

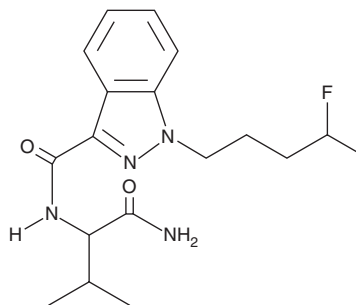
# PRODUCT INFORMATION



## AB-PINACA N-(4-fluoropentyl) isomer

Item No. 15441

**CAS Registry No.:** 2748160-98-5  
**Formal Name:** N-[1-(aminocarbonyl)-2-methylpropyl]-1-(4-fluoropentyl)-1H-indazole-3-carboxamide  
**Synonyms:** AB-PINACA N-(4-fluoropentyl) analog, 4-fluoro AB-PINACA  
**MF:** C<sub>18</sub>H<sub>25</sub>FN<sub>4</sub>O<sub>2</sub>  
**FW:** 348.4  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 209, 302 nm  
**Supplied as:** A solution in methanol  
**Storage:** -20°C  
**Stability:** ≥4 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.<sup>1</sup> Structurally, this designer drug is developed on an indazole base, which distinguishes it from the many JWH compounds having an indolyl base.<sup>1,2</sup> AB-PINACA N-(4-fluoropentyl) isomer is a derivative of AB-PINACA featuring a fluorine atom added to the pentyl group. The physiological and toxicological properties of this compound have not been determined. This product is intended for research and forensic applications.

### References

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).
2. Huffman, J.W., Zengin, G., Wu, M.J., *et al.* Structure-activity relationships for 1-alkyl-3-(1-naphthoyl) indoles at the cannabinoid CB<sub>1</sub> and CB<sub>2</sub> receptors: Steric and electronic effects of naphthoyl substituents. New highly selective CB<sub>2</sub> receptor agonists. *Bioorg. Med. Chem.* **13(1)**, 89-112 (2005).

#### 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