

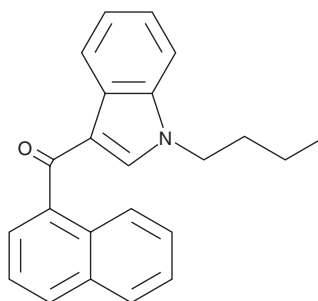
PRODUCT INFORMATION



JWH 073

Item No. 10904

CAS Registry No.: 208987-48-8
Formal Name: (1-butyl-1H-indol-3-yl)-1-naphthalenyl-methanone
MF: C₂₃H₂₁NO
FW: 327.4
Purity: ≥97%
UV/Vis.: λ_{max}: 218, 316 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 073 (Item No. 10904) is a mildly selective agonist of the central cannabinoid (CB₁) receptor derived from the aminoalkylindole WIN 55,212-2. The K_i values for binding CB₁ and the peripheral cannabinoid (CB₂) receptor are 8.9 and 38 nM, respectively for a CB₁:CB₂ ratio of 0.23.¹ Its effects on suppression of spontaneous activity, maximum possible antinociceptive effect in the tail-flick assay, and rectal temperature are comparable to those of WIN 55,212-2 when tested in rats.² This product is intended for forensic and research applications.

References

1. Aung, M.M., Griffin, G., Huffman, J.W., *et al.* Influence of the N-1 alkyl chain length of cannabimimetic indoles upon CB₁ and CB₂ receptor binding. *Drug Alcohol Depend.* **60(2)**, 133-140 (2000).
2. 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, 06/16/2023

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