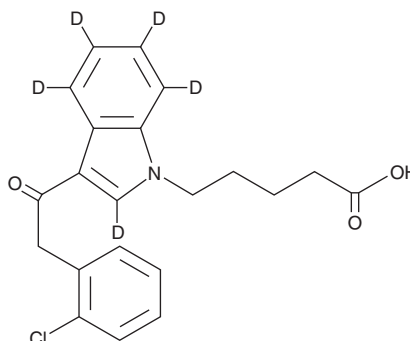


PRODUCT INFORMATION



JWH 203 N-pentanoic acid metabolite-d₅ Item No. 14369

CAS Registry No.: 2749328-26-3
Formal Name: 5-(3-(2-(2-chlorophenyl)acetyl)-1H-indol-1-yl-2,4,5,6,7-d₅)pentanoic acid
MF: C₂₁H₁₅D₅ClNO₃
FW: 374.9
Chemical Purity: ≥98% (JWH 203 N-pentanoic acid metabolite)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d_x); ≤1% d₀
UV/Vis.: λ_{max}: 214, 243, 300 nm
Supplied as: A solution in acetonitrile
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-pentanoic acid metabolite-d₅ (Item No. 14369) is intended for use as an internal standard for the quantification of JWH 203 N-pentanoic acid metabolite (Item No. 14229) 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 is an analgesic chemical from the phenylacetylindole family that acts as a cannabinoid (CB) agonist with K_i values of 8.0 and 7.0 nM at the central (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 CB compounds. Compared to JWH 250, JWH 203 displays slightly more potent binding affinities for the CB₁ and CB₂ CB receptors (JWH 250 K_is = 11 and 33 nM, respectively).¹ JWH 203 N-(5-hydroxypentyl) metabolite is expected to be a metabolite of JWH 203 that would be detectable both in serum and in urine. JWH 203 (Item No. 9000736) is a synthetic cannabinoid (CB) that displays high affinities for both the central CB₁ receptor (K_i = 8.0 nM) and the peripheral CB₂ receptor (K_i = 7.0 nM).¹ JWH 203 N-pentanoic acid metabolite is an expected metabolite of JWH 203, based on the metabolism of similar compounds.² The physiological and toxicological properties of this compound have not been characterized. This product is intended for forensic and research purposes.

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. Moran, C.L., Le, V.H., Chimalakonda, K.C., *et al.* Quantitative measurement of JWH-018 and JWH-073 metabolites excreted in human urine. *Anal. Chem.* **83**(11), 4228-4236 (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, 06/14/2023

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