

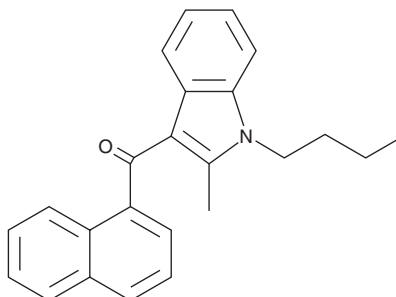
# PRODUCT INFORMATION



**JWH 016**

Item No. 10849

**CAS Registry No.:** 155471-09-3  
**Formal Name:** (1-butyl-2-methyl-1H-indol-3-yl)-1-naphthalenyl-methanone  
**MF:** C<sub>24</sub>H<sub>23</sub>NO  
**FW:** 341.5  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 218, 320 nm  
**Supplied as:** A solution in methanol  
**Storage:** -20°C  
**Stability:** ≥5 years



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

## Description

JWH 016 (Item No. 10849) is an analytical reference standard categorized as a synthetic cannabinoid.<sup>1,2</sup> JWH 016 induces discriminative stimulus effects in rhesus monkeys comparable to those induced by Δ<sup>9</sup>-THC (Item Nos. ISO60157 | 12068).<sup>3</sup> This product is intended for research and forensic 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. 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).
3. Wiley, J.L., Huffman, J.W., Balster, R.L., *et al.* Pharmacological specificity of the discriminative stimulus effects of Δ<sup>9</sup>-tetrahydrocannabinol in rhesus monkeys. *Drug Alcohol Depend.* **40(1)**, 81-86 (1995).

### 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, 07/21/2023

## 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