

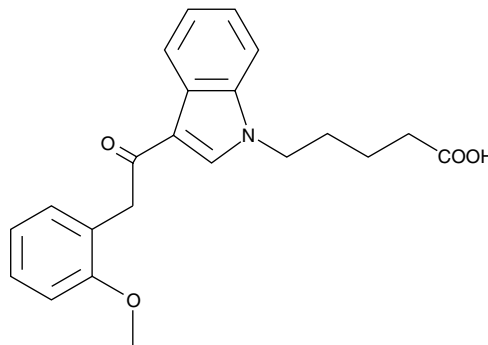
Product Information



JWH 250 N-pentanoic acid metabolite

Item No. 10938

CAS Registry No.: 1379604-65-5
Formal Name: 5-(3-(2-(2-methoxyphenyl)acetyl)-1H-indol-1-yl)pentanoic acid
Synonym: JWH 250 N-(5-carboxypentyl) metabolite
MF: C₂₂H₂₃NO₄
FW: 365.4
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 213, 245, 304 nm



Laboratory Procedures

For long term storage, we suggest that JWH 250 N-pentanoic acid metabolite be stored as supplied at -20°C. It should be stable for at least two years.

JWH 250 N-pentanoic acid metabolite is supplied as a crystalline solid. A stock solution may be made by dissolving the JWH 250 N-pentanoic acid metabolite in the solvent of choice. JWH 250 N-pentanoic acid metabolite is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of JWH 250 N-pentanoic acid metabolite in these solvents is approximately 20 mg/ml.

JWH 250 N-pentanoic acid metabolite is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, JWH 250 N-pentanoic acid metabolite should first be dissolved in DMF and then diluted with the aqueous buffer of choice. JWH 250 N-pentanoic acid metabolite has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

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**, 61-66 (2009).
2. 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**, 1345-1355 (2006).
3. 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).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10938

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog.**

Copyright Cayman Chemical Company, 07/12/2013

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com