

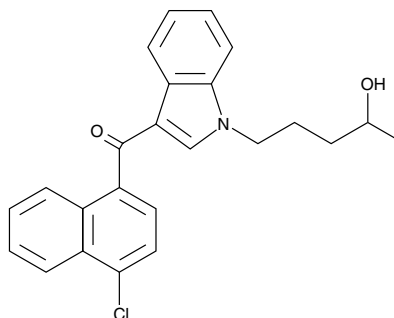
# Product Information



## JWH 398 N-(4-hydroxypentyl) metabolite

Item No. 10943

**CAS Registry No.:** 1537889-06-7  
**Formal Name:** (4-chloronaphthalen-1-yl)(1-(4-hydroxypentyl)-1H-indol-3-yl)methanone  
**MF:** C<sub>24</sub>H<sub>22</sub>ClNO<sub>2</sub>  
**FW:** 391.9  
**Purity:** ≥98%  
**Stability:** ≥1 year at -20°C  
**Supplied as:** A solution in methanol  
**UV/Vis.:** λ<sub>max</sub>: 218, 317 nm



### Laboratory Procedures

For long term storage, we suggest that JWH 398 N-(4-hydroxypentyl) metabolite be stored as supplied at -20°C. It should be stable for at least one year.

JWH 398 N-(4-hydroxypentyl) metabolite is supplied as a solution in methanol. To change the solvent, simply evaporate the methanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of JWH 398 N-(4-hydroxypentyl) metabolite in these solvents is approximately 15, 25, and 20 mg/ml, respectively.

JWH 398 N-(4-hydroxypentyl) metabolite is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

JWH 398 (Item No. 13636) is a synthetic cannabinoid (CB) that activates both central CB<sub>1</sub> and peripheral CB<sub>2</sub> receptors (K<sub>i</sub> = 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-(4-hydroxypentyl) metabolite is an expected phase I metabolite of JWH 398, detectable in serum and urine. While similar hydroxylated phase I metabolites of synthetic CBs retain activity, the physiological properties of this compound have yet to be determined.<sup>4,5</sup> This product is intended for research and forensic applications.

### 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).
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. Brents, L.K., Reichard, E.E., Zimmerman, M., *et al.* Phase I hydroxylated metabolites of the K2 synthetic cannabinoid JWH-018 retain *in vitro* and *in vivo* cannabinoid 1 receptor affinity and activity. *PLoS One* **6(7)**, 1-9 (2011).
5. Brents, L.K., Gallus-Zawala, A., Radomska-Pandya, A., *et al.* Monohydroxylated metabolites of the K2 synthetic cannabinoid JWH-073 retain intermediate to high cannabinoid 1 receptor (CB1R) affinity and exhibit neutral antagonist to partial agonist activity. *Biochem. Pharmacol.* **83**, 952-961 (2012).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/10943](http://www.caymanchem.com/catalog/10943)

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY. NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog**.

Copyright Cayman Chemical Company, 03/19/2014

### Cayman Chemical

#### Mailing address

1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

#### Phone

(800) 364-9897  
(734) 971-3335

#### Fax

(734) 971-3640

#### E-Mail

[custserv@caymanchem.com](mailto:custserv@caymanchem.com)

#### Web

[www.caymanchem.com](http://www.caymanchem.com)