

# PGF<sub>2α</sub> high sensitivity ELISA kit

Most sensitive PGF<sub>2α</sub> ELISA kit available for prostaglandin research.

The PGF<sub>2α</sub> 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 for PGF<sub>2α</sub> commercially available in an enzyme immunoassay. Assay more samples without extraction.

Prostaglandin F<sub>2α</sub> (PGF<sub>2α</sub>) is formed in a variety of cells from PGH<sub>2</sub>, which itself is synthesized from arachidonic acid by the enzyme prostaglandin synthetase. PGF<sub>2α</sub> is often viewed as an antagonist to PGE<sub>2</sub> due to their opposing effects on various tissues. PGF<sub>2α</sub> is a potent bronchoconstrictor and has been implicated in asthma attacks. PGF<sub>2α</sub> 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<sub>2α</sub> have also been associated with preeclampsia.

Citations: 7

[View Online »](#)

## Ordering Information

[Order Online »](#)

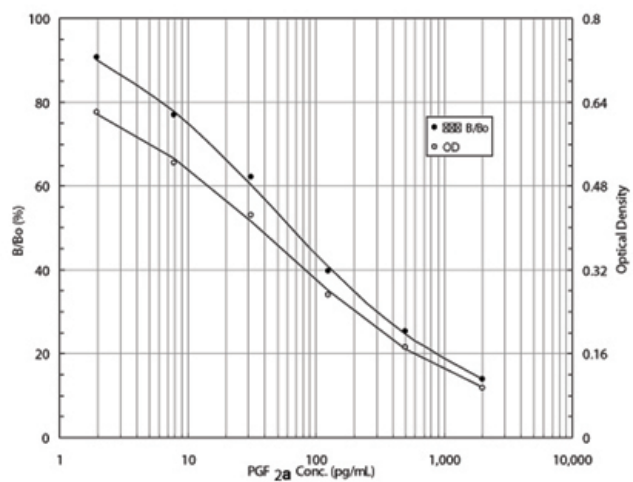
ADI-931-069

5x96 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 F2 $\alpha$

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of PGF<sub>2 $\alpha$</sub>  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 concentrate, 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



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)