

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

Highly sensitive PGE<sub>2</sub> ELISA kit for inflammation and eicosanoid research.

The PGE<sub>2</sub> EIA kit is a colorimetric competitive enzyme immunoassay kit with results in < 3 hours. Absorbance is read at 405 nm. Screen inhibitors of COX II activity by measuring the levels of downstream PGE<sub>2</sub>. Commercially available since 1992, this kit is widely cited in peer-reviewed publications. The non-radioactive ready-to-use liquid color-coded reagents reduce errors.

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: 186

[View Online »](#)

## Ordering Information

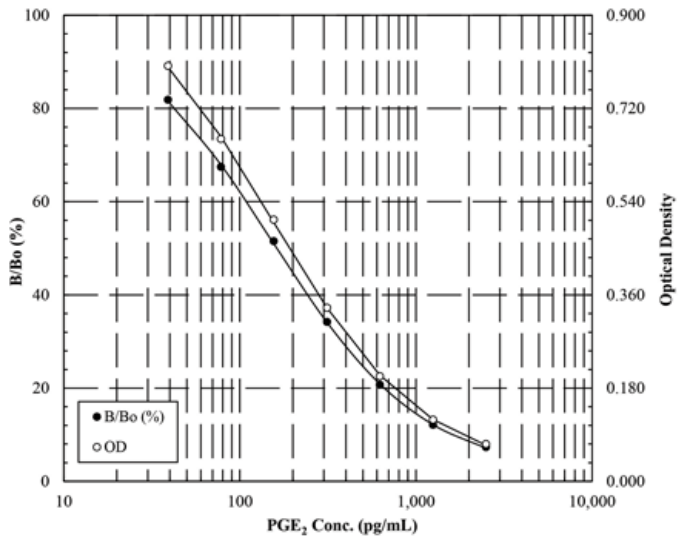
[Order Online »](#)

ADI-900-001	96 wells
-------------	----------

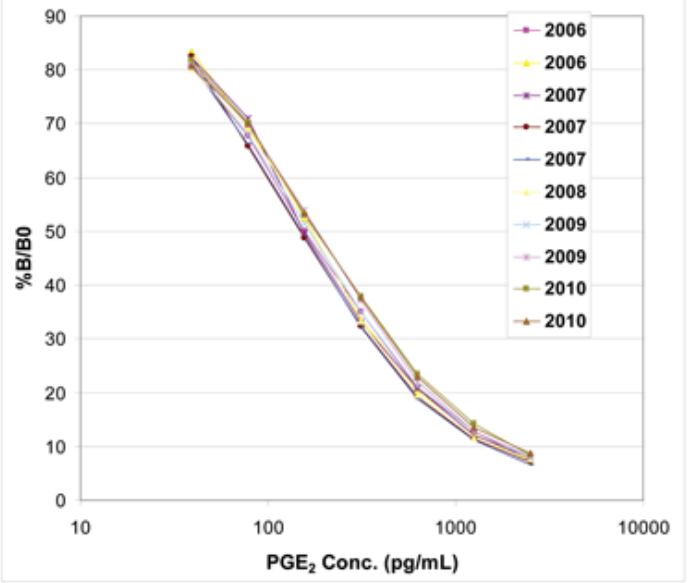
## Manuals, SDS & CofA

[View Online »](#)

- Highly sensitive measurement, detecting as little as 13.4 pg/ml PGE<sub>2</sub>
- Higher throughput format with results in <3 hrs for up to 37 samples in duplicate
- Widely cited in peer reviewed literature
- Ready-to-use liquid color-coded reagents reduce errors
- Reproducible results day-after-day and lot-to-lot



ADI-900-001 PGE<sub>2</sub> Lot-to-lot Consistency



## Handling & Storage

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

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Prostaglandin E2

**Application** Colorimetric detection, ELISA

**Application Notes** For the quantitative determination of PGE<sub>2</sub> in culture supernatants, serum, saliva, urine, and whole blood from any species. Cited sample types include cerebral spinal fluid, dialysate, gingival crevicular fluid, peritoneal exudate fluid and peritoneal exudate cell supernatant, plasma, and tissue.

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

**Contents** GxM IgG Microtiter plate, Conjugate, Antibody, Assay buffer, Wash buffer concentrate, Standard, pNpp Substrate, Stop solution

**Sensitivity** 13.4 pg/ml (range 39.1 – 2,500 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)