

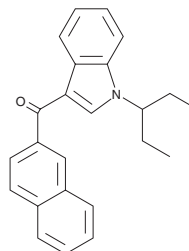
# PRODUCT INFORMATION



## JWH 018 2'-naphthyl-N-(1-ethylpropyl) isomer

Item No. 11586

**CAS Registry No.:** 1869959-37-4  
**Formal Name:** [1-(1-ethylpropyl)-1H-indol-3-yl]-  
2-naphthalenyl-methanone  
**MF:** C<sub>24</sub>H<sub>23</sub>NO  
**FW:** 341.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 216, 253, 325 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥5 years



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

### Description

JWH 018 is a potent synthetic cannabinoid (CB) which has frequently been found in herbal blends.<sup>1-3</sup> JWH 018 2'-naphthyl-N-(1-ethylpropyl) isomer is an isomer of JWH 018, differing by having the naphthyl group attached at the 2, rather than 1, position, as well as an ethylpropyl group in place of the pentyl chain. The biological and toxicological properties of this compound have not been characterized. This product is intended for research and forensic applications.

### References

1. 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).
2. Lindigkeit, R., Boehme, A., Eiserloh, I., *et al.* Spice: A never ending story? *Forensic Sci. Int.* **191(1-3)**, 58-63 (2009).
3. 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).

#### 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, 07/18/2023

#### 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