

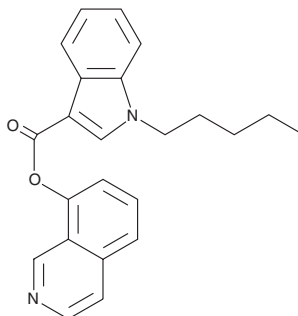
PRODUCT INFORMATION



PB-22 8-hydroxyisoquinoline isomer

Item No. 14558

CAS Registry No.: 1798021-82-5
Formal Name: 1-pentyl-1H-indole-3-carboxylic acid, 8-isoquinolinyl ester
MF: C₂₃H₂₂N₂O₂
FW: 358.4
Purity: ≥98%
UV/Vis.: λ_{max}: 216, 293 nm
Supplied as: A 10 mg/ml solution in acetonitrile
Storage: -20°C
Stability: ≥4 years



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

Description

PB-22 (Item No. 14096) is a synthetic cannabinoid that is structurally analogous to the potent aminoalkylindoles like JWH 018 (Item No. 10900). It differs from JWH 018 by having 8-hydroxyisoquinoline replacing the naphthalene group of JWH 018. PB-22 8-hydroxyisoquinoline isomer is a structural isomer of PB-22. The physiological and toxicological properties of this compound have not been characterized. This product is intended for forensic and research purposes.

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, 12/15/2025

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