

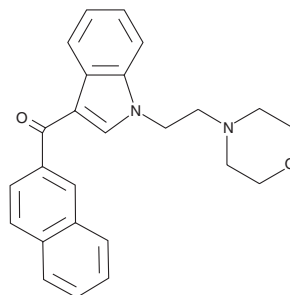
# PRODUCT INFORMATION



## JWH 200 2'-naphthyl isomer

Item No. 9000897

**CAS Registry No.:** 133438-66-1  
**Formal Name:** [1-[2-(4-morpholinyl)ethyl]-1H-indol-3-yl]-2-naphthalenyl-methanone  
**MF:** C<sub>25</sub>H<sub>24</sub>N<sub>2</sub>O<sub>2</sub>  
**FW:** 384.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 215, 253, 326 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 is a 3-indolyl-1-naphthylmethane that acts as a cannabinoid (CB) receptor agonist, binding to the CB<sub>1</sub> receptor with high affinity (IC<sub>50</sub> = 42 nM).<sup>1</sup> This synthetic cannabinoid has been identified as an adulterant of herbal blends.<sup>2</sup> JWH 200 2'-naphthyl isomer differs structurally from JWH 200 by having the naphthyl group attached at the 2' position. The biological and toxicological properties of this compound have not been characterized. This product is intended for forensic purposes.

### References

- 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).
- Kikura-Hanajiri, R., Uchiyama, N., and Goda, Y. Survey of current trends in the abuse of psychotropic substances and plants in Japan. *Leg. Med. (Tokyo)* **13(3)**, 109-115 (2011).

#### 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, 07/10/2023

#### 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