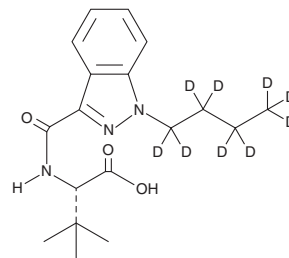


PRODUCT INFORMATION



MDMB-BUTINACA butanoic acid metabolite-d₉ Item No. 42707

Formal Name: N-[(1-butyl-d₉-1H-indazol-3-yl)carbonyl]-3-methyl-L-valine
Synonyms: ADB-BUTINACA 3,3-dimethylbutanoic acid metabolite-d₉,
MDMB-BINACA 3,3-dimethylbutanoic acid metabolite-d₉,
MDMB-BUTINACA 3,3-dimethylbutanoic acid metabolite-d₉
MF: C₁₈H₁₆D₉N₃O₃
FW: 340.5
Chemical Purity: ≥98% (MDMB-BUTINACA butanoic acid metabolite)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₉); ≤1% d₀
Supplied as: A 10 mg/ml solution in acetonitrile
Storage: -20°C
Stability: ≥1 year



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

Description

MDMB-BUTINACA butanoic acid metabolite-d₉ (Item No. 42707) is intended for use as an internal standard for the quantification of MDMB-BUTINACA butanoic acid metabolite (Item No. 28939) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

MDMB-BUTINACA butanoic acid metabolite is categorized as a synthetic cannabinoid metabolite.¹ It is a metabolite of ADB-BUTINACA (Item No. 29350). This product is intended for research and forensic applications.

Reference

1. Kronstrand, R., Norman, C., Vikingsson, S., et al. The metabolism of the synthetic cannabinoids ADB-BUTINACA and ADB-4en-PINACA and their detection in forensic toxicology casework and infused papers seized in prisons. *Drug Test. Anal.* **14**(4), 634-652 (2021).

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, 09/18/2025

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