

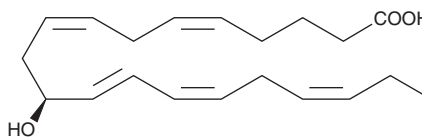
# PRODUCT INFORMATION



## 11(S)-HEPE

Item No. 32510

**CAS Registry No.:** 156473-26-6  
**Formal Name:** 11S-hydroxy-5Z,8Z,12E,14Z,17Z-eicosapentaenoic acid  
**MF:** C<sub>20</sub>H<sub>30</sub>O<sub>3</sub>  
**FW:** 318.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 236 nm  
**Supplied as:** A solution in ethanol  
**Storage:** -20°C  
**Stability:** ≥2 years  
**Special Conditions:** Oxygen and light sensitive



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

### Laboratory Procedures

11(S)-HEPE 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. 11(S)-HEPE 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 11(S)-HEPE is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of 11(S)-HEPE in PBS, pH 7.2 is approximately 0.8 mg/ml. For greater aqueous solubility, 11(S)-HEPE can be directly dissolved in 0.1 M Na<sub>2</sub>CO<sub>3</sub> (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

11(S)-HEPE is a monohydroxy fatty acid derived from EPA, theoretically by the action of COX. Although its arachidonate derived congener 11(S)-HETE has been isolated from cultured neutrophils, it is not known if similar incubations with EPA can produce 11(S)-HEPE.<sup>1</sup>

### Reference

1. Myers, R.F. and Siegel, M.I. Differential effects of anti-inflammatory drugs in lipoxygenase and cyclooxygenase activities of neutrophils from a reverse passive arthus reaction. *Biochem. Biophys. Res. Commun.* **112**, 586-594 (1983).

#### 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, 01/19/2024

#### 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