

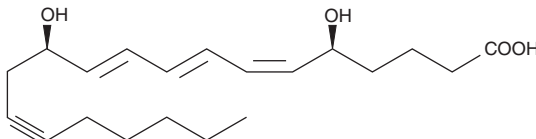
Product Information



14,15-dehydro Leukotriene B₄

Catalog No. 20150

CAS Registry No.: 114616-11-4
Formal Name: 5S,12R-dihydroxy-6Z,8E,10E-eicosatrien-14-ynoic acid
Synonym: 14,15-dehydro LTB₄
MF: C₂₀H₃₀O₄
FW: 334.5
Purity: ≥97%
Stability: ≥1 year at -80°C
Supplied as: A solution in ethanol



Laboratory Procedures

For long term storage, we suggest that 14,15-dehydro leukotriene B₄ (14,15-dehydro LTB₄) be stored as supplied at -80°C. It should be stable for at least one year.

14,15-dehydro LTB₄ 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. 14,15-dehydro LTB₄ is miscible in these solvents.

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-dehydro LTB₄ is needed, the ethanol can be evaporated under a stream of nitrogen and the neat oil dissolved directly in the buffer of choice. 14,15-dehydro LTB₄ is soluble in PBS, pH 7.2, at a concentration of 1 mg/ml. Be certain that your buffers are free of oxygen, transition metal ions, and redox active compounds. We do not recommend storing the aqueous solution for more than one day.

LTB₄ is a dihydroxy fatty acid derived from arachidonic acid through the 5-lipoxygenase pathway.^{1,2} It promotes a number of leukocyte functions including aggregation, stimulation of ion fluxes, enhancement of lysosomal enzyme release, superoxide anion production, chemotaxis, and chemokinesis.^{3,4} At least two LTB₄ receptors, termed BLT₁ and BLT₂, have been identified. 14,15-dehydro LTB₄ is a LTB₄ receptor antagonist that has a higher binding affinity for BLT₁, demonstrating a K_i value of 27 nM, compared to BLT₂, which has a K_i value of 473 nM.⁴ 14,15-dehydro LTB₄ inhibits LTB₄-induced release of lysozymes from rat polymorphonuclear leukocytes with an IC₅₀ value of 1 μM.⁵

References

1. Rådmark, O., Malmsten, C., Samuelsson, B., *et al.* Leukotriene A: Stereochemistry and enzymatic conversion to leukotriene B. *Biochem. Biophys. Res. Commun.* **92**, 954-961 (1980).
2. McGee, J. and Fitzpatrick, F. Enzymatic hydration of leukotriene A₄. *J. Biol. Chem.* **260**, 12832-12837 (1985).
3. Ford-Hutchinson, A.W. Leukotriene B₄ in inflammation. *Crit. Rev. Immunol.* **10**, 1-12 (1990).
4. McMillan, R.M. and Foster, S.J. Leukotriene B₄ and inflammatory disease. *Agents Actions* **24**, 114-119 (1988).
5. Wang, S., Gustafson, E., Pang, L., *et al.* A novel hepatointestinal leukotriene B₄ receptor. *J. Biol. Chem.* **275**, 40686-40694 (2000).
6. Shimazaki, T., Kobayashi, Y., Sato, F., *et al.* Some newly synthesized leukotriene B₄ analogs inhibit LTB₄-induced lysozyme release from rat polymorphonuclear leukocytes. *Prostaglandins* **39**, 459-467 (1990).

Related Product

Leukotriene B₄ - Cat. No. 20110 • U-75302 - Cat. No. 70705 • LY255283 - Cat. No. 70715 • BLT₁ Receptor Monoclonal Antibody - Cat. No. 120111 • BLT₁ Receptor Polyclonal Antibody - Cat. No. 120114

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 under separate cover to the MSDS supervisor at 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, 11/17/2008

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