

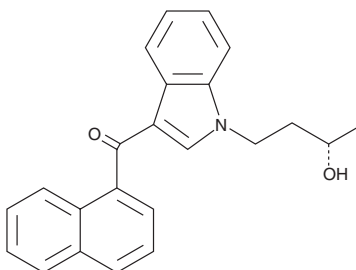
PRODUCT INFORMATION



(S)-(+)-JWH 073 N-(3-hydroxybutyl) metabolite

Item No. 10898

CAS Registry No.: 2748589-83-3
Formal Name: [1-[(3S)-3-hydroxybutyl]-1H-indol-3-yl]-1-naphthalenyl-methanone
MF: C₂₃H₂₁NO₂
FW: 343.4
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 247, 315 nm
Supplied as: A solution in methanol
Storage: -20°C
Stability: ≥2 years



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

Description

(S)-(+)-JWH 073 N-(3-hydroxybutyl) metabolite (Item No. 10898) is an analytical reference standard categorized as a synthetic cannabinoid metabolite.¹ It is a metabolite of JWH 073 (Item No. 10904). This product is intended for research and forensic applications.

Reference

1. Chimalakonda, K.C., Moran, C.L., Kennedy, P.D., *et al.* Solid-phase extraction and quantitative measurement of omega and omega-1 metabolites of JWH-018 and JWH-073 in human urine. *Anal. Chem.* **83**(16), 6381-6388 (2011).

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, 07/21/2023

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