

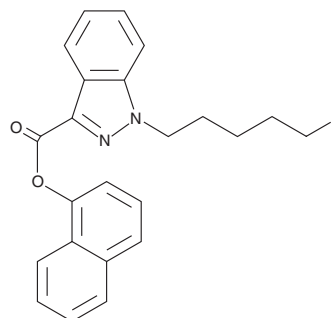
# PRODUCT INFORMATION



## 5-fluoro SDB-005

Item No. 15353

**CAS Registry No.:** 2185863-14-1  
**Formal Name:** 1-(5-fluoropentyl)-1H-indazole-3-carboxylic acid, 1-naphthalenyl ester  
**MF:** C<sub>23</sub>H<sub>21</sub>FN<sub>2</sub>O<sub>2</sub>  
**FW:** 376.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 221, 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

5-fluoro SDB-005 is an analog of SDB-005, a synthetic cannabinoid (CB) that activates the central CB<sub>1</sub> receptor (EC<sub>50</sub> = 21 nM) preferentially over the peripheral CB<sub>2</sub> receptor (EC<sub>50</sub> = 140 nM).<sup>1</sup> While the properties of this compound are not known, the addition of fluorine to the terminal carbon of the alkyl chain in similar cannabimimetic compounds typically increases affinity for both CB receptors significantly.<sup>2</sup> This product is intended for research and forensic applications.

### References

1. Banister, S.D., Wilkinson, S.M., Longworth, M., *et al.* The synthesis and pharmacological evaluation of adamantane-derived indoles: Cannabimimetic drugs of abuse. *ACS Chem. Neurosci.* **4**(7), 1081-1092 (2013).
2. Makriyannis, A. and Deng, H. Cannabimimetic indole derivatives. *University of Connecticut. WO01/28557A1*, (2001).

#### 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, 06/27/2023

#### 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