

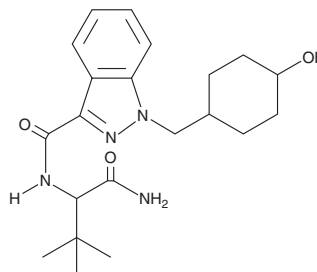
PRODUCT INFORMATION



MAB-CHMINACA metabolite M1

Item No. 43619

CAS Registry No.: 2460433-30-9
Formal Name: N-[1-(aminocarbonyl)-2,2-dimethylpropyl]-1-[(4-hydroxycyclohexyl)methyl]-1H-indazole-3-carboxamide
MF: $C_{21}H_{30}N_4O_3$
FW: 386.5
Purity: $\geq 98\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years



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

Description

MAB-CHMINACA metabolite M1 (Item No. 43619) is an analytical reference standard categorized as a synthetic cannabinoid metabolite.¹ MAB-CHMINACA metabolite M1 is a metabolite of MAB-CHMINACA (Item Nos. 22047 | 16616). This product is intended for research and forensic applications.

Reference

1. Gaunitz, F., Kieliba, T., Thevis, M., *et al.* Solid-phase extraction-liquid chromatography-tandem mass spectrometry method for the qualitative analysis of 61 synthetic cannabinoid metabolites in urine. *Drug Test. Anal.* **12**(1), 27-40 (2019).

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, 06/25/2025

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