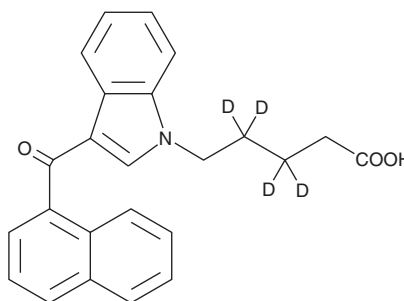


PRODUCT INFORMATION



JWH 018 N-pentanoic acid metabolite-d₄ Item No. 9000867

CAS Registry No.: 1320363-49-2
Formal Name: 3-(1-naphthalenylcarbonyl)-1H-Indole-1-pentanoic-β,β,γ,γ-d₄ acid
MF: C₂₄H₁₇D₄NO₃
FW: 375.5
Chemical Purity: ≥98% (JWH 018 N-pentanoic acid metabolite)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₄); ≤1% d₀
UV/Vis.: λ_{max}: 219, 247, 316 nm
Supplied as: A solution in methanol
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

JWH 018 N-pentanoic acid metabolite-d₄ (Item No. 9000867) is intended for use as an internal standard for the quantification of JWH 018 N-pentanoic acid metabolite (Item No. 9000856) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

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, 07/10/2023

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