

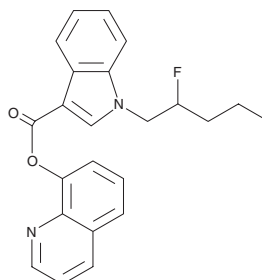
PRODUCT INFORMATION



5-fluoro PB-22 N-(2-fluoropentyl) isomer

Item No. 14546

CAS Registry No.: 2365471-10-7
Formal Name: 1-(2-fluoropentyl)-1H-indole-3-carboxylic acid, 8-quinolinyl ester
MF: $C_{23}H_{21}FN_2O_2$
FW: 376.4
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 215, 230, 293 nm
Supplied as: A 1 mg/ml solution in acetonitrile
Storage: -20°C
Stability: ≥ 4 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 derivative of PB-22 (Item No. 14096), a synthetic cannabinoid. 5-fluoro PB-22 N-(2-fluoropentyl) isomer differs from 5-fluoro PB-22 by having the fluorine atom at the 2 position, rather than the terminal 5 position, of the alkyl group. The physiological and toxicological properties of this compound have not been evaluated. 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, 12/08/2025

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