

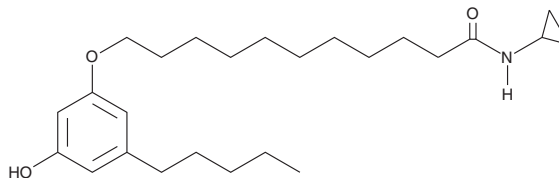
PRODUCT INFORMATION



CB-25

Item No. 10010116

CAS Registry No.: 869376-63-6
Formal Name: N-cyclopropyl-11-(3-hydroxy-5-pentylphenoxy)-undecanamide
MF: C₂₅H₄₁NO₃
FW: 403.6
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

CB-25 is supplied as a crystalline solid. A stock solution may be made by dissolving the CB-25 in the solvent of choice. CB-25 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of CB-25 in ethanol is approximately 50 mg/ml and approximately 30 mg/ml in DMSO and DMF.

CB-25 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CB-25 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. CB-25 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-25 is a stable analog of Δ^9 -tetrahydrocannabinol (THC) and anandamide (AEA). It exhibits high affinity for the central cannabinoid (CB₁) and peripheral cannabinoid (CB₂) receptors with K_i values of 5.2 and 13 nM, respectively. CB-25 behaves as an inverse agonist for the CB₁ receptor as assessed in a cyclic AMP (cAMP) functional assay.{14815,14816}

References

1. Brizzi, A., Brizzi, V., Cascio, M.G., *et al.* Design, synthesis, and binding studies of new potent ligands of cannabinoid receptors. *J. Med. Chem.* **48(23)**, 7343-7350 (2005).
2. Cascio, M.G., Bisogno, T., Palazzo, E., *et al.* *In vitro* and *in vivo* pharmacology of synthetic olivetol- or resorcinol-derived cannabinoid receptor ligands. *Br. J. Pharmacol.* **149(4)**, 431-440 (2006).

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, 01/25/2024

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