

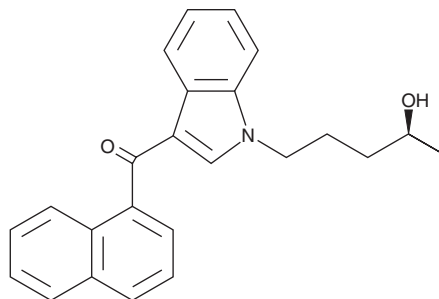
PRODUCT INFORMATION



(S)-(+)-JWH 018 N-(4-hydroxypentyl) metabolite

Item No. 10920

CAS Registry No.: 1454924-73-2
Formal Name: (S)-1-(4-hydroxypentyl)-1H-indol-3-yl(naphthalen-1-yl)methanone
MF: C₂₄H₂₃NO₂
FW: 357.5
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 247, 315 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

(S)-(+)-JWH 018 N-(4-hydroxypentyl) metabolite (Item No. 10920) is a purified enantiomer of JWH 018 N-(4-hydroxypentyl) metabolite (Item No. 10796). Its biological properties have not been examined. This product is intended for forensic purposes.

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, 07/15/2019

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