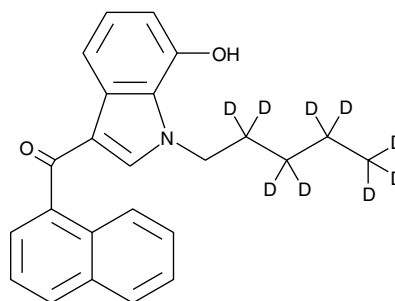


Product Information



JWH 018 7-hydroxyindole metabolite-d₉ Item No. 10715

Formal Name: (7-hydroxy-1-pentyl-1H-indol-3-yl)naphthalen-1-yl)-methanone-2,2',3,3',4,4',5,5,5-d₉
MF: C₂₄H₁₄D₉NO₂
FW: 366.5
Chemical Purity: ≥98%
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₉); ≤1% d₀
Stability: ≥2 years at -20°C
Supplied as: A solution in methanol
UV/Vis.: λ_{max}: 218, 340 nm



Laboratory Procedures

JWH 018 7-hydroxyindole metabolite-d₉ contains nine deuterium atoms at the 2, 2', 3, 3', 4, 4', 5, 5, and 5 positions. It is intended for use as an internal standard for the quantification of JWH 018 7-hydroxyindole metabolite-d₉ by GC- or LC-mass spectrometry (MS). For long term storage, we suggest that JWH 018 7-hydroxyindole metabolite-d₉ be stored as supplied at -20°C. It should be stable for at least two years.

JWH 018 7-hydroxyindole metabolite-d₉ 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 ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of JWH 018 7-hydroxyindole metabolite-d₉ in these solvents is approximately 10 mg/ml.

JWH 018 7-hydroxyindole metabolite-d₉ is used as an internal standard for the quantification of JWH 018 7-hydroxyindole metabolite by stable isotope dilution 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 018 7-hydroxyindole metabolite is a minor monohydroxylated urinary metabolite of JWH 018, a WIN 55,212-2 derivative that is a mildly selective agonist of the peripheral cannabinoid receptor.¹

Reference

1. Sobolevsky, T., Prasolov, I., and Rodchenkov, G. Detection of JWH-018 metabolites in smoking mixture post-administration urine. *Forensic Sci. Int.* **200**, 141-147 (2010).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10715

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY; NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog**.

Copyright Cayman Chemical Company, 08/03/2012