

# 15(S)-HETE ELISA kit

The 15(S)-HETE ELISA kit is a colorimetric competitive enzyme immunoassay kit

The 15(S)-HETE ELISA kit is a colorimetric competitive enzyme immunoassay kit with results overnight + 1 hour. Absorbance is read at 405 nm.

HETEs are byproducts generated by the metabolism of arachidonic acid by lipoxygenases. 12(S)-HETE is the stereospecific hydroxy product from the reduction of 12(S)-hydroperoxy tetraenoic eicosatetraenoic acid [12(S)-HpETE], which itself is a 12-lipoxygenase metabolite of arachidonic acid. 12(S)-HETE is chemotactic and chemokinetic for polymorphonuclear leukocytes and vascular smooth muscle cells. It also acts as a second messenger in angiotensin-II induced aldosterone production. 15-hydroxyeicosatetraenoic acid [15(S)-HETE] is the major hydroxy derivative of arachidonic acid when acted upon by 15-lipoxygenase. It is also the primary monohydroxy acid synthesized by the lipoxygenase activity of Cyclooxygenase-1 upon linoleic acid. 15(S)-HETE acts as a paracrine regulator of smooth muscle and lung neutrophil recruitment, and elevated levels of 15(S)-HETE are associated with asthma, rhinitis, chronic paranasal sinusitis and rheumatoid arthritis.

Citations: 23

[View Online »](#)

## Ordering Information

[Order Online »](#)

ADI-901-051

5x96 wells

## Manuals, SDS & CofA

[View Online »](#)

- Sensitive - very low detection limit
- Easy - few reagent preparation steps save time in the lab
- Trusted - highly cited in peer-reviewed publications
- Reliable - reproducible results day-after-day and lot-after-lot
- 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

**Alternative Name** 15(S)-hydroperoxy tetraenoic eicosatetraenoic acid

**Application** Colorimetric detection, ELISA

**Application Notes** For the quantitative determination of 15(S)-HETE in culture supernatants, plasma, serum, and urine from any species.

**Assay Time** Overnight + 1 hour

**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** 69.21 pg/ml (range 78.1 – 20,000 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-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)