

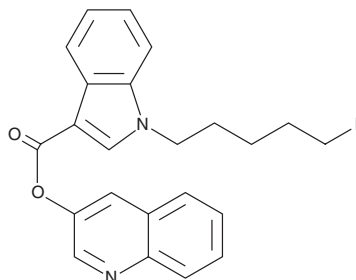
PRODUCT INFORMATION



5-fluoro PB-22 3-hydroxyquinoline isomer

Item No. 14511

CAS Registry No.: 2365471-26-5
Formal Name: 1-(5-fluoropentyl)-1H-indole-3-carboxylic acid, 3-quinolinyl ester
MF: C₂₃H₂₁FN₂O₂
FW: 376.4
Purity: ≥95%
UV/Vis.: λ_{max}: 207, 233, 294 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

5-fluoro PB-22 (Item No. 14095) is a synthetic cannabinoid that is an analog of potent 5-fluoropentyl JWH-type cannabimimetics. 5-fluoro PB-22 3-hydroxyquinoline isomer differs from 5-fluoro PB-22 by having the ester linkage inserted at the three, rather than at the eight, position of the quinoline moiety. The physiological and pharmacological properties of this isomer have not been investigated. This product is intended for forensic and research purposes.

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, 06/19/2023

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