

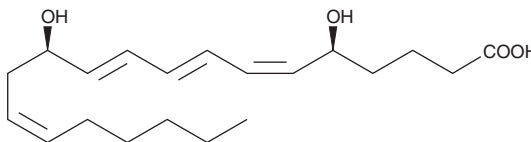
# PRODUCT INFORMATION



## Leukotriene B<sub>4</sub>

Item No. 20110

CAS Registry No.: 71160-24-2  
Formal Name: 5S,12R-dihydroxy-6Z,8E,10E,14Z-eicosatetraenoic acid  
Synonym: LTB<sub>4</sub>  
MF: C<sub>20</sub>H<sub>32</sub>O<sub>4</sub>  
FW: 336.5  
Purity: ≥95%\*  
UV/Vis.: λ<sub>max</sub>: 270 nm  
Supplied as: A solution in ethanol  
Storage: -20°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

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 or dimethyl formamide purged with an inert gas can be used. 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. If an organic solvent-free solution of LTB<sub>4</sub> is needed, the ethanol can be evaporated under a stream of nitrogen and the neat oil dissolved in the buffer of choice. 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. Also, ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. 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>1-3</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. In subnanomolar ranges (3.9 x 10<sup>-10</sup> M), LTB<sub>4</sub> causes chemotaxis and chemokinesis in human polymorphonuclear leukocytes.<sup>4</sup> At higher concentrations, (1.0 x 10<sup>-7</sup> M), LTB<sub>4</sub> leads to neutrophil aggregation and degranulation as well as superoxide anion production.<sup>4,5</sup>

### 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. Ford-Hutchinson, A.W., Bray, M.A., Doig, M.V., *et al.* Leukotriene B, a potent chemokinetic and aggregating substance released from polymorphonuclear leukocytes. *Nature* **286**, 264-265 (1980).
3. McGee, J. and Fitzpatrick, F. Enzymatic hydration of leukotriene A<sub>4</sub>. *J. Biol. Chem.* **260**, 12832-12837 (1985).
4. Ford-Hutchinson, A.W. Leukotriene B<sub>4</sub> in inflammation. *Crit. Rev. Immunol.* **10**, 1-12 (1990).
5. McMillan, R.M. and Foster, S.J. Leukotriene B<sub>4</sub> and inflammatory disease. *Agents Actions* **24**, 114-119 (1988).

\*All cysteinyl leukotrienes may contain a small amount of the 11-*trans* isomer. The purity for all such leukotrienes excludes the 1-4% *trans* isomer which may be present

#### 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, 06/15/2022

#### 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