

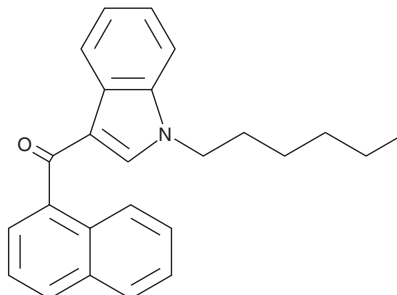
PRODUCT INFORMATION



JWH 019 (exempt preparation)

Item No. 15664

CAS Registry No.: 209414-08-4
Formal Name: (1-hexyl-1H-indol-3-yl)-1-naphthalenyl-methanone
MF: C₂₅H₂₅NO
FW: 355.5
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 247, 315 nm
Supplied as: A 1 mg/ml solution in methanol
Storage: -20°C
Stability: ≥3 years



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

Description

JWH 019 is a cannabimimetic indole that shows a high-affinity for both the central cannabinoid (CB₁) and the peripheral cannabinoid (CB₂) receptors. The affinities of JWH 019 for both CB₁ (K_i = 9.8 nM) and CB₂ (K_i = 5.6 nM) are superior to those of the traditional cannabinoid, Δ⁹-THC (Item No. 12068) (K_i values of 40.7 and 36.4 nM for CB₁ and CB₂, respectively).¹ JWH 019 (exempt preparation) (Item No. 15664) is provided as a DEA exempt preparation. This product is intended for research and forensic 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, 02/14/2025

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