

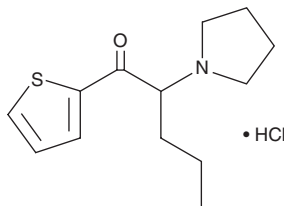
PRODUCT INFORMATION



α -Pyrrolidinopentiothiophenone (hydrochloride)

Item No. 14182

CAS Registry No.: 2748622-52-6
Formal Name: 2-(1-pyrrolidinyl)-1-(2-thienyl)-1-pentanone, monohydrochloride
Synonym: α -PVT
MF: C₁₃H₁₉NOS • HCl
FW: 273.8
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 268, 296 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

N-Pyrrolidinyl-cathinones, like pyrovalerone (Item No. 10817), are designer drugs which act as stimulants. Pyrovalerone is an inhibitor of the transporters for certain monoamine neurotransmitters, including dopamine and norepinephrine, preventing their uptake.^{1,2} α -Pyrrolidinopentiothiophenone is an analog of pyrovalerone in which the benzene ring has been replaced with a thiophene ring. Like pyrovalerone, this compound is expected to be a psychoactive stimulant. Its neurotoxic potential is unknown. 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, 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