

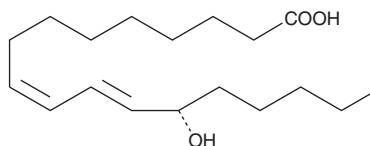
# PRODUCT INFORMATION



## 13(S)-HODE

Item No. 38610

CAS Registry No.: 29623-28-7  
Formal Name: 13S-hydroxy-9Z,11E-octadecadienoic acid  
MF: C<sub>18</sub>H<sub>32</sub>O<sub>3</sub>  
FW: 296.5  
Purity: ≥98%  
UV/Vis.: λ<sub>max</sub>: 234 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

13(S)-HODE 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. The solubility of 13(S)-HODE in these solvents is approximately 50 mg/ml.

13(S)-HODE is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanol solution of 13(S)-HODE should be diluted with the aqueous buffer of choice. The solubility of 13(S)-HODE in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

13(S)-HODE is produced by incubation of linoleic acid with plant and mammalian lipoxygenases. It has been shown to inhibit the adhesion of tumor cells to the endothelium at concentrations around 1 μM,<sup>1,2</sup> and down regulates the expression of the IRGpIIb/IIIa receptor.<sup>3</sup>

### References

1. Buchanan, M.R., Haas, T.A., Lagarde, M., *et al.* 13-Hydroxyoctadecadienoic acid is the vessel wall chemorepellant factor, LOX. *J. Biol. Chem.* **260(30)**, 16056-16059 (1985).
2. Honn, K.V., Nelson, K.K., Renaud, C., *et al.* Fatty acid modulation of tumor cell adhesion to microvessel endothelium and experimental metastasis. *Prostaglandins* **44(5)**, 413-429 (1992).
3. Grossi, I.M., Fitzgerald, L.A., Umbarger, L.A., *et al.* Bidirectional control of membrane expression and/or activation of the tumor cell IRGpIIb/IIIa receptor and tumor cell adhesion by lipoxygenase products of arachidonic acid and linoleic acid. *Cancer Res.* **49(4)**, 1029-1037 (1989).

#### 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, 01/17/2024

#### 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