

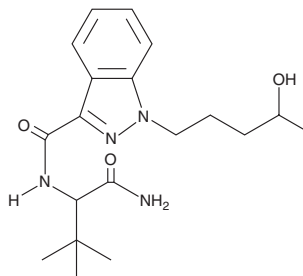
PRODUCT INFORMATION



ADB-PINACA N-(4-hydroxypentyl) metabolite

Item No. 15053

CAS Registry No.: 2748155-91-9
Formal Name: N-[1-(aminocarbonyl)-2,2-dimethylpropyl]-1-(4-hydroxypentyl)-1H-indazole-3-carboxamide
MF: C₁₉H₂₈N₄O₃
FW: 360.5
Purity: ≥98%
UV/Vis.: λ_{max}: 303 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

ADB-PINACA (Item No. 14763) is an analog of AKB48 (Item No. ISO00060), a synthetic cannabinoid also known as APINACA. ADB-PINACA N-(4-hydroxypentyl) metabolite is an expected phase 1 metabolite of ADB-PINACA, based on the known metabolism of similar compounds.^{1,2} The physiological properties of this compound have not been evaluated. This product is intended for forensic and research applications.

References

1. Zhang, Q., Ma, P., Cole, R.B., *et al.* Identification of *in vitro* metabolites of JWH-015, an aminoalkylindole agonist for the peripheral cannabinoid receptor (CB₂) by HPLC-MS/MS. *Anal. Bioanal. Chem.* **386**(5), 1345-1355 (2006).
2. Wintermeyer, A., Möller, I., Thevis, M., *et al.* *In vitro* phase I metabolism of the synthetic cannabinimimetic JWH-018. *Anal. Bioanal. Chem.* **398**(5), 2141-53 (2010).

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