

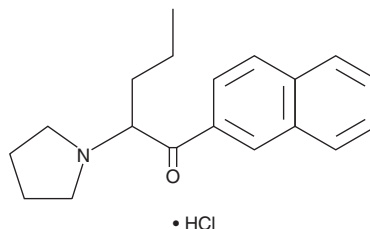
PRODUCT INFORMATION



Naphyrone (hydrochloride) (exempt preparation)

Item No. 16056

CAS Registry No.: 850352-11-3
Formal Name: 1-(2-naphthalenyl)-2-(1-pyrrolidinyl)-
1-pentanone, monohydrochloride
Synonyms: β -Naphyrone, Naphpyrovalerone,
NRG-1, O-2482
MF: $C_{19}H_{23}NO \cdot HCl$
FW: 317.9
Purity: $\geq 97\%$
UV/Vis.: λ_{max} : 253, 295 nm
Supplied as: A solution in methanol
Storage: $-20^{\circ}C$
Stability: ≥ 3 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

Pyrovalerone is a psychoactive compound with stimulatory effects. It is a controlled drug that has been found as a contaminant in products sold as bath salts.¹ Naphyrone (hydrochloride) is an analog of pyrovalerone that is characterized by the substitution of the methylphenyl group of pyrovalerone with a naphthyl group. Naphyrone potently inhibits dopamine, serotonin, and norepinephrine transporters (IC_{50} = 20, 33, and 136 nM, respectively).² Naphyrone, like pyrovalerone, has been detected in bath salt-type products. This compound is intended for forensic and research purposes.

References

1. Kikura-Hanajiri, R., Uchiyama, N., and Goda, Y. Survey of current trends in the abuse of psychotropic substances and plants in Japan. *Leg. Med. (Tokyo)* **13**(3), 109-15 (2011).
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, 01/15/2025

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