

PGF_{2α} high sensitivity ELISA kit

Most sensitive PGF_{2α} ELISA kit available for prostaglandin research.

The PGF_{2α} 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_{2α} commercially available in an enzyme immunoassay. Assay more samples without extraction.

Prostaglandin F_{2α} (PGF_{2α}) is formed in a variety of cells from PGH₂, which itself is synthesized from arachidonic acid by the enzyme prostaglandin synthetase. PGF_{2α} is often viewed as an antagonist to PGE₂ due to their opposing effects on various tissues. PGF_{2α} is a potent bronchoconstrictor and has been implicated in asthma attacks. PGF_{2α} 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_{2α} have also been associated with preeclampsia.

Citations: 8

[View Online »](#)

Ordering Information

[Order Online »](#)

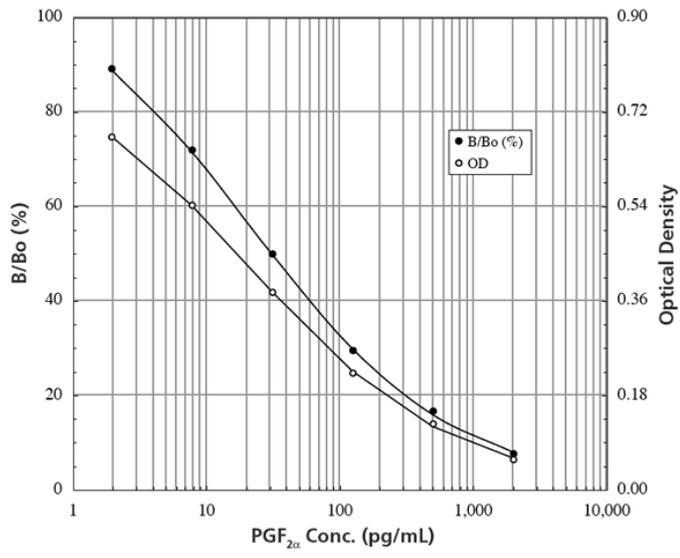
ADI-930-069

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 F2 α

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of PGF_{2 α} 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



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-fr@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