

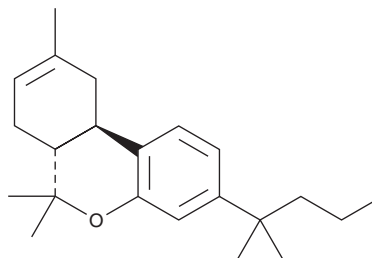
# PRODUCT INFORMATION



## JWH 133

Item No. 10005428

**CAS Registry No.:** 259869-55-1  
**Formal Name:** 3-(1,1-dimethylbutyl)-6aR,7,10,10aR-tetrahydro-6,6,9-trimethyl-6H-dibenzo[b,d]pyran  
**MF:** C<sub>22</sub>H<sub>32</sub>O  
**FW:** 312.5  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 278, 285 nm  
**Supplied as:** A solution in methyl acetate  
**Storage:** -20°C  
**Stability:** ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Description

JWH 133 is a synthetic cannabinoid (CB) that is over 200-fold selective for the peripheral CB<sub>2</sub> receptor (K<sub>i</sub> = 3.4 nM) over the central CB<sub>1</sub> receptor (K<sub>i</sub> = 677 nM).<sup>1</sup> It inhibits the growth of tumors in mice, increasing apoptosis and impairing angiogenesis.<sup>2,3</sup> JWH 133 also suppresses inflammation in experimental colitis and pancreatitis in mice as well as *in vitro* models.<sup>4-6</sup>

### References

1. Huffman, J.W., Little, J., Yu, S., et al. *Bioorg. Med. Chem.* **7**, 2905-14 (1999).
2. Casanova, M.L., Blázquez, C., Martínez-Palacio, J., et al. *J. Clin. Invest.* **111**(1), 43-50 (2003).
3. Blázquez, C., Casanova, M.L., Planas, A., et al. *FASEB J.* **17**, 529-531 (2003).
4. Kimball, E.S., Schneider, C.R., Wallace, N.H., et al. *Am. J. Physiol. Gastrointest. Liver Physiol.* **291**(2), G364-G371 (2006).
5. Michler, T., Storr, M., Kramer, J., et al. *Am. J. Physiol. Gastrointest. Liver Physiol.* **304**(2), G181-G192 (2013).
6. Ramirez, S.H., Haskó, J., Skuba, A., et al. *J. Neurosci.* **32**(12), 4004-4016 (2012).

#### 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, 07/11/2023

#### 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