

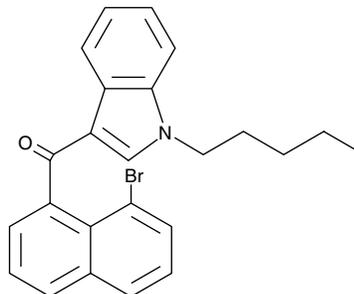
# Product Information



## 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%  
**Stability:** ≥1 year at -20°C  
**Supplied as:** A solution in methanol  
**UV/Vis.:** λ<sub>max</sub>: 224, 315 nm



### Laboratory Procedures

For long term storage, we suggest that JWH 424 be stored as supplied at -20°C. It should be stable for at least one year.

JWH 424 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 424 in ethanol is approximately 30 mg/ml and approximately 5 mg/ml in DMSO and DMF.

JWH 424 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.

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**, 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**, 832-837 (2009).
3. Lindigkeit, R., Boehme, A., Eiserloh, I., *et al.* Spice: A never ending story? *Forensic Sci. Int.* **191**, 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**, 61-66 (2009).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/9001203](http://www.caymanchem.com/catalog/9001203)

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY; NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog.**

Copyright Cayman Chemical Company, 11/09/2012

### Cayman Chemical

#### Mailing address

1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

#### Phone

(800) 364-9897  
(734) 971-3335

#### Fax

(734) 971-3640

#### E-Mail

[custserv@caymanchem.com](mailto:custserv@caymanchem.com)

#### Web

[www.caymanchem.com](http://www.caymanchem.com)