

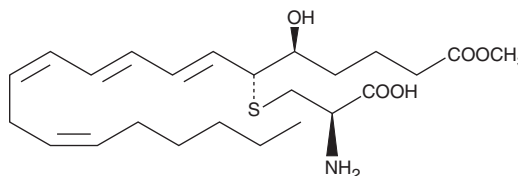
PRODUCT INFORMATION



Leukotriene E₄ methyl ester

Item No. 10007166

CAS Registry No.: 89461-65-4
Formal Name: 5S-hydroxy-6R-(S-cysteinyl)-7E,9E,11Z,14Z-eicosatetraenoic acid, methyl ester
Synonym: LTE₄ methyl ester
MF: C₂₄H₃₉NO₅S
FW: 453.6
Purity: ≥97%
UV/Vis.: λ_{max}: 281 nm
Supplied as: A solution in ethanol
Storage: -80°C
Stability: ≥1 year
Special Conditions: Light-sensitive



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

Laboratory Procedures

Leukotriene E₄ (LTE₄) methyl ester is supplied as a solution in ethanol. To change the solvent, simply 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 LTE₄ methyl ester in these solvents is approximately 50 mg/ml.

LTE₄ methyl ester is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of LTE₄ methyl ester should be diluted with the aqueous buffer of choice. The solubility of LTE₄ methyl ester in PBS (pH 7.2) is approximately 0.1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

LTE₄ is produced by the action of a dipeptidase on LTD₄ leaving only the cysteinyl group still attached to the fatty acid backbone.¹ It is one of the constituents of slow-reacting substance of anaphylaxis (SRS-A).² LTE₄ is considerably less active (8 to 12-fold) than LTC₄ in the biological activities characteristic of cysteinyl LTs.^{1,3} Unlike LTC₄ and LTD₄, LTE₄ accumulates in both plasma and urine. Therefore, urinary excretion of LTE₄ is most often used as an indicator of asthma.⁴⁻⁶ LTE₄ methyl ester is a more lipid soluble form of LTE₄. The biological activity of LTE₄ methyl ester has not been reported.

References

1. Bernström, K. and Hammarström, S. Metabolism of leukotriene D by porcine kidney. *J. Biol. Chem.* **256(18)**, 9579-9582 (1981).
2. Samuelsson, B. Leukotrienes: Mediators of immediate hypersensitivity reactions and inflammation. *Science* **220(4597)**, 568-575 (1983).
3. Lefer, A.M. Leukotrienes as mediators of ischemia and shock. *Biochem. Pharmacol.* **35(2)**, 123-127 (1986).
4. Kumlin, M., Stensvad, F., Larsson, L., *et al.* Validation and application of a new simple strategy for measurements of urinary leukotriene E₄ in humans. *Clin. Exp. Allergy* **25(5)**, 467-479 (1995).
5. Drazen, J.M., O'Brien, J., Sparrow, D., *et al.* Recovery of leukotriene E₄ from the urine of patients with airway obstruction. *Am. Rev. Respir. Dis.* **146(1)**, 104-108 (1992).
6. Kumlin, M., Dahlén, B., Björck, T., *et al.* Urinary excretion of leukotriene E₄ and 11-dehydro-thromboxane B₂ in response to bronchial provocations with allergen, aspirin, leukotriene D₄, and histamine in asthmatics. *Am. Rev. Respir. Dis.* **146(1)**, 96-103 (1992).

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, 12/02/2024

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