

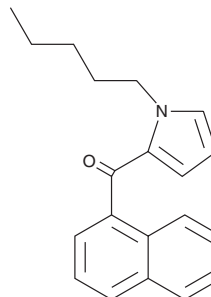
PRODUCT INFORMATION



JWH 030 2-naphthoyl isomer

Item No. 11632

CAS Registry No.: 2365471-38-9
Formal Name: naphthalen-1-yl(1-pentyl-1H-pyrrol-2-yl)methanone
MF: C₂₀H₂₁NO
FW: 291.4
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 303 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 030 is a naphthoyl pyrrole cannabinimetic with specificity for the central cannabinoid (CB₁) receptor (EC₅₀ = 30.5 nM for rat and K_i = 87 nM for mouse) over the peripheral CB₂ receptor (EC₅₀ = 552 nM for human CB₂).^{1,2} JWH 030 2-naphthoyl isomer differs from JWH 030 by having the naphthoyl group attached to the 2 position, rather than the 3 position, of the pyrrole group. The affinity of this isomer for CB receptors and its effects *in vivo* have not been evaluated. This product is intended for forensic and research applications.

References

1. Tarzia, G., Duranti, A., Tontini, A., *et al.* Synthesis and structure-activity relationships of a series of pyrrole cannabinoid receptor agonists. *Bioorg. Med. Chem.* **11(18)**, 3965-3973 (2003).
2. Lainton, J.A.H., Huffman, J.W., Martin, B.R., *et al.* 1-Alkyl-3-(1-naphthoyl)pyrroles: A new class of cannabinoid. *Tetrahedron Lett.* **36(9)**, 1401-1404 (1995).

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, 07/18/2023

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