

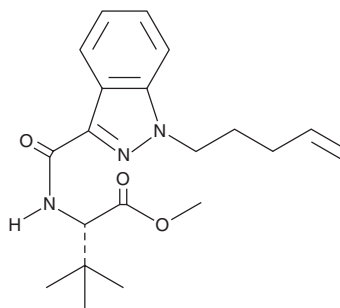
# PRODUCT INFORMATION



## MDMB-4en-PINACA

Item No. 26097

**CAS Registry No.:** 2504100-70-1  
**Formal Name:** 3-methyl-N-[[1-(4-penten-1-yl)-1H-indazol-3-yl]carbonyl]-L-valine, methyl ester  
**Synonyms:** MDMB-PENINACA,  
MDMB-PINACA N1-pentyl-4-en isomer  
**MF:** C<sub>20</sub>H<sub>27</sub>N<sub>3</sub>O<sub>3</sub>  
**FW:** 357.5  
**Purity:** ≥98%  
**Supplied as:** A neat solid  
**Storage:** -20°C  
**Stability:** ≥7 years



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

### Description

MDMB-4en-PINACA (Item No. 26097) is an analytical reference material categorized as a synthetic cannabinoid.<sup>1</sup> MDMB-4en-PINACA substitutes for Δ<sup>9</sup>-THC (Item Nos. ISO60157 | 9003771 | 12068) in a drug discrimination test and induces hypothermia in mice.<sup>2</sup> It has been associated with overdose and found in samples seized by law enforcement.<sup>3,4</sup> MDMB-4en-PINACA is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards for reference materials.

### References

1. Cannaert, A., Sparkes, E., Pike, E., *et al.* Synthesis and *in vitro* cannabinoid receptor 1 activity of recently detected synthetic cannabinoids 4F-MDMB-BICA, 5F-MPP-PICA, MMB-4en-PICA, CUMYL-CBMICA, ADB-BINACA, APP-BINACA, 4F-MDMB-BINACA, MDMB-4en-PINACA, A-CHMINACA, 5F-AB-P7AICA, 5F-MDMB-P7AICA, and 5F-AP7AICA. *ACS Chem. Neurosci.* **11**(24), 4434-4446 (2020).
2. Marusich, J.A., Gamage, T.F., Zhang, Y., *et al.* In vitro and in vivo pharmacology of nine novel synthetic cannabinoid receptor agonists. *Pharmacol. Biochem. Behav.* **220**, 173467 (2022).
3. Manini, A.F., Krotulski, A.J., Schimmel, J., *et al.* Respiratory failure in confirmed synthetic cannabinoid overdose. *Clin. Toxicol. (Phila)* **60**(4), 524-526 (2022).
4. Rodrigues, T.B., Souza, M.P., de Melo Barbosa, L., *et al.* Synthetic cannabinoid receptor agonists profile in infused papers seized in Brazilian prisons. *Forensic Toxicol.* **40**(1), 119-124 (2022).

#### 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, 04/03/2024

#### 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