

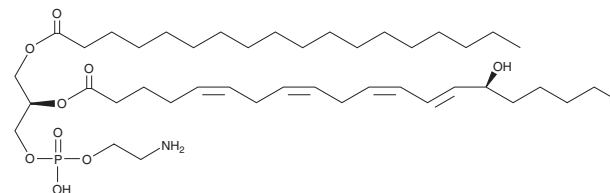
PRODUCT INFORMATION



1-Stearoyl-2-15(S)-HETE-*sn*-glycero-3-PE

Item No. 21139

CAS Registry No.: 947381-58-0
Formal Name: 15S-hydroxy-5Z,8Z,11Z,13E-eicosatetraenoic acid, (1R)-2-[[[(2-aminoethoxy)hydroxyphosphinyl]oxy]-1-[[[(1-oxooctadecyl)oxy]methyl]ethyl ester
Synonyms: 15(S)-HETE-SAPE, 15(S)-Hydroxyeicosatetraenoic Acid-SAPE, 1-Stearoyl-2-15(S)-HETE-*sn*-glycero-3-Phosphatidylethanolamine
MF: C₄₃H₇₈NO₉P
FW: 784.1
Purity: ≥98%
UV/Vis.: λ_{max}: 237 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

1-Stearoyl-2-15(S)-HETE-*sn*-glycero-3-PE is supplied as a solution in ethanol. 1-Stearoyl-2-15(S)-HETE-*sn*-glycero-3-PE is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 1-stearoyl-2-15(S)-HETE-*sn*-glycero-3-PE should be diluted with the aqueous buffer of choice. 1-Stearoyl-2-15(S)-HETE-*sn*-glycero-3-PE has a solubility of approximately 0.3 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method.

Description

1-Stearoyl-2-15(S)-HETE-*sn*-glycero-3-PE is a phospholipid that contains stearic acid (Item No. 10011298) at the *sn*-1 position and 15(S)-HETE (Item No. 34720) at the *sn*-2 position. It is formed in human peripheral monocytes activated by the calcium ionophore A23187 (Item No. 11016) by direct oxidation of 1-stearoyl-2-arachidonoyl-*sn*-glycero-3-PE (SAPE; Item No. 25871) by 15-LO.^{1,2} Phosphoethanolamine (PE) HETEs (PE-HETEs), including 1-stearoyl-2-15(S)-HETE-*sn*-glycero-3-PE, are the main source of esterified HETE in ionophore-activated monocytes.

References

1. Maskrey, B.H., Bermúdez-Fajardo, A., Morgan, A.H., *et al.* Activated platelets and monocytes generate four hydroxyphosphatidylethanolamines via lipoxygenase. *J. Biol. Chem.* **282**(28), 20151-20163 (2007).
2. Morgan, A.H., Hammond, V.J., Morgan, L., *et al.* Quantitative assays for esterified oxylipins generated by immune cells. *Nat. Protoc.* **5**(12), 1919-1931 (2010).

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, 05/05/2022

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