

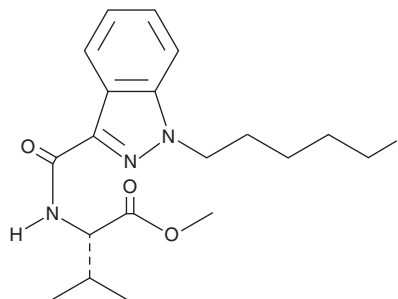
PRODUCT INFORMATION



5-fluoro AMB

Item No. 15489

CAS Registry No.: 1801552-03-3
Formal Name: N-[[1-(5-fluoropentyl)-1H-indazol-3-yl] carbonyl]-L-valine, methyl ester
Synonyms: 5-fluoro AMB-PINACA, 5-fluoro AMP, 5-fluoro MMB-PINACA
MF: C₁₉H₂₆FN₃O₃
FW: 363.4
Purity: ≥98%
UV/Vis.: λ_{max}: 209, 301 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

AB-PINACA (Item No. 14038) is a synthetic cannabinoid recently identified in illegal herbal products.¹ 5-fluoro AMB is an analog of AB-PINACA characterized by the replacement of a primary amine with a methoxy group and the addition of an alkyl-terminal fluorine atom. The physiological and toxicological properties of this compound have not been determined. This product is intended for research and forensic applications.

Reference

1. Uchiyama, N., Matsuda, S., Wakana, D., *et al.* New cannabimimetic indazole derivatives, N-(1-amino-3-methyl-1-oxobutan-2-yl)-1-pentyl-1H-indazole-3-carboxamide (AB-PINACA) and N-(1-amino-3-methyl-1-oxobutan-2-yl)-1-(4-fluorobenzyl)-1H-indazole-3-carboxamide (AB-FUBINACA) identified as designer drugs in illegal products. *Forensic Toxicol.* **31**(1), 93-100 (2013).

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, 06/28/2023

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