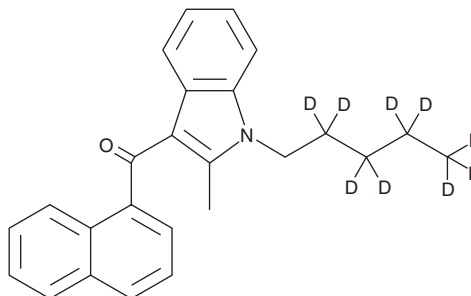


PRODUCT INFORMATION



JWH 007-d₉
Item No. 10486

CAS Registry No.: 1651833-48-5
Formal Name: (2-methyl-1-pentyl-1H-indol-3-yl)-1-naphthalenyl-2,2,3,3,4,4,5,5,5-d₉-methanone
MF: C₂₅H₁₆D₉NO
FW: 364.5
Chemical Purity: ≥98% (JWH 007)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₉); ≤1% d₀
UV/Vis.: λ_{max}: 218, 319 nm
Supplied as: A 1 mg/ml solution in methyl acetate
Storage: -20°C
Stability: ≥4 years



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

Description

JWH 007-d₉ (Item No. 10486) is intended for use as an internal standard for the quantification of JWH 007 (Item No. 10266) 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 007 is a potent cannabinoid (CB) receptor agonist that avidly binds to both CB₁ and CB₂ (K_i = 9.5 and 2.9 nM, respectively).^{1,2} This compares favorably with the binding of Δ⁹-THC, which binds CB₁ and CB₂ with K_i values of 41 and 36 nM, respectively.² Similarly, JWH 007 performs comparably to Δ⁹-THC in murine studies on spontaneous activity, antinociception, hypothermia, and catalepsy.¹ This product is intended for forensic and research purposes.

References

1. Huffman, J.W., Dai, D., Martin, B.R., *et al.* Design, synthesis and pharmacology of cannabimimetic indoles. *Bioorg. Med. Chem. Lett.* **4**(4), 563-566 (1994).
2. 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**(1), 89-112 (2005).

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, 11/19/2025

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