

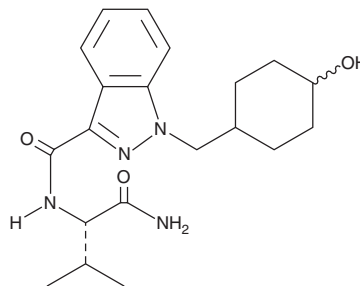
PRODUCT INFORMATION



AB-CHMINACA metabolite M1A

Item No. 16387

CAS Registry No.: 2130879-83-1
Formal Name: N-[(1S)-1-(aminocarbonyl)-2-methylpropyl]-1-[(4-hydroxycyclohexyl)methyl]-1H-indazole-3-carboxamide
MF: C₂₀H₂₈N₄O₃
FW: 372.5
Purity: ≥98% (mixture of diastereomers)
UV/Vis.: λ_{max}: 210, 303 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

AB-CHMINACA metabolite M1A is supplied as a crystalline solid. A stock solution may be made by dissolving the AB-CHMINACA metabolite M1A in the solvent of choice. AB-CHMINACA metabolite M1A is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of AB-CHMINACA metabolite M1A in ethanol and DMSO is approximately 30 mg/ml and approximately 50 mg/ml in DMF.

AB-CHMINACA metabolite M1A is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AB-CHMINACA metabolite M1A should first be dissolved in DMF and then diluted with the aqueous buffer of choice. AB-CHMINACA metabolite M1A has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

AB-CHMINACA (Item No. 15434) is a synthetic cannabinoid (CB) that is an analog of AB-FUBINACA (Item No. 14039), a potent agonist of the central CB₁ receptor (K_i = 0.9 nM).^{1,2} AB-CHMINACA metabolite M1A is a metabolite produced from AB-CHMINACA that can be detected in the urine.³ 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.
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)* (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, 08/03/2022

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