

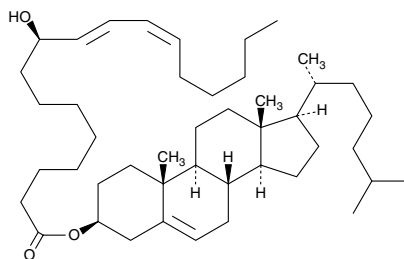
# Product Information



## 9(R)-HODE cholesteryl ester

Item No. 38406

**Formal Name:** 9R-hydroxy-10E,12Z-octadecadienoic acid, cholesteryl ester  
**MF:** C<sub>45</sub>H<sub>76</sub>O<sub>3</sub>  
**FW:** 665.1  
**Purity:** ≥98%  
**Stability:** ≥1 year at -20°C  
**Supplied as:** A solution in ethanol  
**UV/Vis.:** λ<sub>max</sub>: 234 nm ε: 23,000



### Laboratory Procedures

For long term storage, we suggest that 9(R)-HODE cholesteryl ester be stored as supplied at -20°C. It should be stable for at least one year.

9(R)-HODE cholesteryl ester is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of 9(R)-HODE cholesteryl ester in these solvents is approximately 50 mg/ml.

9(R)-HODE is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 9(R)-HODE should be diluted with the aqueous buffer of choice. 9(R)-HODE has a solubility of approximately 10 mg/ml in a 1:10 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

9(R)-HODE cholesteryl ester was originally extracted from atherosclerotic lesions.<sup>1</sup> It remains uncertain whether the oxidized fatty acid portion of the molecule results from enzymatic lipoxygenation or from random lipid peroxidation.<sup>2</sup> 9(R)-HODE cholesteryl ester can be used as a standard for analysis of chiral HODE cholesteryl esters.

### References

1. Brooks, C.J.W., Harland, W.A., Steel, G., et al. Lipids of human atheroma: Isolation of hydroxyoctadecadienoic acids from advanced aortal lesions. *Biochim. Biophys. Acta* 202, 563-566 (1970).
2. Belkner, J., Wiesner, R., Kühn, H., et al. The oxygenation of cholesterol esters by the reticulocyte lipoxygenase. *FEBS Lett.* 279, 110-114 (1991).

### Related Products

(±)9-HODE cholesteryl ester - Item No. 38401 • 9(R)-HODE - Item No. 38405 • 9(S)-HODE cholesteryl ester - Item No. 38411

**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, 04/07/2011

### 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