

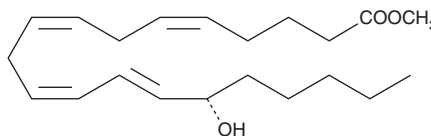
PRODUCT INFORMATION



15(S)-HETE methyl ester

Item No. 21327

CAS Registry No.: 70946-44-0
Formal Name: 15S-hydroxy-5Z,8Z,11Z,13E-eicosatetraenoic acid, methyl ester
Synonym: 15(S)-Hydroxyeicosatetraenoic Acid methyl ester
MF: C₂₁H₃₄O₃
FW: 334.5
Purity: ≥98%
UV/Vis.: λ_{max}: 236 nm
Supplied as: A 100 µg/ml solution in methanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

15(S)-HETE methyl ester is supplied as a solution in methanol. To change the solvent, simply evaporate the methanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. 15(S)-HETE methyl ester 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 15(S)-HETE methyl ester is needed, it can be prepared by evaporating the methanol and directly dissolving the neat oil in aqueous buffers. The solubility of 15(S)-HETE methyl ester in PBS (pH 7.2) is approximately 0.1 mg/ml. For maximum solubility in aqueous buffers, the methanolic solution of 15(S)-HETE methyl ester should be diluted with the aqueous buffer of choice. 15(S)-HETE methyl ester has a solubility of 0.2 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

15(S)-HETE methyl ester is a synthetic derivative of 15(S)-HETE (Item No. 34720), a major arachidonic acid metabolite from the 15-lipoxygenase pathway. Methyl esters of lipids are commonly used in formulations of nutritional supplements.

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, 11/12/2025

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