

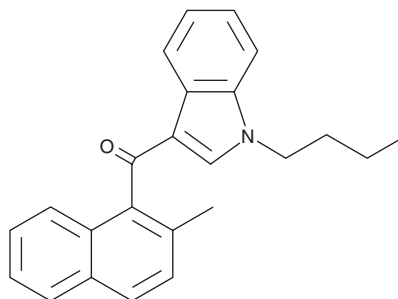
PRODUCT INFORMATION



JWH 073 2-methylnaphthyl analog

Item No. 10848

CAS Registry No.: 1427325-61-8
Formal Name: (1-butyl-1H-indol-3-yl)
(2-methylnaphthalen-1-yl)methanone
MF: C₂₄H₂₃NO
FW: 341.5
Purity: ≥98%
UV/Vis.: λ_{max}: 223, 249, 315 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.

Description

JWH 073 is a synthetic cannabinoid (CB) that activates both the CB₁ and the CB₂ receptor (K_i = 8.9 and 38 nM, respectively).¹ This cannabimimetic compound has been identified in herbal blends.²⁻⁴ JWH 073 2-methylnaphthyl analog (Item No. 10848) differs structurally from JWH 073 by having a methyl group added at the 2 position of the naphthyl moiety. The biological and toxicological properties of this compound have not been characterized. This compound 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. Kikura-Hanajiri, R., Uchiyama, N., and Goda, Y. Survey of current trends in the abuse of psychotropic substances and plants in Japan. *Leg. Med. (Tokyo)* **13(3)**, 109-115 (2011).
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., Ogata, J., *et al.* Chemical analysis of synthetic cannabinoids as designer drugs in herbal products. *Forensic Sci. Int.* **198(1-3)**, 31-38 (2010).

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, 06/16/2023

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