

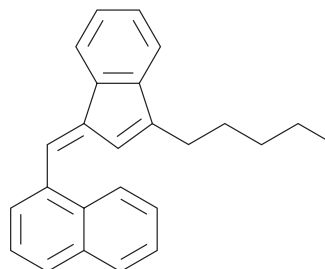
# PRODUCT INFORMATION



**JWH 176**

Item No. 11617

**CAS Registry No.:** 619294-62-1  
**Formal Name:** 1-[(E)-(3-pentyl-1H-inden-1-ylidene)methyl]-naphthalene  
**MF:** C<sub>25</sub>H<sub>24</sub>  
**FW:** 324.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 210, 348 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 176 is a cannabimimetic indene hydrocarbon with high affinity for the central cannabinoid (CB<sub>1</sub>) receptor (K<sub>i</sub> = 26 nM).<sup>1,2</sup> It was originally prepared to examine the roles of hydrogen bonding involving the carbonyl group and aromatic stacking in CB<sub>1</sub> receptor docking studies.<sup>2</sup> This product is intended for forensic and research applications.

## References

1. Huffman, J.W. and Padgett, L.W. Recent developments in the medicinal chemistry of cannabimimetic indoles, pyrroles and indenenes. *Curr. Med. Chem.* **12**, 1395-1411 (2005).
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**, 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, 07/18/2023

## 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