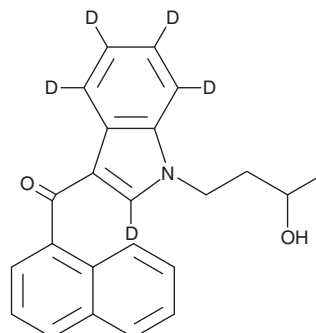


PRODUCT INFORMATION



(±)-JWH 073 N-(3-hydroxybutyl) metabolite-d₅ Item No. 10927

CAS Registry No.: 1413427-47-0
Formal Name: (1-(3-hydroxybutyl)-1H-indol-3-yl)(naphthalen-1-yl)-methanone-2,4,5,6,7-d₅
MF: C₂₃H₁₆D₅NO₂
FW: 348.5
Chemical Purity: ≥98% ((±)-JWH 073 N-(3-hydroxybutyl) metabolite)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀
UV/Vis.: λ_{max}: 218, 246, 315 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 073 N-(3-hydroxybutyl) metabolite-d₅ is intended for use as an internal standard for the quantification of (±)-JWH 073 N-(3-hydroxybutyl) metabolite by GC- or LC-mass spectrometry (MS). JWH 073 is a potent cannabimimetic that has been identified as an adulterant in smoking mixtures. JWH 073 N-(3-hydroxybutyl) metabolite is expected to be produced by hepatic metabolism of JWH 073.^{1,2} The biological actions of this metabolite are unknown.

References

1. Wintermeyer, A., Möller, I., Thevis, M., *et al.* In vitro phase I metabolism of the synthetic cannabimimetic JWH-018. *Anal. Bioanal. Chem.* **398**(5), 2141-2153 (2010).
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).

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, 06/13/2023

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