

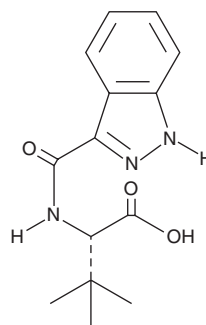
PRODUCT INFORMATION



MDMB-INACA 3,3-dimethylbutanoic acid metabolite

Item No. 38151

CAS Registry No.: 1189552-82-6
Formal Name: N-(1H-indazol-3-ylcarbonyl)-3-methyl-L-valine
Synonym: ADB-INACA 3,3-dimethylbutanoic acid metabolite
MF: C₁₄H₁₇N₃O₃
FW: 275.3
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Description

MDMB-INACA 3,3-dimethylbutanoic acid metabolite (Item No. 38151) is analytical reference standard categorized as a synthetic cannabinoid metabolite.¹ MDMB-INACA 3,3-dimethylbutanoic acid metabolite is a metabolite of 4-fluoro MDMB-BUTINACA (Item Nos. 28186 | 26645). It is also a potential metabolite of MDMB-INACA (Item No. 27248) and ADB-INACA (Item No. 37729) based on the published metabolism of 4-fluoro MDMB-BUTINACA. This product is intended for research and forensic applications.

Reference

1. Leong, H.S., Watanabe, S., Kuzhiumparambil, U., *et al.* Monitoring metabolism of synthetic cannabinoid 4F-MDMB-BINACA via high-resolution mass spectrometry assessed in cultured hepatoma cell line, fungus, liver microsomes and confirmed using urine samples. *Forensic Toxicol.* **39**, 198–212 (2021).

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, 04/11/2025

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