

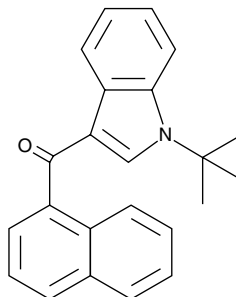
Product Information



JWH 073 N-(1,1-dimethylethyl) isomer

Item No. 9001013

Formal Name:	(1-(tert-butyl)-1H-indol-3-yl) (naphthalen-1-yl)methanone
MF:	C ₂₃ H ₂₁ NO
FW:	327.4
Purity:	≥98%
Stability:	≥2 years at -20°C
Supplied as:	A crystalline solid
UV/Vis.:	λ _{max} : 219, 314 nm



Laboratory Procedures

For long term storage, we suggest that JWH 073 N-(1,1-dimethylethyl) isomer be stored as supplied at -20°C. It should be stable for at least two years.

JWH 073 N-(1,1-dimethylethyl) isomer is supplied as a crystalline solid. A stock solution may be made by dissolving the JWH 073 N-(1,1-dimethylethyl) isomer in the solvent of choice. JWH 073 N-(1,1-dimethylethyl) isomer is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of JWH 073 N-(1,1-dimethylethyl) isomer in ethanol and DMF is approximately 20 mg/ml, and in DMSO it is approximately 10 mg/ml.

JWH 073 N-(1,1-dimethylethyl) isomer is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

JWH 073 is a synthetic cannabinoid (CB) that activates both the central cannabinoid (CB₁) and peripheral cannabinoid (CB₂) receptors (K_i = 8.9 and 38 nM, respectively).¹ This cannabimimetic compound has been found in herbal blends.²⁻⁴ JWH 073 N-(1,1-dimethylethyl) isomer differs structurally from JWH 073 by having a 1,1-dimethylethyl group, rather than a butyl chain, extending from the indole group. The biological activities of this compound have not been characterized. 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**, 133-140 (2000).
2. Kikura-Hanajiri, R., Uchiyama, N., and Goda, Y. Survey of current trends in the abuse of psychotropic substances and plants in Japan. *Leg. Med. (Tokyo)* **13(3)**, 109-15 (2011).
3. Lindigkeit, R., Boehme, A., Eiserloh, I., *et al.* Spice: A never ending story? *Forensic Sci. Int.* **191**, 58-63 (2009).
4. Uchiyama, N., Kikura-Hanajiri, R., Ogata, J., *et al.* Chemical analysis of synthetic cannabinoids as designer drugs in herbal products. *Forensic Sci. Int.* **198**, 31-38 (2010).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/9001013

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL 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 Material 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, 08/02/2012

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