

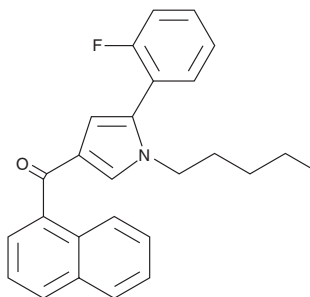
PRODUCT INFORMATION



JWH 307

Item No. 10797

CAS Registry No.: 914458-26-7
Formal Name: [5-(2-fluorophenyl)-1-pentyl-1H-pyrrol-3-yl]-1-naphthalenyl-methanone
MF: C₂₆H₂₄FNO
FW: 385.5
Purity: ≥96%
UV/Vis.: λ_{max}: 217, 250 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 307 is a (1-naphthoyl)pyrrole cannabinimimetic that potently activates both central cannabinoid (CB₁) and peripheral cannabinoid (CB₂) receptors (K_i values of 7.7 and 3.3 nM, respectively).¹ Its physiological, neurological, and toxicological properties have not been evaluated. This compound 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