

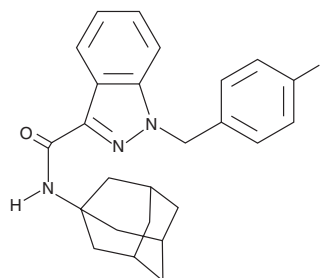
PRODUCT INFORMATION



AKB48 N-(4-fluorobenzyl) analog

Item No. 15937

CAS Registry No.: 2180933-90-6
Formal Name: 1-[(4-fluorophenyl)methyl]-N-tricyclo[3.3.1.1^{3,7}]dec-1-yl-1H-indazole-3-carboxamide
Synonyms: AFB-48, AFUBINACA, FUB-AKB48, FUB-APINACA
MF: C₂₅H₂₆FN₃O
FW: 403.5
Purity: ≥98%
UV/Vis.: λ_{max}: 210, 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

AKB48 (Item No. ISO00060) is a pentyl indazole with structural similarity to the synthetic cannabinoids (CBs) JWH 018 adamantyl carboxamide (Item No. 9001193) and STS-135 (Item No. 11564), which have been sold for recreational use.¹ Generally, quinolones with adamantyl-carboxamide moieties display high affinity for the peripheral cannabinoid (CB₂) receptor but greatly reduced affinity for the central cannabinoid (CB₁) receptor.² AKB48 N-(4-fluorobenzyl) analog is an AKB48 derivative in which the pentyl chain has been replaced with a 4-fluorobenzyl group in order to mimic a structural feature of the potent CB₁ modulator AB-FUBINACA (Item No. 14039). The physiological properties of this compound are not known. AKB48 N-(4-fluorobenzyl) analog is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

References

1. AKB48 (APINACA) and 5F-AKB48 (5F-APINACA), *Drug Enforcement Administration* (2013).
2. 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**(15), 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, 06/29/2023

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