

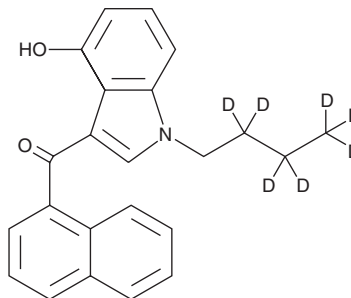
# PRODUCT INFORMATION



## JWH 073 4-hydroxyindole metabolite-d<sub>7</sub> Item No. 10717

**CAS Registry No.:** 2748464-25-5  
**Formal Name:** (1-butyl-(2,2,3,3,4,4,4-d<sub>7</sub>)-4-hydroxy-1H-indol-3-yl)(naphthalen-1-yl)-methanone  
**MF:** C<sub>23</sub>H<sub>14</sub>D<sub>7</sub>NO<sub>2</sub>  
**FW:** 350.5  
**Chemical Purity:** ≥98% (JWH 073 4-hydroxyindole metabolite)

**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>7</sub>); ≤1% d<sub>0</sub>  
**UV/Vis.:** λ<sub>max</sub>: 218, 355 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 073 4-hydroxyindole metabolite-d<sub>7</sub> (Item No. 10717) is intended for use as an internal standard for the quantification of XX (Item No. XX) 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 073 is a mildly selective agonist of the central cannabinoid (CB<sub>1</sub>) receptor derived from the aminoalkylindole WIN 55,212-2. The K<sub>i</sub> values for binding CB<sub>1</sub> and the peripheral cannabinoid (CB<sub>2</sub>) receptor are 8.9 and 38 nM, respectively for a CB<sub>1</sub>:CB<sub>2</sub> ratio of 0.23.<sup>1</sup> JWH 073 is one of several synthetic CBs which have been included in smoking mixtures. JWH 073 4-hydroxyindole metabolite is expected to be a urinary metabolite of JWH 073 based on the metabolism of the closely-related JWH 015 and JWH 018.<sup>2,3</sup>

### References

1. Aung, M.M., Griffin, G., Huffman, J.W., *et al.* Influence of the N-1 alkyl chain length of cannabimimetic indoles upon CB<sub>1</sub> and CB<sub>2</sub> receptor binding. *Drug Alcohol Depend.* **60(2)**, 133-140 (2000).
2. Zhang, Q., Ma, P., Cole, R.B., *et al.* Identification of in vitro metabolites of JWH-015, an aminoalkylindole agonist for the peripheral cannabinoid receptor (CB<sub>2</sub>) by HPLC-MS/MS. *Anal. Bioanal. Chem.* **386(5)**, 1345-1355 (2006).
3. 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