

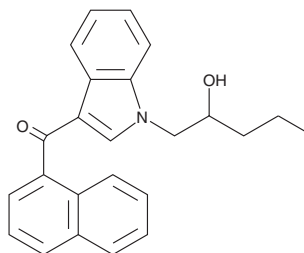
PRODUCT INFORMATION



(±)-JWH 018 N-(2-hydroxypentyl) metabolite

Item No. 11119

CAS Registry No.: 1936529-84-8
Formal Name: [1-(2-hydroxypentyl)-1H-indol-3-yl]-
1-naphthalenyl-methanone
MF: C₂₄H₂₃NO₂
FW: 357.5
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 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 018 N-(2-hydroxypentyl) metabolite is a potential metabolite of JWH 018. Monohydroxylation of the alkyl side chain of JWH 018 occurs through cytochrome P450 action, resulting in metabolites, like (±)-JWH 018 N-(2-hydroxypentyl) metabolite, that may be detectable in serum and, subsequently, in urine.¹⁻³ This product is intended for forensic purposes.

References

1. 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).
2. Sobolevsky, T., Prasolov, I., and Rodchenkov, G. Detection of JWH-018 metabolites in smoking mixture post-administration urine. *Forensic Sci. Int.* **200**, 141-147 (2010).
3. Moran, C.L., Le, V.H., Chimalakonda, K.C., *et al.* Quantitative measurement of JWH-018 and JWH-073 metabolites excreted in human urine. *Anal. Chem.* **83**(11), 4228-4236 (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, 07/07/2023

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