

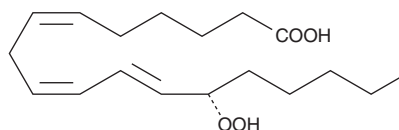
# PRODUCT INFORMATION



## 13(S)-HpOTrE( $\gamma$ )

Item No. 45210

**CAS Registry No.:** 121107-97-9  
**Formal Name:** 13S-hydroperoxy-6Z,9Z,11E-octadecatrienoic acid  
**MF:** C<sub>18</sub>H<sub>30</sub>O<sub>4</sub>  
**FW:** 310.4  
**Purity:**  $\geq$ 98%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 236 nm  
**Supplied as:** A solution in ethanol  
**Storage:** -80°C  
**Stability:**  $\geq$ 2 years



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

### Laboratory Procedures

13(S)-HpOTrE( $\gamma$ ) is supplied as a solution in ethanol. To change the solvent, 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 13(S)-HpOTrE( $\gamma$ ) in these solvents is approximately 50 mg/ml.

Further dilutions of the stock 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. If an organic solvent-free solution of 13(S)-HpOTrE( $\gamma$ ) is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of 13(S)-HpOTrE( $\gamma$ ) in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

13(S)-HpOTrE( $\gamma$ ) is a monohydroxy polyunsaturated fatty acids produced by the action of soybean lipoxygenase-1 (LO-1) on  $\gamma$ -linolenic acid.<sup>1</sup> Further action of soybean LO-1 converts 13(S)-HpOTrE( $\gamma$ ) to all four isomers of 6,13-DiHOTrE.<sup>2</sup> At concentrations greater than 100  $\mu$ M, 13(S)-HpOTrE( $\gamma$ ) inhibits the activity of soybean LO-1.<sup>3</sup>

### References

1. Funk, M.O., Isaac, R., and Porter, N.A. Preparation and purification of lipid hydroperoxides from arachidonic and  $\gamma$ -linolenic acids. *Lipids* **11**, 113-116 (1976).
2. Kim, M.R. and Sok, D.-E. Formation of 6,13-dihydroxyoctadecatrienoic acid isomers from  $\gamma$ -linolenic acid. *Biochem. Biophys. Res. Commun.* **159**, 1154-1160 (1989).
3. Kim, M.R. and Sok, D.-E. Irreversible inhibition of soybean lipoxygenase-1 by hydroperoxy acids as substrates. *Arch. Biochem. Biophys.* **288**, 270-275 (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, 03/07/2024

#### 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