

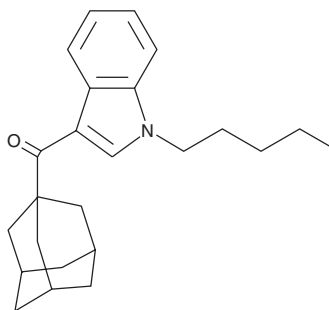
PRODUCT INFORMATION



JWH 018 adamantyl analog

Item No. 9000799

CAS Registry No.: 1345973-49-0
Formal Name: (1-pentyl-1H-indol-3-yl)tricyclo[3.3.1.1^{3,7}]dec-1-yl-methanone
Synonym: AB-001
MF: C₂₄H₃₁NO
FW: 349.5
Purity: ≥97%
UV/Vis.: λ_{max}: 214, 247, 305 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 adamantyl analog is a mildly selective agonist of the peripheral cannabinoid receptor, where the naphthalene ring is substituted with an adamantyl group.¹ Adamantyl substitutions provide structural rigidity and bulk. Though no biological activity has been reported for JWH 018 adamantyl analog, Δ⁸-THC analogs with adamantyl substituted for the carbon side chain demonstrate improved affinity and selectivity of central cannabinoid and peripheral cannabinoid receptor binding.² This synthetic cannabinoid has been identified in a herbal product.³ This product is intended for research and forensic applications.

References

1. Sobolevsky, T., Prasolov, I., and Rodchenkov, G. Detection of JWH-018 metabolites in smoking mixture post-administration urine. *Forensic Sci. Int.* **200(1-3)**, 141-147 (2010).
2. Lu, D., Meng, Z., Thakur, G.A., et al. Adamantyl cannabinoids: A novel class of cannabinergic ligands. *J. Med. Chem.* **48(14)**, 4576-4585 (2005).
3. Kneisel, S., Westphal, F., Bisel, P., et al. Identification and structural characterization of the synthetic cannabinoid 3-(1-adamantoyl)-1-pentylindole as an additive in 'herbal incense'. *J. Mass. Spectrom.* **47(2)**, 195-200 (2012).

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, 07/07/2023

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