

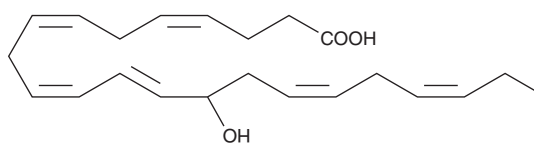
PRODUCT INFORMATION



(±)14-HDHA

Item No. 33550

CAS Registry No.: 87042-40-8
Formal Name: (±)14-hydroxy-4Z,7Z,10Z,12E,16Z,19Z-docosahexaenoic acid
Synonyms: 14-hydroxy Docosahexaenoic Acid, (±)14-HDoHE
MF: C₂₂H₃₂O₃
FW: 344.5
Purity: 98%
UV/Vis.: λ_{max}: 237 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

(±)14-HDHA 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. (±)14-HDHA 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 (±)14-HDHA is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of (±)14-HDHA in PBS (pH 7.2) is approximately 0.8 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

(±)14-HDHA is an autoxidation product of docosahexaenoic acid (DHA) *in vitro*.^{1,2} It is also produced from incubations of DHA in rat liver, brain, and intestinal microsomes.³⁻⁵ DHA is metabolized to 14(S)-HDHA by human platelets along with 11(S)-HDHA.⁵⁻⁷ 14(S)-HDHA is also produced by salmon gills upon stimulation with calcium ionophore.⁸ 14(S)-HDHA is an inhibitor of U-46619-induced human platelet aggregation and rabbit and rat aortic smooth muscle contraction with IC₅₀ values of about 70, 3.6, and 5.3 μM, respectively.^{7,9} (±)14-HDHA is a potential marker of oxidative stress in brain and retina where DHA is an abundant polyunsaturated fatty acid.

References

1. VanRollins, M. and Murphy, R.C. *J. Lipid Res.* **25(5)**, 507-517 (1984).
2. Reynaud, D., Thickitt, C.P., and Pace-Asciak, C.R. *Anal. Biochem.* **214(1)**, 165-170 (1993).
3. VanRollins, M., Baker, R.C., Sprecher, H., *et al. J. Biol. Chem.* **259(9)**, 5776-5783 (1984).
4. Yamane, M., Abe, A., and Yamane, S. *J. Chromatogr.* **652(2)**, 123-136 (1994).
5. Kim, H.Y., Karanian, J.W., Shingu, T., *et al. Prostaglandins* **40(5)**, 473-491 (1990).
6. Aveldaño, M.I. and Sprecher, H. *J. Biol. Chem.* **258(15)**, 9339-9343 (1983).
7. Lagarde, M., Croset, M., Guichardant, M., *et al. Adv. Exp. Med. Bio.* **192**, 327-335 (1985).
8. Bell, J.G., Dick, J.R., and Sargent, J.R. *Lipids* **28**, 819-826 (1993).
9. Karanian, J.W., Kim, H.Y., and Salem, N., Jr. *J. Pharmacol. Exp. Ther.* **270(3)**, 1105-1109 (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, 07/09/2018

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