PRODUCT INFORMATION



5-Fluoropentyl-3-pyridinoylindole (hydrochloride)

Item No. 14295

CAS Registry No.: Formal Name:	2365471-19-6 [1-(5-fluoropentyl)-1H-indol-3-yl]-3- pyridinyl-methanone, monohydrochloride		
MF:	$C_{19}H_{19}FN_2O \bullet HCI$		
FW:	346.8		
Purity:	≥98%		
UV/Vis.:	λ _{max} : 212, 260, 323 nm		• HCI
Supplied as:	A crystalline solid		1101
Storage:	-20°C	N.	
Stability:	≥5 years		
Information represents the product energiantions. Datch energia analytical results are previded on each partitionts of energy			

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

Description

5-Fluoropentyl-3-pyridinoylindole is a synthetic cannabinoid (CB) which shares the aminopentylindole structure of potent CBs like JWH 018 (Item No. 10900). This compound has a novel pyridinoyl group replacing the more common naphthoyl group found in WIN 55,212-2 and related JWH compounds. It also contains an alkyl-terminal fluoro group, which commonly enhances binding to both the central CB1 and peripheral CB₂ receptor.¹ The physiological and toxicological properties of this compound are not known, although it has been detected in herbal mixtures.² This product is intended for forensic and research applications.

References

- 1. Makriyannis, A. and Deng, H. Cannabimimetic indole derivatives. University of Connecticut. US7241799B2, (2001).
- 2. Uchiyama, N., Kawamura, M., Kikura-Hanajiri, R., et al. URB-754: A new class of designer drug and 12 synthetic cannabinoids detected in illegal products. Forensic Sci. Int. 227(1-3), 21-32 (2013).

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.

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, 06/16/2023

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