

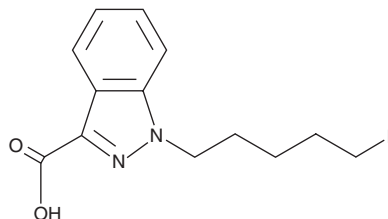
# PRODUCT INFORMATION



## 5-fluoro AB-PINACA 3-carboxyindazole metabolite

Item No. 27733

**CAS Registry No.:** 1535166-43-8  
**Formal Name:** 1-(5-fluoropentyl)-1H-indazole-3-carboxylic acid  
**MF:** C<sub>13</sub>H<sub>15</sub>FN<sub>2</sub>O<sub>2</sub>  
**FW:** 250.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 300 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Description

5-fluoro AB-PINACA 3-carboxyindazole metabolite (Item No. 27733) is an analytical reference standard that is structurally similar to known synthetic cannabinoid metabolites. 5-fluoro AB-PINACA 3-carboxyindazole metabolite is a precursor in the synthesis of 5-fluoro AB-PINACA (Item No. 14755), 5-fluoro ADB-PINACA, (S)-5-fluoro ADB (Item Nos. 25012 | 23631), 5-fluoro AMB (Item No. 15489), and 5-fluoro CUMYL-PINACA.<sup>1-3</sup> This product is intended for research and forensic applications.

### References

1. Banister, S.D., Moir, M., Stuart, J., *et al.* Pharmacology of indole and indazole synthetic cannabinoid designer drugs AB-FUBINACA, ADB-FUBINACA, AB-PINACA, ADB-PINACA, 5F-AB-PINACA, 5F-ADB-PINACA, ADBICA, and 5F-ADBICA. *ACS Chem. Neurosci.* **6(9)**, 1546-1559 (2015).
2. Banister, S.D., Longworth, M., Kevin, R., *et al.* Pharmacology of valinate and tert-leucinate synthetic cannabinoids 5F-AMBICA, 5F-AMB, 5F-ADB, AMB-FUBINACA, MDMB-FUBINACA, MDMB-CHMICA, and their analogues. *ACS Chem. Neurosci.* **7(9)**, 1241-1254 (2016).
3. Banister, S.D., Adams, A., Kevin, R.C., *et al.* Synthesis and pharmacology of new psychoactive substance 5F-CUMYL-P7AICA, a scaffold-hopping analog of synthetic cannabinoid receptor agonists 5F-CUMYL-PICA and 5F-CUMYL-PINACA. *Drug Test Anal.* **11(2)**, 279-291 (2019).

#### 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, 05/22/2024

#### 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