6-keto-PGF1α ELISA kit

Highly sensitive ELISA, detecting as low as 1.40 pg/ml in just 3 hours.

The 6-keto-PGF $_{1\alpha}$ EIA kit is a colorimetric competitive enzyme immunoassay kit with results in < 3 hours. Absorbance is read at 405 nm. This is a highly specific kit with low cross-reactivity to structurally related molecules. The 6-keto-PGF $_{1\alpha}$ EIA kit has been fully validated in-house in a wide variety of sample matrices.

6-keto-PGF $_{1\alpha}$ is a stable hydrolysis product of prostacyclin (PGI $_2$). Due to the short half-life of PGI $_2$ in buffer (2-3 minutes) the production of prostacyclin is typically monitored by measurement of the 6-keto-PGF $_{1\alpha}$ metabolite and its derivative in urine, 2,3-dinor-6-keto-PGF $_{1\alpha}$. PGI $_2$ has autocoid functions that include platelet aggregation, inhibition of vasoconstrictor responses to pressor hormones, platelet anti- and disaggregatory properties, and reproductive functions. The effects are induced by receptor-mediated elevation of cAMP. Prostacyclin has clinical applications in atherosclerosis and protection from myocardial ischemia.

Citations: 26

View Online »

Ordering Information

Order Online »

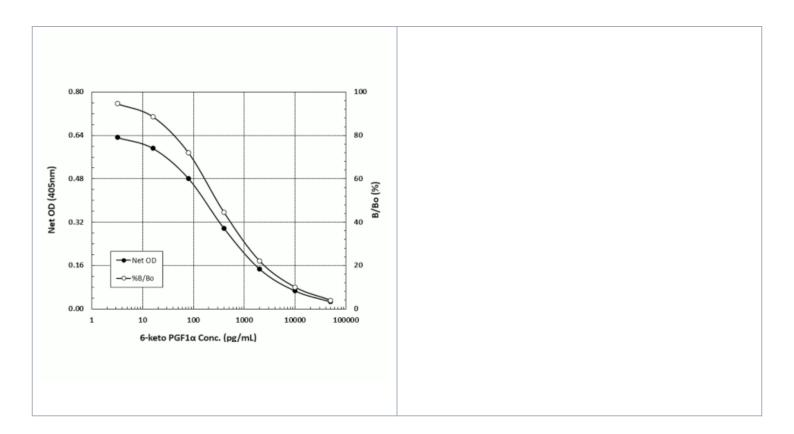
ADI-900-004

96 wells

Manuals, SDS & CofA

View Online »

- Fast results from up to 38 samples in duplicate in just 3 hours
- Highly specific with low crossreactivity with structurally related molecules
- Reliable and fully validated inhouse with a variety of matrices
- Widely published in peer-reviewed journals



Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 6-keto Prostaglandin F1α

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of 6-keto-PGF $_{1\alpha}$ in

culture supernatants, serum, saliva, and urine from any

species.

Assay Time <3 hours

Compatibility This product is compatible with the Absorbance 96 Plate

Reader.

Contents DxS IgG Microtiter plate, Conjugate, Antibody, Assay

buffer, Wash buffer concentrate, Standard, pNpp

Substrate, Stop solution

Sensitivity 1.40 pg/ml (range 3.2 – 50,000 pg/ml)

Species Reactivity Species independent

Wavelength 405 nm

Last modified: May 29, 2024

