

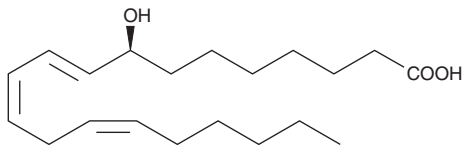
PRODUCT INFORMATION



8(S)-HETrE

Item No. 36360

CAS Registry No.: 889573-69-7
Formal Name: 8S-hydroxy-9E,11Z,14Z-eicosatrienoic acid
MF: C₂₀H₃₄O₃
FW: 322.5
Purity: ≥98%
UV/Vis.: λ_{max}: 236 nm
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years
Special Conditions: Oxygen and light sensitive



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

8(S)-HETrE 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 8(S)-HETrE in these solvents is miscible.

8(S)-HETrE is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 8(S)-HETrE should be diluted with the aqueous buffer of choice. The solubility of 8(S)-HETrE in PBS (pH 7.2) and 0.1 M Na₂CO₃ is approximately 0.8 and 2 mg/ml, respectively. We do not recommend storing the aqueous solution for more than one day.

Description

8(S)-HETrE is a metabolite of the ω-6 fatty acid γ-linolenic acid (GLA; Item No. 90220).¹ It is formed from GLA by 5-lipoxygenase (5-LO) via a dihomo-γ-linolenic acid (Item No. 90230) intermediate. Serum levels of 8(S)-HETrE are decreased in a mouse model of high-fat high-sucrose diet-induced obesity.

Reference

1. Du, Y., Li, D.-X., Lu, D.-Y., *et al.* Lipid metabolism disorders and lipid mediator changes of mice in response to long-term exposure to high-fat and high sucrose diets and ameliorative effects of mulberry leaves. *Food Funct.* **13(8)**, 4576-4591 (2022).

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, 01/16/2024

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