

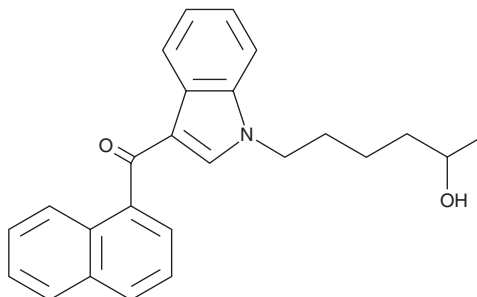
PRODUCT INFORMATION



JWH 019 N-(5-hydroxyhexyl) metabolite

Item No. 14814

CAS Registry No.: 1435934-47-6
Formal Name: [1-(5-hydroxyhexyl)-1H-indol-3-yl]-1-naphthalenyl-methanone
MF: C₂₅H₂₅NO₂
FW: 371.5
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 245, 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 019 (Item No. 13633) is a cannabimimetic indole that shows a high-affinity for both the central cannabinoid (CB₁) (K_i = 9.8 nM) and the peripheral cannabinoid (CB₂) (K_i = 5.6 nM) receptors.¹ JWH 019 N-(5-hydroxyhexyl) metabolite is an expected urinary metabolite of JWH 019 that should be detectable both in serum and urine based on studies of the metabolism of the structurally similar compound, JWH 018.¹⁻⁴ The physiological and toxicological properties of this compound have not been tested. This product is intended for forensic and research applications.

References

1. Aung, M.M., Griffin, G., Huffman, J.W., *et al.* Influence of the N-1 alkyl chain length of cannabimimetic indoles upon CB₁ and CB₂ receptor binding. *Drug Alcohol Depend.* **60(2)**, 133-140 (2000).
2. Teske, J., Weller, J.P., Fieguth, A., *et al.* Sensitive and rapid quantification of the cannabinoid receptor agonist naphthalen-1-yl-(1-pentylindol-3-yl)methanone (JWH-018) in human serum by liquid chromatography-tandem mass spectrometry. *J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.* **878(27)**, 2659-63 (2010).
3. 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).
4. 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).

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