

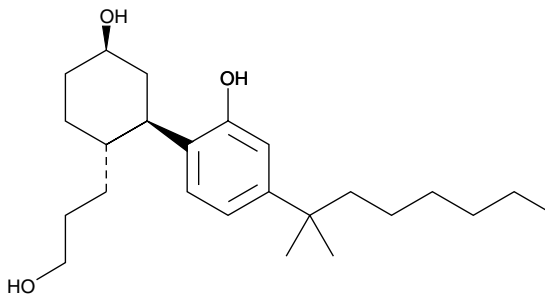
Product Information



(-)-CP 55,940

Item No. 90084

CAS Registry No.: 83002-04-4
Formal Name: 5-(1,1-dimethylheptyl)-2-[(1R,2R,5R)-5-hydroxy-2-(3-hydroxypropyl)cyclohexyl]-phenol
MF: C₂₄H₄₀O₃
FW: 376.6
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 204, 219, 277 nm



Laboratory Procedures

For long term storage, we suggest that (-)-CP 55,940 be stored as supplied at -20°C. It should be stable for at least two years.

(-)-CP 55,940 is supplied as a crystalline solid. A stock solution may be made by dissolving the (-)-CP 55,940 in an organic solvent purged with an inert gas. (-)-CP 55,940 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of (-)-CP 55,940 in these solvents is approximately 30 mg/ml.

(-)-CP 55,940 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, (-)-CP 55,940 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. (-)-CP 55,940 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

(-)-CP 55,940 is a potent, non-selective cannabinoid (CB) receptor agonist with K_i values of 0.58 and 0.69 nM for human recombinant CB₁ and CB₂, respectively.² It is a commonly used reference molecule in CB research and is considerably more potent than Δ⁹-THC in both behavioral tests and in receptor binding assays.¹

References

1. Pertwee, R.G. Pharmacology of cannabinoid receptor ligands. *Current Medicinal Chemistry* **6**, 635-664 (1999).
2. Wiley, J.L., Barrett, R.L., Lowe, J., *et al.* Discriminative stimulus effects of (-)-CP 55,940 and structurally dissimilar cannabinoids in rats. *Neuropharmacology* **34(6)**, 669-676 (1995).

Related Products

(+)-CP 47,497 - Item No. 10910 • (±)-CP 47,497 - Item No. 10913 • (-)-CP 47, 497 - Item No. 10917 • (±)3-*epi* CP 47,497-C8-homolog - Item No. 10918 • (±)-*epi* CP 47, 497 - Item No. 10919 • 2-Arachidonoyl Glycerol - Item No. 62160 • Arachidonoyl Ethanolamide - Item No. 90050 • Cannabidiol - Item No. 90080 • HU-210 - Item No. 90082 • HU-210 (solution) - Item No. 90083 • HU-331 - Item No. 10005673 • JWH 015 - Item No. 10009018

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, 06/20/2011