

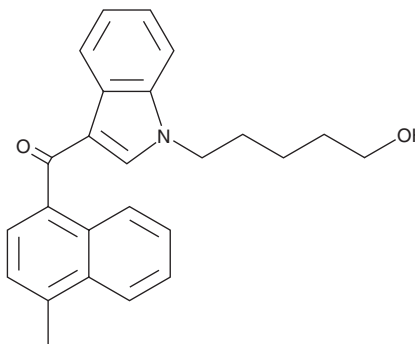
# PRODUCT INFORMATION



## JWH 122 N-(5-hydroxypentyl) metabolite

Item No. 10925

**CAS Registry No.:** 1379604-68-8  
**Formal Name:** [1-(5-hydroxypentyl)-1H-indol-3-yl](4-methyl-1-naphthalenyl)-methanone  
**Synonym:** MAM2201 N-(5-hydroxypentyl) metabolite  
**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, 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) of the naphthoylindole class that is structurally related to JWH 018 and JWH 073. It displays high-affinities for both CB<sub>1</sub> (K<sub>i</sub> = 0.69 nM) and CB<sub>2</sub> (K<sub>i</sub> = 1.2 nM) receptors.<sup>1,2</sup> JWH 122 N-(5-hydroxypentyl) metabolite is a metabolite of JWH 122 (Item No. 10591) that is characterized by monohydroxylation of the N-alkyl chain. Currently, there are no reports of its biological activity.

### References

- Ernst, L., Schiebel, H.M., Theuring, C., *et al.* Identification and characterization of JWH-122 used as new ingredient in "spice-like" herbal incenses. *Forensic Sci. Int.* **208(1-3)**, e31-e35 (2011).
- 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).

#### 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, 06/19/2023

#### 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](http://WWW.CAYMANCHEM.COM)