

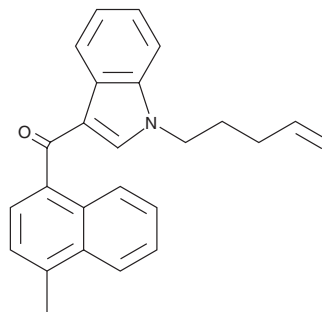
PRODUCT INFORMATION



JWH 122 N-(4-pentenyl) analog

Item No. 11611

CAS Registry No.: 1445577-68-3
Formal Name: (4-methylnaphthalen-1-yl)(1-(pent-4-en-1-yl)-1H-indol-3-yl)methanone
Synonyms: JWH 022 4-methylnaphthyl analog, MAM2201 N-(4-pentenyl) analog
MF: C₂₅H₂₃NO
FW: 353.5
Purity: ≥98%
UV/Vis.: λ_{max}: 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) that displays high affinities for both the central CB₁ receptor (K_i = 0.69 nM) and the peripheral CB₂ receptor (K_i = 1.2 nM).^{1,2} JWH 122 N-(4-pentenyl) analog is structurally related to JWH 122, differing only by the presence of a terminal double bond on the acyl chain. This compound is also structurally related to the cannabimimetic MAM2201. Analogs like this may be impurities or degradants that are found with, and serve as forensic markers for, the parent compounds. The physiological and toxicological properties of this analog are not known. This product is intended for forensic and research 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 CB₁ and CB₂ receptors: steric and electronic effects of naphthoyl substituents. New highly selective CB₂ receptor agonists. *Bioorg. Med. Chem.* **13**, 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 CB₁ cannabinoid receptor. *Bioorg. Med. Chem.* **11**, 539-549 (2003).

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, 07/18/2023

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