

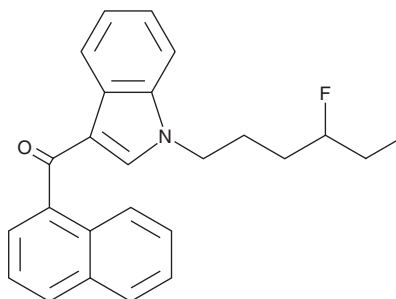
PRODUCT INFORMATION



JWH 019 N-(4-fluorohexyl) isomer

Item No. 15444

CAS Registry No.: 2365471-36-7
Formal Name: [1-(4-fluorohexyl)-1H-indol-3-yl]-1-naphthalenyl-methanone
Synonym: JWH 019 N-(4-fluorohexyl) analog
MF: C₂₅H₂₄FNO
FW: 373.5
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 247, 314 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 a cannabimimetic aminoalkylindole that shows a high-affinity for both the central cannabinoid (CB1) and the peripheral cannabinoid (CB2) receptors.¹ JWH 019 N-(4-fluorohexyl) isomer is an isomer of JWH 019 that has a fluorine atom at the four position of the alkyl chain. The physiological and toxicological properties of this compound are not known. This product is intended for forensic and research applications.

Reference

1. Aung, M.M., Griffin, G., Huffman, J.W., *et al.* Influence of the N-1 alkyl chain length of cannabimimetic indoles upon CB₁ and CB₂ receptor binding. *Drug Alcohol Depend.* **60(2)**, 133-140 (2000).

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