

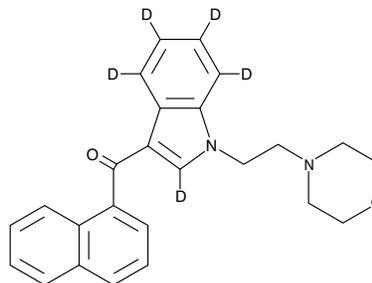
Product Information



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
Stability: ≥2 years at -20°C
Supplied as: A solution in acetonitrile
UV/Vis.: λ_{max}: 219, 247, 314 nm



Laboratory Procedures

JWH 200-d₅ (exempt preparation) contains five deuterium atoms at the 2', 4', 5', 6', and 7' positions. It is intended for use as an internal standard for the quantification of JWH 200 (exempt preparation) by GC- or LC-mass spectrometry (MS). For long term storage, we suggest that JWH 200-d₅ (exempt preparation) be stored as supplied at -20°C. It should be stable for at least two years.

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 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 20 mg/ml.

JWH 200-d₅ (exempt preparation) is used as an internal standard for the quantification of JWH 200 (exempt preparation) by stable isotope dilution 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 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. J. Med. Chem.* **38**, 3094-3105 (1995).
2. Huffman, J.W., Mabon, R., Wu, M.-J., *et al. Bioorg. Med. Chem.* **11**, 539-549 (2003).
3. Compton, D.R., Gold, L.H., Ward, S.J., *et al. J. Pharmacol. Exp. Ther.* **263**(3), 1118-1126 (1992).
4. Pacheco, M., Childers, S.R., Arnold, R., *et al. J. Pharmacol. Exp. Ther.* **257**(1), 170-183 (1991).
5. Bell, M.R., D'Ambra, T.E., Kumar, V., *et al. J. Med. Chem.* **34**, 1099-1110 (1991).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10682

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, 04/14/2015

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

Web

www.caymanchem.com