

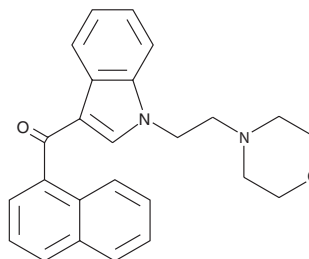
PRODUCT INFORMATION



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: C₂₅H₂₄N₂O₂
FW: 384.5
Purity: ≥98%
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.

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 potentially 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, 01/16/2025

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