

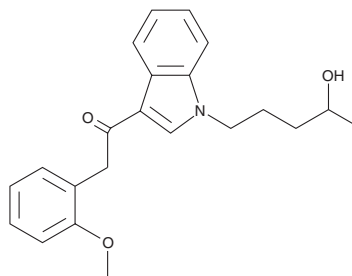
PRODUCT INFORMATION



JWH 250 N-(4-hydroxypentyl) metabolite

Item No. 10939

CAS Registry No.: 1427521-38-7
Formal Name: 1-(1-(4-hydroxypentyl)-1H-indol-3-yl)-2-(2-methoxyphenyl)ethanone
MF: C₂₂H₂₅NO₃
FW: 351.4
Purity: ≥98%
UV/Vis.: λ_{max}: 212, 246, 304 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 250 is a synthetic cannabinoid (CB) which potently activates both CB₁ and CB₂ receptors (K_i values of 11 and 33 nM, respectively).¹ It has been identified in herbal blends.² JWH 250 N-(4-hydroxypentyl) metabolite is expected to be a cytochrome P450 phase I metabolite of JWH 250, detectable both in serum and urine.³⁻⁵ This product is intended for research and forensic applications.

References

1. Huffman, J.W., Szklennik, P.V., Almond, A., *et al.* 1-Pentyl-3-phenylacetylindoles, a new class of cannabimimetic indoles. *Bioorg. Med. Chem. Lett.* **15**(18), 4110-4113 (2005).
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. Wintermeyer, A., Möller, I., Thevis, M., *et al.* In vitro phase I metabolism of the synthetic cannabimimetic JWH-018. *Anal. Bioanal. Chem.* **398**(5), 2141-2153 (2010).
4. Zhang, Q., Ma, P., Cole, R.B., *et al.* Identification of in vitro metabolites of JWH-015, an aminoalkylindole agonist for the peripheral cannabinoid receptor (CB₂) by HPLC-MS/MS. *Anal. Bioanal. Chem.* **386**(5), 1345-1355 (2006).
5. 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, 06/22/2023

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