

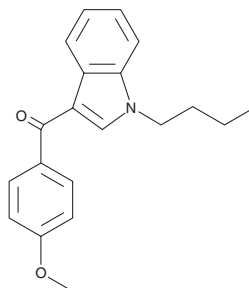
PRODUCT INFORMATION



RCS-4-C4 homolog

Item No. 10798

CAS Registry No.: 1345966-77-9
Formal Name: (4-methoxyphenyl)(1-butyl-1H-indol-3-yl)-methanone
Synonyms: BTM-4, E-4, OBT-199, SR-19
MF: C₂₀H₂₁NO₂
FW: 307.4
Purity: ≥98%
UV/Vis.: λ_{max}: 214, 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 (Item No. 10645) is a cannabimimetic compound similar to JWH 018, a potent cannabinoid receptor agonist. RCS-4, like JWH 018, has been identified in herbal blends. RCS-4-C4 homolog is identical to RCS-4 except the N-1 alkyl chain length has been shortened from C5 to C4. The biological activities of RCS-4-C4 homolog have not been reported.

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, 06/15/2023

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