

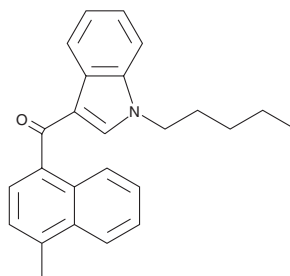
PRODUCT INFORMATION



JWH 122 (exempt preparation)

Item No. 15667

CAS Registry No.: 619294-47-2
Formal Name: (4-methyl-1-naphthalenyl)(1-pentyl-1H-indol-3-yl)-methanone
MF: C₂₅H₂₅NO
FW: 355.5
Purity: ≥98%
UV/Vis.: λ_{max}: 221, 247, 316 nm
Supplied as: A 1 mg/ml solution in methanol
Storage: -20°C
Stability: ≥5 years



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

Description

JWH 122 (exempt preparation) (Item No. 15667) is an analytical reference standard categorized as a synthetic cannabinoid.¹ JWH 122 has been found in Spice-like herbal blends.² JWH 122 is regulated as a Schedule I compound in the United States. JWH 122 (exempt preparation) (Item No. 15667) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

References

1. Cannaert, A., Storme, J., Franz, F., *et al.* Detection and activity profiling of synthetic cannabinoids and their metabolites with a newly developed bioassay. *Anal. Chem.* **88**(23), 11476-11485 (2016).
2. Ernst, L., Schiebel, H.M., Theuring, C., *et al.* Identification and characterization of JWH-122 used as new ingredient in "spice-like" herbal incenses. *Forensic Sci. Int.* **208**(1-3), e31-e35 (2011).

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, 06/29/2026

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