

PRODUCT INFORMATION



JWH 210-d₉
Item No. 10510

CAS Registry No.: 1651833-52-1
Formal Name: (4-ethyl-1-naphthalenyl)(1-pentyl-1H-indol-3-yl)-methanone-d₉

MF: C₂₆H₁₈D₉NO
FW: 378.6

Chemical Purity: ≥98% JWH 210

Deuterium

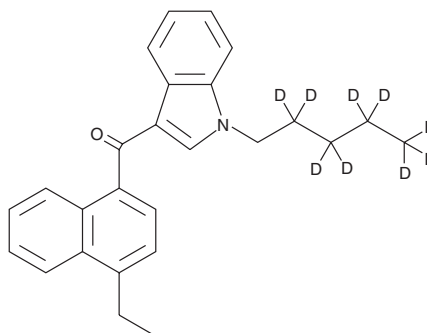
Incorporation: ≥99% deuterated forms (d₁-d₉);
≤1% d₀

UV/Vis.: λ_{max}: 222, 247, 316 nm

Supplied as: A solution in methanol

Storage: -20°C

Stability: As supplied, 1 year from the QC date provided on the Certificate of Analysis, when stored properly



Description

JWH 210-d₉ is intended for use as an internal standard for the quantification of JWH 210 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 210 is a potent cannabimimetic alkylindole, binding the central cannabinoid and peripheral cannabinoid receptors with K_i values of 0.46 and 0.69 nM, respectively.¹ The effects of JWH 210 in whole cells or organisms have not been evaluated.

References

1. 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).

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, 07/20/2016

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