

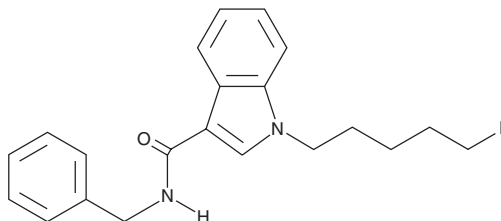
PRODUCT INFORMATION



5-fluoro SDB-006

Item No. 15390

CAS Registry No.: 1776086-02-2
Formal Name: 1-(5-fluoropentyl)-N-(phenylmethyl)-
1H-indole-3-carboxamide
MF: $C_{21}H_{23}FN_2O$
FW: 338.4
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 218, 291 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

SDB-006 (Item No. 15156) is an analog of the cannabimimetic indole JWH 018 adamantyl carboxamide (Item No. 9001193) in which the adamantane cage has been replaced with a phenyl ring.¹ It binds the central cannabinoid (CB_1) and peripheral cannabinoid (CB_2) receptors with EC_{50} values of 19 and 134 nM, respectively.¹ 5-fluoro SDB-006 is an analog of SDB-006 with a fluorine atom added to the terminal carbon of the alkyl chain. While the properties of this compound are not known, this modification of similar cannabimimetic compounds typically increases affinity for both CB receptors significantly.² This product is intended for forensic and research 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