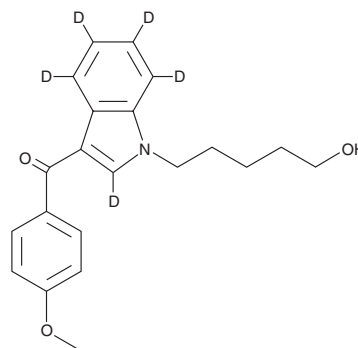


# PRODUCT INFORMATION



## RCS-4 N-(5-hydroxypentyl) metabolite-d<sub>5</sub> Item No. 14364

**CAS Registry No.:** 2703972-65-8  
**Formal Name:** (1-(5-hydroxypentyl)-1H-indol-3-yl-2,4,5,6,7-d<sub>5</sub>)  
(4-methoxyphenyl)methanone  
**MF:** C<sub>21</sub>H<sub>18</sub>D<sub>5</sub>NO<sub>3</sub>  
**FW:** 342.4  
**Chemical Purity:** ≥98% (RCS-4 N-(5-hydroxypentyl) metabolite)  
**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>5</sub>); ≤1% d<sub>0</sub>  
**UV/Vis.:** λ<sub>max</sub>: 214, 265, 313 nm  
**Supplied as:** A solution in acetonitrile  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Description

RCS-4 N-(5-hydroxypentyl) metabolite-d<sub>5</sub> is intended for use as an internal standard for the quantification of RCS-4 N-(5-hydroxypentyl) 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).

RCS-4 (Item No. 10645) is structurally similar to certain JWH compounds, synthetic cannabinoids used in herbal mixtures meant to mimic cannabis. RCS-4 itself has been identified in herbal mixtures.<sup>1</sup> RCS-4 N-(5-hydroxypentyl) metabolite (Item No. 10935) is a potential metabolite of RCS-4. A similar JWH metabolite has been detected in human urine following JWH smoking.<sup>2</sup> This metabolite is almost completely glucuroconjugated in urine samples, requiring enzymolysis of the sample before analysis.<sup>2</sup>

### References

1. Kikura-Hanajiri, R., Uchiyama, N., and Goda, Y. Survey of current trends in the abuse of psychotropic substances and plants in Japan. *Leg. Med. (Tokyo)* **13(3)**, 109-15 (2011).
2. 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, 06/16/2023

#### 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