

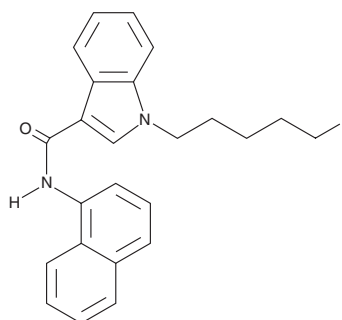
PRODUCT INFORMATION



5-fluoro NNEI

Item No. 14147

CAS Registry No.: 1445580-60-8
Formal Name: 1-(5-fluoropentyl)-N-1-naphthalenyl-1H-indole-3-carboxamide
Synonym: 5-fluoro MN-24
MF: C₂₄H₂₃FN₂O
FW: 374.5
Purity: ≥98%
UV/Vis.: λ_{max}: 221, 299 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 (Item No. 10900) is a potent synthetic cannabinoid (CB) which has been identified in smoking mixtures.^{1,2} NNEI (Item No. 15001) is an analog of JWH 018 which differs by having an amide linker inserted between the naphthalene and ketone groups. 5-fluoro NNEI is a derivative of NNEI in which a fluorine atom has been added to the terminal carbon of the pentyl chain. This modification of other synthetic CBs typically increases the compound's affinity for both CB receptors significantly.³ 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. Moran, C.L., Le, V.H., Chimalakonda, K.C., *et al.* Quantitative measurement of JWH-018 and JWH-073 metabolites excreted in human urine. *Anal. Chem.* **83(11)**, 4228-4236 (2011).
3. Makriyannis, A. and Deng, H. Cannabimimetic indole derivatives. *University of Connecticut WO01/28557A1* (2001).

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, 07/25/2023

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