

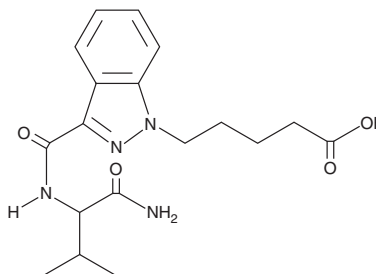
# PRODUCT INFORMATION



## AB-PINACA N-pentanoic acid metabolite

Item No. 15051

**CAS Registry No.:** 1879029-93-2  
**Formal Name:** 3-[[[1-(aminocarbonyl)-2-methylpropyl]amino]carbonyl]-1H-indazole-1-pentanoic acid  
**MF:** C<sub>18</sub>H<sub>24</sub>N<sub>4</sub>O<sub>4</sub>  
**FW:** 360.4  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 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

AB-PINACA (Item No. 14038) is a synthetic cannabinoid (CB) that has been identified in illegal herbal products.<sup>1</sup> AB-PINACA pentanoic acid metabolite is an expected compound produced by phase I metabolism of AB-PINACA *in vivo*, based on the metabolism of comparable alkylindole CBs.<sup>2,3</sup> The physiological and toxicological properties of this compound have not been determined. This product is intended for research and forensic applications.

### References

1. Uchiyama, N., Matsuda, S., Wakana, D., *et al.* New cannabimimetic indazole derivatives, N-(1-amino-3-methyl-1-oxobutan-2-yl)-1-pentyl-1H-indazole-3-carboxamide (AB-PINACA) and N-(1-amino-3-methyl-1-oxobutan-2-yl)-1-(4-fluorobenzyl)-1H-indazole-3-carboxamide (AB-FUBINACA) identified as designer drugs in illegal products. *Forensic Toxicol.* **31**(1), 93-100 (2013).
2. Sobolevsky, T., Prasolov, I., and Rodchenkov, G. Detection of JWH-018 metabolites in smoking mixture post-administration urine. *Forensic Sci. Int.* **200**(1-3), 141-147 (2010).
3. Chimalakonda, K.C., Moran, C.L., Kennedy, P.D., *et al.* Solid-phase extraction and quantitative measurement of omega and omega-1 metabolites of JWH-018 and JWH-073 in human urine. *Anal. Chem.* **83**(16), 6381-6388 (2011).

#### 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, 06/26/2023

#### 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](http://WWW.CAYMANCHEM.COM)