

PRODUCT INFORMATION



MDMB-4en-PINACA butanoic acid metabolite-d₄ Item No. 37592

Formal Name: (S)-3,3-dimethyl-2-(1-(pent-4-en-1-yl)-1H-indazole-3-carboxamido-d₄)butanoic acid

Synonyms: MDMB-PENINACA butanoic acid metabolite-d₄, MDMB-4en-PINACA 3,3-Dimethylbutanoic acid-d₄, MDMB-PINACA N1-pentyl-4-en isomer butanoic acid metabolite-d₄

MF: C₁₉H₂₁D₄N₃O₃

FW: 347.5

Chemical Purity: ≥98% (MDMB-4en-PINACA butanoic acid metabolite)

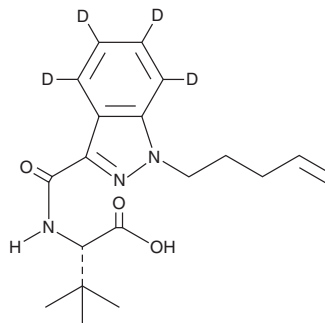
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₄); ≤1% d₀

UV/Vis.: λ_{max}: 301 nm

Supplied as: A solution in acetonitrile

Storage: -20°C

Stability: ≥4 years



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

Description

MDMB-4en-PINACA butanoic acid metabolite-d₄ (Item No. 37592) is intended for use as an internal standard for the quantification of MDMB-4en-PINACA butanoic acid metabolite (Item No. 28740) 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-4en-PINACA butanoic acid metabolite is categorized as a synthetic cannabinoid metabolite.¹ It is a metabolite of MDMB-4en-PINACA (Item Nos. 30761 | 26097). This product is intended for research and forensic applications.

Reference

1. Ozturk, Y.E. and Yeter, O. In vitro phase I metabolism of the recently emerged synthetic MDMB-4en-PINACA and its detection in human urine samples. *J. Anal. Toxicol.* **44**(9), 976-984 (2020).

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, 07/10/2023

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