

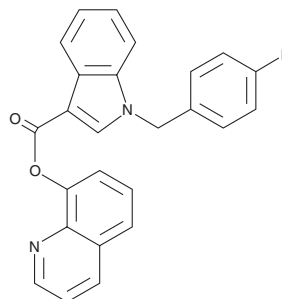
PRODUCT INFORMATION



FUB-PB-22

Item No. 14949

CAS Registry No.: 1800098-36-5
Formal Name: 1-[(4-fluorophenyl)methyl]-1H-indole-3-carboxylic acid, 8-quinoliny ester
MF: C₂₅H₁₇FN₂O₂
FW: 396.4
Purity: ≥98%
UV/Vis.: λ_{max}: 230, 293 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

FUB-PB-22 (Item No. 14949) is an analytical reference standard categorized as a synthetic cannabinoid.¹ FUB-PB-22 has been found in samples seized by law enforcement.² This product is intended for research and forensic applications

References

1. Heww, C., Schoeder, C.T., Pillaiyar, T., *et al.* Pharmacological evaluation of synthetic cannabinoids identified as constituents of spice. *Forensic Toxicol.* **34**, 329-343 (2016).
2. Dronova, M., Smolianitski, E., and Lev, O. Electrooxidation of new synthetic cannabinoids: Voltammetric determination of drugs in seized street samples and artificial saliva. *Anal. Chem.* **88(8)**, 4487-4494 (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, 06/26/2023

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