

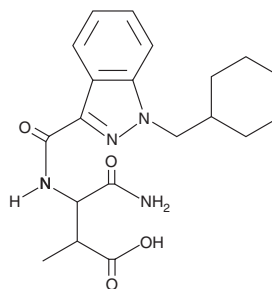
# PRODUCT INFORMATION



## AB-CHMINACA metabolite M6

Item No. 16453

**CAS Registry No.:** 2748161-80-8  
**Formal Name:** 4-amino-3-[[[1-(cyclohexylmethyl)-1H-indazol-3-yl]carbonyl]amino]-2-methyl-4-oxo-butanoic acid  
**MF:** C<sub>20</sub>H<sub>26</sub>N<sub>4</sub>O<sub>4</sub>  
**FW:** 386.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 210, 305 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 a cyclohexyl-substituted derivative of the indazole-based synthetic cannabinoid, AB-FUBINACA (Item No. 14039).<sup>1,2</sup> AB-CHMINACA metabolite M6 is a potential metabolite in which the carboxamide-linked aminocarbonyl-2-methylpropyl side chain associated with both AB-CHMINACA and AB-FUBINACA features an additional carboxyl group. 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-1*H*-indazole-3-carboxamide (AB-PINACA) and *N*-(1-amino-3-methyl-1-oxobutan-2-yl)-1-(4-fluorobenzyl)-1*H*-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).

#### 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