

LTB4 receptor (BLT1) polyclonal antibody

Leukotrienes are metabolic products of arachidonic acid formed by the action of 5-lipoxygenase (5-LO), with the principal products being leukotrienes C4 (LTC4) and B4 (LTB4). LTB4 is a dihydroxy-leukotriene produced in a variety of cells including eosinophils, neutrophils, platelets, and vascular cells, and is a potent leukocyte chemoattractant known to regulate leukocyte activation, cytokine secretion, inflammatory gene transcription, and IgE synthesis. As such, LTB4 has been implicated in a number of disease pathologies including atherosclerosis, asthma, psoriasis, and inflammatory bowel disease. The many actions of LTB4 are associated with two identified mammalian receptors, BLT1 and BLT2. BLT1 is a high-affinity LTB4 receptor expressed on peripheral leukocytes, while BLT2 is a more ubiquitously expressed low affinity LTB4 receptor whose function is not well understood. The BLT1 receptor is transiently expressed by antigen-primed T-cells in humans, and its expression is known to play a critical role in collagen-induced arthritis in mice.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

[View Online »](#)

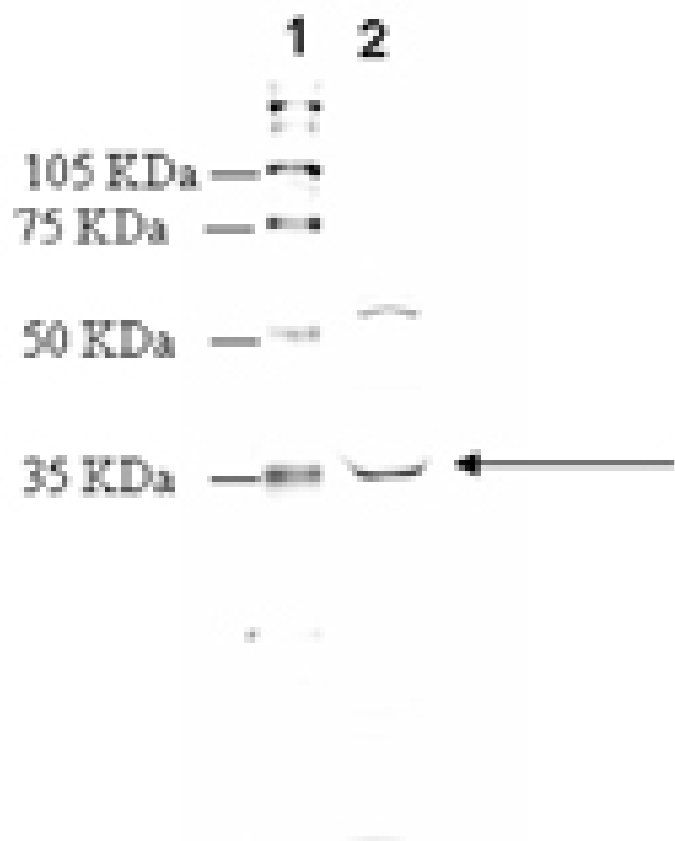
Ordering Information

[Order Online »](#)

ADI-905-793-100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis: MW marker (1) and 50µg rat heart membrane extract (2) probed with LTB4 Receptor (BLT1) pAb at 4.9µg/ml.

Handling & Storage

Handling Keep on ice at all times.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Leukotriene B4 Receptor 1, LTB4-R1

Application WB

Application Notes Predicted MW of ~38kDa. Higher MW species (~50kDa) also observed by Western blot which reflect post-translational modification of the receptor.

Crossreactivity Predicted species reactivity based on epitope sequence homology: mouse.

Formulation Liquid. In PBS containing 50% glycerol and 0.01% sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to the sequence near the N-terminus of rat LTB4 Receptor (BLT1).

Purity Detail Peptide affinity purified.

Species Reactivity Rat

UniProt ID Q9R0Q2

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#)



ENZO LIFE SCIENCES,
INC.
Last modified: May 29, 2024
Phone: +1 949 946 9120
usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info@enzolifesciences.com
be@enzolifesciences.com

France
Phone: +33 472 440 655
info@enzolifesciences.com
fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info@enzolifesciences.com
de@enzolifesciences.com

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