

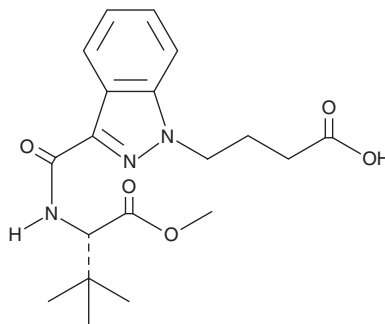
PRODUCT INFORMATION



4-fluoro MDMB-BUTINACA N-butanoic acid metabolite

Item No. 28941

Formal Name: (S)-4-(3-((1-methoxy-3,3-dimethyl-1-oxobutan-2-yl)carbamoyl)-1H-indazol-1-yl)butanoic acid
Synonym: 4-fluoro MDMB-BINACA N-butanoic acid metabolite
MF: C₁₉H₂₅N₃O₅
FW: 375.4
Purity: ≥98%
Supplied as: A solution in methyl acetate
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 metabolite (Item No. 28941) is an analytical reference standard that is structurally similar to known synthetic cannabinoids. 4-fluoro MDMB-BUTINACA N-butanoic acid metabolite is a potential metabolite of MDMB-BUTINACA (Item No. 28739) and 4-fluoro MDMB-BUTINACA (Item Nos. 28186 | 26645) based on the published metabolism of ADB-PINACA (Item Nos. ISO60212 | ISO00150) and 5-fluoro ADB-PINACA (Item No. 14764), respectively.¹ At the time 4-fluoro MDMB-BUTINACA N-butanoic acid metabolite (Item No. 28941) 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. 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, 07/26/2023

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