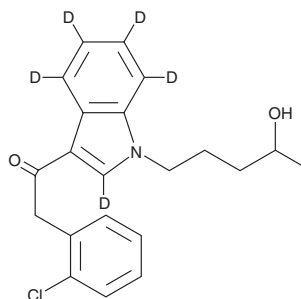


PRODUCT INFORMATION



JWH 203 N-(4-hydroxypentyl) metabolite-d₅ Item No. 14367

CAS Registry No.: 2747915-98-4
Formal Name: 2-(2-chlorophenyl)-1-(1-(4-hydroxypentyl)-1H-indol-3-yl)-2,4,5,6,7-d₅ethan-1-one
MF: C₂₁H₁₇D₅ClNO₂
FW: 360.9
Chemical Purity: ≥98% (JWH 203 N-(4-hydroxypentyl) metabolite)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀
UV/Vis.: λ_{max}: 212, 245, 304 nm
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: ≥4 years



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

Description

JWH 203 N-(4-hydroxypentyl) metabolite-d₅ (Item No. 14367) is intended for use as an internal standard for the quantification of JWH 203 N-(4-hydroxypentyl) metabolite (Item No. 14227) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

JWH 203 (Item No. 9000736) is an analgesic chemical from the phenylacetylindole family that acts as an agonist with K_i values of 8.0 and 7.0 nM at the central cannabinoid (CB₁) and peripheral (CB₂) receptors, respectively.¹ Similar to the related 2'-methoxy compound JWH 250 (Item No. 13634), JWH 203 has a phenylacetyl group in place of the naphthoyl ring used in most aminoalkylindole cannabinoid compounds. Compared to JWH 250, JWH 203 displays slightly more potent binding affinities for the CB₁ and CB₂ receptors (JWH 250 K_s = 11 and 33 nM, respectively).¹ JWH 203 N-(4-hydroxypentyl) metabolite is expected to be a metabolite of JWH 203 that would be detectable both in serum and in urine. While similar hydroxylated phase I metabolites of synthetic CBs retain activity, the physiological properties of this compound have yet to be determined.^{2,3} 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. Brents, L.K., Reichard, E.E., Zimmerman, M., *et al.* Phase I hydroxylated metabolites of the K2 synthetic cannabinoid JWH-018 retain *in vitro* and *in vivo* cannabinoid 1 receptor affinity and activity. *PLoS One* **6**(7), e21917 (2011).
3. Brents, L.K., Gallus-Zawala, A., Radomska-Pandya, A., *et al.* Monohydroxylated metabolites of the K2 synthetic cannabinoid JWH-073 retain intermediate to high cannabinoid 1 receptor (CB1R) affinity and exhibit neutral antagonist to partial agonist activity. *Biochem. Pharmacol.* **83**(7), 952-961 (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, 06/14/2023

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