

# LTB<sub>4</sub> ELISA kit

The LTB<sub>4</sub> ELISA kit is a colorimetric competitive enzyme immunoassay kit

The LTB<sub>4</sub> 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 B<sub>4</sub> (LTB<sub>4</sub>) stimulates leukocyte functions including lysosomal enzyme release, adhesion, and aggregation of polymorphonuclear leukocytes. LTB<sub>4</sub> has been implicated as a potent mediator of inflammatory diseases and immunoregulation. The cysteinyl leukotrienes (LTC<sub>4</sub>, LTD<sub>4</sub>, and LTE<sub>4</sub>) are formed by addition of cysteine derivatives to LTA<sub>4</sub>. The conversion of LTA<sub>4</sub> to LTC<sub>4</sub> by the enzyme LTC<sub>4</sub> 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: 55

[View Online »](#)

## Ordering Information

[Order Online »](#)

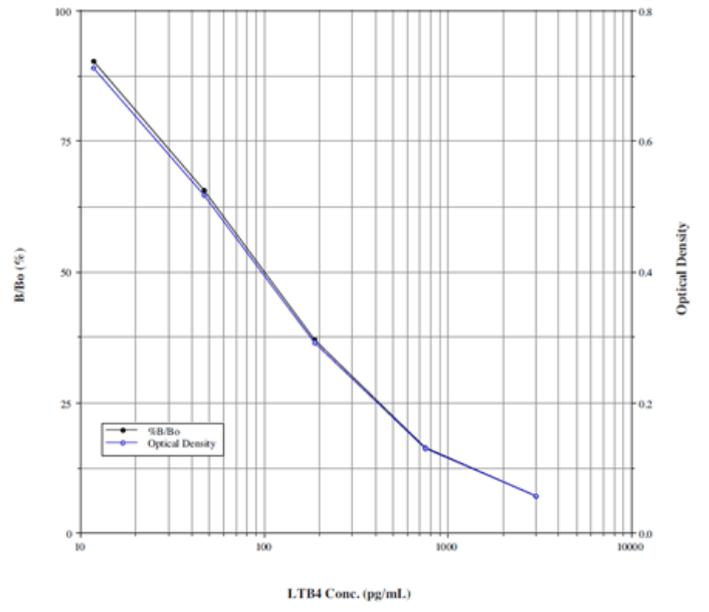
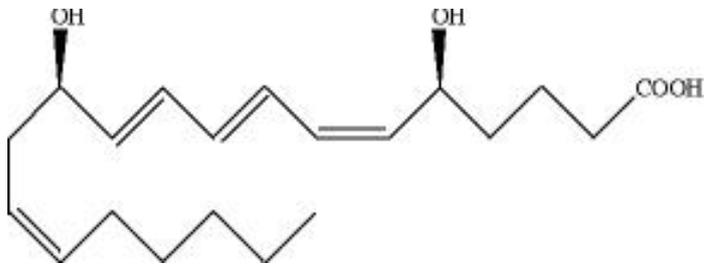
ADI-901-068

5x96 wells

## Manuals, SDS & CofA

[View Online »](#)

- Ultrasensitive measurement of LTB<sub>4</sub>, 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
- Save with bulk kit package!



## 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 LTB<sub>4</sub> 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 Microtiter plate, Conjugate, Antibody, Assay buffer concentrate, 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

**Last modified: May 29, 2024**



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)