

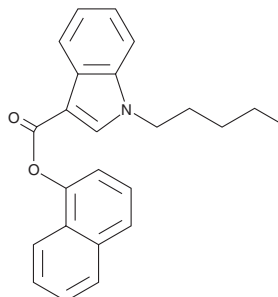
PRODUCT INFORMATION



CBL-018

Item No. 16443

CAS Registry No.: 2365471-54-9
Formal Name: 1-pentyl-1H-indole-3-carboxylic acid,
1-naphthalenyl ester
MF: C₂₄H₂₃NO₂
FW: 357.5
Purity: 357.5
UV/Vis.: λ_{max}: 220, 293 nm
Supplied as: A solution in acetonitrile
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

JWH 018 (Item No. 10900) is a mildly selective agonist of the peripheral cannabinoid (CB₂) receptor with K_i values of 9.0 and 2.94 nM for the central cannabinoid (CB₁) and CB₂ receptors, respectively.¹ CBL-018 is an analog of JWH 018 where the core indole base and the naphthalene group is linked by a carboxylate ester. The physiological and toxicological properties of this compound have not been tested. This product is intended for forensic and research applications.

Reference

1. Aung, M.M., Griffin, G., Huffman, J.W., *et al.* Influence of the N-1 alkyl chain length of cannabimimetic indoles upon CB₁ and CB₂ receptor binding. *Drug Alcohol Depend.* **60(2)**, 133-140 (2000).

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, 06/30/2023

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