

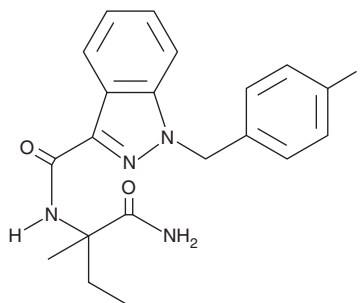
PRODUCT INFORMATION



AB-FUBINACA isomer 2

Item No. 9001529

CAS Registry No.: 2365471-64-1
Formal Name: N-[1-(aminocarbonyl)-1-methylpropyl]-1-[(4-fluorophenyl)methyl]-1H-indazole-3-carboxamide
MF: C₂₀H₂₁FN₄O₂
FW: 368.4
Purity: ≥98%
UV/Vis.: λ_{max}: 209, 301 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

AB-FUBINACA (Item No. 14039) is an adamantylindole-based synthetic cannabinoid (CB) with 10-fold greater affinity for the central CB₁ receptor (K_i = 0.9 nM) than that of JWH 018 (Item No. 10900).^{1,2} AB-FUBINACA was first synthesized by Pfizer as a potent CB₁ receptor modulator for potential therapeutic use, but recently was identified along with AB-PINACA (Item No. 14038) in illegal herbal products.^{1,2} AB-FUBINACA isomer 2 is a regioisomer of AB-FUBINACA. The physiological and toxicological properties of this compound have not been determined. This product is intended for research and forensic applications.

References

1. Uchiyama, N., Matsuda, S., Wakana, D., *et al.* New cannabimimetic indazole derivatives, N-(1-amino-3-methyl-1-oxobutan-2-yl)-1-pentyl-1H-indazole-3-carboxamide (AB-PINACA) and N-(1-amino-3-methyl-1-oxobutan-2-yl)-1-(4-fluorobenzyl)-1H-indazole-3-carboxamide (AB-FUBINACA) identified as designer drugs in illegal products. *Forensic Toxicol.* **31(1)**, 93-100 (2013).
2. Buchler, I.P., Hayes, M.J., Hedge, S.G., *et al.* Indazole derivatives. WO 2009/106982 A1 (2009), 1-283, PCT/IB2009/000432.

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, 07/12/2023

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