

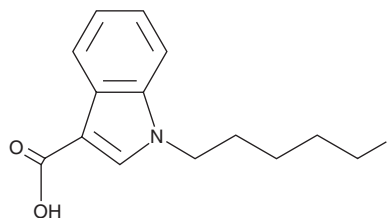
# PRODUCT INFORMATION



## 5-fluoro PB-22 3-carboxyindole metabolite

Item No. 14381

CAS Registry No.: 1432794-98-3  
Formal Name: 1-(5-fluoropentyl)-1H-indole-3-carboxylic acid  
MF:  $C_{14}H_{16}FNO_2$   
FW: 249.3  
Purity:  $\geq 98\%$   
Supplied as: A neat solid  
Storage:  $-20^{\circ}C$   
Stability:  $\geq 6$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Description

5-fluoro PB-22 3-carboxyindole metabolite (Item No. 14381) is an analytical reference material categorized as a synthetic cannabinoid metabolite.<sup>1</sup> It is a metabolite of 5-fluoro PB-22 (Item Nos. 16048 | 14095).<sup>2</sup> This product is intended for research and forensic applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards for reference materials.

### References

1. Davies, B.B., Bayard, C., Larson, S.J., *et al.* Retrospective analysis of synthetic cannabinoid metabolites in urine of individuals suspected of driving impaired. *J. Anal. Toxicol.* **40**(2), 89-96 (2016).
2. Wohlfarth, A., Gandhi, A.S., Pang, S., *et al.* Metabolism of synthetic cannabinoids PB-22 and its 5-fluoro analog, 5F-PB-22, by human hepatocyte incubation and high-resolution mass spectrometry. *Anal. Bioanal. Chem.* **406**(6), 1763-1780 (2014).

#### 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, 08/25/2022

#### 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