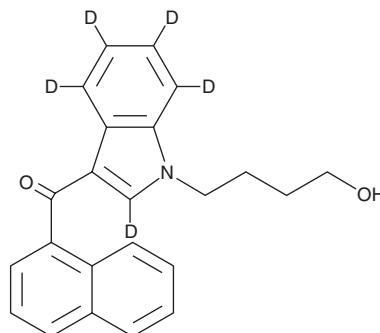


PRODUCT INFORMATION



JWH 073 N-(4-hydroxybutyl) metabolite-d₅ Item No. 10934

CAS Registry No.: 2484976-96-5
Formal Name: (1-(4-hydroxybutyl)-1H-indol-3-yl-2',4',5',6',7'-d₅)
(naphthalen-1-yl)-methanone
MF: C₂₃H₁₆D₅NO₂
FW: 348.5
Chemical Purity: ≥98% (JWH 073 N-(4-hydroxybutyl) metabolite)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀
UV/Vis.: λ_{max}: 218, 247, 314 nm
Supplied as: A 1 mg/ml 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 073 N-(4-hydroxybutyl) metabolite-d₅ (Item No. 10934) is intended for use as an internal standard for the quantification of JWH 073 N-(4-hydroxybutyl) metabolite (Item No. 9000865) 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 073 is a mildly selective agonist of the central cannabinoid (CB₁) receptor derived from the aminoalkylindole WIN 55,212-2. The K_i values for binding CB₁ and the peripheral cannabinoid (CB₂) receptor are 8.9 and 38 nM, respectively for a CB₁:CB₂ ratio of 0.23.¹ JWH 073 is one of several synthetic CBs which have been included in smoking mixtures. JWH 073 N-(4-hydroxybutyl) metabolite is expected to be a major urinary metabolite of JWH 073, characterized by monohydroxylation of the N-alkyl chain, based on the metabolism of the closely-related JWH 015 and JWH 018.^{2,3} In urine samples, this metabolite of JWH 018 is almost completely glucuronidated.³

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. 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).
3. 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, 11/20/2025

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