

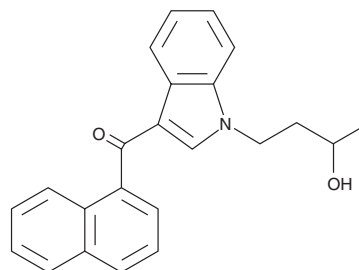
PRODUCT INFORMATION



(±)-JWH 073 N-(3-hydroxybutyl) metabolite

Item No. 10795

CAS Registry No.: 1320363-48-1
Formal Name: (1-(3-hydroxybutyl)-1H-indol-3-yl)
(naphthalen-1-yl)-methanone
MF: C₂₃H₂₁NO₂
FW: 343.4
Purity: ≥98%
UV/Vis.: λ_{max}: 219, 247, 316 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

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 during the metabolism of JWH 073 by human liver microsomes *in vitro*, generated by the oxidation of the aminoalkyl chain.^{1,2} The biological actions of this metabolite are unknown. This product is intended for research and forensic applications.

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, 06/15/2023

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