

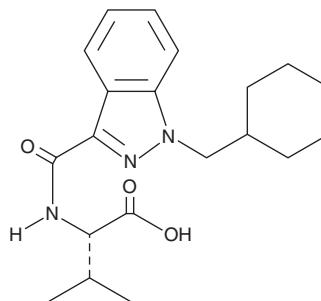
# PRODUCT INFORMATION



## AB-CHMINACA metabolite M2

Item No. 16389

**CAS Registry No.:** 1185887-51-7  
**Formal Name:** N-[[1-(cyclohexylmethyl)-1H-indazol-3-yl]carbonyl]-L-valine  
**MF:** C<sub>20</sub>H<sub>27</sub>N<sub>3</sub>O<sub>3</sub>  
**FW:** 357.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 210, 303 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-CHMINACA (Item No. 15434) is an indazole-based synthetic cannabinoid (CB) that is structurally related to AB-FUBINACA (Item No. 14039), a high affinity ligand of the central CB<sub>1</sub> receptor (K<sub>i</sub> = 0.9 nM).<sup>1,2</sup> AB-CHMINACA metabolite M2 is a metabolite of AB-CHMINACA that can be detected in the urine.<sup>3</sup> The physiological and toxicological properties of this compound are not known. This product is intended for forensic and research 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. Buchler, I.P., Hayes, M.J., Hedge, S.G., *et al.* Indazole derivatives. *Pfizer Inc.* **WO2009/106982A1** (2009).
3. Wurita, A., Hasegawa, K., Minakata, K., *et al.* Identification and quantification of metabolites of AB-CHMINACA in a urine specimen of an abuser. *Leg. Med. (Tokyo)* **19**, (2015).

#### 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, 06/30/2023

#### 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