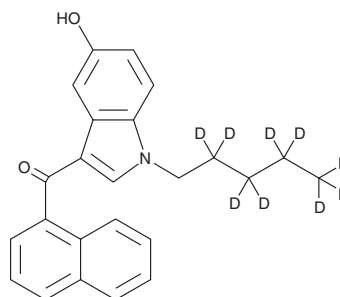


PRODUCT INFORMATION



JWH 018 5-hydroxyindole metabolite-d₉ Item No. 10713

CAS Registry No.: 2748533-34-6
Formal Name: (5-hydroxy-1-pentyl-2,2,3,3,4,4,5,5,5-d₉-1H-indol-3-yl)(naphthalen-1-yl)-methanone
MF: C₂₄H₁₄D₉NO₂
FW: 366.5
Chemical Purity: ≥98% (JWH 018 5-hydroxyindole metabolite)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₉); ≤1% d₀
UV/Vis.: λ_{max}: 219, 256, 280, 324 nm
Supplied as: A solution in methanol
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

JWH 018 5-hydroxyindole metabolite-d₉ is intended for use as an internal standard for the quantification of JWH 018 5-hydroxyindole metabolite (Item No. 9000852) 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 018 5-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 DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of JWH 018 5-hydroxyindole metabolite-d₉ in these solvents is approximately 30 mg/ml.

Description

JWH 018 5-hydroxyindole metabolite is a major monohydroxylated urinary metabolite of JWH 018, a WIN 55,212-2 derivative that is a mildly selective agonist of the peripheral cannabinoid receptor. In urine samples, this metabolite is almost completely glucuronidated.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(1-3)**, 141-147 (2010).

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, 07/18/2023

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