

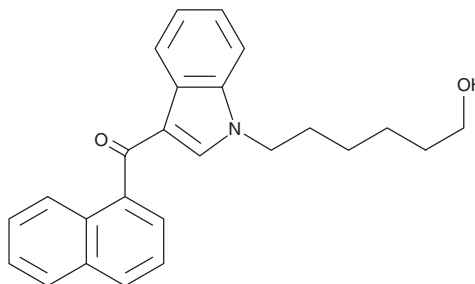
# PRODUCT INFORMATION



## JWH 019 N-(6-hydroxyhexyl) metabolite

Item No. 9000765

**CAS Registry No.:** 1435934-29-4  
**Formal Name:** [1-(6-hydroxyhexyl)-1H-indol-3-yl]-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>: 218, 247, 316 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

Cannabimimetic indoles, including certain JWH compounds, have been identified in herbal blends. JWH 018, a potent agonist of both the central cannabinoid (CB<sub>1</sub>) receptor and peripheral cannabinoid (CB<sub>2</sub>) receptor, has been found in many herbal blends and has been regulated in the United States and other countries. Studies indicate that JWH 018 is rapidly metabolized by the liver and that certain metabolites can be identified in the urine.<sup>1-3</sup> JWH 019 is nearly identical to JWH 018 in structure and activity at both CB receptors.<sup>4</sup> JWH 019 N-(6-hydroxyhexyl) metabolite is an expected metabolite of JWH 019, detectable both in serum and urine. Its biological activities have not been evaluated. This product is intended for research and forensic applications.

### References

1. 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).
2. 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).
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. Aung, M.M., Griffin, G., Huffman, J.W., *et al.* Influence of the N-1 alkyl chain length of cannabimimetic indoles upon CB<sub>1</sub> and CB<sub>2</sub> receptor binding. *Drug Alcohol Depend.* **60(2)**, 133-140 (2000).

#### 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, 03/08/2024

#### 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