

PRODUCT INFORMATION



4(S),11(R)-DiHDoHE

Item No. 9001290

Formal Name: 4S,11R-dihydroxydocosa-5Z,7E,9E,13Z,16Z,19Z-hexaenoic acid

MF: C₂₂H₃₂O₄

FW: 360.5

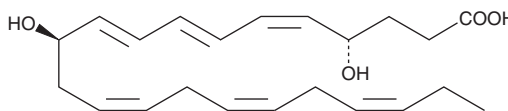
Purity: ≥95%

UV/Vis.: λ_{max}: 270 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

4(S),11(R)-DiHDoHE 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. 4(S),11(R)-DiHDoHE 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 4(S),11(R)-DiHDoHE is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of 4(S),11(R)-DiHDoHE in PBS, pH 7.2, is approximately 0.5 mg/ml. For greater aqueous solubility, 4(S),11(R)-DiHDoHE can be directly dissolved in 0.1 M Na₂CO₃ (solubility of 1 mg/ml) and then diluted with PBS (pH 7.2) to achieve the desired concentration. We do not recommend storing the aqueous solution for more than one day.

Description

Docosahexaenoic acid (DHA; Item No. 90310) is an ω-3 fatty acid that is abundant in the brain and the retina and is known to be important in early development.^{1,2} Recently, various DHA metabolites, including 7(R)-Maresin 1 (Item No. 10878) and 10(S),17(S)-DiHDoHE (Item No. 10008128), have been shown to block infiltration of neutrophils during an acute inflammatory response.^{3,4} 4(S),11(R)-DiHDoHE is a dihydroxy-DHA derivative. The biological actions of this compound have not yet been determined.

References

1. Su, H.-M. Mechanisms of n-3 fatty acid-mediated development and maintenance of learning memory performance. *J. Nutr. Biochem.* **21**(5), 364-373 (2010).
2. Wu, T.C. and Chen, P.H. Health consequences of nutrition in childhood and early infancy. *Pediatr. Neonatol.* **50**(4), 135-142 (2009).
3. Serhan, C.N., Yang, R., Martinod, K., *et al.* Maresins: Novel macrophage mediators with potent antiinflammatory and proresolving actions. *J. Exp. Med.* **206**(1), 15-23 (2009).
4. Serhan, C.N., Gotlinger, K., Hong, S., *et al.* Anti-inflammatory actions of neuroprotectin D1/protectin D1 and its natural stereoisomers: Assignments of dihydroxy-containing docosatrienes. *J. Immunol.* **176**(3), 1848-1859 (2006).

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, 11/13/2019

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