

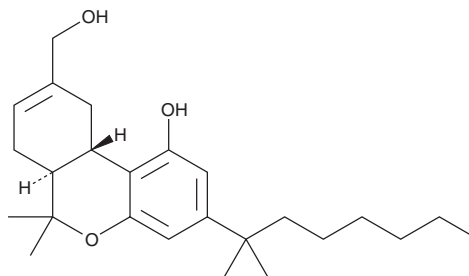
# Product Information



## HU-211

Item No. 10006350

**CAS Registry No.:** 112924-45-5  
**Formal Name:** 3-(1,1-dimethylheptyl)-6aS,7,10,10aS-tetrahydro-1-hydroxy-6,6-dimethyl-6H-dibenzo[b,d]pyran-9-methanol  
**Synonym:** Dexanabinol  
**MF:** C<sub>25</sub>H<sub>38</sub>O<sub>3</sub>  
**FW:** 386.6  
**Purity:** ≥98%  
**Stability:** ≥2 years at -20°C  
**Supplied as:** A crystalline solid



### Laboratory Procedures

For long term storage, we suggest that HU-211 be stored as supplied at -20°C. It should be stable for at least two years.

HU-211 is supplied as a crystalline solid. A stock solution may be made by dissolving the HU-211 in an organic solvent purged with an inert gas. HU-211 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of HU-211 in these solvents is approximately 30 mg/ml.

HU-211 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, HU-211 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. HU-211 has a solubility of approximately 0.25 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

The endocannabinoids present a rich system of central cannabinoid (CB<sub>1</sub>), peripheral cannabinoid (CB<sub>2</sub>), and non-CB receptor-mediated pharmacology that has stimulated research in many fields including memory, weight loss and appetite, neurodegeneration, tumor surveillance, analgesia, and inflammation.<sup>1,2</sup> HU-211 is a synthetic terpene-based cannabinoid devoid of CB<sub>1</sub> and CB<sub>2</sub> agonist activity, thus lacking the psychomotor responses characteristic of Δ<sup>9</sup>-THC. HU-211 does exhibit the neuroprotective, antioxidant properties of other related compounds like cannabidiol.<sup>3</sup> HU-211 also has anti-inflammatory properties derived through inhibition of NF-κB and the resulting decreases in cytokines such as TNFα and interleukin-6.

### References

1. Pertwee, R.G. Pharmacology of cannabinoid receptor ligands. *Current Medicinal Chemistry* **6**, 635-664 (1999).
2. Martin, B.R., Mechoulam, R., and Razdan, R.K. Discovery and characterization of endogenous cannabinoids. *Life Sci* **65**, 573-595 (1999).
3. Juttler, E., Potrovita, I., Tarabin, V., *et al.* The cannabinoid dexanabinol is inhibitor of the nuclear factor-kappa B (NF-κB). *Neuropharmacology* **47**, 580-592 (2004).

### Related Products

Cannabidiol (DEA Schedule I Regulated Compound) - Item No. 90080 • Cannabidiol (solution) - Item No. 90081 • HU-210 (DEA Schedule I Regulated Compound) - Item No. 90082 • HU-308 - Item No. 90086 • CAY10429 - Item No. 10004259

### 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, 03/07/2011