

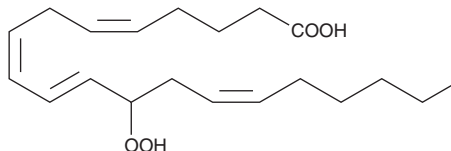
# PRODUCT INFORMATION



## (±)12-HpETE

Item No. 10138

**CAS Registry No.:** 71030-35-8  
**Formal Name:** 12-hydroperoxy-5Z,8Z,10E,14Z-  
eicosatetraenoic acid  
**MF:** C<sub>20</sub>H<sub>32</sub>O<sub>4</sub>  
**FW:** 336.5  
**Purity:** ≥96%  
**UV/Vis.:** λ<sub>max</sub>: 237 nm  
**Supplied as:** A solution in ethanol  
**Storage:** -80°C  
**Stability:** ≥6 months



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

### Laboratory Procedures

(±)12-HpETE 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. (±)12-HpETE is miscible in these solvents.

(±)12-HpETE is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of (±)12-HpETE should be diluted with the aqueous buffer of choice. The solubility of (±)12-HpETE in PBS (pH 7.2) and 0.1 M Na<sub>2</sub>CO<sub>3</sub> is approximately 0.5 and 2 mg/ml, respectively. We do not recommend storing the aqueous solution for more than one day.

### Description

(±)12-HpETE is one of the six monohydroperoxy fatty acids produced by the non-enzymatic oxidation of arachidonic acid (Item No. 90010) and consists of an equal mixture of the R and S isomers. Reduction of the hydroperoxide yields the more stable hydroxyl fatty acid (±)12-HETE (Item No. 34550). The biological activity of (±)12-HpETE has not been documented.

#### 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, 11/13/2023

#### 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