

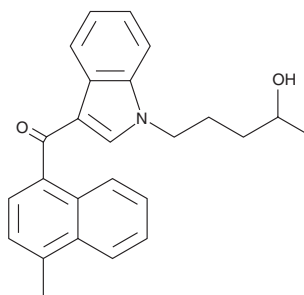
# PRODUCT INFORMATION



## JWH 122 N-(4-hydroxypentyl) metabolite

Item No. 11784

**CAS Registry No.:** 1537889-07-8  
**Formal Name:** [1-(4-hydroxypentyl)-1H-indol-3-yl]  
(4-methyl-1-naphthalenyl)-methanone  
**MF:** C<sub>25</sub>H<sub>25</sub>NO<sub>2</sub>  
**FW:** 371.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 222, 246, 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 122 is a synthetic cannabinoid (CB) that displays high affinities for both the central CB<sub>1</sub> receptor (K<sub>i</sub> = 0.69 nM) and the peripheral CB<sub>2</sub> receptor (K<sub>i</sub> = 1.2 nM).<sup>1,2</sup> JWH 122 N-(4-hydroxypentyl) metabolite is an expected phase I metabolite of JWH 122, detectable in serum and urine. While similar hydroxylated phase I metabolites of synthetic cannabinoids retain activity, the physiological properties of this compound have yet to be determined.<sup>3,4</sup> This product is intended for research and forensic applications.

### References

1. Huffman, J.W., Zengin, G., Wu, M.J., *et al.* Structure-activity relationships for 1-alkyl-3-(1-naphthoyl) indoles at the cannabinoid CB<sub>1</sub> and CB<sub>2</sub> receptors: Steric and electronic effects of naphthoyl substituents. New highly selective CB<sub>2</sub> receptor agonists. *Bioorg. Med. Chem.* **13(1)**, 89-112 (2005).
2. Huffman, J.W., Mabon, R., Wu, M.J., *et al.* 3-indolyl-1-naphthylmethanes: New cannabimimetic indoles provide evidence for aromatic stacking interactions with the CB<sub>1</sub> cannabinoid receptor. *Bioorgan. Med. Chem.* **11(4)**, 539-549 (2003).
3. 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)**, e21917 (2011).
4. 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(7)**, 952-961 (2012).

#### 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, 07/24/2023

#### 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