

PRODUCT INFORMATION



5-fluoro-3,5-AB-PFUPPYCA

Item No. 17181

CAS Registry No.: 2365471-73-2
Formal Name: N-[(1S)-1-(aminocarbonyl)-2-methylpropyl]-1-(5-fluoropentyl)-3-(4-fluorophenyl)-1H-pyrazole-5-carboxamide

Synonyms: 5-fluoro AB-FUPPYCA, 5-fluoro AB-FUPYCA, AZ-037

MF: $C_{20}H_{26}F_2N_4O_2$

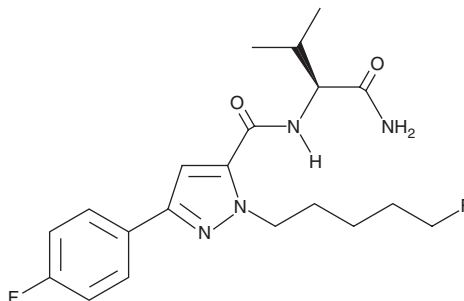
FW: 392.4

Purity: $\geq 98\%$

Supplied as: A neat solid

Storage: -20°C

Stability: ≥ 5 years



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

Description

5-fluoro-3,5-AB-PFUPPYCA (Item No. 17181) is an analytical reference material that is structurally categorized as a synthetic cannabinoid. The structure of highly substituted pyrazoles identified as cannabimimetic designer drugs has been reported.¹ The physiological and toxicological properties of this compound are not known. 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.

Reference

1. Girreser, U., Rösner, P., and Vasilev, A. Structure elucidation of the designer drug N-(1-amino-3,3-dimethyl-1-oxobutan-2-yl)-1-(5-fluoropentyl)-3-(4-fluorophenyl)-pyrazole-5-carboxamide and the relevance of predicted ^{13}C NMR shifts - a case study. *Drug Test Anal.* (2015).

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, 08/12/2024

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