

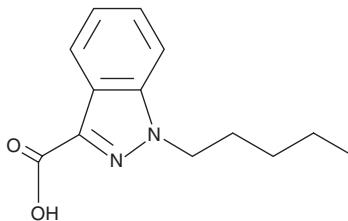
# PRODUCT INFORMATION



## AB-PINACA 3-carboxyindazole metabolite

Item No. 27732

CAS Registry No.: 1283576-17-9  
Formal Name: 1-pentyl-1H-indazole-3-carboxylic acid  
MF:  $C_{13}H_{16}N_2O_2$   
FW: 232.3  
Purity:  $\geq 98\%$   
UV/Vis.:  $\lambda_{\max}$ : 300 nm  
Supplied as: A crystalline solid  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 4$  years



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

### Description

AB-PINACA 3-carboxyindazole metabolite (Item No. 27732) is an analytical reference standard categorized as a synthetic cannabinoid metabolite.<sup>1</sup> AB-PINACA 3-carboxyindazole metabolite is a metabolite of APINAC (Item No. 18589). This product is intended for research and forensic applications.

### Reference

1. Hwang, J., Hwang, J., Ganganna, B., *et al.* Metabolic and pharmacokinetic characterization of a new synthetic cannabinoid APINAC in rats. *Forensic Toxicol.* **36**, 88-101 (2017).

#### 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, 05/20/2024

#### 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