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 $\rm B_4$ (LTB $_4$) stimulates leukocyte functions including lysosomal enzyme release, adhesion, and aggregation of polymorphonuclear leukocytes. LTB $_4$ has been implicated as a potent mediator of inflammatory diseases and immunoregulation. The cysteinyl leukotrienes (LTC $_4$, LTD $_4$, and LTE $_4$) are formed by addition of cysteine derivatives to LTA $_4$. The conversion of LTA $_4$ to LTC $_4$ by the enzyme LTC $_4$ 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: 54

View Online »

Ordering Information

Order Online »

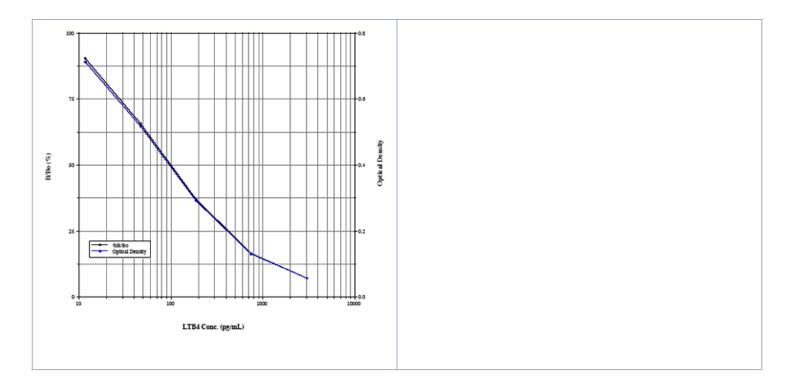
ADI-900-068

96 wells

Manuals, SDS & CofA

View Online »

- Ultrasensitive measurement of LTB₄, detecting as little as 5.63 pg/ml
- Widely cited in numerous peerreviewed 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 dayafter-day and lot-after-lot
- Fully quantitative results that surpass semi-quantitaive 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 LTB $_{\!\scriptscriptstyle A}$ 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

Last modified: May 29, 2024

