

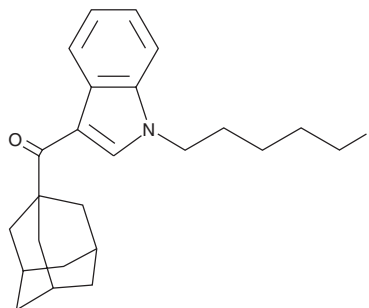
PRODUCT INFORMATION



5-fluoro JWH 018 adamantyl analog

Item No. 14432

CAS Registry No.: 1364933-62-9
Formal Name: [1-(5-fluoropentyl)-1H-indol-3-yl]tricyclo[3.3.1.1^{3,7}]dec-1-yl-methanone
Synonym: 5-fluoro-AB-001, AM2201 adamantyl analog
MF: C₂₄H₃₀FNO
FW: 367.5
Purity: ≥98%
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 is an analog of JWH 018, a mildly selective agonist of the peripheral cannabinoid (CB) 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, Δ⁸-THC analogs with adamantyl substituted for the carbon side chain demonstrate improved affinity and selectivity of central CB₁ and peripheral CB₂ binding.² 5-fluoro JWH 018 adamantyl analog is a derivative of JWH 018 adamantyl (Item No. 9000799), having a fluorine atom at the terminal carbon of the pentyl chain. The physiological and toxicological properties of this compound are not known. This product is intended for forensic and research 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).

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, 06/19/2023

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