

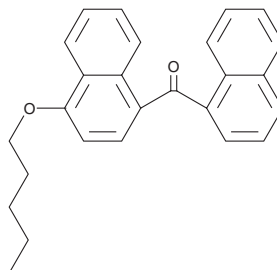
Product Information



CB-13

Item No. 10010398

CAS Registry No.: 432047-72-8
Formal Name: 1-naphthalenyl[4-(pentyllox)-1-naphthalenyl]-methanone
Synonym: CRA-13
MF: C₂₆H₂₄O₂
FW: 368.5
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that CB-13 be stored as supplied at -20°C. It should be stable for at least two years.

CB-13 is supplied as a crystalline solid. A stock solution may be made by dissolving the CB-13 in an organic solvent purged with an inert gas. CB-13 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of CB-13 in these solvents is approximately 0.2, 5, and 20 mg/ml, respectively.

If aqueous stock solutions are required for biological experiments, they can best be prepared by diluting the organic solvent into aqueous buffers or isotonic saline. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

CB-13 is a dual agonist of the, central cannabinoid (CB₁) (IC₅₀ = 15 nM) and peripheral cannabinoid (CB₂) (IC₅₀ = 98 nM) receptors.¹ In rats, oral CB-13 (3 mg/kg) potently blocks CB₁-dependent neuropathic mechanical hyperalgesia, shows limited brain penetration, and exhibits good oral bioavailability.¹ CB-13 does not possess genotoxic potential, as indicated by negative results in both chromosome aberration and reverse mutation assays.¹ The effects of CB-13 on other aspects of CB signaling have not been assessed.

Reference

1. Dziadulewicz, E.K., Bevan, S.J., Brain, C.T., *et al.* Naphthalen-1-yl-(4-pentylloxynaphthalen-1-yl)methanone: A potent, orally bioavailable human CB₁/CB₂ dual agonist with antihyperalgesic properties and restricted central nervous system penetration. *J. Med. Chem.* **50**, 3851-3856 (2007).

Related Products

JWH 018 - Cat. No. 13169 • JWH 073 - Cat. No. 13170 • JWH 200 - Cat. No. 13171 • (±)-CP 55,940 - Cat. No. 13241 • (±)-CP 47,497 - Cat. No. 16851 • 2-Arachidonoyl Glycerol - Cat. No. 62160 • 2-Arachidonoyl Glycerol ether - Cat. No. 62165 • Arachidonoyl Ethanolamide - Cat. No. 90050 • Arachidonoyl-2'-Fluoroethylamide - Cat. No. 90054 • HU-210 (DEA Schedule I Regulated Compound) - Cat. No. 90082 • (-)-CP 55,940 - Cat. No. 90084 • Arachidonoyl Cyclopropylamide - Cat. No. 91053 • Arachidonoyl 2'-Chloroethylamide - Cat. No. 91054 • JWH 015 - Cat. No. 90009018 • WIN 55212-2 (mesylate) - Cat. No. 1009023 • O-2545 - Cat. No. 10009195 • L-759,633 - Cat. No. 10009280

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog.**

Copyright Cayman Chemical Company, 02/21/2011