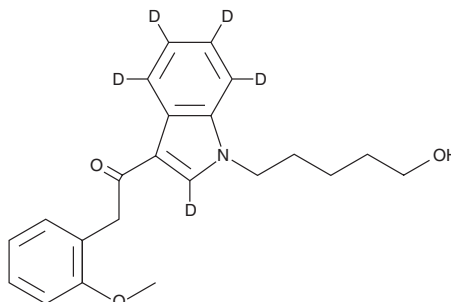


PRODUCT INFORMATION



JWH 250 N-(5-hydroxypentyl) metabolite-d₅ Item No. 11474

CAS Registry No.: 2748212-89-5
Formal Name: 1-[1-(5-hydroxypentyl)-1H-indol-3-yl-2,4,5,6,7-d₅]-2-(2-methoxyphenyl)-ethanone
MF: C₂₂H₂₀D₅NO₃
FW: 356.5
Chemical Purity: ≥98% JWH 250 N-(5-hydroxypentyl) metabolite
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀
UV/Vis.: λ_{max}: 212, 245, 303 nm
Supplied as: A solution in methanol
Storage: -20°C
Stability: ≥1 year



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

Laboratory Procedures

JWH 250 N-(5-hydroxypentyl) metabolite-d₅ is intended for use as an internal standard for the quantification of JWH 250 N-(5-hydroxypentyl) metabolite (Item No. 9000767) 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 250 N-(5-hydroxypentyl) metabolite-d₅ is supplied as a solution in methanol. To change the solvent, simply evaporate the methanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of JWH 250 N-(5-hydroxypentyl) metabolite-d₅ in these solvents is approximately 14 and 20 mg/ml, respectively.

Description

Cannabimimetic indoles, including certain “JWH” compounds, have been identified in herbal blends.¹ Hydroxylated and glucuronidated metabolites of two closely-related compounds, JWH 015 and JWH 018, have been identified from *in vitro* liver microsomal metabolism and from urine, respectively.²⁻⁴ JWH 250 is a cannabimimetic indole that is structurally- and functionally-related to JWH 015 and JWH 018. JWH 250 N-(5-hydroxypentyl) metabolite is expected to be a metabolite of JWH 250 that would be detectable both in serum and in urine.

References

1. Uchiyama, N., Kikura-Hanajiri, R., Kawahara, N., *et al. Forensic Toxicol.* **27**, 61-66 (2009).
2. Zhang, Q., Ma, P., Cole, R.B., *et al. Anal. Bioanal. Chem.* **386**, 1345-1355 (2006).
3. Sobolevsky, T., Prasolov, I., and Rodchenkov, G. *Forensic Sci. Int.* **200**, 141-147 (2010).
4. Moran, C.L., Le, V.H., Chimalakonda, K.C., *et al. 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, 05/23/2022

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