

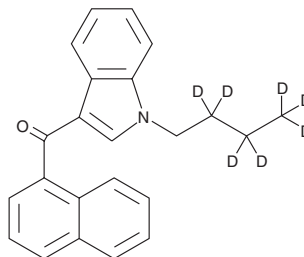
Product Information



JWH 073-d₇ (exempt preparation)

Item No. 9000868

CAS Registry No.: 1415744-43-2
Formal Name: [1-(butyl-2,2,3,3,4,4,4-d₇)-1H-indol-3-yl]-1-naphthalenyl-methanone
MF: C₂₃H₁₄D₇NO
FW: 334.5
Chemical Purity: ≥98% JWH 073-d₇ (exempt preparation)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₇); ≤1% d₀
Stability: ≥2 years at -20°C
Supplied as: A solution in methanol
UV/Vis.: λ_{max}: 218, 246, 315 nm



Laboratory Procedures

JWH 073-d₇ (exempt preparation) contains seven deuterium atoms at the 2, 2', 3, 3', 4, 4, and 4' positions. It is intended for use as an internal standard for the quantification of JWH 073 (exempt preparation) by GC- or LC-mass spectrometry (MS). For long term storage, we suggest that JWH 073-d₇ (exempt preparation) be stored as supplied at -20°C. It should be stable for at least two years.

JWH 073-d₇ (exempt preparation) is supplied as a solution in methanol. To change the solvent, simply evaporate the methanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of JWH 073-d₇ (exempt preparation) in these solvents is approximately 10 mg/ml.

JWH 073-d₇ (exempt preparation) is used as an internal standard for the quantification of JWH 073 (exempt preparation) by stable isotope dilution MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated *versus* unlabeled).

Description

JWH 073 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

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 and Alcohol Dependence* **60**, 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 LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 06/30/2015

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com