

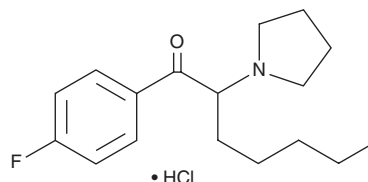
PRODUCT INFORMATION



4-fluoro PV8 (hydrochloride)

Item No. 15356

CAS Registry No.: 2748485-01-8
Formal Name: 1-(4-fluorophenyl)-2-(pyrrolidin-1-yl)heptan-1-one, monohydrochloride
Synonyms: 4-fluoro α -PHPP, *p*-fluoro α -PHPP, *p*-fluoro-PV8, *para*-fluoro α -PHPP, *para*-fluoro-PV8
MF: C₁₇H₂₄FNO • HCl
FW: 313.8
Purity: \geq 98%
UV/Vis.: λ_{max} : 254 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

PV8 (Item No. 14762) is an analog of the regulated psychoactive compound α -pyrrolidinopentiophenone. 4-fluoro PV8 is a *para*-fluoro analog of PV8. 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, 06/20/2023

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