

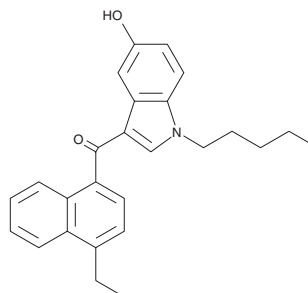
PRODUCT INFORMATION



JWH 210 5-hydroxyindole metabolite

Item No. 9000771

CAS Registry No.: 1427325-81-2
Formal Name: (4-ethyl-1-naphthalenyl)(5-hydroxy-1-pentyl-1H-indol-3-yl)-methanone
MF: C₂₆H₂₇NO₂
FW: 385.5
Purity: ≥98%
UV/Vis.: λ_{max}: 221, 256, 282, 323 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 210 is a potent cannabimimetic alkylindole that has been identified in extracts from herbal blends. Structurally and functionally similar alkylindoles are rapidly metabolized by the liver and secreted in the urine, with 5-hydroxyindole metabolites detected both in liver secretions and in urine.¹⁻³ JWH 210 5-hydroxyindole metabolite is an expected metabolite of JWH 210, detectable both in serum and urine. Its biological activities have yet to be determined. 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).

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