

PRODUCT INFORMATION



CH-PIATA N-pentanoic acid metabolite

Item No. 37603

Formal Name: 5-(3-(2-(cyclohexylamino)-2-oxoethyl)-1H-indol-1-yl)pentanoic acid

Synonyms: CH-PIACA N-pentanoic acid metabolite, CHX-PIACA N-pentanoic acid metabolite, CHX-PIATA N-pentanoic acid metabolite, Cyclohexyl-PIATA N-pentanoic acid metabolite

MF: C₂₁H₂₈N₂O₃

FW: 356.5

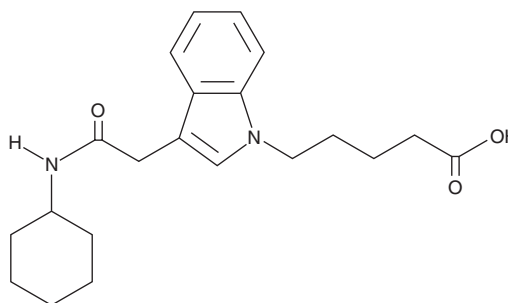
Purity: ≥97%

UV/Vis.: λ_{max}: 225 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥4 years



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

Description

CH-PIATA N-pentanoic acid metabolite (Item No. 37603) is an analytical reference standard that is structurally similar to known synthetic cannabinoids. CH-PIATA N-pentanoic acid metabolite is a potential metabolite of CH-PIATA (Item No. 36558) based on the published metabolism of 5-fluoro ABICA (Item No. 9001555).¹ At the time CH-PIATA N-pentanoic acid metabolite (Item No. 37603) was made available for purchase, specific metabolism data had not been published. Contact us if updated information on this molecule is now available. This product is intended for research and forensic applications.

Reference

1. 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, 04/11/2025

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