

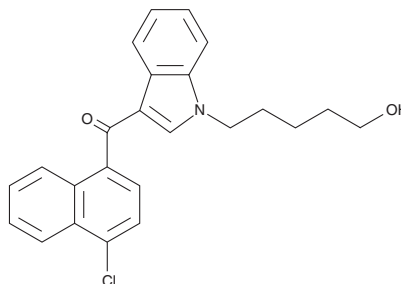
# PRODUCT INFORMATION



## JWH 398 N-(5-hydroxypentyl) metabolite

Item No. 9000770

**CAS Registry No.:** 1379604-69-9  
**Formal Name:** (4-chloro-1-naphthalenyl)[1-(5-hydroxypentyl)-1H-indol-3-yl]-methanone  
**MF:** C<sub>24</sub>H<sub>22</sub>ClNO<sub>2</sub>  
**FW:** 391.9  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 219, 315 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 is a synthetic cannabinoid (CB) that activates both CB<sub>1</sub> and CB<sub>2</sub> receptors (K<sub>i</sub> values of 2.3 and 2.8 nM, respectively).<sup>1</sup> It has been reported to be an adulterant of herbal products.<sup>2,3</sup> JWH 398 N-(5-hydroxypentyl) metabolite is a potential metabolite of JWH 398. Metabolism of structurally similar compounds leads to monohydroxylation of the N-alkyl chain, detectable in urine.<sup>4</sup> This product is intended for research and forensic applications.

### 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)**, 109-115 (2011).
3. Dresen, S., Ferreirós, N., Pütz, M., et al. Monitoring of herbal mixtures potentially containing synthetic cannabinoids as psychoactive compounds. *J. Mass Spectrom.* **45(10)**, 1186-94 (2010).
4. Sobolevsky, T., Prasolov, I., and Rodchenkov, G. Detection of JWH-018 metabolites in smoking mixture post-administration urine. *Forensic Sci. Int.* **200(1-3)**, 141-147 (2010).

#### 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