

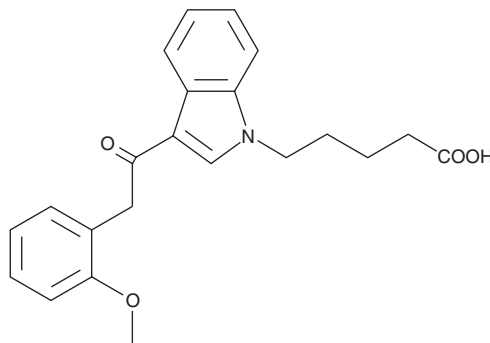
PRODUCT INFORMATION



JWH 250 N-pentanoic acid metabolite

Item No. 10938

CAS Registry No.: 1379604-65-5
Formal Name: 3-[2-(2-methoxyphenyl)acetyl]-1H-indole-1-pentanoic acid
Synonym: JWH 250 N-(5-carboxypentyl) metabolite
MF: C₂₂H₂₃NO₄
FW: 365.4
Purity: ≥98%
UV/Vis.: λ_{max}: 213, 245, 304 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.¹ Metabolites of two closely-related compounds, JWH 015 and JWH 018, have been identified from *in vitro* liver microsomal metabolism and from urine, respectively.^{2,3} JWH 250 is a cannabimimetic indole that is structurally- and functionally-related to JWH 015 and JWH 018. JWH 250 N-pentanoic acid metabolite is expected to be a metabolite of JWH 250 that would be detectable both in serum and in urine.

References

1. Uchiyama, N., Kikura-Hanajiri, R., Kawahara, N., *et al.* Identification of a cannabimimetic indole as a designer drug in a herbal product. *Forensic Toxicol.* **27(2)**, 61-66 (2009).
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. Zhang, Q., Ma, P., Cole, R.B., *et al.* Identification of *in vitro* metabolites of JWH-015, an aminoalkylindole agonist for the peripheral cannabinoid receptor (CB₂) by HPLC-MS/MS. *Anal. Bioanal. Chem.* **386(5)**, 1345-1355 (2006).

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, 06/22/2023

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