

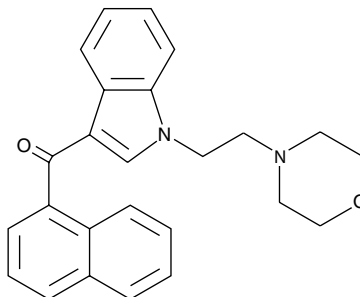
# Product Information



## JWH 200 (exempt preparation)

Item No. 13171

**CAS Registry No.:** 103610-04-4  
**Formal Name:** [1-[2-(4-morpholinyl)ethyl]-1H-indol-3-yl]-1-naphthalenyl-methanone  
**MF:** C<sub>25</sub>H<sub>24</sub>N<sub>2</sub>O<sub>2</sub>  
**FW:** 384.5  
**Purity:** ≥98%  
**Stability:** ≥2 years at -20°C  
**Supplied as:** A solution in acetonitrile



### Laboratory Procedures

For long term storage, we suggest that JWH 200 (exempt preparation) be stored as supplied at -20°C. It should be stable for at least two years.

JWH 200 (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 DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of JWH 200 (exempt preparation) in these solvents is approximately 20 mg/ml.

JWH 200 (exempt preparation) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, JWH 200 (exempt preparation) should first be dissolved in DMF and then diluted with the aqueous buffer of choice. JWH 200 (exempt preparation) has a solubility of approximately 0.1 mg/ml in a 1:6 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

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 murine 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**, 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. *Bioorg. Med. Chem.* **11**, 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**, 1099-1110 (1991).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/13171](http://www.caymanchem.com/catalog/13171)

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog**.

Copyright Cayman Chemical Company, 01/15/2014

### Cayman Chemical

#### Mailing address

1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

#### Phone

(800) 364-9897  
(734) 971-3335

#### Fax

(734) 971-3640

#### E-Mail

[custserv@caymanchem.com](mailto:custserv@caymanchem.com)

#### Web

[www.caymanchem.com](http://www.caymanchem.com)