

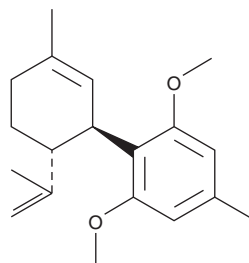
# Product Information



## O-1918

Item No. 10004914

**CAS Registry No.:** 536697-79-7  
**Formal Name:** 1,3-dimethoxy-5-methyl-2-[(1R,6R)-3-methyl-6-(1-methylethenyl)-2-cyclohexen-1-yl]-benzene  
**MF:** C<sub>19</sub>H<sub>26</sub>O<sub>2</sub>  
**FW:** 286.4  
**Purity:** ≥98%  
**Stability:** ≥1 year at -20°C  
**Supplied as:** A solution in methyl acetate



### Laboratory Procedures

For long term storage, we suggest that O-1918 be stored as supplied at -20°C. It should be stable for at least one year.

O-1918 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 O-1918 in these solvents is approximately 30 mg/ml.

O-1918 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of O-1918 should be diluted with the aqueous buffer of choice. O-1918 has a solubility of approximately 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.

Abnormal cannabidiol is a synthetic regioisomer of cannabidiol that fails to elicit either central cannabinoid (CB<sub>1</sub>) or peripheral cannabinoid (CB<sub>2</sub>) responsiveness and is without psychotropic activity. It induces endothelium-dependent vasodilation *via* a CB<sub>1</sub>/CB<sub>2</sub>/nitric oxide-independent mechanism.<sup>1</sup> O-1918 is a cannabidiol analog that acts as a selective antagonist of abnormal cannabidiol at the non-CB<sub>1</sub>/CB<sub>2</sub> endothelial receptor. It does not bind to CB<sub>1</sub> or CB<sub>2</sub> receptors at concentrations up to 30 μM and inhibits the vasorelaxant effects of abnormal cannabidiol *in vitro* and in whole animals.<sup>2</sup> It also blocks the abnormal cannabidiol-induced activation of the phosphatidylinositol 3-kinase/Akt pathway in human umbilical vein endothelial cells.<sup>2</sup>

### References

1. Járai, Z., Wagner, J.A., Varga, K., *et al.* Cannabinoid-induced mesenteric vasodilation through an endothelial site distinct from CB<sub>1</sub> or CB<sub>2</sub> receptors. *Proc. Natl. Acad. Sci. USA* **96(24)**, 14136-14141 (1999).
2. Offertáler, L., Mo, F.-M., Bátka, S., *et al.* Selective ligands and cellular effectors of a G protein-coupled endothelial cannabinoid receptor. *Mol. Pharmacol.* **63(3)**, 699-705 (2003).

### Related Products

Cannabidiol (DEA Schedule I Regulated Compound) - Item No. 90080 • Cannabidiol (solution) - Item No. 90081 • CAY10429 - Item No. 10004259 • N-Arachidonoyl L-Serine - Item No. 10005455 • O-1602 - Item No. 10006803 • O-1821 - Item No. 10006804

### 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