

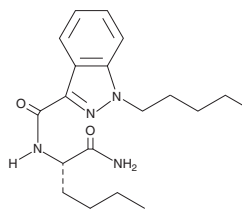
# PRODUCT INFORMATION



## ADB-PINACA isomer 3

Item No. 9001527

**Formal Name:** (S)-N-(1-amino-1-oxohexan-2-yl)-1-pentyl-1H-indazole-3-carboxamide  
**MF:** C<sub>19</sub>H<sub>28</sub>N<sub>4</sub>O<sub>2</sub>  
**FW:** 344.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 209, 302 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥5 years



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

### Description

ADB-PINACA (Item No. 14763) is an analog of AKB48 (Item No. ISO00060), a synthetic cannabinoid recently identified in herbal products.<sup>1</sup> ADB-PINACA isomer 3 is a regioisomer of ADB-PINACA. 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., Kawamura, M., Kikura-Hanajiri, R., *et al.* Identification of two new-type synthetic cannabinoids, N-(1-adamantyl)-1-pentyl-1H-indole-3-carboxamide (APICA) and N-(1-adamantyl)-1-pentyl-1H-indazole-3-carboxamide (APINACA), and detection of five synthetic cannabinoids, AM-1220, AM-2233, AM-1241, CB-13 (CRA-13), and AM-1248, as designer drugs in illegal products. *Forensic Toxicol.* **30**, 114-125 (2012).

#### 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, 07/12/2023

#### 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