

# Cysteinyl Leukotriene ELISA kit

Highly sensitive ELISA kit enabling detection  
of LTC<sub>4</sub>/D<sub>4</sub>/E<sub>4</sub> in 4 hours.

The Cysteinyl Leukotriene ELISA kit is a colorimetric competitive enzyme immunoassay kit with results in 4 hours. Absorbance is read at 405 nm. Safe non-radioactive measurement of LTC<sub>4</sub>/D<sub>4</sub>/E<sub>4</sub>. Ready-to-use liquid color-coded reagents reduce errors.

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: 24

[View Online »](#)

## Ordering Information

[Order Online »](#)

ADI-901-070

5x96 wells

## Manuals, SDS & CofA

[View Online »](#)

- Sensitive measurement of LTC<sub>4</sub>/D<sub>4</sub>/E<sub>4</sub>, detecting as little as 26.6 pg/ml
- High throughput format with results in 4 hours for up to 38 samples in duplicate
- 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 standard and conjugate at -20°C.

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Application** Colorimetric detection, ELISA

**Application Notes** For the quantitative determination of Cysteinyl Leukotrienes (LTC<sub>4</sub>/D<sub>4</sub>/E<sub>4</sub>) in culture supernatants and urine from any species.

**Assay Time** 4 hours

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

**Contents** GxM Microtiter plate, Conjugate, Antibody, Assay buffer concentrate, Wash buffer concentrate, Standard, pNpp Substrate, Stop solution

**Sensitivity** 26.6 pg/ml (range 78.1 – 2,500 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-  
eu@enzolifesciences.com](mailto:info-eu@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)