

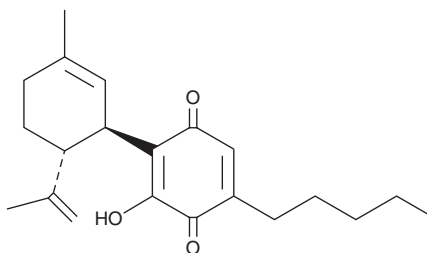
Product Information



HU-331

Item No. 10005673

CAS Registry No.: 137252-25-6
Formal Name: 3-hydroxy-2-[(1R,6R)-3-methyl-6-(1-methylethenyl)-2-cyclohexen-1-yl]-5-pentyl-2,5-cyclohexadiene-1,4-dione
MF: C₂₁H₂₈O₃
FW: 328.5
Purity: ≥95%
Stability: ≥1 year at -20°C
Supplied as: A solution in methyl acetate



Laboratory Procedures

For long term storage, we suggest that HU-331 be stored as supplied at -20°C. It should be stable for at least one year. HU-331 is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of HU-331 in these solvents is at least 20 mg/ml.

HU-331 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of HU-331 should be diluted with the aqueous buffer of choice. HU-331 has a solubility of 0.5 mg/ml in a 1:1 solution of ethanol: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 (CB₁), peripheral cannabinoid (CB₂), 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.^{1,2} HU-331 is a hydroxylquinone cannabidiol analog that exhibits potent antineoplastic activity on a variety of human cancer cell lines.³ It effectively inhibits the growth of human Raji and Jurkat lymphoma cells *in vitro* with an EC₅₀ of approximately 0.2 µg/ml and an EC₈₀ of 1.56 µg/ml. HU-331 also inhibits the growth of HT-29 colon carcinoma cells, which have been inoculated into nude mice, by more than 50% at a dose of 5 mg/kg.³

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. Kogan, N.M., Rabinowitz, R., Levi, P., *et al.* Synthesis and antitumor activity of quinonoid derivatives of cannabinoids. *J. Med. Chem.* **47**, 3800-3806 (2004).

Related Products

(+)-CP 47,497 - Item No. 10910 • (±)-CP 47,497 - Item No. 10913 • (-)-CP 47,497 - Item No. 10917 • Cannabidiol - Item No. 90080 • HU-210 - Item No. 90082 • HU-308 - Item No. 90086 • HU-211 - Item No. 10006350

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, 05/05/2011