

Leukotriene B4

Prostanoid receptor ligand

Activator of PPAR α . Stimulates leukocyte functions including chemokinesis, chemotaxis, lysosomal enzyme release, adhesion, stimulation of ion fluxes, superoxyde anion production and C3b receptor expression. Induces vascular permeability, NK cell activity, and bronchoconstriction. UVmax: 270 nm (50,000).

Citations: 13

[View Online »](#)

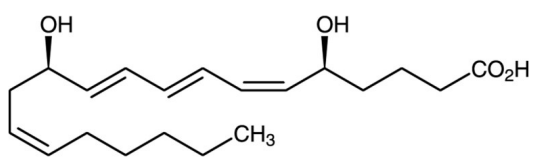
Ordering Information

[Order Online »](#)

BML-LB004-1000	1mg
BML-LB004-0050	50µg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -80°C.
Handling	Open ampule and aliquot into suitable tubes or vials. Store at -80°C or colder. For experiments with cells or tissues, aliquots of ethanol solutions may be added directly to aqueous media. For HPLC, aliquots of ethanol solutions should be evaporated to dryness under a stream of nitrogen and HPLC mobile phase added to the desired concentration.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	LTB4, 5,12-Dihydroxy-[S-[R*,S*-(E,Z,Z,E)]]-6,8,10,14-eicosatetraenoic acid
Appearance	Colorless liquid.
CAS	71160-24-2
Couple Target	Leukotriene receptor, PPAR, Prostanoid receptor
Couple Type	Activator, Ligand
Formula	C ₂₀ H ₃₂ O ₄
Formulation	Oil dissolved in ethanol.
MI	14: 5455
MW	336.5
Purity	≥97% (HPLC)
Source	Synthetic.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)