

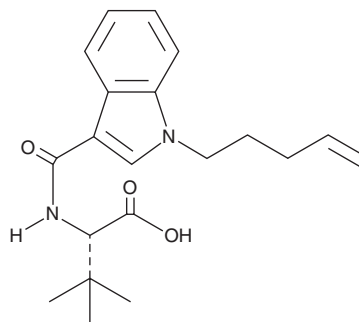
# PRODUCT INFORMATION



## MDMB-4en-PICA butanoic acid metabolite

Item No. 28943

**Formal Name:** (S)-3,3-dimethyl-2-(1-(pent-4-en-1-yl)-1H-indole-3-carboxamido)butanoic acid  
**Synonym:** MDMB-4en-PICA 3,3-dimethylbutanoic acid metabolite  
**MF:** C<sub>20</sub>H<sub>26</sub>N<sub>2</sub>O<sub>3</sub>  
**FW:** 342.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 218, 290 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

MDMB-4en-PICA butanoic acid metabolite (Item No. 28943) is an analytical reference standard that is structurally similar to known synthetic cannabinoids. MDMB-4en-PICA butanoic acid metabolite is a potential metabolite of MDMB-4en-PICA (Item No. 28541) based on the published metabolism of MDMB-PICA.<sup>1</sup> At the time MDMB-4en-PICA butanoic acid metabolite (Item No. 28943) 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. Franz, F., Jechle, H., Wilde, M., et al. Structure-metabolism relationships of valine and *tert*-leucine-derived synthetic cannabinoid receptor agonists: A systematic comparison of the in vitro phase I metabolism using pooled human liver microsomes and high-resolution mass spectrometry. *Forensic Toxicol.* **37**, 316-329 (2019).

#### 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, 07/25/2023

#### 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](http://WWW.CAYMANCHEM.COM)