

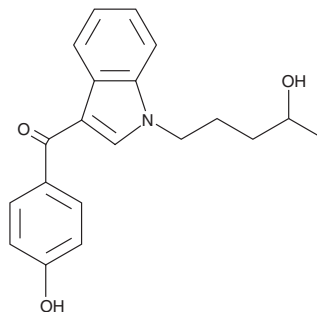
PRODUCT INFORMATION



RCS-4 M9 metabolite

Item No. 11977

CAS Registry No.: 1630022-96-6
Formal Name: [1-(4-hydroxypentyl)-1H-indol-3-yl]
(4-hydroxyphenyl)-methanone
Synonym: RCS-4 N-(4-hydroxypentyl)phenol metabolite
MF: C₂₀H₂₁NO₃
FW: 323.4
Purity: ≥98%
UV/Vis.: λ_{max}: 214, 262, 320 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

RCS-4 is a synthetic cannabinoid (CB) which has been detected in herbal mixtures.¹ RCS-4 M9 metabolite (Item No. 11977) is an expected metabolite of RCS-4 (Item No. 10645), characterized by hydroxylation on both the alkyl and phenyl groups. Such dihydroxylated metabolites have been shown to result during the metabolism of similar synthetic CBs both *in vitro*, with liver microsomes, and *in vivo*.²⁻⁴ This product is 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), 109-115 (2011).
2. Wintermeyer, A., Möller, I., Thevis, M., et al. In vitro phase I metabolism of the synthetic cannabimimetic JWH-018. *Anal. Bioanal. Chem.* **398**(5), 2141-2153 (2010).
3. 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 (CB₂) by HPLC-MS/MS. *Anal. Bioanal. Chem.* **386**(5), 1345-1355 (2006).
4. 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, 07/24/2023

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