

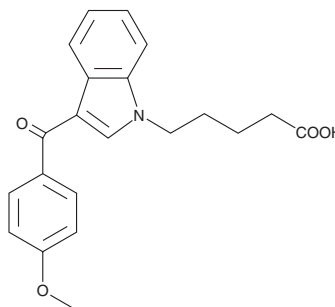
PRODUCT INFORMATION



RCS-4 N-(5-carboxypentyl) metabolite

Item No. 10937

CAS Registry No.: 1427521-39-8
Formal Name: 5-(3-(4-methoxybenzoyl)-1H-indol-1-yl)pentanoic acid
Synonym: RCS-4 N-pentanoic acid metabolite
MF: C₂₁H₂₁NO₄
FW: 351.4
Purity: ≥98%
UV/Vis.: λ_{max}: 213, 263, 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 which is structurally similar to JWH 018. Like JWH 018, RCS-4 has been detected in herbal products.¹ RCS-4 N-(5-carboxypentyl) metabolite is an expected metabolite of RCS-4. Carboxylation of the N-alkyl chain of JWH 018 occurs during *in vivo* metabolism, resulting in a 5-carboxypentyl metabolite that is detectable in the urine.²

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, 06/19/2023

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