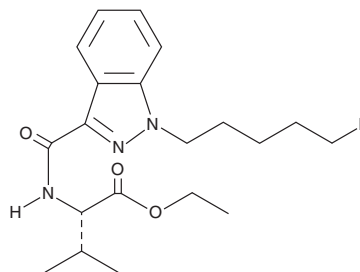


PRODUCT INFORMATION

5-fluoro AEB

Item No. 17735

CAS Registry No.: 2365471-51-6
Formal Name: N-[[1-(5-fluoropentyl)-1H-indazol-3-yl]carbonyl]-L-valine, ethyl ester
Synonym: 5-fluoro EMB-PINACA
MF: C₂₀H₂₈FN₃O₃
FW: 377.5
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥2 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 identified in illegal herbal products.¹ Variations of AB-PINACA have been synthesized including derivatives that include a fluorine atom added to the terminal carbon of the pentyl group. 5-fluoro AEB is an analog of AB-PINACA characterized by the replacement of the carboxamide with ethyl acetate and the addition of an alkyl-terminal fluorine atom. The physiological and toxicological properties of this compound have not been determined. 5-fluoro AEB 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.

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, 01/17/2023

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