

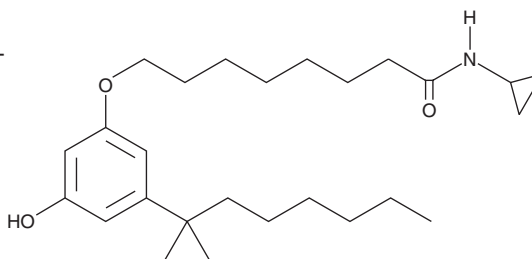
PRODUCT INFORMATION



CB-86

Item No. 13289

CAS Registry No.: 1150586-64-3
Formal Name: N-cyclopropyl-8-[3-(1,1-dimethylheptyl)-5-hydroxyphenoxy]-octanamide
MF: C₂₆H₄₃NO₃
FW: 417.6
Purity: ≥98%
UV/Vis.: λ_{max}: 206, 275 nm
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

CB-86 is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of CB-86 in these solvents is approximately 30 mg/ml.

CB-86 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of CB-86 should be diluted with the aqueous buffer of choice. CB-86 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CB-86 is a resorcinol-anandamide hybrid compound that acts as a partial agonist for the central cannabinoid (CB₁) receptor and a neutral antagonist for the peripheral cannabinoid (CB₂) receptor with K_i values of 5.6 and 7.9 nM, respectively.¹ Its demonstrated potency is similar to that of WIN 55,212-2 (K_i = 21 and 2.1 nM for CB₁ and CB₂, respectively).¹ CB-86 (1 mg/kg) exhibits antinociceptive effects in mice treated with formalin.¹

Reference

1. Brizzi, A., Brizzi, V., Cascio, M.G., *et al.* New resorcinol-anandamide “hybrids” as potent cannabinoid receptor ligands endowed with antinociceptive activity *in vivo*. *J. Med. Chem.* **52(8)**, 2506-2514 (2009).

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, 02/26/2024

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