

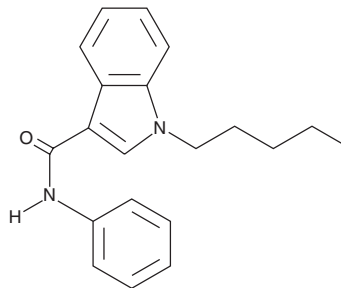
PRODUCT INFORMATION



SDB-006 N-phenyl analog

Item No. 16044

CAS Registry No.: 1430634-87-9
Formal Name: 1-pentyl-N-phenyl-1H-indole-3-carboxamide
Synonyms: phenyl-PICA, SDB-005
MF: C₂₀H₂₂N₂O
FW: 306.4
Purity: ≥98%
UV/Vis.: λ_{max}: 216, 259, 300 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 N-phenyl analog is a cannabimimetic indole that binds the central cannabinoid (CB₁) and peripheral cannabinoid (CB₂) receptors with EC₅₀ values of 21 and 140 nM, respectively.¹ It was discovered during research of the related compound JWH 018 adamantyl carboxamide (Item No. 9001193), which has been sold illicitly in herbal blends.¹ This product is intended for research and forensic applications.

Reference

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).

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, 06/29/2023

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