

PRODUCT INFORMATION



MDMB-PICA-d₉

Item No. 39946

Formal Name: 3-methyl-N-[(1-pentyl-2,2,3,3,4,4,5,5,5-d₉-1H-indol-3-yl)carbonyl]-L-valine, methyl ester

MF: C₂₁H₂₁D₉N₂O₃

FW: 367.5

Chemical Purity: ≥98% (MDMB-PICA)

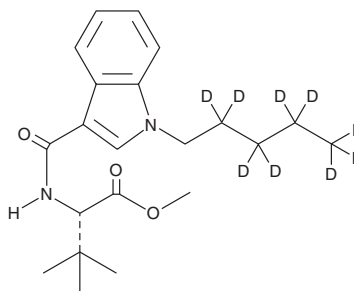
Deuterium

Incorporation: ≥99% deuterated forms (d₁-d₉); ≤1% d₀

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

MDMB-PICA-d₉ (Item No. 39946) is intended for use as an internal standard for the quantification of MDMB-PICA (Item No. 39729) 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-PICA is categorized as a synthetic cannabinoid.^{1,2} This product is intended for research and forensic applications.

References

1. Banister, S.D., Longworth, M., Kevin, R., *et al.* Pharmacology of valinate and tert-leucinate synthetic cannabinoids 5F-AMBICA, 5F-AMB, 5F-ADB, AMB-FUBINACA, MDMB-FUBINACA, MDMB-CHMICA, and their analogues. *ACS Chem. Neurosci.* **7**(9), 1241-1254 (2016).
2. Gioé-Gallo, C., Ortigueira, S., Brea, J., *et al.* Pharmacological insights emerging from the characterization of a large collection of synthetic cannabinoid receptor agonists designer drugs. *Biomed. Pharmacother.* **164**, 114934 (2023)

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, 07/12/2024

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