

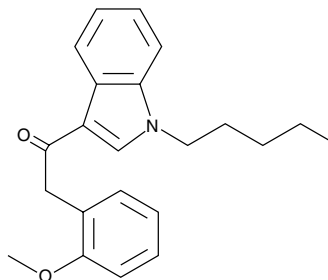
Product Information



JWH 250

Item No. 13634

CAS Registry No.: 864445-43-2
Formal Name: 1-(1-pentyl-1H-indol-3-yl)-2-(2-methoxyphenyl)-ethanone
MF: C₂₂H₂₅NO₂
FW: 335.2
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that JWH 250 be stored as supplied at -20°C. It should be stable for at least two years. JWH 250 is supplied as a crystalline solid. A stock solution may be made by dissolving the JWH 250 in the solvent of choice. JWH 250 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 250 is approximately 2 mg/ml in ethanol and approximately 20 mg/ml in DMSO and DMF.

If aqueous stock solutions are required for biological experiments, they can best be prepared by diluting the organic solvent into aqueous buffers or isotonic saline. Ensure that 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 250 is a cannabimimetic indole that shows a high-affinity for both the central cannabinoid (CB₁) and the peripheral cannabinoid (CB₂) receptors. The affinities of JWH 250 for both CB₁ (K_i = 11 nM) and CB₂ (K_i = 33 nM) receptors are comparable to those of the traditional CB, Δ⁹-tetrahydrocannabinol (Δ⁹-THC; K_i = 40.7 and 36.4 nM, respectively).¹

Reference

1. Huffman, J.W., Szklennik, P.V., Almond, A., *et al.* 1-Pentyl-3-phenylacetylindoles, a new class of cannabimimetic indoles. *Bioorg. Medicinal Chem. Letters* **15**, 4110-4113 (2005).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/13634

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, 12/09/2011

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