

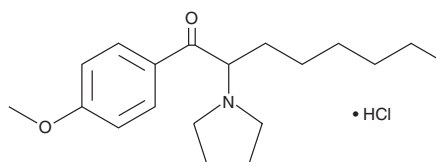
PRODUCT INFORMATION



4-methoxy PV9 (hydrochloride)

Item No. 15064

CAS Registry No.: 2748590-60-3
Formal Name: 1-(4-methoxyphenyl)-2-(pyrrolidin-1-yl) octan-1-one, monohydrochloride
Synonyms: MeO- α -POP, 4-MeO PV9, 4-methoxy α -POP, *para*-methoxy PV9
MF: C₁₉H₂₉NO₂ • HCl
FW: 339.9
Purity: ≥98%
UV/Vis.: λ_{max} : 225, 291 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

PV9 (Item No. 15062) is an analog of α -pyrrolidinopentiophenone (Item No. 9001083) in which the pentyl chain is elongated by three carbons. 4-methoxy PV9 differs from PV9 by having a methoxy group at the four, or *para*, position of the phenyl group. The physiological and toxicological properties of this compound are not known. This product is intended for forensic and research applications.

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