

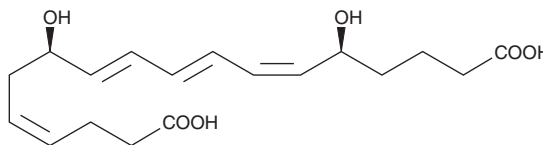
PRODUCT INFORMATION



18-carboxy dinor Leukotriene B₄

Item No. 20170

CAS Registry No.: 102674-12-4
Formal Name: 7R,14S-dihydroxy-4Z,8E,10E,12Z-octadecatetraenedioic acid
Synonym: 18-carboxy dinor LTB₄
MF: C₁₈H₂₆O₆
FW: 338.4
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

18-carboxy dinor LTB₄ 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 18-carboxy dinor LTB₄ 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 18-carboxy dinor LTB₄ is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of 18-carboxy dinor LTB₄ in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

18-carboxy dinor LTB₄ is a b-oxidation metabolite of LTB₄.¹ In the liver, LTB₄ is rapidly metabolized to 20-carboxy LTB₄, which then undergoes b-oxidation to 18-carboxy dinor LTB₄.²

References

1. Harper, T.W., Garrity, M.J., and Murphy, R.C. Metabolism of leukotriene B₄ in isolated rat hepatocytes. Identification of a novel 18-carboxy-19,20-dinor leukotriene B₄ metabolite. *J. Biol. Chem.* **261**, 5414-5418 (1986).
2. Shirley, M.A. and Murphy, R.C. Metabolism of leukotriene B₄ in isolated rat hepatocytes. Involvement of 2,4-dienoyl-coenzyme a reductase in leukotriene B₄ metabolism. *J. Biol. Chem.* **265**, 16288-16295 (1990).

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, 04/26/2022

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