

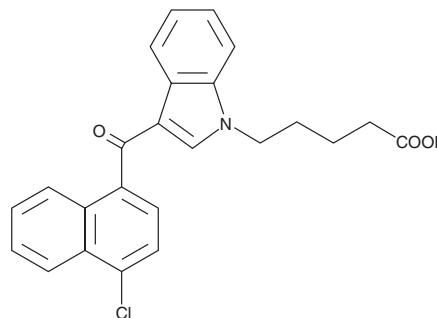
PRODUCT INFORMATION



JWH 398 N-pentanoic acid metabolite

Item No. 10942

CAS Registry No.: 1537889-10-3
Formal Name: 3-[[4-chloro-1-naphthalenyl]carbonyl]-1H-indole-1-pentanoic acid
Synonym: JWH 398 N-(5-carboxypentyl) metabolite
MF: C₂₄H₂₀ClNO₃
FW: 405.9
Purity: ≥98%
UV/Vis.: λ_{max}: 217, 317 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

JWH 398 (Item No. 13636) is a synthetic cannabinoid (CB) that acts as an agonist at both the central CB₁ receptor and the peripheral CB₂ receptor (K_is = 2.3 and 2.8 nM, respectively).¹ It has been detected as an additive of herbal mixtures.² JWH 398 N-pentanoic acid metabolite is an expected phase I metabolite of JWH 398. The biological and toxicological properties of this compound have not been characterized. This product is intended for forensic and research purposes.

References

1. Huffman, J.W. Cannabimimetic indoles, pyrroles, and indenes: Structure-activity relationships and receptor interactions. *The cannabinoid receptors*. Reggio, P.H., editor, *Humana Press* (2009).
2. 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).

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, 06/22/2023

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