

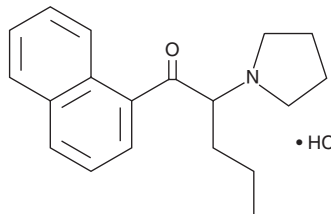
PRODUCT INFORMATION



Naphyrone 1-naphthyl isomer (hydrochloride)

Item No. 11240

CAS Registry No.: 1349245-31-3
Formal Name: 1-(naphthalen-1-yl)-2-(pyrrolidin-1-yl)pentan-1-one, monohydrochloride
Synonym: α -Naphyrone
MF: C₁₉H₂₃NO • HCl
FW: 317.9
Purity: \geq 95%
UV/Vis.: λ_{max} : 212, 242, 323 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

Naphyrone (Item No. 10517) is an analog of the psychoactive compound pyrovalerone, characterized by the substitution of the methylphenyl group of pyrovalerone with 2-naphthyl. Like pyrovalerone, naphyrone has been detected in products marketed as party pills or bath salts.¹ Naphyrone 1-naphthyl isomer (hydrochloride) (Item No. 11240) is a structural isomer of naphyrone, having the naphthyl group attached at the 1, rather than 2, position. The physiological and toxicological properties of this compound have not been reported. This product is intended for forensic and research applications.

Reference

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)**, 109-115 (2011).

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, 07/07/2023

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