

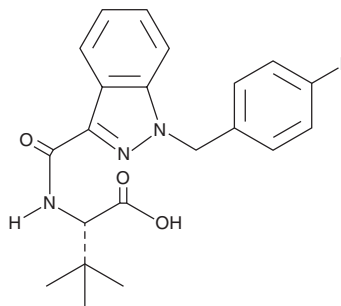
PRODUCT INFORMATION



MDMB-FUBINACA metabolite M1

Item No. 20061

CAS Registry No.: 2693397-47-4
Formal Name: N-[[1-[(4-fluorophenyl)methyl]-1H-indazol-3-yl]carbonyl]-3-methyl-L-valine
Synonym: MDMB-FUBINACA 3,3-dimethylbutanoic acid
MF: C₂₁H₂₂FN₃O₃
FW: 383.4
Purity: ≥98%
UV/Vis.: λ_{max}: 302 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-FUBINACA metabolite M1 (Item No. 20061) is an analytical reference standard that is structurally classified as a synthetic cannabinoid. It is an expected metabolite of the indazole-based synthetic cannabinoid MDMB-FUBINACA (Item No. 16966) and also the presumptive metabolite of ADB-FUBINACA (Item No. 14292). The physiological and toxicological properties of this compound have not been determined. This product is intended for research and forensic applications.

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, 06/20/2023

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