

Product Information



(+)-CP 55,940

Item No. 13608

Formal Name: 2-((1S,2S,5S)-5-hydroxy-2-(3-hydroxypropyl)cyclohexyl)-5-(2-methyloctan-2-yl)phenol

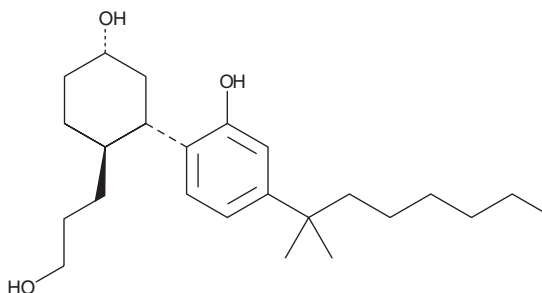
MF: C₂₄H₄₀O₃

FW: 376.6

Purity: ≥98%

Stability: ≥2 years at -20°C

Supplied as: A crystalline solid



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 the solvent of choice. (+)-CP 55,940 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. 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 bicyclic mimetic of Δ^9 -THC that binds both cannabinoid receptors and has potent analgesic properties.¹⁻³ The isomer (-)-CP 55,940, which structurally resembles the A and C ring of Δ^9 -THC, is a potent analgesic.² (+)-CP 55,940 is an enantiomer purified from the (±)-CP 55,940 racemic mixture. The functional characteristics of this isomer have not been studied. However, there is no significant difference between the racemate and (-) isomer in analgesic activity in four of five tests, suggesting that (+)-CP 55,940 is biologically active.²

References

1. Devane, W.A., Dysarz, F.A., III, Johnson, M.R., *et al.* Determination and characterization of a cannabinoid receptor in rat brain. *Mol. Pharmacol.* **34**, 605-613 (1988).
2. Howlett, A.C., Johnson, M.R., Melvin, L.S., *et al.* Nonclassical cannabinoid analgetics inhibit adenylate cyclase: Development of a cannabinoid receptor model. *Mol. Pharmacol.* **33**, 297-302 (1987).
3. Pertwee, R.G. Pharmacology of cannabinoid receptor ligands. *Curr. Med. Chem.* **6**, 635-664 (1999).

Related Products

(±)3-*epi* CP 47,497-C8-homolog - Item No. 10918 • (±)-*epi* CP 47, 497 - Item No. 10919 • JWH 018 - Item No. 13169 • JWH 073 - Item No. 13170 • JWH 200 - Item No. 13171 • (±)-CP 47,497-C8-homolog - Item No. 13216 • (-)-CP 47,497 - Item No. 13218 • (+)-CP 47,497 - Item No. 13219 • (±)-CP 55,940 - Item No. 13241 • JWH 019 - Item No. 13633 • JWH 250 - Item No. 13634 • (±)-CP 47,497 - Item No. 16851 • HU-210 (DEA Schedule I Regulated Compound) - Item No. 90082 • (-)-CP 55,940 - Item No. 90084 • HU-308 - Item No. 90086 • HU-331 - Item No. 10005673 • JWH 015 - Item No. 10009018 • WIN 55212-2 (mesylate) - Item No. 10009023

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, 05/06/2011