

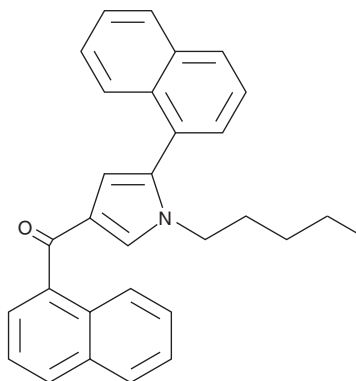
PRODUCT INFORMATION



JWH 309

Item No. 10830

CAS Registry No.: 914458-42-7
Formal Name: 1-naphthalenyl[5-(1-naphthalenyl)-1-pentyl-1H-pyrrol-3-yl]-methanone
MF: C₃₀H₂₇NO
FW: 417.6
Purity: ≥98%
UV/Vis.: λ_{max}: 220, 258 nm
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: ≥4 years



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

Description

JWH 309 is a synthetic cannabinoid (CB) that displays high affinities for both the central CB₁ receptor (K_i = 41 nM) and the peripheral CB₂ receptor (K_i = 49 nM).¹ The physiological and toxicological properties of this compound have not been examined. This product is intended for forensic and research applications.

Reference

1. 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 CB₁ and CB₂ 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