

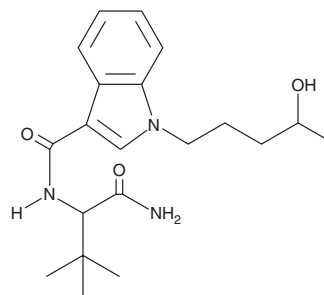
# PRODUCT INFORMATION



## ADBICA N-(4-hydroxypentyl) metabolite

Item No. 15240

**CAS Registry No.:** 2460433-26-3  
**Formal Name:** N-[1-(aminocarbonyl)-2,2-dimethylpropyl]-1-(4-hydroxypentyl)-1H-indole-3-carboxamide  
**Synonym:** ADB-PICA N-(4-hydroxypentyl) metabolite  
**MF:** C<sub>20</sub>H<sub>29</sub>N<sub>3</sub>O<sub>3</sub>  
**FW:** 359.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 218, 290 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

ADBICA (Item No. 14293) is a synthetic cannabinoid (CB) that has recently been identified in herbal blends.<sup>1</sup> Structurally, it is a modified aminoalkylindole like many cannabimimetics with pronounced activity at both CB receptors.<sup>2</sup> ADBICA N-(4-hydroxypentyl) metabolite is a potential phase 1 metabolite of ADBICA, based on the known metabolism of other aminoalkylindoles.<sup>3,4</sup> The physiological and toxicological properties of this compound have not been determined. This product is intended for forensic and research applications.

### References

1. Uchiyama, N., Matsuda, S., Kawamura, M., *et al.* Two new-type cannabimimetic quinolinyll carboxylates, QUPIC and QUCHIC, two new cannabimimetic carboxamide derivatives, ADB-FUBINACA and ADBICA, and five synthetic cannabinoids detected with a thiophene derivative α-PVT and an opioid receptor agonist AH-7921 identified in illegal products. *Forensic Toxicol.* **31(2)**, 223-240 (2013).
2. Aung, M.M., Griffin, G., Huffman, J.W., *et al.* Influence of the N-1 alkyl chain length of cannabimimetic indoles upon CB<sub>1</sub> and CB<sub>2</sub> receptor binding. *Drug Alcohol Depend.* **60(2)**, 133-140 (2000).
3. Wintermeyer, A., Möller, I., Thevis, M., *et al.* *In vitro* phase I metabolism of the synthetic cannabimimetic JWH-018. *Anal. Bioanal. Chem.* **398(5)**, 2141-53 (2010).
4. 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