

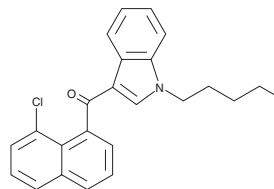
PRODUCT INFORMATION



JWH 398 8-chloronaphthyl isomer

Item No. 9001028

CAS Registry No.: 1366067-88-0
Formal Name: (8-chloronaphthalen-1-yl)(1-pentyl-1H-indol-3-yl) methanone
MF: C₂₄H₂₂NO
FW: 375.9
Purity: ≥98%
UV/Vis.: λ_{max}: 223, 311 nm
Supplied as: A solution in methanol
Storage: -20°C
Stability: ≥4 years



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

Description

JWH 398 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 8-chloronaphthyl isomer differs structurally from JWH 398 by having chlorine positioned on the naphthyl rings at the 8, rather than the 4, position. The biological and toxicological properties of this compound have not been characterized. This product is intended for forensic purposes.

References

- 1 Huffman, J.W. The cannabinoid receptors. Humana, New York (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, 06/20/2023

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