

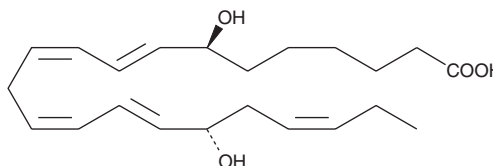
PRODUCT INFORMATION



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: 7,17-dihydroxy-8E,10Z,13Z,15E,19Z-docosapentaenoic acid
Synonyms: 7(S),17(S)-hydroxy DPA, Resolvin D5 n-3 DPA
MF: C₂₂H₃₄O₄
FW: 362.5
Purity: ≥98%
UV/Vis.: λ_{max}: 244 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

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

Description

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**(3), 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**(7146), 869-874 (2007).

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, 09/11/2024

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