

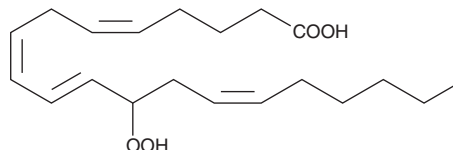
# PRODUCT INFORMATION



## (±)12-HpETE

Item No. 10138

CAS Registry No.: 71030-35-8  
Formal Name: 12-hydroperoxy-5Z,8Z,10E,14Z-eicosatetraenoic acid  
MF:  $C_{20}H_{32}O_4$   
FW: 336.5  
Purity:  $\geq 96\%$   
Stability:  $\geq 6$  months at  $-80^{\circ}C$   
Supplied as: A solution in ethanol  
UV/Vis.:  $\lambda_{max}$ : 237 nm  $\epsilon$ : 27,000



### Laboratory Procedures

For long term storage, we suggest that (±)12-HpETE be stored as supplied at  $-80^{\circ}C$ . It should be stable for at least six months.

(±)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.

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-HpETE is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of (±)12-HpETE in PBS (pH 7.2) is approximately 0.5 mg/ml. For greater aqueous solubility, (±)12-HpETE can be directly dissolved in 0.1 M  $Na_2CO_3$  (2 mg/ml) and then diluted with PBS (pH 7.2) to achieve the desired concentration or pH. Store aqueous solutions of (±)12-HpETE on ice and use within 12 hours of preparation.

### 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, 08/13/2015

#### 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