

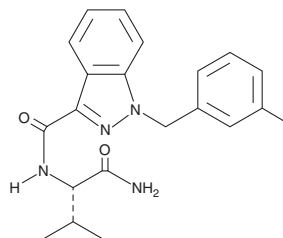
# PRODUCT INFORMATION



## AB-FUBINACA 3-fluorobenzyl isomer

Item No. 9001523

**CAS Registry No.:** 1185282-19-2  
**Formal Name:** N-[(1S)-1-(aminocarbonyl)-2-methylpropyl]-1-[(3-fluorophenyl)methyl]-1H-indazole-3-carboxamide  
**MF:** C<sub>20</sub>H<sub>21</sub>FN<sub>4</sub>O<sub>2</sub>  
**FW:** 368.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 209, 302 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 indazole-based synthetic cannabinoid (CB) that potently binds the central CB<sub>1</sub> receptor (K<sub>i</sub> = 0.9 nM).<sup>1-2</sup> This compound has been identified along with AB-PINACA (Item No. 14038) in illegal herbal products.<sup>1-2</sup> AB-FUBINACA 3-fluorobenzyl isomer is an isomer of AB-FUBINACA that differs by having the fluoro group at the three, rather than four, position on the benzyl group. This compound binds CB<sub>1</sub> with a K<sub>i</sub> value of 51.1 nM.<sup>2</sup> 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, 07/11/2023

### 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