

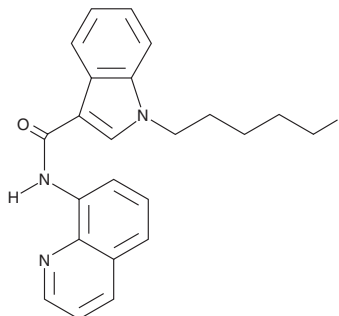
PRODUCT INFORMATION



AM2201 8-quinolinyl carboxamide

Item No. 14892

CAS Registry No.: 2365471-42-5
Formal Name: 1-(5-fluoropentyl)-N-8-quinolinyl-1H-indole-3-carboxamide
MF: C₂₃H₂₂FN₃O
FW: 375.4
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 244, 307, 330 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

Several cannabimimetic quinolinyl carboxylates (PB-22 (Item No. 14096) and BB-22 (Item No. 14099)) and cannabimimetic carboxamide derivatives (ADB-FUBINACA (Item No. 14292) and ADBICA (Item No. 14293)) have been identified as designer drugs in illegal products.¹ AM 2201 8-quinolinyl carboxamide is a derivative of these compounds where 8-quinolinol replaces the naphthalene group of AM2201 (Item No. 10707), resulting in a chemical structure similar to that of 5-fluoro PB-22 (Item No. 14095). The biological actions of AM 2201 8-quinolinyl carboximide are unknown. This product is intended for forensic and research applications.

Reference

1. Uchiyama, N., Matsuda, S., Kawamura, M., *et al.* Two new-type cannabimimetic quinolinyl carboxylates, QUPIC and QUCHIC, two new cannabimimetic carboxamide derivatives, ADB-FUBINACA and ADBICA, and five synthetic cannabinoids detected with a thiophene derivative α -PVT and an opioid receptor agonist AH-7921 identified in illegal products. *Forensic Toxicol.* **31(2)**, 223-240 (2013).

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, 06/22/2023

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