# **Product Information**



## **Oxy-16**

Item No. 9000395

CAS Registry No.: 596-94-1

Formal Name: cholest-5-ene-3\beta,20,22R-triol Synonym: 20(R),22(R)-dihydroxy Cholesterol

MF:  $C_{27}H_{46}O_3$ 418.7 FW: **Purity:** ≥95%

Stability: ≥2 years at -20°C A crystalline solid Supplied as:

### **Laboratory Procedures**

For long term storage, we suggest that oxy-16 be stored as supplied at -20°C. It should be stable for at least two years. Oxy-16 is supplied as a crystalline solid. A stock solution may be made by dissolving the oxy-16 in the solvent of choice. Oxy-16 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of oxy-16 in ethanol is approximately 10 mg/ml and approximately 1.5 mg/ml in DMSO

If aqueous stock solutions are required for biological experiments, they can best be prepared by dissolving the compound in aqueous buffers or isotonic saline. 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.

Naturally occurring oxysterols are products of cholesterol oxidation that can stimulate the hedgehog (Hh) signaling pathway in target cells associated with cardiovascular disease and with bone formation. Depending on the target cell, activation of Hh signaling can modulate inflammatory responses to additional atherogenic factors such as lesion-producing macrophages or enable osteoblast differentiation.<sup>2,3</sup> Oxy-16 is a synthetic oxysterol compound that may function as an antagonist of hedgehog activity. Published data regarding its efficacy remains forthcoming.

### References

- 1. Dwyer, J.R., Sever, N., Carlson, M., et al. Oxysterols are novel activators of the hedgehog signaling pathway in pluripotent mesenchymal cells. J. Biol. Chem. 282(12), 8959-8968 (2007).
- Björkhem, I. and Diczfalusy, U. Oxysterols: Friends, foes, or just fellow passengers? Arterioscler. Thromb. Vasc. Biol. 22, 734-742 (2002).
- 3. Amantea, C.M., Kim, W.-K., Meliton, V., et al. Oxysterol-induced osteogenic differentiation of marrow stromal cells is regulated by Dkk-1 inhibitable and PI3-kinase mediated signaling. J. Cell. Biochem. 105(2), 424-436 (2008).

#### **Related Products**

For a list of related products please visit: www.caymanchem.com/catalog/9000395

WARNING: This product is for laboratory research only: not for administration to humans. Not for human or veterinary DIAGNOSTIC OR THERAPEUTIC USE.

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

ayman 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 ose, 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

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 y 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, 06/15/2012

### Cayman Chemical

### **Mailing address**

1180 E. Ellsworth Road Ann Arbor, MI 48108 USA

### **Phone**

(800) 364-9897 (734) 971-3335

(734) 971-3640

custserv@caymanchem.com

www.caymanchem.com