

Cysteinyl Leukotriene ELISA kit

Highly sensitive ELISA kit enabling detection
of LTC₄/D₄/E₄ 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 LTC₄/D₄/E₄. Ready-to-use liquid color-coded reagents reduce errors.

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: 24

[View Online »](#)

Ordering Information

[Order Online »](#)

ADI-901-070

5x96 wells

Manuals, SDS & CofA

[View Online »](#)

- Sensitive measurement of LTC₄/D₄/E₄, 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 (LTC₄/D₄/E₄) 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

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

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

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

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