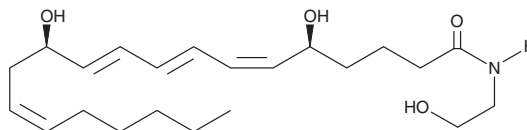


PRODUCT INFORMATION



Leukotriene B₄ Ethanolamide Item No. 20112

CAS Registry No.: 877459-63-7
Formal Name: N-(2-hydroxyethyl)-5S,12R-dihydroxy-6Z,8E,10E,14Z-eicosatetraenamide
Synonym: LTB₄-EA
MF: C₂₂H₃₇NO₄
FW: 379.5
Purity: ≥97%
UV/Vis.: λ_{max}: 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

LTB₄-EA 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 LTB₄-EA in these solvents is approximately 100 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 LTB₄-EA is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of LTB₄-EA in PBS (pH 7.2) is approximately 12 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

The effects of LTB₄ are mediated by two known receptors, BLTR₁ and BLTR₂.^{1,2} LTB₄ is a high affinity ligand for BLTR₁, and many of its pro-inflammatory effects are believed to be transduced through this receptor. BLTR₂ is more inegmatic, in that LTB₄ is not a high-affinity ligand, nor is it clear that BLTR₂ activation promotes inflammation. LTB₄-EA is a theoretical 5-lipoxygenase metabolite of the endocannabinoid anandamide. In CHO cells transfected with human BLTR₁, LTB₄-EA was a potent antagonist with about three times greater affinity for the receptor (K_i=1.22 nM versus 3.88 nM) than native LTB₄ itself. LTB₄-EA also antagonized the LTB₄-induced contractions of guinea pig lung parenchyma with an EC₅₀ of 10 nM.³ LTB₄-EA thus represents a potential endogenous anti-inflammatory compound functioning as a natural antagonist of BLTR1.

References

1. Yokomizo, T., Izumi, T., Chang, K., *et al.* A G-protein-coupled receptor for leukotriene B₄ that mediates chemotaxis. *Nature* **387(6633)**, 620-624 (1997).
2. Yokomizo, T., Kato, K., Terawaki, K., *et al.* A second leukotriene B₄ receptor, BLT2: A new therapeutic target in inflammation and immunological disorders. *J. Exp. Med.* **193(3)**, 421-431 (2000).
3. Pharmacological characterization of LTB₄ ethanolamide: Interaction with leukotriene (BLT) and vanilloid (TRPV1) receptors, (2003), 121 in 13th Annual ICRS Cannabinoid Symposium.

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, 11/15/2021

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