

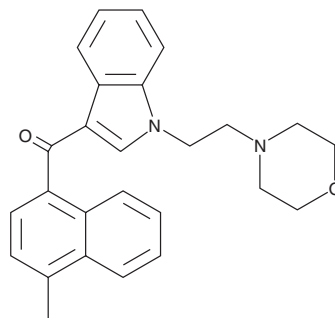
# PRODUCT INFORMATION



## JWH 193

Item No. 13980

**CAS Registry No.:** 133438-58-1  
**Formal Name:** (4-methyl-1-naphthalenyl)  
[1-[2-(4-morpholinyl)ethyl]-1H-  
indol-3-yl]-methanone  
**MF:** C<sub>26</sub>H<sub>26</sub>N<sub>2</sub>O<sub>2</sub>  
**FW:** 398.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 222, 313 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 193 (Item No. 13980) is a synthetic cannabinoid (CB) which binds the central CB<sub>1</sub> receptor with high affinity (K<sub>i</sub> = 6 nM).<sup>1</sup> This aminoalkylindole shares structural features with the antinociceptive CB<sub>1</sub> agonists pravadoline (Item No. 10006973) and WIN 55,212-2.<sup>1</sup> The physiological and toxicological properties of this compound have not been investigated. This product is intended for research and forensic applications.

### Reference

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

#### 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