

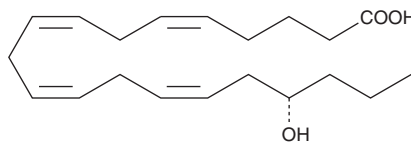
# PRODUCT INFORMATION



## 17(S)-HETE

Item No. 10011305

**CAS Registry No.:** 183509-25-3  
**Formal Name:** 17S-hydroxy-5Z,8Z,11Z,14Z-eicosatetraenoic acid  
**Synonym:** 17(S)-Hydroxyeicosatetraenoic Acid  
**MF:** C<sub>20</sub>H<sub>32</sub>O<sub>3</sub>  
**FW:** 320.5  
**Purity:** ≥98%  
**Supplied as:** A solution in ethanol  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

17(S)-HETE 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. 17(S)-HETE is miscible in organic solvents such as DMSO and dimethyl formamide. 17(S)-HETE is soluble in 0.1 M Na<sub>2</sub>CO<sub>3</sub> at a concentration of approximately 2 mg/ml

17(S)-HETE is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 17(S)-HETE should be diluted with the aqueous buffer of choice. The solubility of 17(S)-HETE in PBS (pH 7.2) is approximately 0.8 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Electrolyte and fluid transport in the kidney are regulated in part by arachidonic acid and its metabolites. 17-HETE is a cytochrome P450 (CYP450) metabolite of arachidonic acid that has stereospecific effects on sodium transport in the kidney. 17(S)-HETE inhibits proximal tubule ATPase activity by as much as 70% at a concentration of 2 μM.<sup>1</sup>

### Reference

1. Carroll, M.A., Balazy, M., Margiotta, P., *et al.* Cytochrome P-450-dependent HETEs: Profile of biological activity and stimulation by vasoactive peptides. *Am. J. Physiol.* **271**(4 Pt 2), R863-R869 (1996).

#### 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, 09/30/2022

#### 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