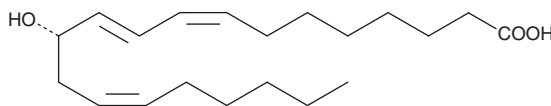


# PRODUCT INFORMATION



## 12(S)-HETrE Item No. 18483

CAS Registry No.: 72710-10-2  
Formal Name: 12S-hydroxy-8Z,10E,14Z-eicosatrienoic acid  
MF:  $C_{20}H_{34}O_3$   
FW: 322.5  
Purity:  $\geq 98\%$   
UV/Vis.:  $\lambda_{\max}$ : 235 nm  
Supplied as: A 100  $\mu\text{g/ml}$  solution in ethanol  
Storage:  $-20^\circ\text{C}$   
Stability:  $\geq 2$  years



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

### Laboratory Procedures

12(S)-HETrE 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 12(S)-HETrE in these solvents is miscible.

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 12(S)-HETrE is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of 12(S)-HETrE in PBS, pH 7.2, is approximately 0.8 mg/ml. For greater aqueous solubility, 12(S)-HETrE can be directly dissolved in 0.1 M  $\text{Na}_2\text{CO}_3$  (solubility of 2 mg/ml) and then diluted with PBS (pH 7.2) to achieve the desired concentration or pH. We do not recommend storing the aqueous solution for more than one day.

### Description

12(S)-HETrE is produced by 12-lipoxygenase oxidation of dihomo- $\gamma$ -linolenic acid (DGLA).<sup>1</sup> 12(S)-HETrE is reported to inhibit agonist-mediated platelet activation ( $\text{IC}_{50} = 40 \mu\text{M}$ ),  $\alpha$  granule secretion, integrin  $\alpha\text{IIb}\beta 3$  activation, Rap1 activation, and thrombin-induced clot retraction *in vitro*.<sup>1</sup>

### Reference

1. Ikei, K.N., Yeung, J., Apopa, P.L., *et al.* Investigations of human platelet-type 12-lipoxygenase: Role of lipoxygenase products in platelet activation. *J. Lipid Res.* **53**(12), 2546-2559 (2012).

#### 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, 02/05/2026

#### 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