

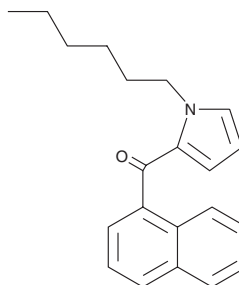
# PRODUCT INFORMATION



## JWH 031 2'-isomer

Item No. 11631

CAS Registry No.: 1797106-18-3  
Formal Name: (1-hexyl-1H-pyrrol-2-yl)-1-naphthalenyl-methanone  
MF: C<sub>21</sub>H<sub>23</sub>NO  
FW: 305.4  
Purity: ≥98%  
Stability: ≥1 year at -20°C  
Supplied as: A solution in methanol  
UV/Vis.: λ<sub>max</sub>: 218, 302 nm



### Laboratory Procedures

For long term storage, we suggest that JWH 031 2'-isomer be stored as supplied at -20°C. It should be stable for at least one year.

JWH 031 2'-isomer is supplied as a solution in methanol. To change the solvent, simply evaporate the methanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of JWH 031 2'-isomer in ethanol is approximately 20 mg/ml and approximately 10 mg/ml in DMSO and DMF.

JWH 031 2'-isomer is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

### Description

JWH 031 (Item No. 10824) is a synthetic cannabinoid (CB) with a relatively low binding affinity for the central CB<sub>1</sub> receptor (K<sub>i</sub> = 399 nM) that is efficacious in reducing spontaneous activity and increasing antinociception in mice.<sup>1</sup> JWH 031 2'-isomer differs from JWH 031 by having the naphthyl group at the 2' position, rather than the 5' position, of the pyrrole group. The physiological properties of this compound are unknown. This product is intended for forensic and research purposes.

### Reference

1. Wiley, J.L., Compton, D.R., Dai, D., *et al.* Structure-activity relationships of indole- and pyrrole-derived cannabinoids. *J. Pharmacol. Exp. Ther.* **285**(3), 995-1004 (1998).

#### 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, 09/11/2015

#### 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