

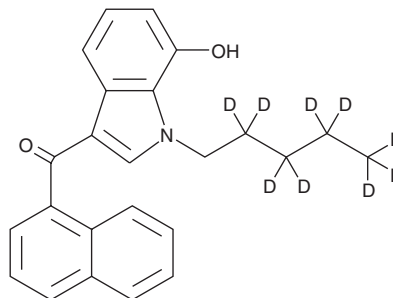
# PRODUCT INFORMATION



## JWH 018 7-hydroxyindole metabolite-d<sub>9</sub> Item No. 10715

**CAS Registry No.:** 2749349-17-3  
**Formal Name:** (7-hydroxy-1-pentyl-2,2,3,3,4,4,5,5,5-d<sub>9</sub>-1H-indol-3-yl)-1-naphthalenyl-methanone  
**MF:** C<sub>24</sub>H<sub>14</sub>D<sub>9</sub>NO<sub>2</sub>  
**FW:** 366.5  
**Chemical Purity:** ≥98% (JWH 018 7-hydroxyindole metabolite)

**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>9</sub>); ≤1% d<sub>0</sub>  
**UV/Vis.:** λ<sub>max</sub>: 218, 340 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.

### Description

JWH 018 7-hydroxyindole metabolite-d<sub>9</sub> (Item No. 10715) is intended for use as an internal standard for the quantification of JWH 018 7-hydroxyindole metabolite 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 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.<sup>1</sup>

### 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, 06/13/2023

### 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