

15(S)-HETE ELISA kit

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

The 15(S)-HETE EIA 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

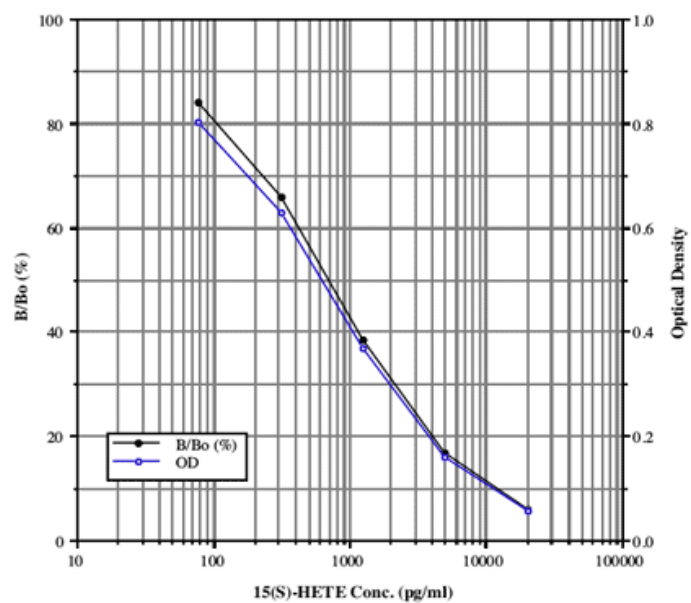
[Order Online »](#)

ADI-900-051	96 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



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 IgG Microtiter plate, Conjugate, Antibody, Assay buffer, 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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)