

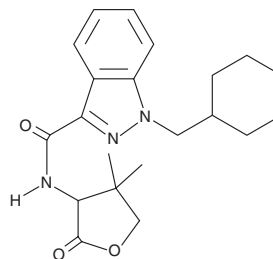
# PRODUCT INFORMATION



## MAB-CHMINACA metabolite M10

Item No. 17817

**CAS Registry No.:** 2748153-59-3  
**Formal Name:** 1-(cyclohexylmethyl)-N-(tetrahydro-4,4-dimethyl-2-oxo-3-furanyl)-1H-indazole-3-carboxamide  
**Synonym:** ADB-CHMINACA metabolite M10  
**MF:** C<sub>21</sub>H<sub>27</sub>N<sub>3</sub>O<sub>3</sub>  
**FW:** 369.5  
**Purity:** ≥98%  
**Supplied as:** A neat solid  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Description

MAB-CHMINACA (Item No. 16616) is an indazole-based synthetic cannabinoid (CB) structurally similar to AB-CHMINACA (Item No. 15434), both of which have been found in herbal blend products and are classified as Schedule I substances by the US Drug Enforcement Administration.<sup>1,2</sup> MAB-CHMINACA is a potent agonist of the CB<sub>1</sub> receptor with a reported K<sub>i</sub> value of 0.289 nM.<sup>3</sup> MAB-CHMINACA metabolite M10 is a potential metabolite of MAB-CHMINACA with two sites of metabolism (the terminal amide has been hydrolyzed to a carboxylic acid, and the isopropyl group has been hydroxylated). This compound readily cyclizes to the lactone. The physiological and toxicological properties of this compound are not known. This product is intended for forensic and research applications.

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

### References

1. Hasegawa, K., Wurita, A., Minakata, K., *et al.* Postmortem distribution of MAB-CHMINACA in body fluids and solid tissues of a human cadaver. *Forensic Toxicol.* **33**, 380-387 (2015).
2. Wurita, A., Hasegawa, K., Minakata, K., *et al.* Identification and quantitation of 5-fluoro-ADB-PINACA and MAB-CHMINACA in dubious herbal products. *Forensic Toxicol.* **33**, 213-220 (2015).
3. Buchler, I.P., Hayes, M.J., Hedge, S.G., *et al.* Indazole derivatives. WO 2009/106982 A1 (2009), 1-283, PCT/IB2009/000432.

#### 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, 08/04/2022

#### 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