

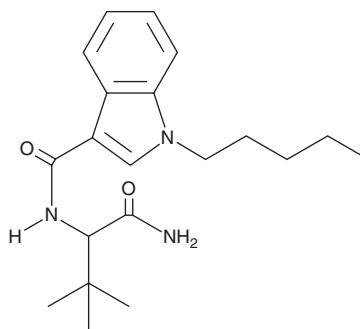
# PRODUCT INFORMATION



## ADBICA

Item No. 14293

**CAS Registry No.:** 1445583-48-1  
**Formal Name:** N-(1-amino-3,3-dimethyl-1-oxobutan-2-yl)-1-pentyl-1H-indole-3-carboxamide  
**Synonym:** ADB-PICA  
**MF:** C<sub>20</sub>H<sub>29</sub>N<sub>3</sub>O<sub>2</sub>  
**FW:** 343.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 218, 290 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

ADBICA is a synthetic cannabinoid (CB) that has recently been identified in herbal blends.<sup>1</sup> Its name refers to its structure, which has a 1-amino-3,3-dimethyl-1-oxobutan-2-yl group linked to a 1-pentyl-1H-indole-3-carboxamide base at the amide group. The base is similar to that of the aminoalkylindoles which potentially activate both CB receptors.<sup>2</sup> The physiological and toxicological properties of this compound have not been determined. This product is intended for forensic and research applications.

### References

1. Uchiyama, N., Matsuda, S., Kawamura, M., *et al.* Two new-type cannabimimetic quinoliny 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).
2. Aung, M.M., Griffin, G., Huffman, J.W., *et al.* Influence of the N-1 alkyl chain length of cannabimimetic indoles upon CB<sub>1</sub> and CB<sub>2</sub> receptor binding. *Drug Alcohol Depend.* **60(2)**, 133-140 (2000).

#### 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, 06/16/2023

#### 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