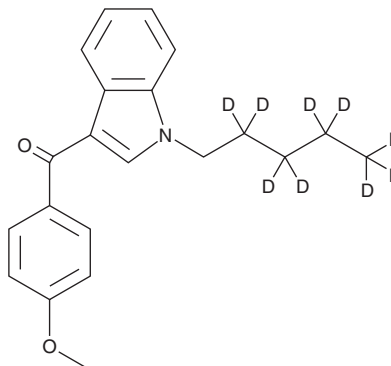


PRODUCT INFORMATION



RCS-4-d₉
Item No. 10513

CAS Registry No.: 2749316-95-6
Formal Name: (4-methoxyphenyl)(1-pentyl-d₉-1H-indol-3-yl)methanone
Synonyms: BTM-4-d₉, E-4-d₉, OBT-199-d₉, SR-19-d₉
MF: C₂₁H₁₄D₉NO₂
FW: 330.5



Chemical Purity: ≥98% (RCS-4)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₉); ≤1% d₀
UV/Vis.: λ_{max}: 213, 265, 321 nm
Supplied as: A solution in methanol
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-d₉ (Item No. 10513) is intended for use as an internal standard for the quantification of RCS-4 (Item Nos. 10645 | 15666) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

RCS-4 is structurally similar to known synthetic cannabinoids. It has been found in several herbal incense products.¹ RCS-4-d₉ 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)**, 109-115 (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, 05/20/2022

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