

# PGE<sub>2</sub> FPIA kit

The PGE<sub>2</sub> FPIA kit is a homogeneous mix-and-read fluorescence polarization enzyme immunoassay kit

The PGE<sub>2</sub> FPIA kit is a homogeneous mix-and-read fluorescence polarization enzyme immunoassay kit with. Results are obtained within just 30 minutes and the generated signal is stable for up to 20 hours. The PGE<sub>2</sub> FPIA kit is ideal for high throughput screening (HTS) applications. A convenient homogeneous mix-and-read assay that is amendable to miniaturization and automation, this kit has a robust Z factor >0.8.

Prostaglandin E<sub>2</sub> (PGE<sub>2</sub>) is an extensively studied prostaglandin owing to its predominance in inflammation, cancer, atherosclerosis, autoimmune disease, and sepsis. Oxidation of arachidonic acid by prostaglandin synthases (COX-1 and COX-2) produces prostaglandin H<sub>2</sub> (PGH<sub>2</sub>), which is further metabolized by PGE synthases into its major product, PGE<sub>2</sub>. PGE<sub>2</sub> mediates autocrine and paracrine signaling by binding to G-protein coupled receptors (EP1, EP2, EP3, EP4) on the cell surface, functioning to modulate phospholipase C and adenylate cyclase activity. PGE<sub>2</sub> has been of great interest as a therapeutic target, either by modulation of its synthesis by COX inhibitors (NSAIDS) or by modulation of its receptors by downregulation or binding antagonists. PGE<sub>2</sub> production in a variety of tissues has been shown to modulate numerous physiological processes including natriuresis in the kidney, smooth muscle elasticity in the vasculature, and the inflammatory response to damaged tissues by monocytes and macrophages.

Citations: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

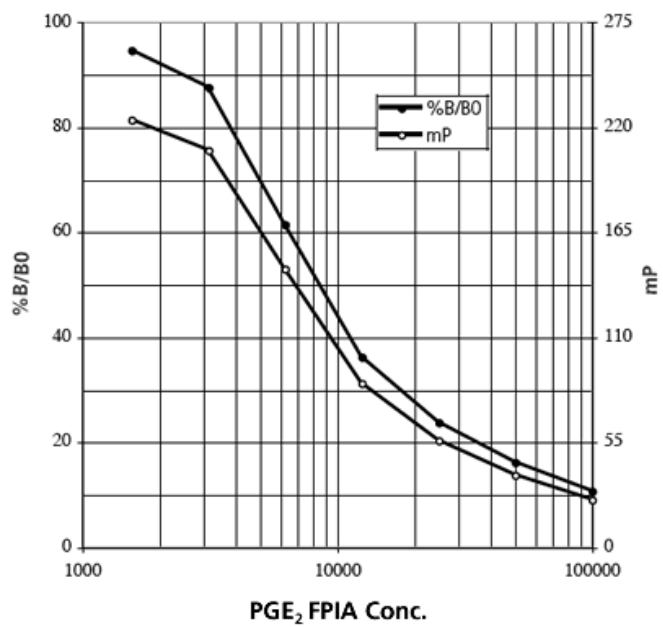
ADI-920-001

100 tests

## Manuals, SDS & CofA

[View Online »](#)

- Higher throughput - homogeneous mix-and-read assay
- Flexible - signal generated can be read after 30 minutes or several hours
- Robust - Z' Factor >0.8



## Handling & Storage

**Use/Stability** Store all components at 4°C, except standard and conjugate concentrate at -20°C.

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Prostaglandin E2

**Application** Fluorescence microscopy

**Application Notes** For the quantitative determination of PGE<sub>2</sub> in culture supernatants from any species.

**Assay Time** 30 minutes

**Compatibility** This product is compatible with the [Absorbance 96 Plate Reader](#).

**Contents** Conjugate concentrate, Antibody, Assay buffer concentrate, Standard

**Sensitivity** 684 pg/ml (range 1,562 – 100,000 pg/ml)

**Species Reactivity** Species independent



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-  
fr@enzolifesciences.com](mailto:info-fr@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)