

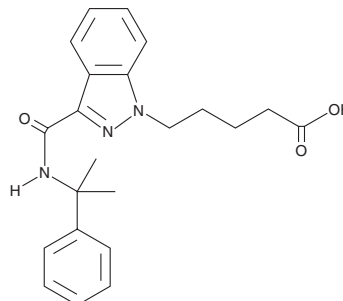
PRODUCT INFORMATION



5-fluoro CUMYL-PINACA N-pentanoic acid metabolite

Item No. 31286

Formal Name: 5-(3-((2-phenylpropan-2-yl)carbamoyl)-1H-indazol-1-yl)pentanoic acid
MF: C₂₂H₂₅N₃O₃
FW: 379.5
Purity: ≥98%
UV/Vis.: λ_{max}: 302 nm
Supplied as: A 10 mg/ml solution in acetonitrile
Storage: -20°C
Stability: ≥2 years



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

Description

5-fluoro CUMYL-PINACA N-pentanoic acid metabolite (Item No. 31286) is an analytical reference standard that is structurally similar to known synthetic cannabinoids. 5-fluoro CUMYL-PINACA N-pentanoic acid metabolite is a metabolite of 5-fluoro CUMYL-PINACA.¹ It is also a potential metabolite of CUMYL-PINACA based on the published metabolism of 5-fluoro CUMYL-PINACA.^{1,2} This product is intended for research and forensic applications.

References

1. Staeheli, S.N., Poetzsch, M., Veloso, V.P., *et al.* In vitro metabolism of the synthetic cannabinoids CUMYL-PINACA, 5F-CUMYL-PINACA, CUMYL-4CN-BINACA, 5F-CUMYL-P7AICA and CUMYL-4CN-B7AICA. *Drug Test Anal.* **10**(1), 148-157 (2018).
2. Franz, F., Jechle, H., Wilde, M., *et al.* Structure-metabolism relationships of valine and *tert*-leucine-derived synthetic cannabinoid receptor agonists: A systematic comparison of the in vitro phase I metabolism using pooled human liver microsomes and high-resolution mass spectrometry. *Forensic Toxicol.* **37**, 316-329 (2019).

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, 10/27/2025

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