

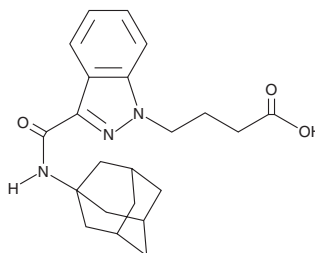
PRODUCT INFORMATION



4-fluoro ABUTINACA N-butanoic acid metabolite

Item No. 30749

Formal Name: 4-(3-(((3s,5s,7s)-adamantan-1-yl)carbamoyl)-1H-indazol-1-yl)butanoic acid
MF: C₂₂H₂₇N₃O₃
FW: 381.5
Purity: ≥98%
UV/Vis.: λ_{max}: 210, 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

4-fluoro ABUTINACA N-butanoic acid metabolite (Item No. 30749) is an analytical reference standard that is structurally similar to known synthetic cannabinoids. 4-fluoro ABUTINACA N-butanoic acid metabolite is a potential metabolite of 4-fluoro ABUTINACA (Item No. 30724) based on the published metabolism of 5-fluoro AKB48 (Item No. 12065).¹ At the time 4-fluoro ABUTINACA N-butanoic acid metabolite was made available for purchase, specific metabolism data had not been published. Contact us if updated information on this molecule is now available. This product is intended for research and forensic applications.

Reference

1. Holm, N.B., Pedersen, A.J., Dalsgaard, P.W., *et al.* Metabolites of 5F-AKB-48, a synthetic cannabinoid receptor agonist, identified in human urine and liver microsomal preparations using liquid chromatography high-resolution mass spectrometry. *Drug Test Anal.* **7**(3), 199-206 (2015).

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, 06/28/2023

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