

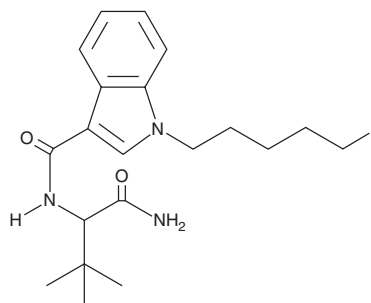
# PRODUCT INFORMATION



## 5-fluoro ADBICA

Item No. 14766

**CAS Registry No.:** 1863065-82-0  
**Formal Name:** N-[1-(aminocarbonyl)-2,2-dimethylpropyl]-1-(5-fluoropentyl)-1H-indole-3-carboxamide  
**Synonym:** 5-fluoro ADB-PICA  
**MF:** C<sub>20</sub>H<sub>28</sub>FN<sub>3</sub>O<sub>2</sub>  
**FW:** 361.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 217, 290 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥5 year



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

### Description

ADBICA is a synthetic cannabinoid that has recently been identified in herbal blends.<sup>1</sup> 5-fluoro ADBICA is a derivative of ADBICA featuring a fluorine atom added to the terminal carbon of the pentyl group. The physiological and toxicological properties of this compound have not been determined. This product is intended for forensic and research applications.

### Reference

1. Uchiyama, N., Matsuda, S., Kawamura, M., *et al.* Two new-type cannabimimetic quinolinyl carboxylates, QUPIC and QUCHIC, two new cannabimimetic carboxamide derivatives, ADB-FUBINACA and ADBICA, and five synthetic cannabinoids detected with a thiophene derivative α-PVT and an opioid receptor agonist AH-7921 identified in illegal products. *Forensic Toxicol.* **31(2)**, 223-240 (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, 06/22/2023

### 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