

# Urinary prostacyclin ELISA kit

The Urinary Prostacyclin ELISA kit is a  
colorimetric competitive enzyme  
immunoassay kit

The Urinary Prostacyclin EIA kit is a colorimetric competitive enzyme immunoassay kit with results in < 3 hours. Absorbance is read at 405 nm. Quantitate  $\text{PGI}_2$  by measuring the stable metabolites in one simple assay.

Prostacyclin ( $\text{PGI}_2$ ) is involved in platelet aggregation, vasoconstriction, and reproductive functions.  $\text{PGI}_2$  has a half life of 60 minutes in plasma but only 2 to 3 minutes in buffer. The production of  $\text{PGI}_2$  is typically monitored by measurement of 6-keto- $\text{PGF}_{1\alpha}$  and 2,3-dinor-6-keto- $\text{PGF}_{1\alpha}$ . 6-keto- $\text{PGF}_{1\alpha}$  is produced by the non-enzymatic hydration of  $\text{PGI}_2$ , and is further metabolized in urine to the 2,3-dinor derivative. A number of pharmaceuticals alter and/or inhibit the synthesis of  $\text{PGI}_2$ . Methods to measure  $\text{PGI}_2$  in urine typically involve HPLC, gas chromatography/mass spectrometry, or radioimmunoassay and enzyme immunoassay for 6-keto- $\text{PGF}_{1\alpha}$ . The conversion of  $\text{PGI}_2$  into 6-keto- $\text{PGF}_{1\alpha}$  and 2,3-dinor-6-keto- $\text{PGF}_{1\alpha}$  is shown below.

Citations: 5

[View Online »](#)

## Ordering Information

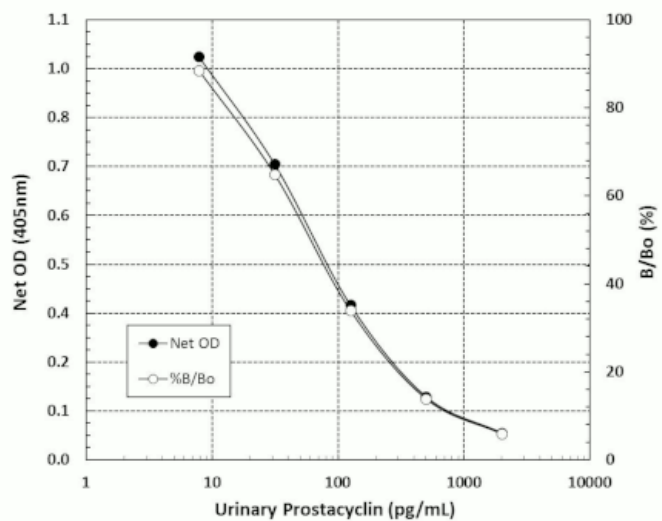
[Order Online »](#)

ADI-901-025	5x96 wells
-------------	------------

## Manuals, SDS & CofA

[View Online »](#)

- Quantitate  $\text{PGI}_2$  - measuring the stable breakdown products 6-keto- $\text{PGF}_{1\alpha}$  and 2,3-dinor-6keto- $\text{PGF}_{1\alpha}$  in one simple assay
- Less stress - much easier than GC/MS with results from up to 39 samples in duplicate in under 3 hours
- Specific - low cross-reactivity with related eicosanoids
- Save with bulk kit package!



## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name PGI2, Prostaglandin I2

Application Colorimetric detection, ELISA

**Application Notes** For the quantitative determination of urinary prostacyclin metabolites in culture supernatants and urine from any species.

Assay Time <3 hours

Compatibility This product is compatible with the [Absorbance 96 Plate Reader](#).

Contents DxS Microtiter plate, Conjugate, Antibody, Assay buffer concentrate, Wash buffer concentrate, Standard, pNpp Substrate, Stop solution

Sensitivity 6.58 pg/ml (range 7.81 – 2,000 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)