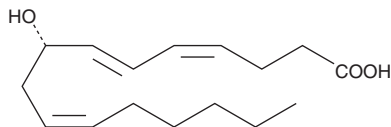


PRODUCT INFORMATION

tetranor-12(S)-HETE

Item No. 10007207

CAS Registry No.: 121842-79-3
Formal Name: 8S-hydroxy-4Z,6E,10Z-hexadecatrienoic acid
Synonym: tetranor-12(S)-Hydroxyeicosatetraenoic Acid
MF: C₁₆H₂₆O₃
FW: 266.4
Purity: ≥98%
UV/Vis.: λ_{max}: 234 nm
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥1 year



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

Laboratory Procedures

tetranor-12(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. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of tetranor-12(S)-HETE in these solvents is at least 25 mg/ml.

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 tetranor-12(S)-HETE is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of tetranor-12(S)-HETE in PBS (pH 7.2) is at least 0.5 mg/ml. For greater aqueous solubility, tetranor-12(S)-HETE can be directly dissolved in 0.1 M Na₂CO₃ (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

12(S)-HETE is a product of arachidonic acid metabolism through the 12-lipoxygenase pathway. It is primarily found in platelets, leukocytes, and to a lesser extent in smooth muscle cells.¹ It enhances tumor cell adhesion to endothelial cells, fibronectin, and the subendothelial matrix.² tetranor-12(S)-HETE is the major β-oxidation product resulting from peroxisomal metabolism of 12(S)-HETE in numerous tissues, and Lewis lung carcinoma cells.^{3,4} No biological function has yet been determined for tetranor-12(S)-HETE. Some data indicate it may play a role in controlling the inflammatory response in injured corneas.⁴ In some diseases (e.g., Zellweger's Syndrome) peroxisomal abnormalities result in the inability of cells to metabolize 12(S)-HETE, which may be responsible for symptoms of the disease. The tetranor derivative of 12(S)-HETE is available as a research tool for the elucidation of the metabolic fate of its parent compound.

References

1. Evans, J.F., Leblanc, Y., Fitzsimmons, B.J., *et al. Biochim. Biophys. Acta* **917**, 406-410 (1987).
2. Grossi, I.M., Fitzgerald, L.A., Umbarger, L.A., *et al. Cancer Res.* **49**, 1029-1037 (1989).
3. Wigren, J., Herbertsson, H., Tollbom, Ö., *et al. J. Lipid Res.* **34**, 625-631 (1993).
4. Hurst, J.S., Bazan, H.P., and Balazy, M. *Exp. Eye Res.* **59**, 97-106 (1994).

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, 05/05/2022

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