

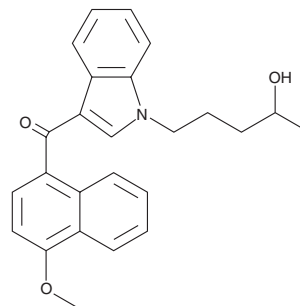
PRODUCT INFORMATION



JWH 081 N-(4-hydroxypentyl) metabolite

Item No. 14348

CAS Registry No.: 2748622-62-8
Formal Name: [1-(4-hydroxypentyl)-1H-indol-3-yl]
(4-methoxy-1-naphthalenyl)-methanone
MF: C₂₅H₂₅NO₃
FW: 387.5
Purity: ≥98%
UV/Vis.: λ_{max}: 212, 313 nm
Supplied as: A solution in acetonitrile
Storage: -20°C
Stability: ≥5 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

JWH 081 (Item No. 10579) is a cannabimimetic indole related to JWH 015 (Item No. 10009018) and JWH 018 (Item No. 10900) that shows preference for the central cannabinoid (CB₁) receptor (K_i = 1.2 nM) over the peripheral cannabinoid (CB₂) receptor (K_i = 12.4 nM).¹⁻² JWH 081 N-(4-hydroxypentyl) metabolite is an expected product of the phase I metabolism of JWH 081, based on studies on the metabolism of the related synthetic cannabinoids JWH 018 and JWH 073 (Item No. 10904).³⁻⁴ The physiological and toxicological properties of this compound are not known. This product is intended for research and forensic purposes.

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**, 133-140 (2000).
2. 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).
3. Wintermeyer, A., Möller, I., Thevis, M., *et al.* *In vitro* phase I metabolism of the synthetic cannabimimetic JWH-018. *Anal. Bioanal. Chem.* **398**, 2141-53 (2010).
4. Chimalakonda, K.C., Moran, C.L., Kennedy, P.D., *et al.* Solid-phase extraction and quantitative measurement of omega and omega-1 metabolites of JWH-018 and JWH-073 in human urine. *Anal. Chem.* **83**(16), 6381-6388 (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, 06/16/2023

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