

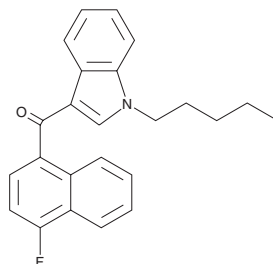
PRODUCT INFORMATION



JWH 412

Item No. 11971

CAS Registry No.: 1364933-59-4
Formal Name: (4-fluoro-1-naphthalenyl)(1-pentyl-1H-indol-3-yl)-methanone
MF: C₂₄H₂₂FNO
FW: 359.4
Purity: ≥98%
UV/Vis.: λ_{max}: 217, 317 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 412 is a positional isomer of AM2201 (Item No. 10707) that was identified in the 'herbal mixture' XoXo.¹ The substitution of hydrogen against fluorine in the 4-position of the naphthyl moiety enhances the binding affinity to the central cannabinoid (CB₁) receptor compared to the unsubstituted JWH 018.¹ JWH 412 demonstrates K_i values of 7.2 and 3.2 nM at CB₁ and peripheral CB₂ receptors, respectively.² This product is intended for research and forensic purposes.

References

1. Moosmann, B., Kneisel, S., Girreser, U., *et al.* Separation and structural characterization of the synthetic cannabinoids JWH-412 and 1-[(5-fluoropentyl)-1H-indol-3yl]-(4-methylnaphthalen-1-yl)methanone using GC-MS, NMR analysis and a flash chromatography system. *Forensic Sci. Int.* **220(1-3)**, e17-e22 (2012).
2. Smith, V.J. Synthesis and pharmacology of N-alkyl-3-(halonaphthoyl)indoles, *Ph.D. diss., Clemson University*, (2008).

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, 07/24/2023

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