

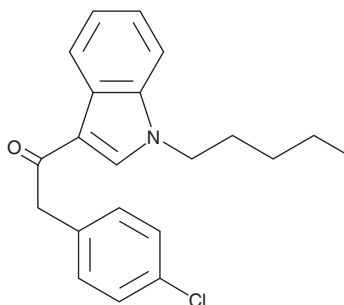
PRODUCT INFORMATION



JWH 203 4-chlorophenyl isomer

Item No. 10867

CAS Registry No.: 864445-58-9
Formal Name: 2-(4-chlorophenyl)-1-(1-pentyl-1H-indol-3-yl)-ethanone
Synonyms: JWH 203 4-chloro isomer, JWH 206
MF: C₂₁H₂₂ClNO
FW: 339.9
Purity: ≥98%
UV/Vis.: λ_{max}: 212, 246, 306 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 4-chlorophenyl isomer (Item No. 10867) is an analytical reference standard categorized as a synthetic cannabinoid.^{1,2} JWH 203 4-chlorophenyl isomer reduces spontaneous locomotor activity, but does not induce antinociception or a decrease body temperature, in mice.² This product is intended for research and forensic applications.

References

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).
2. Wiley, J.L., Marusich, J.A., Martin, B.R., *et al.* 1-Pentyl-3-phenylacetylindoles and JWH 018 share in vivo cannabinoid profiles in mice. *Drug Alcohol Depend.* **123**(1-3), 148-153 (2012).

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, 09/13/2023

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