

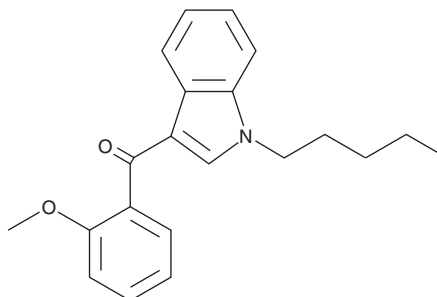
PRODUCT INFORMATION



RCS-4 2-methoxy isomer

Item No. 10865

CAS Registry No.: 1345966-76-8
Formal Name: (2-methoxyphenyl)(1-pentyl-1H-indol-3-yl)-methanone
MF: C₂₁H₂₃NO₂
FW: 321.4
Purity: ≥98%
UV/Vis.: λ_{max}: 214, 252, 313 nm
Supplied as: A solution in methanol
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 (CB) that has been detected as an adulterant of herbal mixtures.¹ RCS-4 2-methoxy isomer is an analog of RCS-4 that differs only in the location of the methoxy group on the phenyl ring. The affinity of RCS-4 2-methoxy isomer for CB receptors, or its physiological effects, have not been documented. However, RCS-4 2-methoxy isomer structurally resembles JWH 250 (Item No. 13634), a cannabimimetic indole that shows a high-affinity for both CB₁ and CB₂ (K_i = 11 and 33 nM, respectively).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. Huffman, J.W., Szklennik, P.V., Almond, A., et al. 1-Pentyl-3-phenylacetylindoles, a new class of cannabimimetic indoles. *Bioorg. Med. Chem. Lett.* **15(18)**, 4110-4113 (2005).

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, 06/16/2023

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