

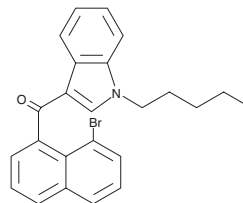
# PRODUCT INFORMATION



## JWH 424

Item No. 9001203

**CAS Registry No.:** 1366068-04-3  
**Formal Name:** (8-bromonaphthalen-1-yl)(1-pentyl-1H-indol-3-yl)methanone  
**MF:** C<sub>24</sub>H<sub>22</sub>BrNO  
**FW:** 420.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 224, 315 nm  
**Supplied as:** A solution in methanol  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Description

JWH 018 is a synthetic cannabinoid (CB) that potently activates the central CB<sub>1</sub> and peripheral CB<sub>2</sub> receptors (K<sub>i</sub> = 9.0 and 2.94 nM, respectively).<sup>1</sup> This cannabimimetic compound has frequently been found in herbal blends. <sup>2-4</sup> JWH 424 is an 8-bromonaphthyl derivative of JWH 018 which shows a reduced selectivity for CB<sub>1</sub> over CB<sub>2</sub> (K<sub>i</sub> = 20.9 and 5.4 nM, respectively). The properties of this compound *in vivo* are not known. This product is intended for forensic and research applications.

### 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. Auwärter, V., Dresen, S., Weinmann, W., *et al.* 'Spice' and other herbal blends: Harmless incense or cannabinoid designer drugs? *J. Mass Spectrom.* **44(5)**, 832-837 (2009).
3. Lindigkeit, R., Boehme, A., Eiserloh, I., *et al.* Spice: A never ending story? *Forensic Sci. Int.* **191(1-3)**, 58-63 (2009).
4. Uchiyama, N., Kikura-Hanajiri, R., Kawahara, N., *et al.* Identification of a cannabimimetic indole as a designer drug in a herbal product. *Forensic Toxicol.* **27(2)**, 61-66 (2009).

#### 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, 07/06/2023

#### 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