

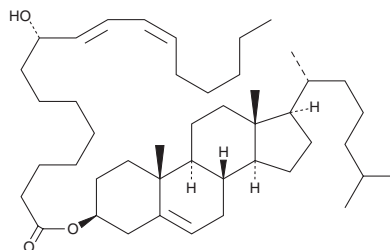
# PRODUCT INFORMATION



## 9(S)-HODE cholesteryl ester

Item No. 38411

**CAS Registry No.:** 143442-54-0  
**Formal Name:** 9S-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%  
**UV/Vis.:** λ<sub>max</sub>: 234 nm ε: 23,000  
**Supplied as:** A solution in ethanol  
**Storage:** -20°C  
**Stability:** ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

9(S)-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 ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of 9(S)-HODE cholesteryl ester in these solvents is approximately 50 mg/ml.

9(S)-HODE cholesteryl ester is sparingly soluble (<20 µg/ml in PBS pH 7.2) in aqueous buffers. Therefore, further dilutions of the organic solvent solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. 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.

### Description

9(S)-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(S)-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(3)**, 563-566 (1970).
2. Belkner, J., Wiesner, R., Kühn, H., *et al.* The oxygenation of cholesterol esters by the reticulocyte lipoxygenase. *FEBS Lett.* **279(1)**, 110-114 (1991).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 01/17/2024

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM