

PRODUCT INFORMATION



Leukotriene B₄-3-aminopropylamide

Item No. 20114

CAS Registry No.: 89596-43-0

Formal Name: N-(3-aminopropyl)-5S,12R-dihydroxy-6Z,8E,10E,14Z-eicosatetraenamide

Synonym: LTB₄-3-aminopropylamide

MF: C₂₃H₄₀N₂O₃

FW: 392.6

Purity: ≥95%

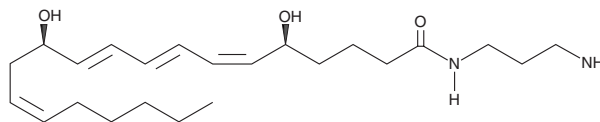
UV/Vis.: λ_{max}: 270 nm

Supplied as: A 100 µg/ml 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

Leukotriene B₄-3-aminopropylamide (LTB₄-3-aminopropylamide) 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. LTB₄-3-aminopropylamide 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 LTB₄-3-aminopropylamide is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of LTB₄-3-aminopropylamide in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

The effects of LTB₄ are mediated by two receptors, BLT₁ and BLT₂.^{1,2} LTB₄-3-aminopropylamide is an analog of LTB₄ that exhibits potent and selective binding to the BLT₁ receptor, having K_i values of 5.1 and 1,227 nM at BLT₁ and BLT₂, respectively.³

References

1. Yokomizo, T., Izumi, T., Chang, K., *et al.* A G-protein-coupled receptor for leukotriene B₄ that mediates chemotaxis. *Nature* **387**, 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**, 421-431 (2000).
3. Wang, S., Gustafson, E., Pang, L., *et al.* A novel hepatointestinal leukotriene B₄ receptor. *J. Biol. Chem.* **275**, 40686-40694 (2000).

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, 12/15/2025

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