

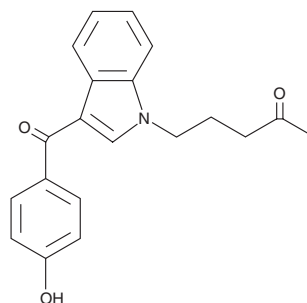
PRODUCT INFORMATION



RCS-4 M11 metabolite

Item No. 11979

CAS Registry No.: 1952336-78-5
Formal Name: 5-[3-(4-hydroxybenzoyl)-1H-indol-1-yl]-2-pentanone
Synonym: RCS-4 N-(4-oxo-pentyl)phenol metabolite
MF: C₂₀H₁₉NO₃
FW: 321.4
Purity: ≥95%
UV/Vis.: λ_{max}: 213, 273, 319 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 (Item No. 10645) is a synthetic cannabinoid which has been detected in herbal mixtures.¹ RCS-4 M11 metabolite is a major urinary metabolite of RCS-4.² It is derived by O-demethylation and oxidation of the N-pentyl chain to a ketone on the parent compound. This metabolite is potentially useful as a metabolic marker for the identification of RCS-4 ingestion and 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)**, 109-15 (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 cannabimimetic, 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, 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