

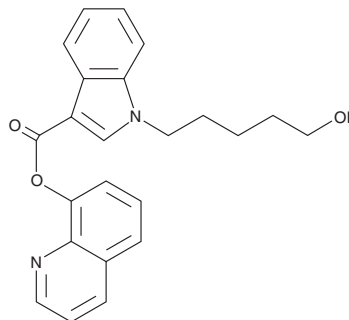
# PRODUCT INFORMATION



## PB-22 N-(5-hydroxypentyl) metabolite

Item No. 14387

**CAS Registry No.:** 1884343-30-9  
**Formal Name:** 1-(5-hydroxypentyl)-1H-indole-3-carboxylic acid, 8-quinolinyl ester  
**MF:** C<sub>23</sub>H<sub>22</sub>N<sub>2</sub>O<sub>3</sub>  
**FW:** 374.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 216, 230, 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

PB-22 N-(5-hydroxypentyl) metabolite is an expected metabolite of PB-22 (Item No. 14096), based on the metabolism of JWH 018 (Item No. 10900).<sup>1</sup> PB-22 is a synthetic cannabinoid that is structurally analogous to the potent aminoalkylindoles, like JWH 018. The physiological and toxicological properties of this compound have not been characterized. This product is intended for forensic and research purposes.

### Reference

1. Moran, C.L., Le, V.H., Chimalakonda, K.C., *et al.* Quantitative measurement of JWH-018 and JWH-073 metabolites excreted in human urine. *Anal. Chem.* **83**(11), 4228-4236 (2011).

#### 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