

PGE₂ high sensitivity ELISA kit

Ultrasensitive PGE₂ ELISA kit for inflammation and eicosanoid research.

The PGE₂ high sensitivity ELISA kit is a colorimetric competitive enzyme immunoassay kit with results overnight + 1 hour.

Prostaglandin E₂ (PGE₂) 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 H₂ (PGH₂), which is further metabolized by PGE synthases into its major product, PGE₂. PGE₂ 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. PGE₂ 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. PGE₂ 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: 60

[View Online »](#)

Ordering Information

[Order Online »](#)

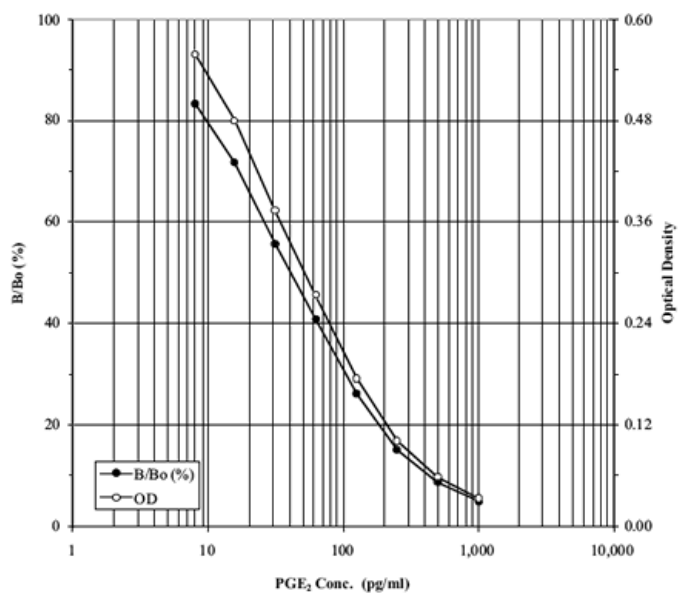
ADI-930-001

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Ultrasensitive measurement, detecting as little as 8.26 pg/ml PGE₂
- Validated in a wide variety of sample matrices
- Reliable and reproducible results lot-after-lot and day-after-day
- Fully quantitative results that surpass semi-quantitative Western blot analysis



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 PGE₂ in culture supernatants, serum, saliva, urine, and whole blood from any species. Cited sample types include plasma.

Assay Time Overnight + 1 hour

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 8.26 pg/ml (range 7.8 – 1,000 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)