

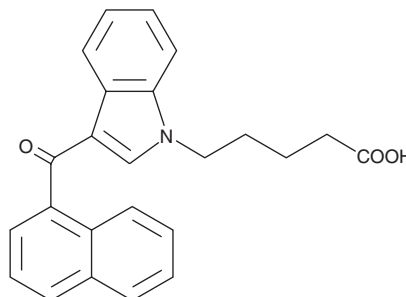
PRODUCT INFORMATION



JWH 018 N-pentanoic acid metabolite

Item No. 9000856

CAS Registry No.: 1254475-87-0
Formal Name: 3-(1-naphthalenylcarbonyl)-1H-indole-1-pentanoic acid
MF: C₂₄H₂₁NO₃
FW: 371.4
Purity: ≥95%
UV/Vis.: λ_{max}: 218, 316 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 is a mildly selective agonist of the peripheral cannabinoid (CB₂) receptor, derived from the aminoalkylindole WIN 55,212-2. The K_i values for binding central cannabinoid (CB₁) and CB₂ receptors are 9.0 and 2.94 nM, respectively, for a CB₁:CB₂ ratio of 3.06.¹ JWH 018 is one of several synthetic CBs which have been included in smoking mixtures. JWH 018 N-pentanoic acid metabolite is a minor urinary metabolite of JWH 018, characterized by carboxylation of the N-alkyl chain.² In urine samples, this metabolite is almost completely glucuronidated.² This product is intended for research and forensic applications.

References

1. 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).
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).

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