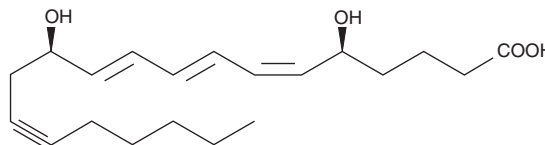


# PRODUCT INFORMATION



## 14,15-dehydro Leukotriene B<sub>4</sub> Item No. 20150

**CAS Registry No.:** 114616-11-4  
**Formal Name:** 5S,12R-dihydroxy-6Z,8E,10E-eicosatrien-14-ynoic acid  
**Synonym:** 14,15-dehydro LTB<sub>4</sub>  
**MF:** C<sub>20</sub>H<sub>30</sub>O<sub>4</sub>  
**FW:** 334.5  
**Purity:** ≥97%  
**UV/Vis.:** λ<sub>max</sub>: 270 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

14,15-dehydro LTB<sub>4</sub> 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 LTB<sub>4</sub> 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 LTB<sub>4</sub> 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 LTB<sub>4</sub> 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.

### Description

LTB<sub>4</sub> is a dihydroxy fatty acid derived from arachidonic acid through the 5-lipoxygenase pathway.<sup>1,2</sup> It promotes a number of leukocyte functions including aggregation, stimulation of ion fluxes, enhancement of lysosomal enzyme release, superoxide anion production, chemotaxis, and chemokinesis.<sup>3,4</sup> At least two LTB<sub>4</sub> receptors, termed BLT<sub>1</sub> and BLT<sub>2</sub>, have been identified. 14,15-dehydro LTB<sub>4</sub> is a LTB<sub>4</sub> receptor antagonist that has a higher binding affinity for BLT<sub>1</sub>, demonstrating a K<sub>i</sub> value of 27 nM, compared to BLT<sub>2</sub>, which has a K<sub>i</sub> value of 473 nM.<sup>4</sup> 14,15-dehydro LTB<sub>4</sub> inhibits LTB<sub>4</sub>-induced release of lysozymes from rat polymorphonuclear leukocytes with an IC<sub>50</sub> value of 1 μM.<sup>5</sup>

### References

1. Rådmark, O., Malmsten, C., Samuelsson, B., *et al. Biochem. Biophys. Res. Commun.* **92**, 954-961 (1980).
2. McGee, J. and Fitzpatrick, F. *J. Biol. Chem.* **260**, 12832-12837 (1985).
3. Ford-Hutchinson, A.W. *Crit. Rev. Immunol.* **10**, 1-12 (1990).
4. McMillan, R.M. and Foster, S.J. *Agents Actions* **24**, 114-119 (1988).
5. Wang, S., Gustafson, E., Pang, L., *et al. J. Biol. Chem.* **275**, 40686-40694 (2000).
6. Shimazaki, T., Kobayashi, Y., Sato, F., *et al. Prostaglandins* **39**, 459-467 (1990).

#### 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, 04/26/2022

#### 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