

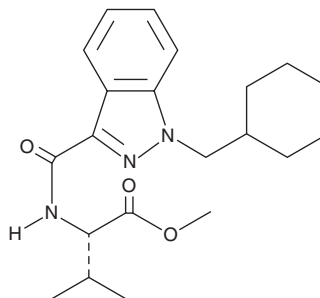
PRODUCT INFORMATION



MA-CHMINACA

Item No. 16421

CAS Registry No.: 1971007-96-1
Formal Name: N-[[1-(cyclohexylmethyl)-1H-indazol-3-yl] carbonyl]-L-valine, methyl ester
Synonyms: AMB-CHMINACA, AMB N-methylcyclohexyl analog, MAB-AB-CHMINACA, MMB-CHMINACA
MF: C₂₁H₂₉N₃O₃
FW: 371.5
Purity: ≥98%
UV/Vis.: λ_{max}: 210, 302 nm
Supplied as: A solution in acetonitrile
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 CB₁ receptor (K_i = 0.9 nM).^{1,2} AB-CHMINACA metabolite M2 is a potential derivative resulting from the replacement of a terminal amide NH₂ with a hydroxyl group. MA-CHMINACA is an analog of AB-CHMINACA metabolite M2 in which the hydroxyl moiety is replaced with a methyl ester group. 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. 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, 06/16/2023

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