

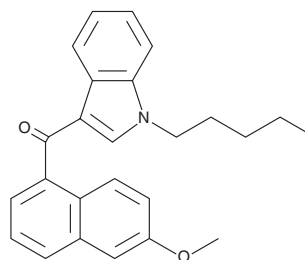
PRODUCT INFORMATION



JWH 081 6-methoxynaphthyl isomer

Item No. 9001047

CAS Registry No.: 824961-41-3
Formal Name: (6-methoxy-1-naphthalenyl)(1-pentyl-1H-indol-3-yl)-methanone
MF: C₂₅H₂₅NO₂
FW: 371.5
Purity: ≥98%
UV/Vis.: λ_{max}: 225, 317 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Description

JWH 081 is a synthetic cannabinoid (CB) that shows a high-affinity for the central cannabinoid (CB₁) receptor with a K_i value of 1.2 nM and ten-fold reduced affinity for the peripheral cannabinoid (CB₂) receptor (K_i = 12.4 nM).^{1,2} JWH 081 6-methoxynaphthyl isomer differs structurally from JWH 081 by having the methoxy group attached to the naphthyl rings at the 6 position, instead of the 4 position. The biological and toxicological properties of this compound have not been characterized. This product is intended for forensic and research purposes.

References

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).
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(1)**, 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, 09/29/2025

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