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 PGI₂ by measuring the stable metabolites in one simple assay.

Prostacyclin (PGI2) 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- $PGF_{1\alpha}$ and 2,3-dinor-6-keto- $PGF_{1\alpha}$. 6-keto- $PGF_{1\alpha}$ 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- $PGF_{1\alpha}$. The conversion of PGI_2 into 6-keto- $PGF_{1\alpha}$ and 2,3-dinor-6-keto- $PGF_{1\alpha}$ is shown below.

Citations: 7

View Online »

Ordering Information

Order Online »

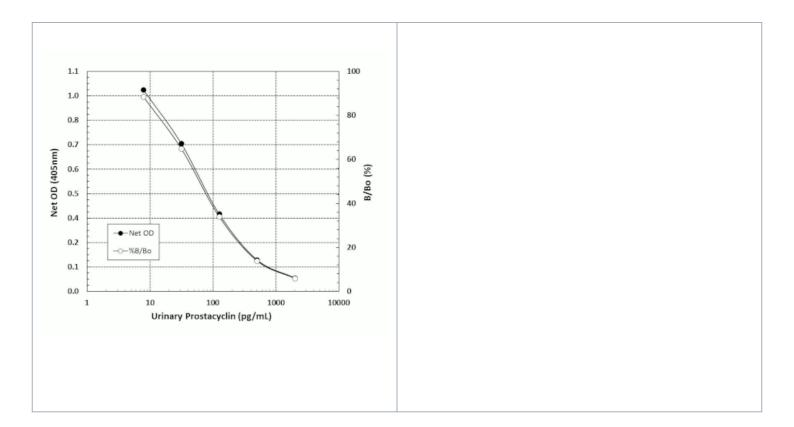
ADI-901-025

5x96 wells

Manuals, SDS & CofA

View Online »

- Quantitate PGI₂ measuring the stable breakdown products 6keto-PGF_{1α} and 2,3-dinor-6keto-PGF_{1α} 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

