

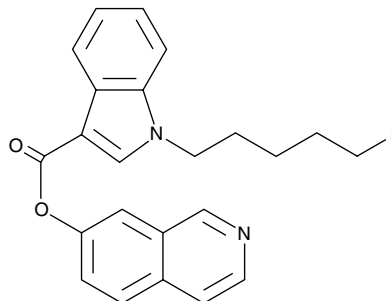
# Product Information



## 5-fluoro PB-22 7-hydroxyisoquinoline isomer

Item No. 14552

**Formal Name:** isoquinolin-7-yl 1-(5-fluoropentyl)-1H-indole-3-carboxylate  
**MF:** C<sub>23</sub>H<sub>21</sub>FN<sub>2</sub>O<sub>2</sub>  
**FW:** 376.4  
**Purity:** ≥98%  
**Stability:** ≥2 years at -20°C  
**Supplied as:** A crystalline solid  
**UV/Vis.:** λ<sub>max</sub>: 217, 294 nm



### Laboratory Procedures

For long term storage, we suggest that 5-fluoro PB-22 7-hydroxyisoquinoline isomer be stored as supplied at -20°C. It should be stable for at least two years.

5-fluoro PB-22 7-hydroxyisoquinoline isomer is supplied as a crystalline solid. A stock solution may be made by dissolving the 5-fluoro PB-22 7-hydroxyisoquinoline isomer in the solvent of choice. 5-fluoro PB-22 7-hydroxyisoquinoline isomer is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of 5-fluoro PB-22 7-hydroxyisoquinoline isomer in these solvents is approximately 10 and 11 mg/ml, respectively.

5-fluoro PB-22 7-hydroxyisoquinoline isomer is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 5-fluoro PB-22 7-hydroxyisoquinoline isomer should first be dissolved in DMF and then diluted with the aqueous buffer of choice. 5-fluoro PB-22 7-hydroxyisoquinoline isomer has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

5-fluoro PB-22 is a synthetic cannabinoid that is an analog of potent 5-fluoropentyl JWH-type cannabimimetics. 5-fluoro PB-22 7-hydroxyisoquinoline isomer differs from 5-fluoro PB-22 by having the quinoline group replaced with an isoquinoline group attached at its seven position. The physiological and pharmacological properties of this isomer have not been investigated. This product is intended for forensic and research purposes.

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/14552](http://www.caymanchem.com/catalog/14552)

### Cayman Chemical

**Mailing address**  
1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

**Phone**  
(800) 364-9897  
(734) 971-3335

**Fax**  
(734) 971-3640

**E-Mail**  
custserv@caymanchem.com

**Web**  
[www.caymanchem.com](http://www.caymanchem.com)

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 02/18/2014