

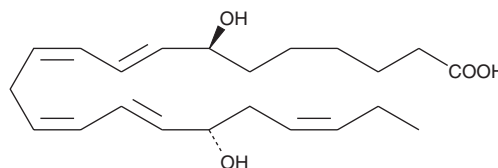
Product Information



7(S),17(S)-dihydroxy-8(E),10(Z),13(Z), 15(E),19(Z)-Docosapentaenoic Acid

Item No. 10546

CAS Registry No.: 887752-13-8
Formal Name: 7S,17S-dihydroxy-8E,10Z,13Z,15E,19Z-docosapentaenoic acid
Synonym: 7(S),17(S)-hydroxy DPA
MF: C₂₂H₃₄O₄
FW: 362.5
Purity: ≥98%
Stability: ≥1 year at -20°C
Supplied as: A solution in ethanol



Laboratory Procedures

For long term storage, we suggest that 7(S),17(S)-dihydroxy-8(E),10(Z),13(Z),15(E),19(Z)-docosapentaenoic acid (7(S),17(S)-hydroxy DPA) be stored as supplied at -20°C. It should be stable for at least one year.

7(S),17(S)-hydroxy DPA 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 7(S),17(S)-hydroxy DPA in these solvents is approximately 100 mg/ml.

If aqueous stock solutions are required for biological experiments, they can best be prepared by diluting the organic solvent into aqueous buffers or isotonic saline. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Novel oxylipins, referred to as docosanoids, have been derived from C22 polyunsaturated fatty acids. 7(S),17(S)-hydroxy DPA is a DPA-derived analog of the 17(S)-dihydroxy series of docosanoids known as protectins. Protectin D1, a DHA-derived dihydroxy fatty acid, exhibits potent anti-inflammatory activities.¹⁻³ Potentially, 7(S),17(S)-hydroxy DPA demonstrates similar properties; however, its biological activity has yet to be determined.

References

1. 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**, 1848-1859 (2006).
2. Ariel, A. and Serhan, C.N. Resolvins and protectins in the termination program of acute inflammation. *Trends Immunol.* **28(4)**, 176-183 (2007).
3. Schwab, J.M., Chiang, N., Arita, M., *et al.* Resolvin E1 and protectin D1 activate inflammation-resolution programmes. *Nature* **447**, 869-875 (2007).

Related Products

Docosapentaenoic Acid - Item No. 90165 • 17-keto-7(Z),10(Z),13(Z),15(E),19(Z)-Docosapentaenoic Acid - Item No. 9000347

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 03/17/2011

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com