

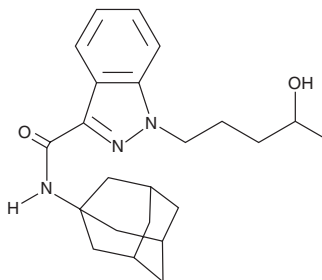
PRODUCT INFORMATION



AKB48 N-(4-hydroxypentyl) metabolite

Item No. 14221

CAS Registry No.: 1843184-41-7
Formal Name: 1-(4-hydroxypentyl)-N-tricyclo[3.3.1.1.3,7]dec-1-yl-1H-indazole-3-carboxamide
Synonym: APINACA N-(4-hydroxypentyl) metabolite
MF: C₂₃H₃₁N₃O₂
FW: 381.5
Purity: ≥98%
UV/Vis.: λ_{max}: 210, 302 nm
Supplied as: A 10 mg/ml solution in methanol
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

AKB48 is a pentyl indazole with structural similarity to JWH 018 adamantyl carboxamide (Item No. 9001193) and STS-135 (Item No. 11564), which are synthetic cannabinoids that may be sold for recreational use.¹ AKB48 N-(4-hydroxypentyl) metabolite is an expected phase 1 metabolite of AKB48, based on the known metabolism of similar compounds.² This product is intended for research and forensic applications.

References

1. Drug Enforcement Administration AKB48 (APINACA) and 5F-AKB48 (5F-APINACA), (2013).
2. Wintermeyer, A., Möller, I., Thevis, M., *et al.* *In vitro* phase I metabolism of the synthetic cannabimimetic JWH-018. *Anal. Bioanal. Chem.* **398**, 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, 12/08/2025

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