

PRODUCT INFORMATION



5-fluoro AKB48

Item No. 12065

CAS Registry No.: 1400742-13-3

Formal Name: N-((3s,5s,7s)-adamantan-1-yl)-1-(5-fluoropentyl)-1H-indazole-3-carboxamide

Synonyms: AKB48 N-(5-fluoropentyl) analog, 5-fluoro APINACA, APINACA 5-fluoropentyl analog

MF: $C_{23}H_{30}FN_3O$

FW: 383.5

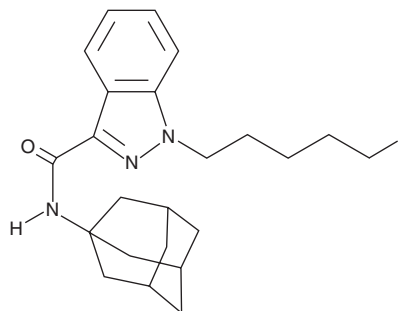
Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 209, 302 nm

Supplied as: A crystalline solid

Storage: $-20^{\circ}C$

Stability: ≥ 2 years



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

Description

AKB48 (Item No. ISO00060) is a pentylindazole with structural similarity to JWH018 adamantyl carboxamide (Item No. 9001193) and STS-135 (Item No. 11564), which are synthetic cannabinoids (CBs) that may be sold for recreational use. 5-fluoro AKB48 differs structurally from AKB48 by having fluorine at the terminal carbon of the pentyl chain. While the physiological properties of this compound are not known, quinolones with adamantyl-carboxamide moieties display high affinity for the peripheral CB_2 receptor but greatly reduced affinity for the central CB_1 receptor.¹ This product is intended for research and forensic applications.

Reference

1. Pasquini, S., De Rosa, M., Pedani, V., *et al.* Investigations on the 4-quinolone-3-carboxylic acid motif. 4. Identification of new potent and selective ligands for the cannabinoid type 2 receptor with diverse substitution patterns and antihyperalgesic effects in mice. *J. Med. Chem.* **54**, 5444-5453 (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, 10/23/2025

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