

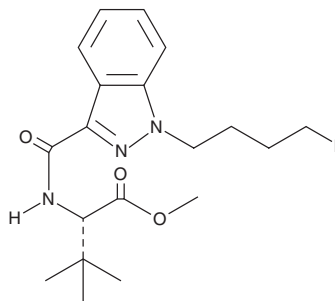
PRODUCT INFORMATION



4-fluoro MDMB-BUTINACA

Item No. 26645

CAS Registry No.: 2390036-46-9
Formal Name: N-[[1-(4-fluorobutyl)-1H-indazol-3-yl] carbonyl]-3-methyl-L-valine, methyl ester
Synonym: 4-fluoro MDMB-BINACA
MF: C₁₉H₂₆FN₃O₃
FW: 363.4
Purity: ≥98%
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

4-fluoro MDMB-BUTINACA (Item No. 26645) is an analytical reference material categorized as a synthetic cannabinoid.¹ 4-fluoro MDMB-BUTINACA has been associated with fatalities.² It is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards for reference materials.

References

1. Sparkes, E., Boyd, R., Chen, S., *et al.* Synthesis and pharmacological evaluation of newly detected synthetic cannabinoid receptor agonists AB-4CN-BUTICA, MMB-4CN-BUTINACA, MDMB-4F-BUTICA, MDMB-4F-BUTINACA and their analogs. *Front. Psychiatry* **13**, 1010501 (2022).
2. Simon, G., Tóth, D., Heckmann, V., *et al.* Simultaneous fatal poisoning of two victims with 4F-MDMB-BINACA and ethanol. *Forensic Toxicol.* (2022).

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, 10/25/2022

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