

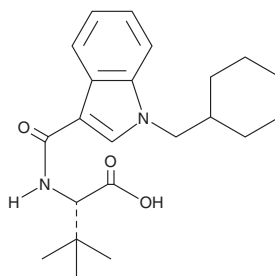
PRODUCT INFORMATION



MDMB-CHMICA metabolite M2

Item No. 19316

CAS Registry No.: 2460730-12-3
Formal Name: (S)-2-(1-(cyclohexylmethyl)-1H-indole-3-carboxamido)-3,3-dimethylbutanoic acid
MF: C₂₂H₃₀N₂O₃
FW: 370.5
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 292, 367 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Description

MDMB-CHMICA metabolite M2 (Item No. 19316) is an analytical reference standard categorized as a synthetic cannabinoid metabolite.¹ It is a metabolite of MDMB-CHMICA (Item No. 16965). This product is intended for research and forensic applications.

Reference

1. Gaunitz, F., Lehmann, S., Thomas, A., *et al.* Post-mortem distribution of the synthetic cannabinoid MDMB-CHMICA and its metabolites in a case of combined drug intoxication. *Int. J. Legal Med.* **132(6)**, 1645-1657 (2018).

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, 06/14/2023

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