

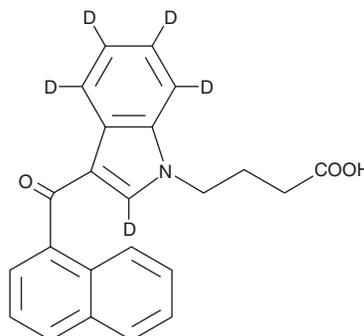
PRODUCT INFORMATION



JWH 073 N-butanoic acid metabolite-d₅

Item No. 9000870

CAS Registry No.: 1320363-50-5
Formal Name: 4-(3-(1-naphthoyl)-1H-indol-1-yl)-2,4,5,6,7-d₅-butanoic acid
MF: C₂₃H₁₄D₅NO₃
FW: 362.4
Chemical Purity: ≥98% (JWH 073 N-butanoic acid metabolite)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀
UV/Vis.: λ_{max}: 218, 246, 314 nm
Supplied as: A solution in methanol
Storage: -20°C
Stability: ≥1 year



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

Description

JWH 073 N-butanoic acid metabolite-d₅ (Item No. 9000870) is intended for use as an internal standard for the quantification of JWH 073 N-butanoic acid metabolite (Item No. 9000866) 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 N-butanoic acid metabolite, characterized by carboxylation of the N-alkyl chain, is expected to be a urinary metabolite of JWH 073 based on the metabolism of the closely-related JWH 015 and JWH 018.^{1,2} This product is intended for research and forensic applications.

References

1. 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).
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