PGE2 CLIA kit

The PGE₂ CLIA kit is a sensitive chemiluminescent competitive enzyme immunoassay kit

The ${\rm PGE}_2$ CLIA kit is a sensitive chemiluminescent competitive enzyme immunoassay kit for the quantitative determination of ${\rm PGE}_2$ in biological fluids from any species.

Prostaglandin E_2 (PGE $_2$) 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 (PGH $_2$), which is further metabolized by PGE synthases into its major product, PGE $_2$. PGE $_2$ 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 $_2$ 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 $_2$ 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: 9

View Online »

Ordering Information

Order Online »

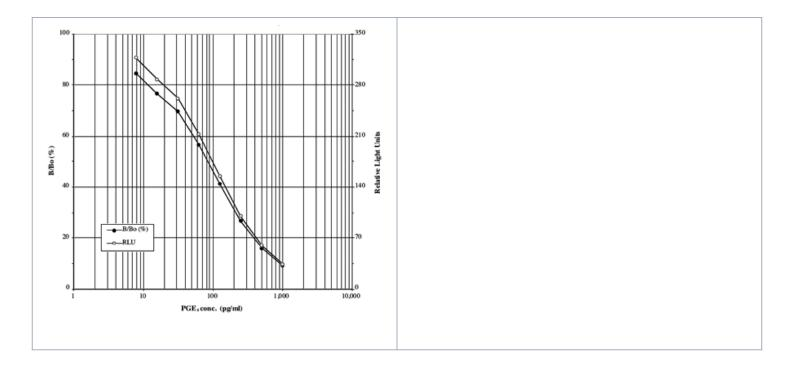
ADI-910-001

96 wells

Manuals, SDS & CofA

View Online »

- Validated in numerous sample matrices
- Results in just 3 hours
- Sensitive to 6.03 pg/ml
- Results from up to 36 samples in duplicate per plate



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 Chemiluminescence

Application Notes For the quantitative determination of PGE₂ in culture

supernatants, serum, saliva, urine, and whole blood from

any species.

Assay Time 3 hours

Compatibility This product is compatible with the Absorbance 96 Plate

Reader.

Contents White GxM IgG Microtiter plate, Conjugate, Antibody,

Assay buffer, Wash buffer concentrate, Standard, CLIA

Substrate

Sensitivity 6.03 pg/ml (range 7.81 – 1,000 pg/ml)

Species Reactivity Species independent

Last modified: May 29, 2024

