

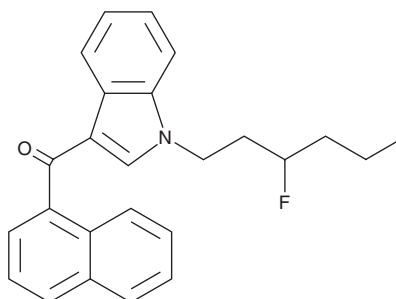
PRODUCT INFORMATION



JWH 019 N-(3-fluorohexyl) isomer

Item No. 15443

CAS Registry No.: 2365471-37-8
Formal Name: [1-(3-fluorohexyl)-1H-indol-3-yl]-1-naphthalenyl-methanone
Synonym: JWH 019 N-(3-fluorohexyl) analog
MF: C₂₅H₂₄FNO
FW: 373.5
Purity: ≥98%
UV/Vis.: λ_{max}: 219, 245, 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 019 (Item No. 13633) is the N-hexyl homolog of the aminoalkylindole cannabimimetic JWH 018 (Item No. 10900). JWH 019 N-(3-fluorohexyl) isomer is a homolog of JWH 019 that has a fluorine atom at the three position of the hexyl chain. The physiological and toxicological properties of this compound are not known. This product is intended for forensic and research applications.

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, 06/28/2023

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