

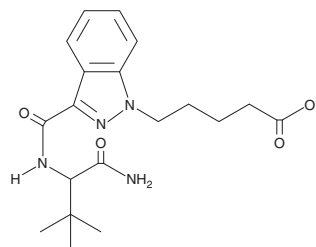
PRODUCT INFORMATION



ADB-PINACA pentanoic acid metabolite

Item No. 15055

CAS Registry No.: 2460433-29-6
Formal Name: 3-[[[1-(aminocarbonyl)-2,2-dimethylpropyl] amino]carbonyl]-1H-indazole-1-pentanoic acid
MF: C₁₉H₂₆N₄O₄
FW: 374.4
Purity: ≥95%
UV/Vis.: λ_{max}: 302 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 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), the pentyl indazole with structural similarity to JWH 018 adamantyl carboxamide (Item No. 9001193) and STS-135 (Item No. 11564), which are synthetic cannabinoids (CBs) with potential to be sold for recreational abuse.¹ ADB-PINACA pentanoic acid metabolite is an expected compound produced by phase I metabolism of ADB-PINACA *in vivo*, based on the metabolism of comparable alkylindole CBs.^{2,3} The physiological and toxicological properties of this compound have not been determined. This product is intended for research and forensic applications.

References

1. Drug Enforcement Administration AKB48 (APINACA) and 5F-AKB48 (5F-APINACA), (2013).
2. Sobolevsky, T., Prasolov, I., and Rodchenkov, G. Detection of JWH-018 metabolites in smoking mixture post-administration urine. *Forensic Sci. Int.* **200**, 141-147 (2010).
3. Chimalakonda, K.C., Moran, C.L., Kennedy, P.D., *et al.* Solid-phase extraction and quantitative measurement of omega and omega-1 metabolites of JWH-018 and JWH-073 in human urine. *Anal. Chem.* **83**(16), 6381-6388 (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, 06/14/2023

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