

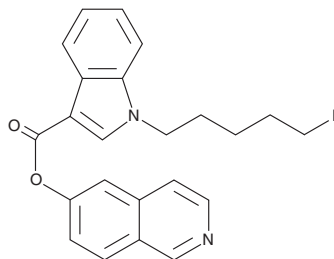
# PRODUCT INFORMATION



## 5-fluoro PB-22 6-hydroxyisoquinoline isomer

Item No. 14551

**CAS Registry No.:** 2365471-44-7  
**Formal Name:** 1-(5-fluoropentyl)-1H-indole-3-carboxylic acid, 6-isoquinolinyl ester  
**MF:** C<sub>23</sub>H<sub>21</sub>FN<sub>2</sub>O<sub>2</sub>  
**FW:** 376.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 219, 293 nm  
**Supplied as:** A 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 synthetic cannabinoid that is an analog of potent 5-fluoropentyl JWH-type cannabimimetics. 5-fluoro PB-22 6-hydroxyisoquinoline isomer differs from 5-fluoro PB-22 by having the quinoline group replaced with an isoquinoline group attached at its six position. 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, 06/20/2023

#### 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