

PRODUCT INFORMATION

ADB-5'Br-PINACA 3,3-dimethylbutanoic acid metabolite

Item No. 42007

Formal Name: (S)-2-(5-bromo-1-pentyl-1H-indazole-3-carboxamido)-3,3-dimethylbutanoic acid

Synonym: ADB-P-5'Br-INACA 3,3-dimethylbutanoic acid metabolite

MF: C₁₉H₂₆BrN₃O₃

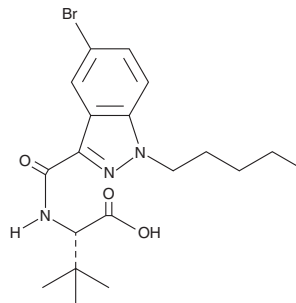
FW: 424.3

Purity: ≥98%

Supplied as: A 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-5'Br-PINACA 3,3-dimethylbutanoic acid metabolite (Item No. 42007) is an analytical reference standard categorized as a synthetic cannabinoid metabolite.¹ ADB-5'Br-PINACA 3,3-dimethylbutanoic acid metabolite is a metabolite of ADB-5'Br-PINACA (Item No. 37758). This product is intended for research and forensic applications.

Reference

1. Rautio, T., Obrist, R., Krebs, L., *et al.* In vitro metabolism study of ADB-P-5Br-INACA and ADB-4en-P-5Br-INACA using human hepatocytes, liver microsomes, and in-house synthesized references. *Drug Test Anal.* (2024).

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, 09/25/2024

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