

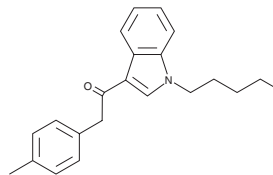
# PRODUCT INFORMATION



## JWH 251 4-methylphenyl isomer

Item No. 9001022

**CAS Registry No.:** 864445-41-0  
**Formal Name:** 2-(4-methylphenyl)-1-(1-pentyl-1H-indol-3-yl)-ethanone  
**MF:** C<sub>22</sub>H<sub>25</sub>NO  
**FW:** 319.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 212, 246, 305 nm  
**Supplied as:** A solution in methanol  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Description

JWH 251 is a cannabimimetic indole that shows a high affinity for the central cannabinoid (CB<sub>1</sub>) receptor with a K<sub>i</sub> value of 29 nM.<sup>1</sup> This compound has been identified as an adulterant in herbal blends.<sup>2</sup> JWH 251 4-methylphenyl isomer differs from JWH 251 by having the methyl group at the 4 position, rather than the 2 position, on the phenyl group. The biological and toxicological properties of this compound have not been characterized. This product is intended for forensic purposes.

### References

- 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).
- 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).

#### 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, 06/20/2023

#### 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