

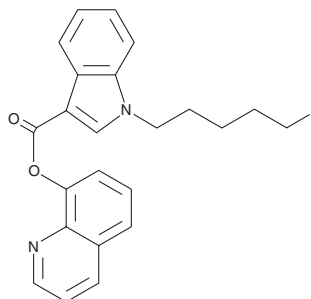
PRODUCT INFORMATION



5-fluoro PB-22 (exempt preparation)

Item No. 16048

CAS Registry No.: 1400742-41-7
Formal Name: 1-(5-fluoropentyl)-8-quinolinyl ester-1H-indole-3-carboxylic acid
Synonym: 5-fluoro QUPIC
MF: C₂₃H₂₁FN₂O₂
FW: 376.4
Purity: ≥98%
Supplied as: A 100 µg/ml solution in acetonitrile
Storage: -20°C
Stability: ≥3 years



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

Description

5-fluoro PB-22 (exempt preparation) (Item No. 16048) is an analytical reference standard categorized as a synthetic cannabinoid.^{1,2} 5-fluoro PB-22 is regulated as a Schedule I compound in the United States. 5-fluoro PB-22 (exempt preparation) (Item No. 16048) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

References

1. Banister, S.D., Stuart, J., Kevin, R.C., *et al.* Effects of bioisosteric fluorine in synthetic cannabinoid designer drugs JWH-018, AM-2201, UR-144, XLR-11, PB-22, 5F-PB-22, APICA, and STS-135. *ACS Chem. Neurosci.* **6(8)**, 1445-1458 (2015).
2. Costain, W.J., Tauskela, J.S., Rasquinha, I., *et al.* Pharmacological characterization of emerging synthetic cannabinoids in HEK293T cells and hippocampal neurons. *Eur. J. Pharmacol.* **786(2016)**, 234-235 (2016).

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, 08/31/2022

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