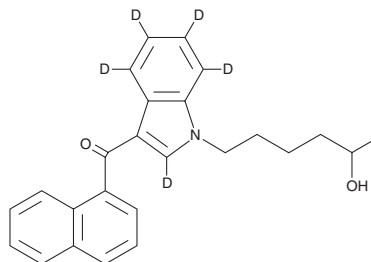


PRODUCT INFORMATION



JWH 019 N-(5-hydroxyhexyl) metabolite-d₅ Item No. 14687

CAS Registry No.: 2747917-84-4
Formal Name: (1-(5-hydroxyhexyl)-1H-indol-3-yl-2,4,5,6,7-d₅)
(naphthalen-1-yl)methanone
MF: C₂₅H₂₀D₅NO₂
FW: 376.5
Chemical Purity: ≥98% (JWH 019 N-(5-hydroxyhexyl) metabolite)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀
UV/Vis.: λ_{max}: 218, 247, 316 nm
Supplied as: A solution in methanol
Storage: -20°C
Stability: ≥4 years



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

Description

JWH 019 N-(5-hydroxyhexyl) metabolite-d₅ (Item No. 14687) is intended for use as an internal standard for the quantification of JWH 019 N-(5-hydroxyhexyl) metabolite (Item No. 14814) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Cannabimimetic indoles, including certain JWH compounds, have been identified in herbal blends. JWH 018, a potent agonist of both the central cannabinoid (CB₁) receptor and peripheral cannabinoid (CB₂) 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.¹⁻³ JWH 019 is nearly identical to JWH 018 in structure and activity at both CB receptors.⁴ JWH 019 N-(5-hydroxyhexyl) metabolite is an expected metabolite of JWH 019, detectable both in serum and urine. Its biological activities have not been evaluated.

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 CB1 and CB2 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, 06/21/2023

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