

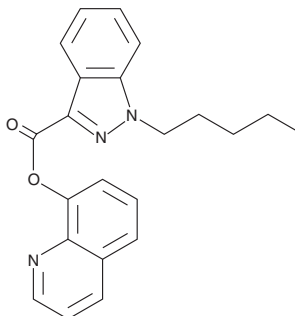
PRODUCT INFORMATION



NPB-22

Item No. 15535

CAS Registry No.: 1445579-61-2
Formal Name: 1-pentyl-1H-indazole-3-carboxylic acid, 8-quinolinyl ester
MF: C₂₂H₂₁N₃O₂
FW: 359.4
Purity: ≥98%
UV/Vis.: λ_{max}: 226, 302 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

The cannabimimetic quinolinyl carboxylate PB-22 (Item No. 14096) is an analog of JWH 018 (Item No. 10900) that has been identified as a designer drug in illegal products.¹ NPB-22 is an indazole-based derivative of PB-22 (Item No. 14096). The physiological and toxicological properties of this compound are not known. 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, 06/28/2023

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