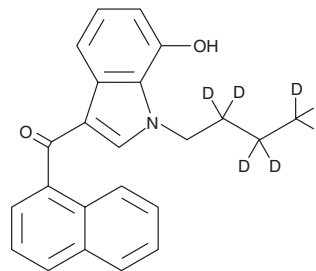


PRODUCT INFORMATION



JWH 073 7-hydroxyindole metabolite-d₇ Item No. 10720

CAS Registry No.: 2749316-88-7
Formal Name: (1-butyl-2,2,3,3,4,4,4-d₇-7-hydroxy-1H-indol-3-yl)-1-naphthalenyl-methanone
MF: C₂₃H₁₄D₇NO₂
FW: 350.5
Purity: ≥99% deuterated forms (d₁-d₇); ≤1% d₀
UV/Vis.: λ_{max}: 218, 338 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.

Laboratory Procedures

JWH 073 7-hydroxyindole metabolite-d₇ is supplied as a solution in methanol. A stock solution may be made by dissolving the JWH 073 7-hydroxyindole metabolite-d₇ in the solvent of choice, which should be purged with an inert gas. JWH 073 7-hydroxyindole metabolite-d₇ is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of JWH 073 7-hydroxyindole metabolite-d₇ in these solvents is approximately 20 mg/ml.

Description

JWH 073 7-hydroxyindole metabolite-d₇ contains seven deuterium atoms at the 2, 2', 3, 3', 4, 4, and 4 positions. It is intended for use as an internal standard for the quantification of JWH 073 7-hydroxyindole metabolite by GC- or LC-mass spectrometry (MS). 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 7-hydroxyindole metabolite is expected to be a urinary metabolite of JWH 073 based on the metabolism of the closely-related JWH 015 and JWH 018.^{2,3}

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**, 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**, 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).

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, 07/18/2023

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