

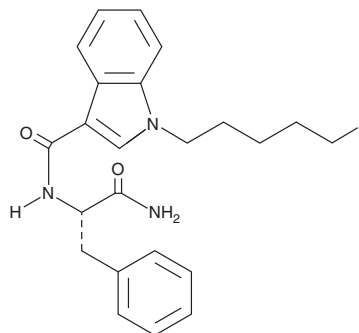
PRODUCT INFORMATION



PX 1

Item No. 16201

CAS Registry No.: 2221100-71-4
Formal Name: N-[(1S)-2-amino-2-oxo-1-(phenylmethyl)ethyl]-1-(5-fluoropentyl)-1H-indole-3-carboxamide
Synonyms: 5-fluoro APP-PICA, SRF-30
MF: C₂₃H₂₆FN₃O₂
FW: 395.5
Purity: ≥98%
UV/Vis.: λ_{max}: 216, 291 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

AM2201 (Item No. 10707) is a potent synthetic cannabinoid (CB) with K_i values of 1.0 and 2.6 nM for the central cannabinoid (CB₁) and peripheral cannabinoid (CB₂) receptors, respectively.¹ PX 1 is a derivative of AM2201 with a PINACA-like upper chain where the isopropyl group has been replaced with a benzyl group. The physiological and toxicological properties of this compound have not been determined. This product is intended for research and forensic applications.

Reference

1. Makriyannis, A. and Deng, H. Cannabimimetic indole derivatives. *University of Connecticut*, WO01/28557A1 (2001).

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, 06/29/2023

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