

PRODUCT INFORMATION



1-Stearoyl-2-15(S)-HETE-*sn*-glycero-3-PE-d₁₁

Item No. 27931

Formal Name: 15S-hydroxy-5Z,8Z,11Z,13E-eicosatetraenoic acid-16,16,17,17,18,18,19,19,20,20,20-d₁₁, (1R)-2-[[[(2-aminoethoxy)hydroxyphosphinyl]oxy]-1-[[[(1-oxooctadecyl)oxy]methyl]ethyl] ester

Synonyms: 15(S)-HETE-SAPE-d₁₁, 15(S)-Hydroxyeicosatetraenoic Acid-SAPE-d₁₁, 1-Stearoyl-2-15(S)-HETE-*sn*-glycero-3-Phosphatidylethanolamine-d₁₁

MF: C₄₃H₆₇D₁₁NO₉P

FW: 795.1

Chemical Purity: ≥98% (1-Stearoyl-2-15(S)-HETE-*sn*-glycero-3-PE)

Deuterium

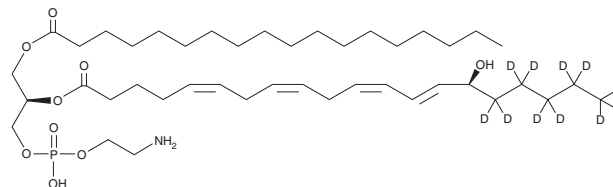
Incorporation: ≥99% deuterated forms (d₁-d₁₁); ≤1% d₀

UV/Vis.: λ_{max}: 235 nm

Supplied as: A 500 µg/ml 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-d₁₁ is intended for use as an internal standard for the quantification of 1-stearoyl-2-15(S)-HETE-*sn*-glycero-3-PE (Item No. 21139) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

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-lipoxygenase (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, 12/19/2025

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