

LTB₄ ELISA kit

Ultrasensitive, well-validated LTB₄ ELISA kit
for inflammation research

The LTB₄ EIA kit is a colorimetric competitive enzyme immunoassay kit with results in 4 hours. Absorbance is read at 405 nm. Assay up to 39 samples in duplicate with a broad dynamic range that is suitable for a wide variety of sample matrices. Document your results to scientists world-wide with this highly cited kit.

Leukotrienes are major products of 5-lipoxygenase metabolism of arachidonic acid. Leukotriene B₄ (LTB₄) stimulates leukocyte functions including lysosomal enzyme release, adhesion, and aggregation of polymorphonuclear leukocytes. LTB₄ has been implicated as a potent mediator of inflammatory diseases and immunoregulation. The cysteinyl leukotrienes (LTC₄, LTD₄, and LTE₄) are formed by addition of cysteine derivatives to LTA₄. The conversion of LTA₄ to LTC₄ by the enzyme LTC₄ synthase limits the rate of cysteinyl leukotriene formation. Cysteinyl leukotrienes are potent lipid mediators in inflammation, inducing smooth muscle contractions and increased capillary permeability.

Citations: 53

[View Online »](#)

Ordering Information

[Order Online »](#)

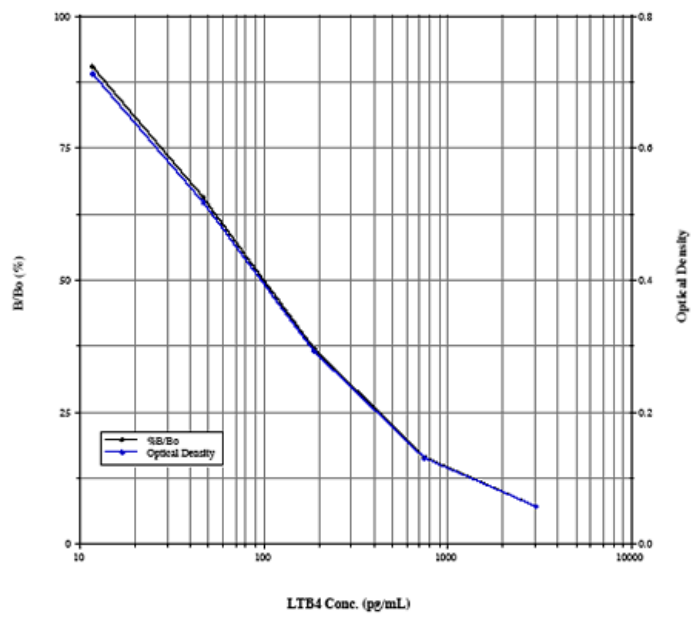
ADI-900-068

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Ultrasensitive measurement of LTB₄, detecting as little as 5.63 pg/ml
- Widely cited in numerous peer-reviewed publications
- Broad dynamic range suitable for a large variety of samples
- High throughput format with results in 4 hours for up to 39 samples in duplicate
- Highly reproducible results day-after-day and lot-after-lot
- Fully quantitative results that surpass semi-quantitative Western blot analysis



Handling & Storage

Use/Stability Store all components at +4°C, except conjugate at -20°C.

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Leukotriene B4

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of LTB₄ in culture supernatants, plasma, saliva, and urine from any species. Cited sample types include bronchioalveolar lavage fluid, cell lysate, crevicular fluid, tissue, serum and whole blood.

Assay Time 4 hours

Compatibility This product is compatible with the [Absorbance 96 Plate Reader](#).

Contents GxR IgG Microtiter plate, Conjugate, Antibody, Assay buffer, Wash buffer concentrate, Standard, pNpp Substrate, Stop solution

Sensitivity 5.63 pg/ml (range 11.7 – 3,000 pg/ml)

Species Reactivity Species independent

Wavelength 405 nm



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-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

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