

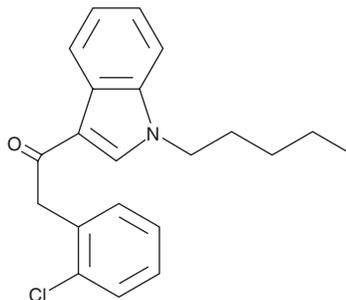
# PRODUCT INFORMATION



**JWH 203**

Item No. 9000736

**CAS Registry No.:** 864445-54-5  
**Formal Name:** 2-(2-chlorophenyl)-1-(1-pentyl-1H-indol-3-yl)-ethanone  
**MF:** C<sub>21</sub>H<sub>22</sub>ClNO  
**FW:** 339.9  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 212, 246, 304 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 203 is an analgesic chemical from the phenylacetylindole family that acts as a cannabinoid (CB) agonist with K<sub>i</sub> values of 8.0 and 7.0 nM at the CB<sub>1</sub> and CB<sub>2</sub> receptors, respectively.<sup>1</sup> Similar to the related 2'-methoxy compound JWH 250, JWH 203 has a phenylacetyl group in place of the naphthoyl ring used in most aminoalkylindole CB compounds. Compared to JWH 250, JWH 203 displays slightly more potent binding affinities for the CB<sub>1</sub> and CB<sub>2</sub> receptors (JWH 250 K<sub>i</sub>s = 11 and 33 nM, respectively).<sup>1</sup> This product is intended for research and forensic applications.

## Reference

1. Huffman, J.W., Szklennik, P.V., Almond, A., *et al.* 1-Pentyl-3-phenylacetylindoles, a new class of cannabimimetic indoles. *Bioorg. Med. Chem. Lett.* **15(18)**, 4110-4113 (2005).

### 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, 07/06/2023

## 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