

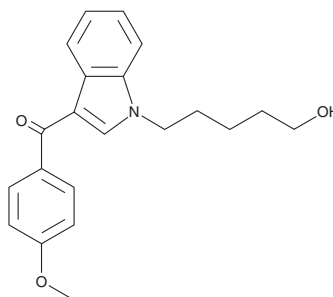
# PRODUCT INFORMATION



## RCS-4 N-(5-hydroxypentyl) metabolite

Item No. 10935

**CAS Registry No.:** 1379604-66-6  
**Formal Name:** (1-(5-hydroxypentyl)-1H-indol-3-yl)  
(4-methoxyphenyl)-methanone  
**MF:** C<sub>21</sub>H<sub>23</sub>NO<sub>3</sub>  
**FW:** 337.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 214, 262, 319 nm  
**Supplied as:** A 10 mg/ml 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 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> This product is intended for research and forensic applications.

### 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)**, 109-115 (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, 11/20/2025

#### 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