## PGF2α ELISA kit

The PGF $_{2\alpha}$  ELISA kit is a colorimetric competitive enzyme immunoassay kit

The PGF $_{2\alpha}$  EIA kit is a colorimetric competitive enzyme immunoassay kit with results in < 3 hours. Absorbance is read at 405 nm. Broad dynamic range is suitable for a wide variety of sample types. Ready-to-use liquid color-coded reagents are safe, non-radioactive, and reduce laboratory errors.

Prostaglandin  $F_{2\alpha}$  (PGF $_{2\alpha}$ ) is formed in a variety of cells from PGH $_2$ , which itself is synthesized from arachidonic acid by the enzyme prostaglandin synthetase. PGF $_{2\alpha}$  is often viewed as an antagonist to PGE $_2$  due to their opposing effects on various tissues. PGF $_{2\alpha}$  is a potent bronchoconstrictor and has been implicated in asthma attacks. PGF $_{2\alpha}$  is also involved in reproductive functions including corpus luteum regulation, uterine contractions, and sperm motility. This has led to its use in terminating pregnancies and inducing labor at term. High levels of PGF $_{2\alpha}$  have also been associated with preeclampsia.

Citations: 35

View Online »

**Ordering Information** 

Order Online »

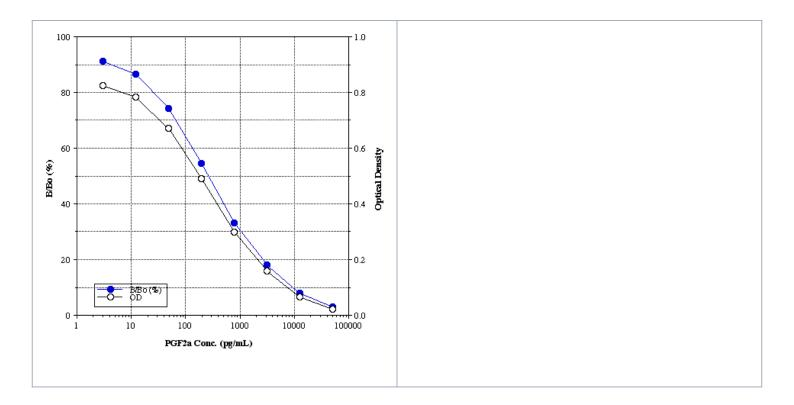
ADI-900-069

96 wells

Manuals, SDS & CofA

**View Online** »

- Broad dynamic range suitable for many sample types
- Stable liquid color-coded reagents reduce errors
- Convenient storage conditions
- Results from up to 36 samples in duplicate in less than 3 hours
- Reproducible results day-after-day and lot-after-lot



## **Handling & Storage**

Long Term Storage +4°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

Alternative Name Prostaglandin F2α

**Application** Colorimetric detection, ELISA

**Application Notes** For the quantitative determination of PGF $_{2\alpha}$  in culture

supernatants, plasma, serum, saliva, and urine from any species. Cited sample types include cell lysate and whole

blood.

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** 6.71 pg/ml (range 3.05 – 50,000 pg/ml)

Species Reactivity Species independent

Wavelength 405 nm

Last modified: May 29, 2024

