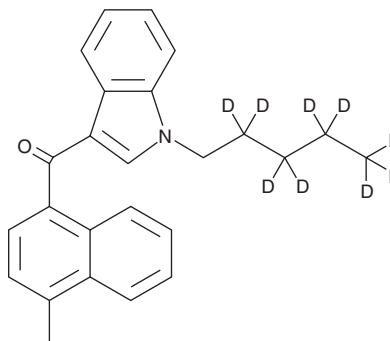


PRODUCT INFORMATION



JWH 122-d₉ (exempt preparation) Item No. 22868

CAS Registry No.: 1651833-50-9
Formal Name: (4-methyl-1-naphthalenyl)(1-pentyl-1H-indol-3-yl)-methanone-2,2',3,3',4,4',5,5,5-d₉
MF: C₂₅H₁₆D₉NO
FW: 364.5
Chemical Purity: ≥98% (JWH 122)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₉); ≤1% d₀
UV/Vis.: λ_{max}: 222, 247, 316 nm
Supplied as: A solution in methanol
Storage: -20°C
Stability: ≥1 year



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

Description

JWH 122-d₉ (exempt preparation) (Item No. 22868) is intended for use as an internal standard for the quantification of JWH 122 (Item No. 10591) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

JWH 122 is a synthetic cannabinoid (CB) that displays high-affinities for both CB₁ (K_i = 0.69 nM) and CB₂ (K_i = 1.2 nM) receptors.^{1,2} JWH 122-d₉ is regulated as a Schedule I compound in the United States. JWH 122-d₉ (exempt preparation) (Item No. 22868) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

References

- Huffman, J.W., Zengin, G., Wu, M.-J., *et al.* Structure-activity relationships for 1-alkyl-3-(1-naphthoyl) indoles at the cannabinoid CB₁ and CB₂ receptors: steric and electronic effects of naphthoyl substituents. New highly selective CB₂ receptor agonists. *Bioorg. Med. Chem.* **13**, 89-112 (2005).
- Huffman, J.W., Mabon, R., Wu, M.-J., *et al.* 3-indolyl-1-naphthylmethanes: New cannabimimetic indoles provide evidence for aromatic stacking interactions with the CB₁ cannabinoid receptor. *Bioorg. Med. Chem.* **11**, 539-549 (2003).

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, 08/24/2017

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