

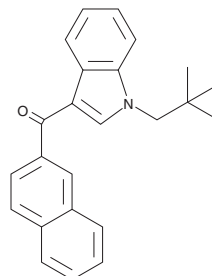
PRODUCT INFORMATION



JWH 018 2'-naphthyl-N-(2,2-dimethylpropyl) isomer

Item No. 9001008

CAS Registry No.: 1869954-38-0
Formal Name: [1-(2,2-dimethylpropyl)-1H-indol-3-yl]-2-naphthalenyl-methanone
MF: C₂₄H₂₃NO
FW: 341.5
Purity: ≥98%
UV/Vis.: λ_{max}: 216, 252, 325 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

JWH 018 (1-pentyl-3-(1-naphthoyl)indole) is a synthetic cannabinoid (CB) that potently activates the central CB₁ and peripheral CB₂ receptors (K_i = 9.0 and 2.94 nM, respectively).¹ This cannabimimetic compound has frequently been found in herbal blends.²⁻⁴ JWH 018 2'-naphthyl-N-(2,2-dimethylpropyl) isomer differs structurally from JWH 018 by having the naphthyl group attached at the 2' position and 2,2-dimethylpropyl in place of the pentyl chain. The biological activities of this compound have not been characterized.

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. Auwärter, V., Dresen, S., Weinmann, W., *et al.* 'Spice' and other herbal blends: Harmless incense or cannabinoid designer drugs? *J. Mass Spectrom.* **4(5)**, 832-837 (2009).
3. Lindigkeit, R., Boehme, A., Eiserloh, I., *et al.* Spice: A never ending story? *Forensic Sci. Int.* **191(1-3)**, 58-63 (2009).
4. Uchiyama, N., Kikura-Hanajiri, R., Kawahara, N., *et al.* Identification of a cannabimimetic indole as a designer drug in a herbal product. *Forensic Toxicol.* **27(2)**, 61-66 (2009).

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, 12/13/2016

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