

PGF_{2α} ELISA kit

The PGF_{2α} ELISA kit is a colorimetric competitive enzyme immunoassay kit

The PGF_{2α} 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_{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: 37

[View Online »](#)

Ordering Information

[Order Online »](#)

ADI-901-069

5x96 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
- 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 Prostaglandin F2 α

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of PGF_{2 α} 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 Microtiter plate, Conjugate, Antibody, Assay buffer concentrate, 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



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-be@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