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, 06/13/2023

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