

PRODUCT INFORMATION



JWH 018 2'-naphthyl-N-(1,1-dimethylpropyl) isomer

Item No. 9001007

CAS Registry No.: 1869951-99-4

Formal Name: [1-(1,1-dimethylpropyl)-1H-indol-3-yl]-2-naphthalenyl-methanone

MF: C₂₄H₂₃NO

FW: 341.5

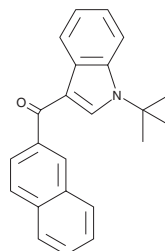
Purity: ≥98%

UV/Vis.: λ_{max}: 216, 252, 325 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: As supplied, 2 years from the QC date provided on the Certificate of Analysis, when stored properly



Description

JWH 018 is a synthetic cannabinoid (CB) that potently activates the central cannabinoid (CB₁) and peripheral cannabinoid (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-(1,1-dimethylpropyl) isomer differs from JWH 018 structurally by having a 1,1-dimethylpropyl chain, rather than a pentyl chain, extending from the indole group and the naphthyl moiety attached at its 2 position. The biological activities of this compound have not been characterized. This product is intended for forensic 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**, 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.* **44**, 832-837 (2009).
3. Lindigkeit, R., Boehme, A., Eiserloh, I., *et al.* Spice: A never ending story? *Forensic Sci. Int.* **191**, 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**, 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