

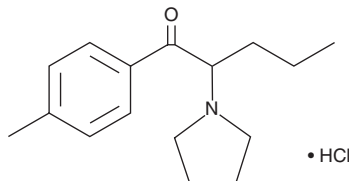
PRODUCT INFORMATION



Pyrovalerone (hydrochloride)

Item No. 10836

CAS Registry No.: 1147-62-2
Formal Name: 1-(4-methylphenyl)-2-(1-pyrrolidinyl)-1-pentanone, monohydrochloride
Synonyms: Centroton, Thymergix, Valerophenone
MF: C₁₆H₂₃NO • HCl
FW: 281.8
Purity: ≥95%
UV/Vis.: λ_{max}: 255 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

Pyrovalerone (hydrochloride) is an inhibitor of the transporters for certain monoamine neurotransmitters, including dopamine and norepinephrine, preventing their uptake.^{1,2} It has been used clinically for the treatment of chronic fatigue, but is used infrequently due to problems with abuse and dependence. Pyrovalerone is a DEA Schedule V controlled substance. Pyrovalerone (hydrochloride) is for use in the forensic analysis of samples that may contain this compound. This product, the hydrochloride salt of pyrovalerone, has superior solubility in aqueous solvents, compared to the free base (Item No. 10817).

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, 06/15/2023

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