

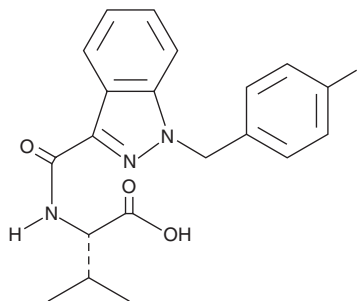
PRODUCT INFORMATION



AB-FUBINACA metabolite 3

Item No. 17761

CAS Registry No.: 1877243-60-1
Formal Name: N-[[1-[(4-fluorophenyl)methyl]-1H-indazol-3-yl]carbonyl]-L-valine
Synonyms: FUB-AMB 3-butanoic Acid metabolite, FUB-AMB 3-methylbutanoic Acid metabolite, MMB-FUBINACA metabolite 1
MF: C₂₀H₂₀FN₃O₃
FW: 369.4
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥3 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

AB-FUBINACA metabolite 3 (Item No. 17761) is an analytical reference material categorized as a synthetic cannabinoid.¹ AB-FUBINACA metabolite 3 is a metabolite of AB-FUBINACA (Item Nos. 14039 | ISO60211) and MMB-FUBINACA (Item Nos. 9001960 | 31090).^{1,2} 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.

References

1. Webb, H.D.J., Finlay, D.B., Chen, S., *et al.* Characterisation of AMB-FUBINACA metabolism and CB1-mediated activity of its acid metabolite. *Forensic Toxicol.* **41(1)**, 114-125 (2023).
2. Castaneto, M.S., Wohlfarth, A., Pang, S., *et al.* Identification of AB-FUBINACA metabolites in human hepatocytes and urine using high-resolution mass spectrometry. *Forensic Toxicol.* **33(2)**, 295-310 (2015).

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 its 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, 01/25/2023

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