

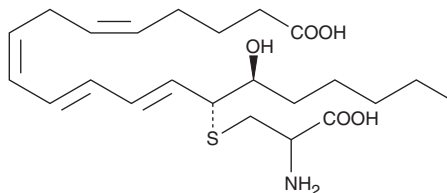
Product Information



14,15-Leukotriene E₄

Catalog No. 10011362

CAS Registry No.:	1000852-57-2
Formal Name:	15S-hydroxy-14R-(S-cysteinyl)-5Z,8Z,10E,12E-eicosatetraenoic acid
Synonyms:	Eoxin E ₄ , EXE ₄ , 14,15-LTE ₄
MF:	C ₂₃ H ₃₇ NO ₅ S
FW:	439.6
Purity:	≥97%
Stability:	≥1 year at -80°C
Supplied as:	A solution in methanol



Laboratory Procedures

For long term storage, we suggest that 14,15-leukotriene E₄ (14,15-LTE₄) be stored as supplied at -80°C. It should be stable for at least one year.

14,15-LTE₄ 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. The solubility of 14,15-LTE₄ 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 14,15-LTE₄ is needed, it can be prepared by evaporating the methanol and directly dissolving the neat oil in aqueous buffers. The solubility of 14,15-LTE₄ in PBS, pH 7.2, is approximately 100 µg/ml. We do not recommend storing the aqueous solution for more than one day.

Leukotrienes (LTs) are a group of acute inflammatory mediators derived from arachidonic acid in leukocytes. The majority of these metabolites are formed through the 5-lipoxygenase (5-LO) pathway.¹ 14,15-LTE₄ is a metabolite of 14,15-LTC₄ and 14,15-LTD₄, an alternate class of LTs synthesized by a pathway involving the dual actions of 15- and 12-LOs on arachidonic acid *via* 15-HpETE and 14,15-LTA₄ intermediates.²⁻⁵ These metabolites are classified as eoxins because they are formed mostly by eosinophils.⁴ Mast cells and nasal polyps can synthesize 14,15-LTC₄ as well, however metabolism to 14,15-LTE₄ in these cells and tissue has not been documented. 14,15-LTE₄ increases vascular permeability of human endothelial cell monolayers with about 10-fold less potency than LTC₄, but approximately 100-fold greater potency than histamine.⁴

References

1. Luo, M., Lee, S., and Brock, T.G. Leukotriene synthesis by epithelial cells. *Histol. Histopathol.* **18**, 587-595 (2003).
2. Yokoyama, C., Shinjo, F., Yoshimoto, T., *et al.* Arachidonate 12-lipoxygenase purified from porcine leukocytes by immunoaffinity chromatography and its reactivity with hydroperoxyeicosatetraenoic acids. *J. Biol. Chem.* **261**, 16714-16721 (1986).
3. Bryant, R.W., Schewe, T., Rapoport, S.M., *et al.* Leukotriene formation by a purified reticulocyte lipoxygenase enzyme. Conversion of arachidonic acid and 15-hydroperoxyeicosatetraenoic acid to 14,15-leukotriene A₄. *J. Biol. Chem.* **260**, 3548-3555 (1985).
4. Feltenmark, S., Gautam, N., Brunnström, Å., *et al.* Eoxins are proinflammatory arachidonic acid metabolites produced *via* the 15-lipoxygenase-1 pathway in human eosinophils and mast cells. *Proc. Natl. Acad. Sci. USA* **105**(2), 680-685 (2008).
5. Sailesh, S., Kumar, Y.V.K., Prasad, M., *et al.* Sheep uterus dual lipoxygenase in the synthesis of 14,15-leukotrienes. *Arch. Biochem. Biophys.* **315**(2), 362-368 (1994).

Related Products

Leukotriene E₄ - Cat. No. 20410 • N-acetyl Leukotriene E₄ - Cat. No. 20420 • Leukotriene E₄ EIA Kit - Cat. No. 520411 • Cysteinyl Leukotriene EIA Kit - Cat. No. 520501 • 14,15-Leukotriene C₄ EIA Kit - Cat. No. 10006748 • Cysteinyl Leukotriene Express EIA Kit - Cat. No. 10009291 • 14,15-Leukotriene C₄ - Cat. No. 10011360

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants **only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 09/08/2009

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com