PGF_{2α} high sensitivity ELISA kit

Most sensitive PGF $_{2\alpha}$ ELISA kit available for prostaglandin research.

The PGF $_{2\alpha}$ high sensitivity EIA kit is a colorimetric competitive enzyme immunoassay kit with results overnight + 3 hours. Absorbance is read at 405 nm. Simply the most sensitive quantification fo PGF $_{2\alpha}$ commercially available in an enzyme immunoassay. Assay more samples without extraction.

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: 6

View Online »

Ordering Information

Order Online »

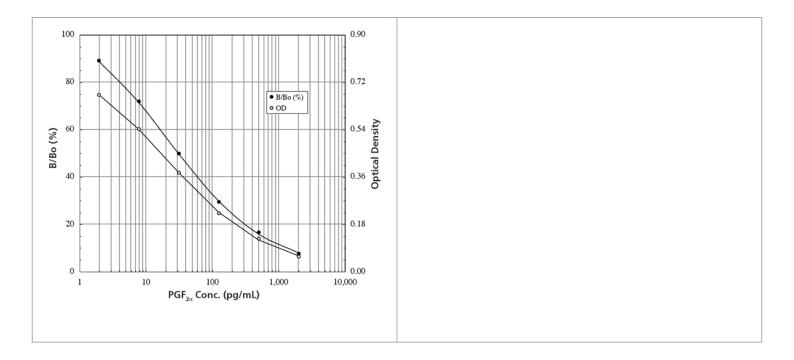
ADI-930-069

96 wells

Manuals, SDS & CofA

View Online »

- Convenient assay most samples without extraction
- Robust fully validated in multiple complex sample matrices
- Results from up to 37 samples in duplicate per 96-well plate



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, milk, plasma, serum, saliva, and urine from

any species.

Assay Time Overnight + 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 0.98 pg/ml (range 1.95 – 2,000 pg/ml)

Species Reactivity Species independent

Wavelength 405 nm

Last modified: May 29, 2024

