

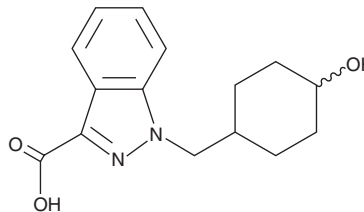
PRODUCT INFORMATION



AB-CHMINACA metabolite M5A

Item No. 16394

CAS Registry No.: 2207957-90-0
Formal Name: 1-((4-hydroxycyclohexyl)methyl)-1H-indazole-3-carboxylic acid
MF: C₁₅H₁₈N₂O₃
FW: 274.3
Purity: ≥98% (mixture of diastereomers)
UV/Vis.: λ_{max}: 210, 299 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 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 M5A is a potential metabolite produced during metabolism of AB-CHMINACA in the liver. 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. *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