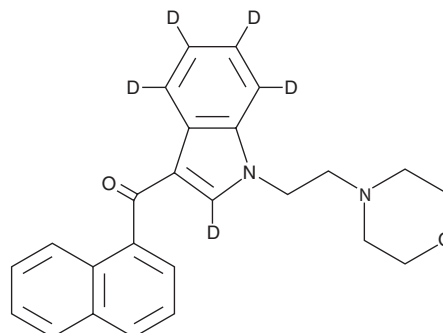


# PRODUCT INFORMATION



**JWH 200-d<sub>5</sub>**  
Item No. 10903

**CAS Registry No.:** 1651833-51-0  
**Formal Name:** [1-[2-(4-morpholinyl)ethyl]-1H-indol-3-yl-2',4',5',6',7'-d<sub>5</sub>]-1-naphthalenyl-methanone  
**MF:** C<sub>25</sub>H<sub>19</sub>D<sub>5</sub>N<sub>2</sub>O<sub>2</sub>  
**FW:** 389.5  
**Chemical Purity:** ≥98% (JWH 200)  
**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>5</sub>); ≤1% d<sub>0</sub>  
**UV/Vis.:** λ<sub>max</sub>: 219, 247, 314 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥5 years



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

## Description

JWH 200-d<sub>5</sub> (Item No. 10903) is intended for use as an internal standard for the quantification of JWH 200 (Item No. 10902) 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 is an aminoalkylindole that acts as a cannabinoid (CB) receptor ligand. It binds to the CB<sub>1</sub> receptor with high-affinity (IC<sub>50</sub> = 7.8-42 nM).<sup>1,2</sup> The effects of JWH 200 in locomotor activity, tail-flick latency, hypothermia, and ring-immobility tests are comparable or superior to Δ<sup>9</sup>-THC or WIN 55,212.<sup>3</sup> It potently inhibits the contraction of electrically-stimulated mouse vas deferens (IC<sub>50</sub> = 3.7-6.0 nM).<sup>4,5</sup>

## 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 CB<sub>1</sub> 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 Δ<sup>9</sup>-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, 06/16/2023

## 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