

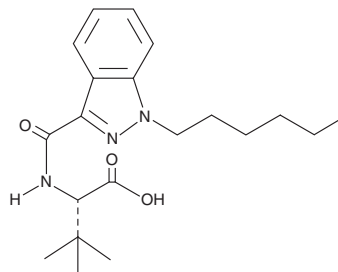
PRODUCT INFORMATION



5-fluoro ADB metabolite 7

Item No. 20178

CAS Registry No.: 2365471-39-0
Formal Name: N-[[1-(5-fluoropentyl)-1H-indazol-3-yl] carbonyl]-3-methyl-L-valine
Synonyms: 5F-ADB-M7,
5-fluoro MDMB-PINACA metabolite 7
MF: C₁₉H₂₆FN₃O₃
FW: 363.4
Purity: ≥98%
UV/Vis.: λ_{max}: 211, 302 nm
Supplied as: A solution in acetonitrile
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 ADB metabolite 7 (Item No. 20178) is an analytical reference standard categorized as a synthetic cannabinoid metabolite.¹ 5-fluoro ADB metabolite 7 is a metabolite of 5-fluoro ADB (Item Nos. 25012 | 23631 | 16603). It has been found in postmortem blood samples from individuals suspected of using 5-fluoro ADB. This product is intended for research and forensic applications.

Reference

1. Seither, J.Z., Reidy, L.J., and Boland, D.M. Identification and quantification of 5-fluoro ADB and the 5-fluoro ADB ester hydrolysis metabolite in postmortem blood samples by LC-MS/MS. *J. Anal. Toxicol.* 44(2), 133-139 (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, 06/21/2023

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