

# PRODUCT INFORMATION



## ADB-FUBINACA

Item No. 14292

**CAS Registry No.:** 1445583-51-6  
**Formal Name:** N-[1-(aminocarbonyl)-2,2-dimethylpropyl]-1-[(4-fluorophenyl)methyl]-1H-indazole-3-carboxamide

**MF:** C<sub>21</sub>H<sub>23</sub>FN<sub>4</sub>O<sub>2</sub>

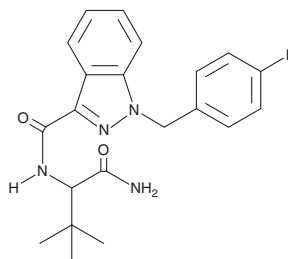
**FW:** 382.4

**Purity:** ≥98%

**Supplied as:** A neat solid

**Storage:** -20°C

**Stability:** As supplied, 3 years from the QC date provided on the Certificate of Analysis, when stored properly



### Description

ADB-FUBINACA is a synthetic cannabinoid 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 (ADB) group linked to a 4-fluorobenzyl-1H-indazole-3-carboxamide (FUBINACA) base at the amide group. The physiological and toxicological properties of this compound have not been determined. This product is intended for forensic and research applications.

This product is a qualified Reference Material (RM) that has been manufactured and tested to meet ISO17025 and Guide 34 guidelines. These materials are tested using validated analytical methods on qualified instrumentation to ensure traceability of measurements. All traceable RMs may be distinguished by their CofAs and can be downloaded using the batch number located on the product label. For a representative CofA please contact our technical support.

### Reference

1. Uchiyama, N., Matsuda, S., Kawamura, M., *et al.* Two new-type cannabimimetic quinolonyl carboxylates, QUPIC and QUCHIC, two new cannabimimetic carboxamide derivatives, ADB-FUBINACA and ADBICA, and five synthetic cannabinoids detected with a thiophene derivative a-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, 07/07/2016

#### 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