

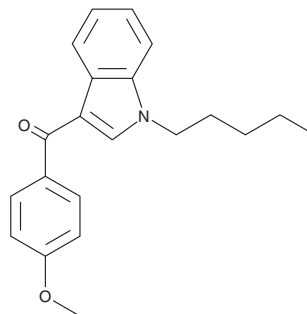
PRODUCT INFORMATION



RCS-4

Item No. 10645

CAS Registry No.: 1345966-78-0
Formal Name: (4-methoxyphenyl)(1-pentyl-1H-indol-3-yl)methanone
Synonyms: BTM-4, E-4, OBT-199, SR-19
MF: C₂₁H₂₃NO₂
FW: 321.4
Purity: ≥98%
UV/Vis.: λ_{max}: 214, 263, 320 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥1 year



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

Description

RCS-4 (Item No. 10645) is an analytical reference standard structurally similar to known synthetic cannabinoids. It has been found in various herbal incense products.¹ RCS-4 is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

Reference

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).

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, 09/15/2017

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