

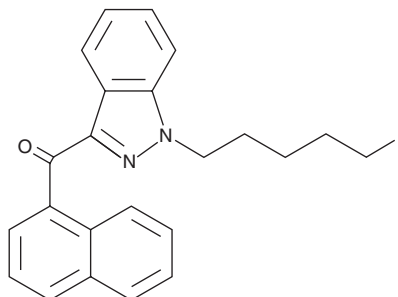
PRODUCT INFORMATION



THJ2201 (exempt preparation)

Item No. 22860

CAS Registry No.: 1801552-01-1
Formal Name: [1-(5-fluoropentyl)-1H-indazol-3-yl]-1-naphthalenyl-methanone
Synonyms: AM2201 indazole analog,
5-fluoropentyl JWH 018 indazole analog,
5-fluoro THJ 018
MF: C₂₃H₂₁FN₂O
FW: 360.4
Purity: ≥98%
UV/Vis.: λ_{max}: 219, 320 nm
Supplied as: A 1 mg/ml solution in methanol
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

THJ2201 (exempt preparation) (Item No. 22860) is an analytical reference standard categorized as a synthetic cannabinoid.¹ THJ2201 is regulated as a Schedule 1 compound in the United States. THJ2201 (exempt preparation) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

Reference

1. Uchiyama, N., Shimokawa, Y., Kawamura, M., *et al.* Chemical analysis of a benzofuran derivative, 2-(2-ethylaminopropyl)benzofuran (2-EAPB), eight synthetic cannabinoids, five cathinone derivatives, and five other designer drugs newly detected in illegal products. *Forensic Toxicol.* **32**(2), 266-281 (2014).

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, 09/13/2017

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