

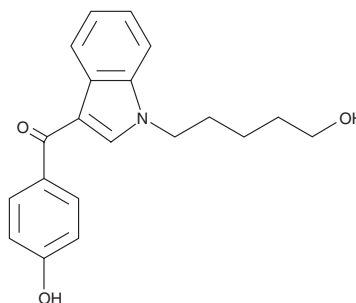
# PRODUCT INFORMATION



## RCS-4 M10 metabolite

Item No. 11978

**CAS Registry No.:** 1952336-77-4  
**Formal Name:** (1-(5-hydroxypentyl)-1H-indol-3-yl)(4-hydroxyphenyl)methanone  
**Synonym:** RCS-4 N-(5-hydroxypentyl)phenol metabolite  
**MF:** C<sub>20</sub>H<sub>21</sub>NO<sub>3</sub>  
**FW:** 323.4  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 214, 272, 319 nm  
**Supplied as:** A crystalline solid  
**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 (Item No. 10645) is a synthetic cannabinoid which has been detected in herbal mixtures.<sup>1</sup> RCS-4 M10 metabolite is a major urinary metabolite of RCS-4, characterized by O-demethylation/monohydroxylation of the N-pentyl chain.<sup>2</sup> Its metabolism has been derived using GC-MS analysis.<sup>2</sup> This metabolite is potentially useful as a metabolic marker for the identification of RCS-4 ingestion and is therefore intended for research and forensic purposes.

### 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. Kavanagh, P., Grigoryev, A., Melnik, A., et al. The identification of the urinary metabolites of 3-(4-methoxybenzoyl)-1-pentylindole (RCS-4), a novel cannabinimetic, by gas chromatography-mass spectrometry. *J. Anal. Toxicol.* **36**, 303-311 (2012).

#### 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](http://WWW.CAYMANCHEM.COM)