12(S)-HETE ELISA kit

Sensitive ELISA kit enabling detection of 12(S)-HETE in culture supernatants and plasma from any species.

The 12(S)-HETE EIA kit is a colorimetric competitive enzyme immunoassay kit with results in 5 hours. Absorbance is read at 405 nm. Fast, accurate, and reliable results are obtained with ready-to-use liquid color-coded reagents that reduce error.

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: 39

View Online »

Ordering Information

Order Online »

ADI-901-050

5x96 wells

Manuals, SDS & CofA

View Online »

- Easy-to-use color-coded reagents and pre-coated plates reduce errors and save time in the lab
- High throughput format, obtain results from up to 39 samples in duplicate in just 5 hours
- Highly sensitive assay, measure as little as 146 pg/ml
- Reliable and 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 12(S)-hydroperoxy tetraenoic eicosatetraenoic acid

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of 12(S)-HETE in culture

supernatants and plasma from any species. Cited sample types include serum, renal interstitial fluid and tissue.

Assay Time 5 hours

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 146 pg/ml (range 195 – 50,000 pg/ml)

Species Reactivity Species independent

Wavelength 405 nm

Last modified: May 29, 2024

