

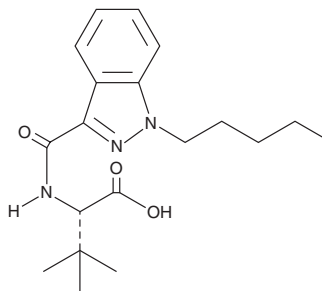
PRODUCT INFORMATION



MDMB-PINACA butanoic acid metabolite

Item No. 46332

CAS Registry No.: 2104034-59-3
Formal Name: 3-methyl-N-[(1-pentyl-1H-indazol-3-yl) carbonyl]-L-valine
Synonyms: ADB-PINACA butanoic acid metabolite, MDMB-PINACA 3,3-dimethylbutanoic acid metabolite
MF: $C_{19}H_{27}N_3O_3$
FW: 345.4
Purity: $\geq 98\%$
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

MDMB-PINACA butanoic acid metabolite (Item No. 46332) is an analytical reference standard categorized as a synthetic cannabinoid metabolite.¹ MDMB-PINACA butanoic acid metabolite is a metabolite of ADB-PINACA (Item Nos. ISO60212 | ISO00150). This product is intended for research and forensic applications.

Reference

1. Longworth, M., Connor, M., Banister, S.D., *et al.* Synthesis and pharmacological profiling of the metabolites of synthetic cannabinoid drugs APICA, STS-135, ADB-PINACA, and 5F-ADB-PINACA. *ACS Chem. Neurosci.* **8**(8), 1673-1680 (2017).

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, 07/30/2026

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