

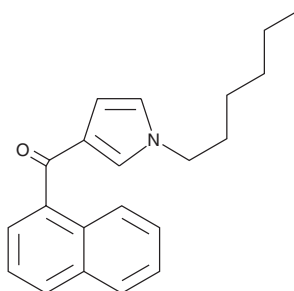
PRODUCT INFORMATION



JWH 031

Item No. 10824

CAS Registry No.: 162934-74-9
Formal Name: (1-hexyl-1H-pyrrol-3-yl)-1-naphthalenyl-methanone
MF: C₂₁H₂₃NO
FW: 305.4
Purity: ≥98%
UV/Vis.: λ_{max}: 217, 281 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 031 is a synthetic cannabinoid (CB) with a relatively low binding affinity for the central CB₁ receptor (K_i = 399 nM).¹ Despite its low affinity for CB₁, this pyrrole-based cannabinoid is efficacious in reducing spontaneous activity and increasing antinociception in mice.¹ This product is intended for forensic and research applications.

Reference

1. Wiley, J.L., Compton, D.R., Dai, D., *et al.* Structure-activity relationships of indole- and pyrrole-derived cannabinoids. *J. Pharmacol. Exp. Ther.* **285(3)**, 995-1004 (1998).

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