

PRODUCT INFORMATION

4-fluoro MDMB-BUTINACA N-(butanoic acid) 3,3-dimethylbutanoic acid metabolite

Item No. 31643

Formal Name: (S)-2-(1-(3-carboxypropyl)-1H-indazole-3-carboxamido)-3,3-dimethylbutanoic acid
Synonyms: 4-fluoro MDMB-BINACA N-(butanoic acid) butanoic acid metabolite, 4-fluoro MDMB-BINACA N-(butanoic acid) ester hydrolysis metabolite, 4-fluoro MDMB-BUTINACA N-(butanoic acid) butanoic acid metabolite

MF: C₁₈H₂₃N₃O₅

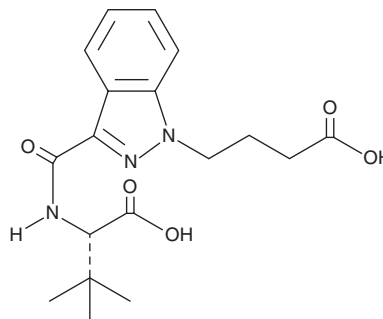
FW: 361.4

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

4-fluoro MDMB-BUTINACA N-(butanoic acid) 3,3-dimethylbutanoic acid metabolite (Item No. 31643) is an analytical reference standard categorized as a synthetic cannabinoid metabolite.¹ 4-fluoro MDMB-BUTINACA N-(butanoic acid) 3,3-dimethylbutanoic acid metabolite is a metabolite of 4-fluoro MDMB-BUTINACA (Item Nos. 28186 | 26645). This product is intended for research and forensic applications.

Reference

1. Haschimi, B., Mogler, L., Halter, S., *et al.* Detection of the recently emerged synthetic cannabinoid 4F-MDMB-BINACA in “legal high” products and human urine specimens. *Drug Test Anal.* **11(9)**, 1377-1386 (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, 06/22/2023

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