

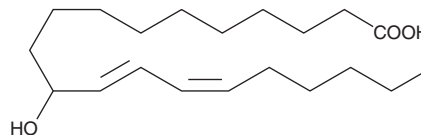
# PRODUCT INFORMATION



## (±)11-HEDE

Item No. 37500

**CAS Registry No.:** 5598-37-8  
**Formal Name:** 11-hydroxy-12E,14Z-eicosadienoic acid  
**Synonym:** 11-hydroxy-12,14-Eicosadienoic Acid  
**MF:** C<sub>20</sub>H<sub>36</sub>O<sub>3</sub>  
**FW:** 324.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 236 nm  
**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

(±)11-HEDE 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 ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of (±)11-HEDE in these solvents is approximately 50 mg/ml.

(±)11-HEDE is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of (±)11-HEDE should be diluted with the aqueous buffer of choice. The solubility of (±)11-HEDE in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

(±)11-HEDE is a racemic mixture of the monohydroxy fatty acids 11(S)-HEDE (Item No. 37510) and 11(R)-HEDE (Item No. 37505). 11-HEDE is formed from eicosadienoic acid by cyclooxygenase (COX) and in macrophages.<sup>1,2</sup>

### References

1. Hemler, M.E., Crawford, C.G., and Lands, W.E.M. Lipoxygenation activity of purified prostaglandin-forming cyclooxygenase. *Biochemistry* **17**(9), 1772-1779 (1978).
2. Rabinovitch, H., Durand, J., Gualde, N., *et al.* Metabolism of polyunsaturated fatty acids by mouse peritoneal macrophages: The lipoxygenase metabolic pathway. *Agents Actions* **11**(6-7), 580-583 (1981).

#### 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, 01/16/2024

#### 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