

# PRODUCT INFORMATION



## JWH 200 analog 1

Item No. 9001951

**Formal Name:** (E)-3,4,4-trimethyl-1-(1-(2-morpholinoethyl)-1H-indol-3-yl)pent-2-en-1-one

**MF:** C<sub>22</sub>H<sub>30</sub>N<sub>2</sub>O<sub>2</sub>

**FW:** 354.5

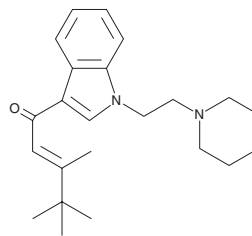
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 213, 254, 321 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 (Item No. 10902) is a synthetic cannabinoid (CB) receptor agonist that binds to the CB<sub>1</sub> receptor with high-affinity (IC<sub>50</sub> = 7.8-42 nM).<sup>1-2</sup> Structurally, JWH 200 features a (4-morpholinyl)ethyl group attached to the indole amide of a 1H-indol-3-yl-(1-naphthyl)-methanone scaffold. JWH 200 analog 1 differs from JWH 200 by having 2,3,3-trimethyl-1-butene in place of the naphthyl group. The physiological and toxicological properties of this compound are not known. This product is intended for forensic and research applications.

### 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. *Bioorg. Med. Chem.* **11(4)**, 539-549 (2003).

#### 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, 07/14/2023

#### 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](http://WWW.CAYMANCHEM.COM)