

PGE₁ ELISA kit

The PGE₁ ELISA kit is a colorimetric competitive enzyme immunoassay kit

The PGE₁ EIA kit is a colorimetric competitive enzyme immunoassay kit with results in < 3 hours. Absorbance is read at 405 nm. Assay up to 38 samples in duplicate per kit with excellent detection sensitivity.

Prostaglandin E₁ (PGE₁) is synthesized from dihomo- γ -linolenic acid (DGLA). PGE₁ has been shown to have a number of biological actions, including vasodilation, proliferation of vascular smooth muscle cells, platelet aggregation and has been shown to have insulin-like actions. Its effects are induced by receptor mediated elevation of cAMP. PGE₁ is the major prostaglandin in semen.

Citations: 9

[View Online »](#)

Ordering Information

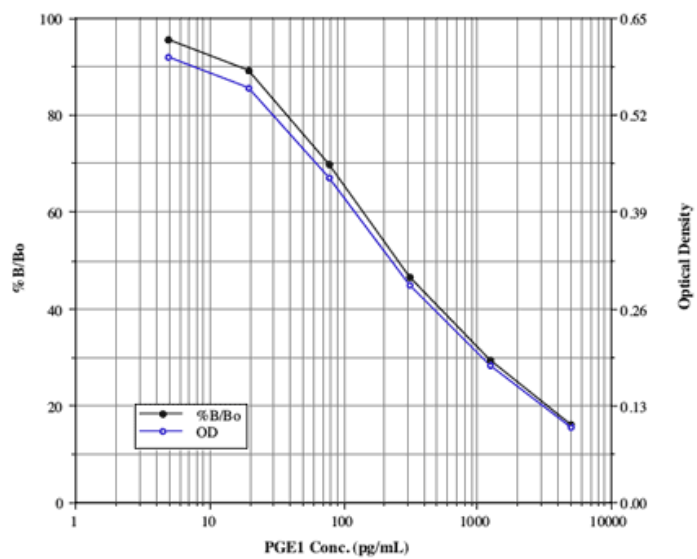
[Order Online »](#)

ADI-900-005	96 wells
-------------	----------

Manuals, SDS & CofA

[View Online »](#)

- Results in from up to 38 samples in duplicate in less than 3 hours
- Validated in multiple same matrices
- Suitable for use with any species
- Reproducible results day-after-day and lot-after-lot



Handling & Storage

Use/Stability Store all components 4°C, except conjugate -20°C.

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Prostaglandin E1

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of PGE₁ in culture supernatants, plasma, serum, saliva, and urine from any species. Cited sample types include simulated interstitial lung fluid.

Assay Time <3 hours

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

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

Sensitivity 5.58 pg/ml (range 4.88 – 5,000 pg/ml)

Species Reactivity Species independent

Wavelength 405 nm

Last modified: May 29, 2024



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)