

# PRODUCT INFORMATION



## 4-fluoro MDMB-BUTINACA butanoic acid metabolite-d<sub>4</sub>

Item No. 37590

**Formal Name:** (S)-2-(1-(4-fluorobutyl)-1H-indazole-3-carboxamido-4,5,6,7-d<sub>4</sub>)-3,3-dimethylbutanoic acid

**Synonyms:** 4-fluoro MDMB-BINACA  
3,3-dimethylbutanoic acid-d<sub>4</sub>,  
4-fluoro MDMB-BINACA ester hydrolysis metabolite-d<sub>4</sub>

**MF:** C<sub>18</sub>H<sub>20</sub>D<sub>4</sub>FN<sub>3</sub>O<sub>3</sub>

**FW:** 353.4

**Chemical Purity:** ≥98% (4-fluoro MDMB-BUTINACA butanoic acid metabolite)

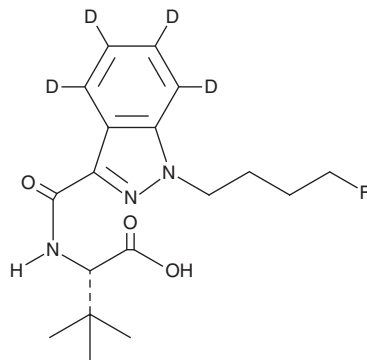
**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>4</sub>); ≤1% d<sub>0</sub>

**UV/Vis.:** λ<sub>max</sub>: 301 nm

**Supplied as:** A 5 mg/ml solution in acetonitrile

**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 butanoic acid metabolite-d<sub>4</sub> (Item No. 37590) is intended for use as an internal standard for the quantification of 4-fluoro MDMB-BUTINACA butanoic acid metabolite (Item No. 28187) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

4-fluoro MDMB-BUTINACA butanoic acid metabolite is categorized as a synthetic cannabinoid metabolite.<sup>1</sup> It is a metabolite of 4-fluoro MDMB-BUTINACA (Item No. 26645).<sup>1</sup> 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, 10/27/2025

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