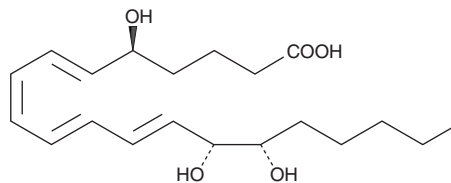


PRODUCT INFORMATION



Lipoxin B₄ Item No. 90420

CAS Registry No.: 98049-69-5
Formal Name: 5S,14R,15S-trihydroxy-6E,8Z,10E,12E-eicosatetraenoic acid
Synonym: LXB₄
MF: C₂₀H₃₂O₅
FW: 352.5
Purity: ≥95%
Supplied as: A solution in ethanol
Storage: -80°C
Stability: ≥1 year
Special Conditions: Oxygen and light sensitive



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

Laboratory Procedures

For long term storage, we suggest that lipoxin B₄ (LXB₄) be stored as supplied at -80°C. It should be stable for at least one year.

LXB₄ 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. It is recommended that this product be stored and handled in an ethanol solution. Lipoxins can isomerize and degrade when put into freeze thaw conditions and/or in solvents such as DMF or DMSO. If diluted with an aqueous buffer, this product should be discarded immediately after use.

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 LXB₄ is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of LXB₄ in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Lipoxin B₄ (LXB₄) is a positional isomer of LXA₄ (Item No. 90410) produced by the metabolism of 15-HETE (Item No. 34700) or 15(S)-HpETE (Item No. 44720) by human leukocytes.^{1,2} At a concentration of 100 nM, LXB₄ inhibits polymorphonuclear leukocyte (PMN) migration stimulated by leukotriene B₄ (LTB₄; Item No. 20110) and inhibits LTB₄-induced adhesion of PMNs with an IC₅₀ value of 0.3 nM.³

References

1. Serhan, C.N., Hamberg, M., Samuelsson, B., *et al.* On the stereochemistry and biosynthesis of lipoxin B. *Proc. Natl. Acad. Sci. USA* 83, 1983-1987 (1986).
2. Serhan, C.N., Hamberg, M., and Samuelsson, B. Lipoxins: Novel series of biologically active compounds formed from arachidonic acid in human leukocytes. *Proc. Natl. Acad. Sci. USA* 81, 5335-5339 (1984).
3. Papayianni, A., Serhan, C.N., and Brady, H.R. Lipoxin A₄ and B₄ inhibit leukotriene-stimulated interactions of human neutrophils and endothelial cells. *J. Immunol.* 156, 2264-2272 (1996).

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, 04/30/2018

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