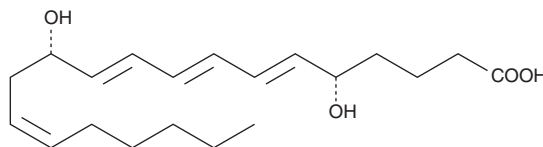


# PRODUCT INFORMATION



## 6-*trans*-12-*epi* Leukotriene B<sub>4</sub> Item No. 35265

**CAS Registry No.:** 71548-19-1  
**Formal Name:** 5S,12S-dihydroxy-6E,8E,10E,14Z-eicosatetraenoic acid  
**Synonym:** 6-*trans*-12-*epi* LTB<sub>4</sub>  
**MF:** C<sub>20</sub>H<sub>32</sub>O<sub>4</sub>  
**FW:** 336.5  
**Purity:** ≥97%  
**UV/Vis.:** λ<sub>max</sub>: 268 nm ε: 50,000 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

6-*trans*-12-*epi* Leukotriene B<sub>4</sub> (6-*trans*-12-*epi* 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. The solubility of 6-*trans*-12-*epi* LTB<sub>4</sub> 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 6-*trans*-12-*epi* LTB<sub>4</sub> is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of 6-*trans*-12-*epi* LTB<sub>4</sub> in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

6-*trans*-12-*epi* LTB<sub>4</sub> is a non-enzymatic hydrolysis product of LTA<sub>4</sub>. It was originally isolated from glycogen-induced rabbit peritoneal polymorphonuclear leukocytes (PMNL).<sup>1,2</sup> 6-*trans*-12-*epi* LTB<sub>4</sub> is weakly chemotactic for PMNL, with approximately 20 times less potency than LTB<sub>4</sub>, and has no effect on the aggregation response.<sup>3</sup>

### References

- Borgeat, P. and Samuelsson, B. Metabolism of arachidonic acid in polymorphonuclear leukocytes. *J. Biol. Chem.* **254**(16), 7865-7869 (1979).
- Borgeat, P. and Samuelsson, B. Arachidonic acid metabolism in polymorphonuclear leukocytes: Unstable intermediate in formation of dihydroxy acids. *Proc. Natl. Acad. Sci. USA* **76**(7), 3213-3217 (1979).
- Lee, T.H., Mencia-Huerta, J.M., Shih, C., *et al.* Characterization and biologic properties of 5,12-dihydroxy derivatives of eicosapentaenoic acid, including leukotriene B<sub>5</sub> and the double lipoxygenase product. *J. Biol. Chem.* **259**(4), 2383-2389 (1984).

#### 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, 01/12/2023

#### 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