

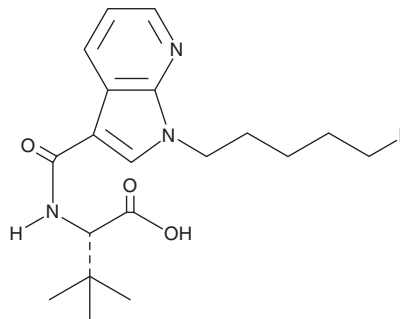
# PRODUCT INFORMATION



## 5-fluoro MDMB-7-PAICA butanoic acid metabolite

Item No. 28693

**CAS Registry No.:** 2712863-53-9  
**Formal Name:** N-[[1-(5-fluoropentyl)-1H-pyrrolo[2,3-b]pyridin-3-yl]carbonyl]-3-methyl-L-valine  
**Synonyms:** 7'-N-5-fluoro ADB butanoic acid metabolite, 5-fluoro MDMB-P7AICA butanoic acid metabolite  
**MF:** C<sub>19</sub>H<sub>26</sub>FN<sub>3</sub>O<sub>3</sub>  
**FW:** 363.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 217, 242, 292 nm  
**Supplied as:** A neat solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Description

5-fluoro MDMB-7-PAICA butanoic acid metabolite (Item No. 28693) is an analytical reference standard categorized as a synthetic cannabinoid metabolite.<sup>1</sup> It is a metabolite of 5-fluoro MDMB-7-PAICA (Item No. 25286). This product is intended for research and forensic applications.

### Reference

1. Richter, L.H.J., Herrmann, J., Andreas, A., et al. Tools for studying the metabolism of new psychoactive substances for toxicological screening purposes - a comparative study using pooled human liver S9, HepaRG cells, and zebrafish larvae. *Toxicol. Lett.* **305**, 73-80 (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, 07/26/2023

#### 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