

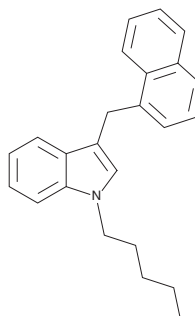
PRODUCT INFORMATION



JWH 175

Item No. 11201

CAS Registry No.: 619294-35-8
Formal Name: 3-(1-naphthalenylmethyl)-1-pentyl-1H-indole
MF: C₂₄H₂₅N
FW: 327.5
Purity: ≥98%
UV/Vis.: λ_{max}: 223, 283 nm
Supplied as: A solution in acetonitrile
Storage: -20°C
Stability: ≥4 years



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

Description

JWH 175 is a synthetic cannabinoid (CB) that potently activates the central CB₁ receptor (K_i = 22 nM).¹ As a naphthylmethylandole, it is structurally related to JWH 018. The properties of this compound at the peripheral CB₂ receptor or *in vivo* have not been characterized. This product is intended for forensic and research purposes.

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, 07/07/2023

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