

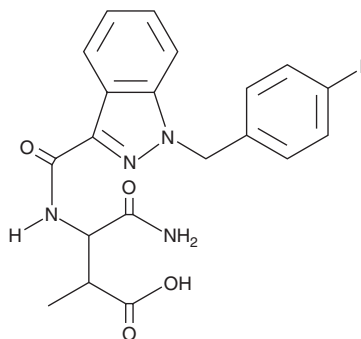
PRODUCT INFORMATION



AB-FUBINACA metabolite 2B

Item No. 17947

CAS Registry No.: 2460433-24-1
Formal Name: 4-amino-3-[[[1-[(4-fluorophenyl)methyl]-1H-indazol-3-yl]carbonyl]amino]-2-methyl-4-oxo-butanoic acid
MF: C₂₀H₁₉FN₄O₄
FW: 398.4
Purity: ≥95%
UV/Vis.: λ_{max}: 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 Nos. ISO60211 | ISO00134 | 14039) is an indazole-based synthetic cannabinoid (CB) that shows high affinity for the CB₁ receptor (K_i = 0.9 nM).^{1,2} AB-FUBINACA metabolite 2B is an expected carboxy metabolite of AB-FUBINACA, based on the major metabolites produced from similar synthetic CBs in humans.^{3,4} AB-FUBINACA metabolite 2B is a single diastereomer, but its relative stereochemistry is unknown. The other diastereomer, AB-FUBINACA metabolite 2A, is sold as Item No. 15529. The physiological and toxicological properties of this compound are not known. This product is intended for forensic and research applications.

References

1. Uchiyama, N., Matsuda, S., Wakana, D., *et al.* New cannabimimetic indazole derivatives, N-(1-amino-3methyl-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. *Pfizer Inc.* **WO2009/106982A1** (2009).
3. Sobolevsky, T., Prasolov, I., and Rodchenkov, G. Detection of JWH-018 metabolites in smoking mixture post-administration urine. *Forensic Sci. Int.* **200(1-3)**, 141-147 (2010).
4. Moran, C.L., Le, V.H., Chimalakonda, K.C., *et al.* Quantitative measurement of JWH-018 and JWH-073 metabolites excreted in human urine. *Anal. Chem.* **83(11)**, 4228-4236 (2011).

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, 07/06/2023

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