

6-keto-PGF_{1α} ELISA kit

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

The 6-keto-PGF_{1α} 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_{1α} EIA kit has been fully validated in-house in a wide variety of sample matrices.

6-keto-PGF_{1α} is a stable hydrolysis product of prostacyclin (PGI₂). Due to the short half-life of PGI₂ in buffer (2-3 minutes) the production of prostacyclin is typically monitored by measurement of the 6-keto-PGF_{1α} metabolite and its derivative in urine, 2,3-dinor-6-keto-PGF_{1α}. PGI₂ has autocooid 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: 24

[View Online »](#)

Ordering Information

[Order Online »](#)

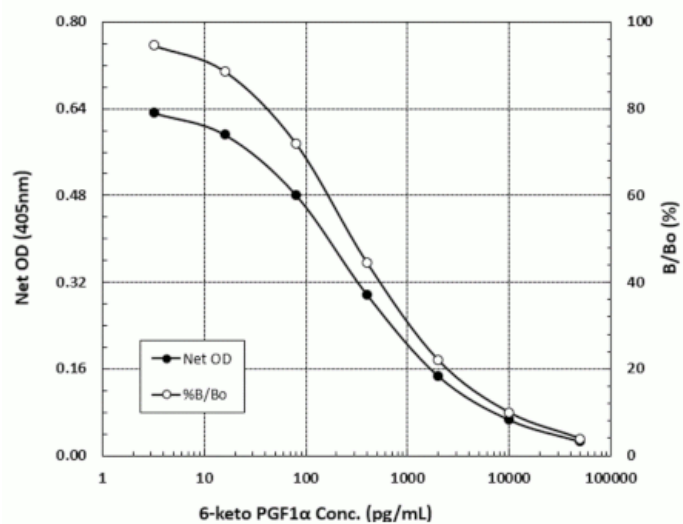
ADI-900-004

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Fast results from up to 38 samples in duplicate in just 3 hours
- Highly specific with low cross-reactivity with structurally related molecules
- Reliable and fully validated in-house 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 F1 α

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of 6-keto-PGF_{1 α} 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



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-
be@enzolifesciences.com](mailto:info-be@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)