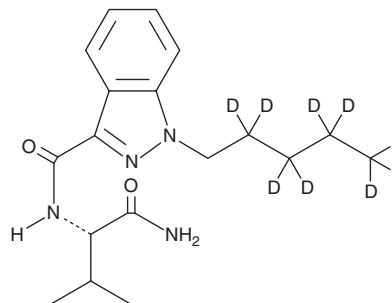


PRODUCT INFORMATION



AB-PINACA-d₉ Item No. 15057

CAS Registry No.: 2484976-98-7
Formal Name: N-[(1S)-1-(aminocarbonyl)-2-methylpropyl]-1-pentyl-2,2,3,3,4,4,5,5,5-d₉-1H-indazole-3-carboxamide
MF: C₁₈H₁₇D₉N₄O₂
FW: 339.5
Chemical Purity: ≥98% (AB-PINACA)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₉); ≤1% d₀
UV/Vis.: λ_{max}: 302 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

AB-PINACA-d₉ (Item No. 15057) is intended for use as an internal standard for the quantification of AB-PINACA (Item Nos. 14038 | ISO60210) 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).

AB-PINACA is categorized as a cannabinoid.¹ AB-PINACA-d₉ is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

Reference

1. Uchiyama, N., Matsuda, S., Wakana, D., *et al.* New cannabimimetic indazole derivatives, N-(1-amino-3-methyl-1-oxobutan-2-yl)-1-pentyl-1H-indazole-3-carboxamide (AB-PINACA) and N-(1-amino-3-methyl-1-oxobutan-2-yl)-1-(4-fluorobenzyl)-1H-indazole-3-carboxamide (AB-FUBINACA) identified as designer drugs in illegal products. *Forensic Toxicol.* **31(1)**, 93-100 (2013).

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, 06/26/2023

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