

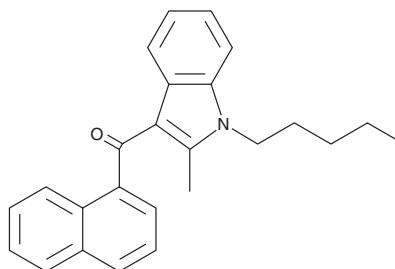
PRODUCT INFORMATION



JWH 007

Item No. 10266

CAS Registry No.: 155471-10-6
Formal Name: (2-methyl-1-pentyl-1H-indol-3-yl)-1-naphthalenyl-methanone
MF: C₂₅H₂₅NO
FW: 355.5
Purity: ≥98%
UV/Vis.: λ_{max}: 219, 320 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.

Laboratory Procedures

JWH 007 is supplied as a solution in methanol. To change the solvent, simply evaporate the methanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of JWH 007 in these solvents is approximately 20 mg/ml.

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.

Description

JWH 007 is a potent cannabinoid (CB) receptor agonist that avidly binds to both CB₁ and CB₂ (K_i = 9.5 and 2.9 nM, respectively).^{1,2} This compares favorably with the binding of Δ⁹-THC, which binds CB₁ and CB₂ with K_i values of 41 and 36 nM, respectively.² Similarly, JWH 007 performs comparably to Δ⁹-THC in mouse studies on spontaneous activity, antinociception, hypothermia, and catalepsy.¹

References

1. Huffman, J.W., Dai, D., Martin, B.R., *et al.* Design, synthesis and pharmacology of cannabimimetic indoles. *Bioorg. Med. Chem. Lett.* **4**(4), 563-566 (1994).
2. Huffman, J.W., Zengin, G., Wu, M.-J., *et al.* Structure-activity relationships for 1-alkyl-3-(1-naphthoyl) indoles at the cannabinoid CB₁ and CB₂ receptors: steric and electronic effects of naphthoyl substituents. New highly selective CB₂ receptor agonists. *Bioorg. Med. Chem.* **13**, 89-112 (2005).

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, 07/18/2023

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