

PRODUCT INFORMATION



9(S)-HpODE

Item No. 48410

CAS Registry No.: 29774-12-7

Formal Name: 9S-hydroperoxy-10E,12Z-octadecadienoic acid

MF: $C_{18}H_{32}O_4$

FW: 312.4

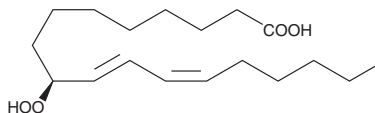
Purity: $\geq 98\%$

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

9(S)-HpODE 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 9(S)-HpODE 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 9(S)-HpODE is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of 9(S)-HpODE in PBS, pH 7.2, is approximately 1 mg/ml. 9(S)-HpODE is highly unstable in aqueous solutions. We recommend that 9(S)-HpODE diluted in aqueous solution be used as soon as possible, preferably within 15 minutes.

Description

9(S)-HpODE is produced by the action of arachidonate 5-lipoxygenase on linoleic acid. It can be further metabolized by potato hydroperoxide dehydratase to colneleic acid.^{1,2}

References

1. Galliard, T. and Phillips, D.R. The enzymic conversion of linoleic acid into 9-(nona-1',3'-dienoxy)non-8-enoic acid, a novel unsaturated ether derivative isolated from homogenates of *Solanum tuberosum* tubers. *Biochem. J.* **129**, 743-753 (1972).
2. Fahlstadius, P. and Hamberg, M. Stereospecific removal of the *pro-R* hydrogen at C-8 of (9S)-hydroperoxy octadecadienoic acid in the biosynthesis of colneleic acid. *J. Chem. Soc. Perkin Trans. 1*, 2027-2030 (1990).

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