

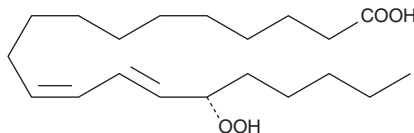
PRODUCT INFORMATION



15(S)-HpEDE

Item No. 47720

CAS Registry No.: 145375-41-3
Formal Name: 15S-hydroperoxy-11Z,13E-eicosadienoic acid
MF: C₂₀H₃₆O₄
FW: 340.5
Purity: ≥95%
UV/Vis.: λ_{max}: 234 nm
Supplied as: A solution in ethanol
Storage: -80°C
Stability: ≥2 years



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

Laboratory Procedures

15(S)-HpEDE 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 DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of 15(S)-HpEDE 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 15(S)-HpEDE is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of 15(S)-HpEDE in PBS, pH 7.2, is approximately 1 mg/ml. Store aqueous solutions of 15(S)-HpEDE on ice and use within 12 hours of preparation. Although the aqueous solutions of 15(S)-HpEDE may be stable for more than 12 hours, we strongly recommend using a fresh preparation each day.

Description

15(S)-HpEDE is a monohydroperoxy polyunsaturated fatty acid produced by the action of 15-lipoxygenase on eicosadienoic acid. Although the biological activities of 15(S)-HpEDE have not been well characterized, it is expected to behave similarly to 15(S)-HpETE (Item No. 44720).

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, 03/11/2024

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