

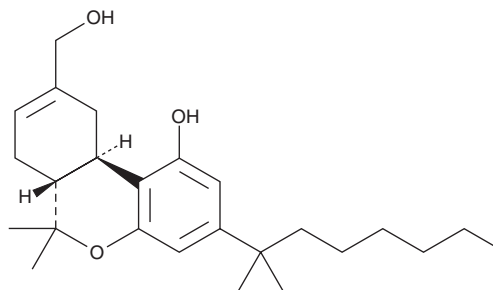
Product Information



HU-210

Item No. 90082 • Lot No. XXXXXX • Size: XX mg

CAS Registry No.: 112830-95-2
Formal Name: 3-(1,1'-dimethylheptyl)-6aR,7,10,10aR-tetrahydro-1-hydroxy-6,6-dimethyl-6H-dibenzo[b,d]pyran-9-methanol
MF: C₂₅H₃₈O₃
FW: 386.6
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
Packaged by: Cayman Chemical Company



Laboratory Procedures

For long term storage, we suggest that HU-210 be stored as supplied at -20°C. It will be stable for at least two years.

HU-210 is supplied as a crystalline solid. A stock solution may be made by dissolving the HU-210 in an organic solvent. HU-210 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of HU-210 in these solvents is at least 50 mg/ml. HU-210 will be stable for at least six months in these solvents if stored at -20°C.

HU-210 is practically insoluble in aqueous solutions (<20 µg/ml), therefore we suggest that further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. 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.

HU-210 is a synthetic agonist analog of Δ⁹-THC, which is the primary psychoactive component of marijuana. It is a potent CB₁ and CB₂ receptor agonist. HU-210 binds to neuroblastoma cell membrane CB₁ receptors with about the same affinity as CP-55940.¹ In murine behavior models (hypothermia, analgesia, hypoactivity, catalepsy) the ED₅₀ value is 5-20 µg/kg.²

References

- Slipetz, D.M., O'Neill, G.P., Gavreau, L., *et al.* Activation of the human peripheral cannabinoid receptor results in inhibition of adenylyl cyclase. *Mol. Pharmacol.* **48**, 352-361 (1995).
- Howlett, A.C., Berglund, B., and Melvin, L.S. Cannabinoid receptor agonists and antagonists. *Current Pharm. Design* **1**, 343-354 (1995).

Related Products

(±)-CP 47,497-C8-homolog-d7 (solution) - Item No. 10686 • (+)-CP 47,497 - Item No. 10910 • (±)-CP 47,497 - Item No. 10913 • (-)-CP 47,497 - Item No. 10917 • Cannabidiol - Item No. 90080 • Cannabidiol (solution) - Item No. 90081 • HU-210 (solution) - Item No. 90083 • CP-55,940 - Item No. 90084 • CB₁ Receptor Polyclonal Antibody - Item No. 101500 • CB₂ Receptor (human) Polyclonal Antiserum - Item No. 101550

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