

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 PGI_2 by measuring the stable metabolites in one simple assay.

Prostacyclin (PGI_2) is involved in platelet aggregation, vasoconstriction, and reproductive functions. PGI_2 has a half life of 60 minutes in plasma but only 2 to 3 minutes in buffer. The production of 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 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 PGI_2 . Methods to measure 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 PGI_2 into 6-keto- $\text{PGF}_{1\alpha}$ and 2,3-dinor-6-keto- $\text{PGF}_{1\alpha}$ is shown below.

Citations: 8

[View Online »](#)

Ordering Information

[Order Online »](#)

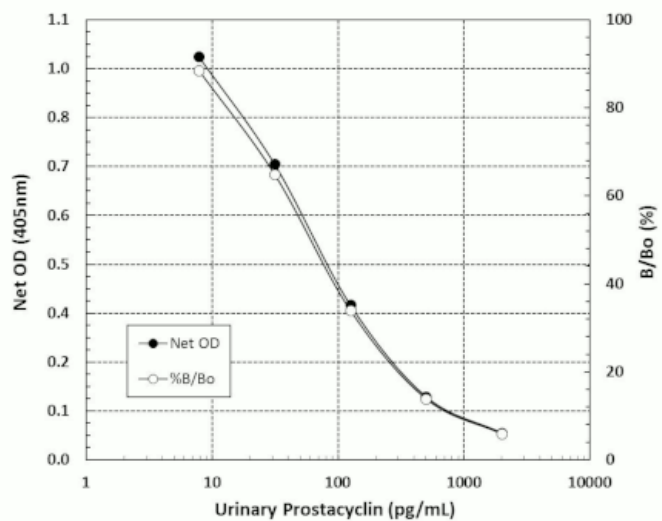
ADI-901-025

5x96 wells

Manuals, SDS & CofA

[View Online »](#)

- Quantitate PGI_2 - measuring the stable breakdown products 6-keto- $\text{PGF}_{1\alpha}$ and 2,3-dinor-6-keto- $\text{PGF}_{1\alpha}$ 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

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com