

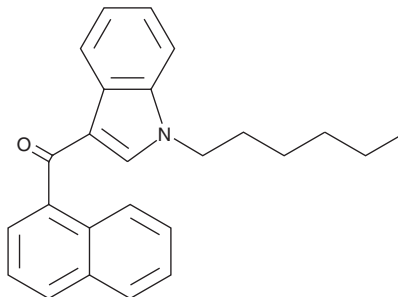
Product Information



JWH 019

Item No. 13633

CAS Registry No.: 209414-08-4
Formal Name: (1-hexyl-1H-indol-3-yl)-1-naphthalenyl-methanone
MF: C₂₅H₂₅NO
FW: 355.5
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that JWH 019 be stored as supplied at -20°C. It should be stable for at least two years.

JWH 019 is supplied as a crystalline solid. A stock solution may be made by dissolving the JWH 019 in the solvent of choice. JWH 019 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of JWH 019 in these solvents is approximately 2.5, 30, and 20 mg/ml, respectively.

If aqueous stock solutions are required for biological experiments, they can best be prepared by diluting the organic solvent into aqueous buffers or isotonic saline. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

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-tetrahydrocannabinol (Δ⁹-THC; K_i values of 40.7 and 36.4 nM, for CB₁ and CB₂, respectively).¹

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 and Alcohol Dependence* **60**, 133-140 (2000).

Related Products

AM694 - Item No. 10567 • JWH 081 - Item No. 10579 • JWH 222 - Item No. 10591 • RCS-8 - Item No. 10636 • KM 233 - Item No. 10640 • RCS-4 - Item No. 10645 • JWH 098 - Item No. 10680 • JWH 302 - Item No. 10722 • JWH 016 - Item No. 10849 • JWH 020 - Item No. 10850 • JWH 018 - Item No. 13169 • JWH 073 - Item No. 13170 • JWH 200 - Item No. 13171 • JWH 250 - Item No. 13634 • JWH 398 - Item No. 13636 • JWH 019 5-hydroxyindole metabolite - Item No. 9000764 • AM630 - Item No. 10006974 • JWH 015 - Item No. 10009018

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY; NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 09/19/2011

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com