

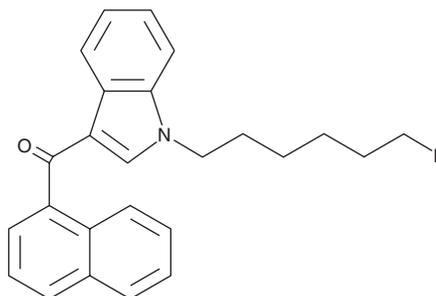
# PRODUCT INFORMATION



## JWH 019 N-(6-fluorohexyl) isomer

Item No. 15446

**CAS Registry No.:** 2365471-82-3  
**Formal Name:** [1-(6-fluorohexyl)-1H-indol-3-yl]-1-naphthalenyl-methanone  
**Synonym:** JWH 019 N-(6-fluorohexyl) analog  
**MF:** C<sub>25</sub>H<sub>24</sub>FNO  
**FW:** 373.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 218, 248, 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

JWH 019 (Item No. 13633) is the N1-hexyl homolog of the cannabimimetic indole JWH 018 (Item No. 10900). It shows a high-affinity for both the central cannabinoid (CB<sub>1</sub>) (K<sub>i</sub> = 9.8 nM) and the peripheral cannabinoid (CB<sub>2</sub>) (K<sub>i</sub> = 5.6 nM) receptors.<sup>1</sup> JWH 019 N-(6-fluorohexyl) isomer differs from JWH 019 by having a fluorine at the 6 position of the hexyl chain. The physiological and toxicological properties of this compound have not been tested. 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 CB<sub>1</sub> and CB<sub>2</sub> 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, 07/29/2022

#### 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