

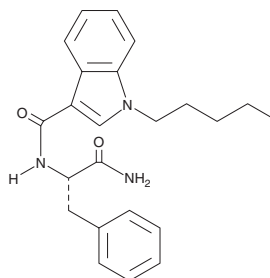
PRODUCT INFORMATION



APP-PICA

Item No. 17613

CAS Registry No.: 2365471-53-8
Formal Name: N-[(1S)-2-amino-2-oxo-1-(phenylmethyl)ethyl]-1-pentyl-1H-indole-3-carboxamide
MF: C₂₃H₂₇N₃O₂
FW: 377.5
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥2 years



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

Description

PX 1 (5-fluoro APP-PICA; Item No. 16201) is a synthetic cannabinoid (CB) that is structurally analogous to AM2201 (Item No. 10707), a potent CB receptor agonist.¹ APP-PICA is an analog of PX 1 that lacks the alkyl-terminal fluorine atom. The physiological and toxicological properties of this compound are not known. This product is intended for forensic and research applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards.

Reference

1. Makriyannis, A. and Deng, H. Cannabimimetic indole derivatives. WO 01/28557 A1 (2001), 1-25, PCT/US00/28832.

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, 08/04/2022

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