

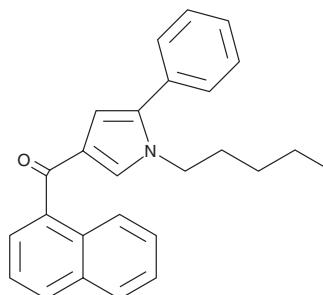
# PRODUCT INFORMATION



## JWH 145

Item No. 10825

**CAS Registry No.:** 914458-19-8  
**Formal Name:** 1-naphthalenyl(1-pentyl-5-phenyl-1H-pyrrol-3-yl)-methanone  
**MF:** C<sub>26</sub>H<sub>25</sub>NO  
**FW:** 367.5  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 217, 254 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

Several aminoalkylindole cannabinoids (CB), including JWH 018 and JWH 073, have been detected in herbal blends.<sup>1,2</sup> JWH 145 (Item No. 10825) is a (1-naphthoyl)pyrrole analog of JWH 018 that potently activates both CB<sub>1</sub> and CB<sub>2</sub> receptors (K<sub>i</sub> values of 14 and 6.4 nM, respectively).<sup>3</sup> Its biological properties have not been evaluated. This compound is intended for forensic and research applications.

### References

1. Kikura-Hanajiri, R., Uchiyama, N., and Goda, Y. Survey of current trends in the abuse of psychotropic substances and plants in Japan. *Leg. Med. (Tokyo)* **13(3)**, 109-115 (2011).
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. Huffman, J.W., Padgett, L.W., Isherwood, M.L., et al. 1-Alkyl-2-aryl-4-(1-naphthoyl)pyrroles: New high affinity ligands for the cannabinoid CB<sub>1</sub> and CB<sub>2</sub> receptors. *Bioorg. Med. Chem. Lett.* **16(20)**, 5432-5435 (2006).

#### 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, 06/15/2023

#### 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