PGE₂ FPIA kit

The PGE₂ FPIA kit is a homogeneous mixand-read fluorescence polarization enzyme immunoassay kit

The PGE₂ 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 ₂ 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_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: 1

View Online »

Ordering Information

Order Online »

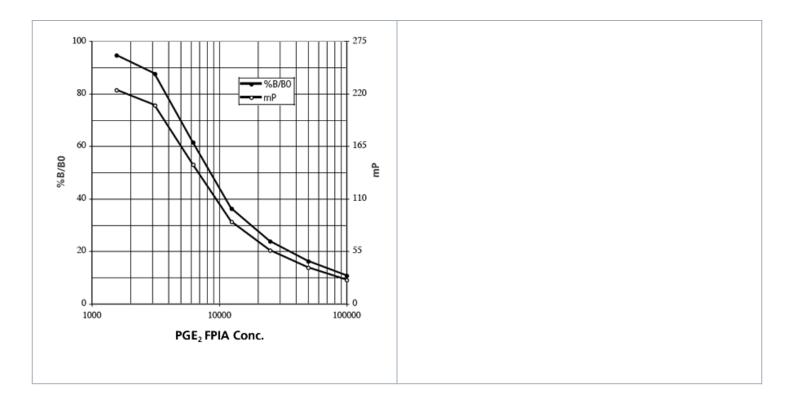
ADI-920-001

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 PGE₂ 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

Last modified: May 29, 2024

