

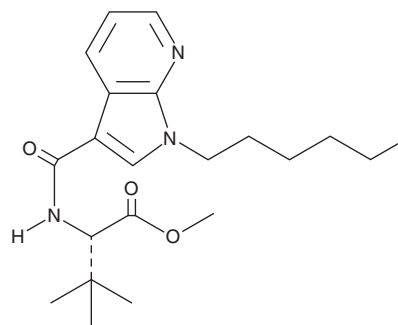
PRODUCT INFORMATION



5-fluoro MDMB-7-PAICA

Item No. 25286

CAS Registry No.: 2377403-81-9
Formal Name: N-[[1-(5-fluoropentyl)-1H-pyrrolo[2,3-b]pyridin-3-yl]carbonyl]-3-methyl-L-valine, methyl ester
Synonyms: 7'N-5-fluoro ADB, 5-fluoro MDMB-P7AICA
MF: C₂₀H₂₈FN₃O₃
FW: 377.5
Purity: ≥98%
UV/Vis.: λ_{max}: 244, 293 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

5-fluoro MDMB-7-PAICA (Item No. 25286) is an analytical reference standard categorized as a synthetic cannabinoid.¹ 5-fluoro MDMB-7-PAICA has been found as an impurity in 4'N 5F-ADB5 samples seized by law enforcement.² This product is intended for research and forensic applications.

References

1. Cannaert, A., Sparkes, E., Pike, E., *et al.* Synthesis and in vitro cannabinoid receptor 1 activity of recently detected synthetic cannabinoids 4F-MDMB-BICA, 5F-MPP-PICA, MMB-4en-PICA, CUMYL-CBMICA, ADB-BINACA, APP-BINACA, 4F-MDMB-BINACA, MDMB-4en-PINACA, A-CHMINACA, 5F-AB-P7AICA, 5F-MDMB-P7AICA, and 5F-AP7AICA. *ACS Chem. Neurosci.* **11(24)**, 4434-4446 (2020).
2. Nacionalni Forenzični Laboratorij. Analytical report - 5F-MDMB-P7AICA (C₂₀H₂₈FN₃O₃) methyl 2-[[1-(5-fluoropentyl)-1H-pyrrolo[2,3-b]pyridin-3-yl]formamido]-3,3-dimethylbutanoate. (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, 01/04/2024

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