

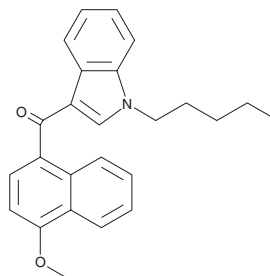
PRODUCT INFORMATION



JWH 081 (exempt preparation)

Item No. 15665

CAS Registry No.: 210179-46-7
Formal Name: (4-methoxy-1-naphthalenyl)(1-pentyl-1H-indol-3-yl)-methanone
MF: C₂₅H₂₅NO₂
FW: 371.5
Purity: ≥98%
UV/Vis.: λ_{max}: 213, 318 nm
Supplied as: A 1 mg/ml solution in methanol
Storage: -20°C
Stability: ≥2 years



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

Description

JWH 081 (exempt preparation) (Item No. 15665) is an analytical reference standard categorized a synthetic cannabinoid.^{1,2} JWH 081 has been found in Spice/K2-type herbal blends and may have neurotoxic properties.³ JWH 081 is regulated as a Schedule I compound in the United States. JWH 081 (exempt preparation) (Item No. 15665) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

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).
3. Cha, H.J., Seong, Y.H., Song, M.J., *et al.* Neurotoxicity of synthetic cannabinoids JWH-081 and JWH-210. *Biomol. Ther. (Seoul)* **23(6)**, 597-603 (2015).

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, 09/20/2017

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