

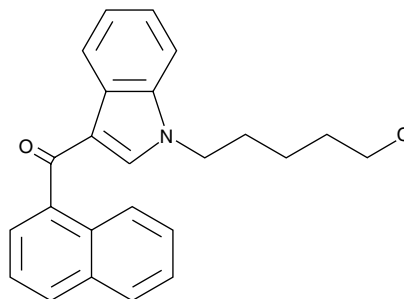
Product Information



JWH 018 N-(5-chloropentyl) analog

Item No. 10521

CAS Registry No.: 1445578-56-2
Formal Name: (1-(5-chloropentyl)-1H-indol-3-yl)
(naphthalen-1-yl)methanone
MF: C₂₄H₂₂ClNO
FW: 375.9
Purity: ≥95%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 218, 246, 314 nm



Laboratory Procedures

For long term storage, we suggest that JWH 018 N-(5-chloropentyl) analog be stored as supplied at -20°C. It should be stable for at least two years.

JWH 018 N-(5-chloropentyl) analog is supplied as a crystalline solid. A stock solution may be made by dissolving the JWH 018 N-(5-chloropentyl) analog in the solvent of choice. JWH 018 N-(5-chloropentyl) analog is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of JWH 018 N-(5-chloropentyl) analog in DMSO is approximately 10 mg/ml and approximately 30 mg/ml in DMF.

JWH 018 N-(5-chloropentyl) analog is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, JWH 018 N-(5-chloropentyl) analog should first be dissolved in DMF and then diluted with the aqueous buffer of choice. JWH 018 N-(5-chloropentyl) analog has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

JWH 018 is a synthetic cannabinoid (CB) that potently activates the central cannabinoid (CB₁) and peripheral cannabinoid (CB₂) receptors (K_i = 9.0 and 2.94 nM, respectively).¹ This cannabimimetic compound has frequently been found in herbal blends.²⁻³ JWH 018 N-(5-chloropentyl) analog differs structurally from JWH 018 by having chlorine added to the 5 position of the pentyl chain. 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. Auwärter, V., Dresen, S., Weinmann, W., *et al.* 'Spice' and other herbal blends: Harmless incense or cannabinoid designer drugs? *J. Mass Spectrom.* **44**, 832-837 (2009).
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., Kawahara, N., *et al.* Identification of a cannabimimetic indole as a designer drug in a herbal product. *Forensic Toxicol.* **27**, 61-66 (2009).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10521

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, 08/06/2013

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