

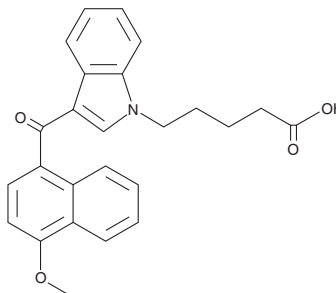
# PRODUCT INFORMATION



## JWH 081 N-pentanoic acid metabolite

Item No. 14353

**CAS Registry No.:** 1537889-08-9  
**Formal Name:** 3-[[4-methoxy-1-naphthalenyl]carbonyl]-1H-indole-1-pentanoic acid  
**MF:** C<sub>25</sub>H<sub>23</sub>NO<sub>4</sub>  
**FW:** 401.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 212, 315 nm  
**Supplied as:** A solution in acetonitrile  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Description

JWH 081 (Item No. 10579) is a cannabimimetic indole related to JWH 015 (Item No. 10009018) and JWH 018 (Item No. 10900) that shows preference for the central cannabinoid (CB<sub>1</sub>) receptor (K<sub>i</sub> = 1.2 nM) over the peripheral cannabinoid (CB<sub>2</sub>) receptor (K<sub>i</sub> = 12.4 nM).<sup>1,2</sup> JWH 081 N-pentanoic acid metabolite is an expected metabolite of JWH 081 that would be detectable both in serum and in urine. The physiological and toxicological properties of this compound are not known. This product is intended for research and forensic purposes.

### References

1. Aung, M.M., Griffin, G., Huffman, J.W., *et al.* Influence of the N-1 alkyl chain length of cannabimimetic indoles upon CB<sub>1</sub> and CB<sub>2</sub> receptor binding. *Drug Alcohol Depend.* **60(2)**, 133-140 (2000).
2. Huffman, J.W., Zengin, G., Wu, M.J., *et al.* Structure-activity relationships for 1-alkyl-3-(1-naphthoyl) indoles at the cannabinoid CB<sub>1</sub> and CB<sub>2</sub> receptors: Steric and electronic effects of naphthoyl substituents. New highly selective CB<sub>2</sub> receptor agonists. *Bioorg. Med. Chem.* **13(1)**, 89-112 (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, 06/13/2023

#### 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