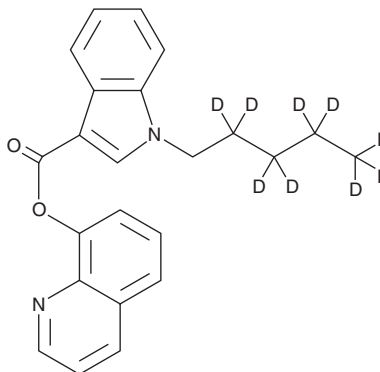


PRODUCT INFORMATION



PB-22-d₉
Item No. 14298

CAS Registry No.: 2484977-00-4
Formal Name: 1-pentyl-2,2,3,3,4,4,5,5,5-d₉-1H-indole-3-carboxylic acid, 8-quinolinyl ester
Synonym: QUPIC-d₉
MF: C₂₃H₁₃D₉N₂O₂
FW: 367.5
Chemical Purity: ≥98% (PB-22)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₉); ≤1% d₀
UV/Vis.: λ_{max}: 203, 217, 230, 294 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

PB-22-d₉ (Item No. 14298) is intended for use as an internal standard for the quantification of PB-22 (Item Nos. ISO00122 | 18203) 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).

PB-22 is categorized as a synthetic cannabinoid.¹ It mimics the effects of Δ⁹-THC (Item Nos. 12068 | ISO60157). PB-22-d₉ is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

Reference

1. Banister, S.D., Stuart, J., Kevin, R.C., *et al.* Effects of bioisosteric fluorine in synthetic cannabinoid designer drugs JWH-018, AM-2201, UR-144, XLR-11, PB-22, 5F-PB-22, APICA, and STS-135. *ACS Chem. Neurosci.* **6**(8), 1445-1458 (2015).

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