

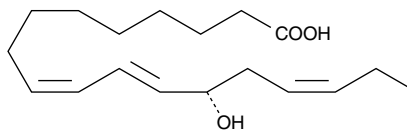
Product Information



13(S)-HOTrE

Item No. 39620

CAS Registry No:	87984-82-5
Formal Name:	13S-hydroxy-9Z,11E,15Z-octadecatrienoic acid
MF:	C ₁₈ H ₃₀ O ₃
FW:	294.4
Purity:	≥98%
Stability:	≥2 years at -20°C
Supplied as:	A solution in ethanol
UV/Vis:	λ _{max} : 234 nm ε: 23,000
Misc:	Oxygen and light sensitive



Laboratory Procedures

For long term storage, we suggest that 13(S)-HOTrE be stored as supplied at -20°C. It should be stable for at least two years.

13(S)-HOTrE is supplied as a solution in ethanol. To change the solvent, evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as dimethyl formamide or DMSO purged with an inert gas can be used. The solubility of 13(S)-HOTrE in these solvents is approximately 50 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 13(S)-HOTrE is needed, the ethanol can be evaporated under a stream of nitrogen and the neat oil dissolved in the buffer of choice. The solubility of 13(S)-HOTrE in PBS (pH 7.2) is approximately 1 mg/ml. More concentrated aqueous solutions of 13(S)-HOTrE can be prepared using concentrated basic buffers (pH > 8.0 and ionic strength ≥ 0.1 M). Add 400 µl of cold buffer (0°C) per mg of 13(S)-HOTrE and vortex vigorously until completely dissolved. Store aqueous solutions of 13(S)-HOTrE on ice and use within twelve hours. We do not recommend storing the aqueous solution for more than one day.

13(S)-HOTrE is the 15-lipoxygenase (15-LO) product of linolenic acid. It has been detected in cell membranes and as the cholesteryl ester associated with the lesions of atherosclerosis, and in the biomembranes of soybeans exposed to 15-LO.^{1,2}

References

1. Belkner, J., Wiesner, R., Kühn, H., *et al.* The oxygenation of cholesterol esters by the reticulocyte lipoxygenase. *FEBS Lett.* **279**, 110-114 (1991).
2. Maccarrone, M., van Aarle, P.G.M., Veldink, G.A., *et al.* In vitro oxygenation of soybean biomembranes by lipoxygenase-2. *Biochim. Biophys. Acta* **1190**, 164-169 (1994).

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

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, 04/08/2010