

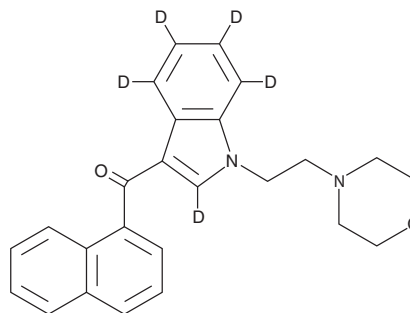
PRODUCT INFORMATION



JWH 200-d₅ (exempt preparation)

Item No. 10682

CAS Registry No.: 1651833-51-0
Formal Name: [1-[2-(4-morpholinyl)ethyl]-1H-indol-3-yl-2',4',5',6',7'-d₅]-1-naphthalenyl-methanone
MF: C₂₅H₁₉D₅N₂O₂
FW: 389.5
Chemical Purity: ≥98% JWH 200-d₅ (exempt preparation)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀
UV/Vis.: λ_{max}: 219, 247, 314 nm
Supplied as: A solution in acetonitrile
Storage: -20°C
Stability: ≥3 years



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

Laboratory Procedures

JWH 200-d₅ (exempt preparation) is intended for use as an internal standard for the quantification of JWH 200 (exempt preparation) (Item No. 13171) 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 200-d₅ (exempt preparation) is supplied as a solution in acetonitrile. To change the solvent, simply evaporate the acetonitrile under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as methyl acetate, methanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of JWH 200-d₅ (exempt preparation) in these solvents is approximately 5, 2, 20, and 25 mg/ml, respectively.

Description

JWH 200 is an aminoalkylindole that acts as a cannabinoid (CB) receptor ligand. It binds to the CB₁ receptor with high-affinity (IC₅₀ = 7.8-42 nM).^{1,2} The effects of JWH 200 in locomotor activity, tail-flick latency, hypothermia, and ring-immobility tests are comparable or superior to Δ⁹-THC or WIN 55,212.³ It potently inhibits the contraction of electrically-stimulated murine vas deferens (IC₅₀ = 3.7-6.0 nM).^{4,5}

References

1. Eissenstat, M.A., Bell, M.R., D'Ambra, T.E., *et al.* Aminoalkylindoles: Structure-activity relationships of novel cannabinoid mimetics. *J. Med. Chem.* **38**(16), 3094-3105 (1995).
2. 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. *Bioorgan. Med. Chem.* **11**(4), 539-549 (2003).
3. Compton, D.R., Gold, L.H., Ward, S.J., *et al.* Aminoalkylindole analogs: Cannabimimetic activity of a class of compounds structurally distinct from Δ⁹-tetrahydrocannabinol. *J. Pharmacol. Exp. Ther.* **263**(3), 1118-1126 (1992).
4. Pacheco, M., Childers, S.R., Arnold, R., *et al.* Aminoalkylindoles: Actions on specific G-protein-linked receptors. *J. Pharmacol. Exp. Ther.* **257**(1), 170-183 (1991).
5. Bell, M.R., D'Ambra, T.E., Kumar, V., *et al.* Antinociceptive (aminoalkyl) indoles. *J. Med. Chem.* **34**(3), 1099-1110 (1991).

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, 01/17/2025

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