

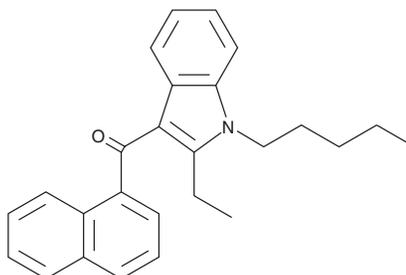
PRODUCT INFORMATION



JWH 116

Item No. 13977

CAS Registry No.: 619294-64-3
Formal Name: (2-ethyl-1-pentyl-1H-indol-3-yl)-
1-naphthalenyl-methanone
MF: C₂₆H₂₇NO
FW: 369.5
Purity: ≥98%
UV/Vis.: λ_{max}: 219, 250, 321 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 116 (Item No. 13977) is a synthetic cannabinoid (CB) receptor ligand from the naphthoylindole family. It is the indole 2-ethyl derivative of the related compound JWH 018 (Item Nos. 10900 and 13169) and has a K_i value of 52 nM for the CB₁ receptor.¹ This product is intended for research and forensic applications.

Reference

1. Huffman, J.W., Mabon, R., Wu, M.J., *et al.* 3-indolyl-1-naphthylmethanes: New cannabimimetic indoles provide evidence for aromatic stacking interactions with the CB₁ cannabinoid receptor. *Bioorgan. Med. Chem.* **11(4)**, 539-549 (2003).

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, 07/14/2023

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