

8-iso-PGF_{2α} ELISA kit

This is a colorimetric competitive immunoassay kit for the quantitative determination of free 8-iso-Prostaglandin F_{2α} in biological fluids such as urine and in culture supernatants.

The measurement of F₂-isoprostanes has emerged as one of the most reliable approaches to assess oxidative stress status *in vivo*, providing an important tool to explore the role of oxidative stress in the pathogenesis of human disease. This is a colorimetric competitive immunoassay kit for the quantitative determination of free 8-iso-Prostaglandin F_{2α} in biological fluids such as urine and in culture supernatants. To measure 8-iso-Prostaglandin F_{2α} levels in serum, plasma or other tissue samples, use the Direct 8-iso-PGF_{2α} EIA Kit.

8-iso-PGF_{2α} is produced *in vivo* by both non-cyclooxygenase and cyclooxygenase dependant mechanisms from arachidonic acid. It has been shown to be a potent vasoconstrictor, a potential mediator of hepatorenal syndrome and a mutagen in 3T3 cells and vascular smooth muscle cells. 8-iso-PGF_{2α} has been shown to circulate in plasma and is excreted in urine.

Prostaglandin F_{2α} (PGF_{2α}) is formed in a variety of cells from PGH₂, which itself is synthesized from arachidonic acid by the enzyme prostaglandin synthetase. PGF_{2α} is often viewed as an antagonist to PGE₂ due to their opposing effects on various tissues. PGF_{2α} is a potent bronchoconstrictor and has been implicated in asthma attacks. PGF_{2α} is also involved in reproductive functions including corpus luteum regulation, uterine contractions, and sperm motility. This has led to its use in terminating pregnancies and inducing labor at term. High levels of PGF_{2α} have also been associated with preeclampsia.

Citations: 35

[View Online »](#)

Ordering Information

[Order Online »](#)

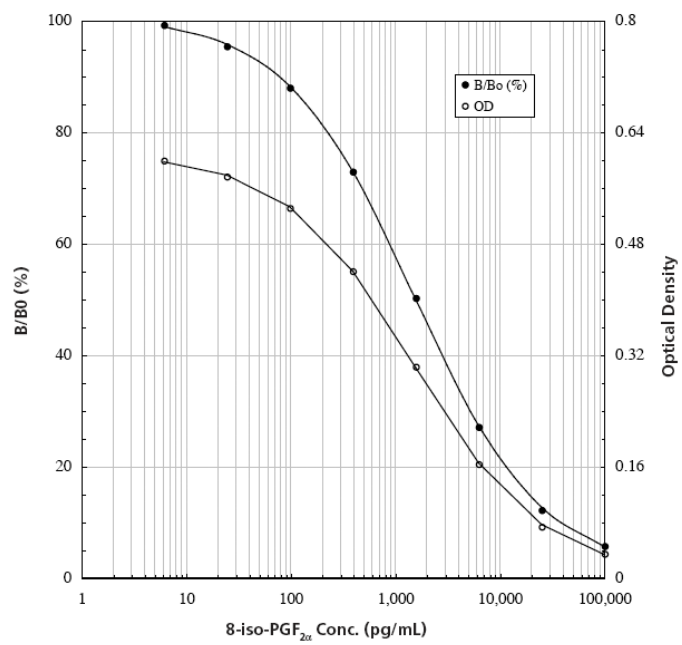
ADI-900-010

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Fast - results from up to 36 samples in duplicate in less than 3 hours
- Ready-to-use - liquid color-coded reagents and pre-coated plates reduce errors
- Trusted - published in peer-reviewed journals
- Reliable - reproducible results day-after-day and lot-after-lot



Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 8-iso-Prostaglandin F2α

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of 8-iso-PGF_{2α} in culture supernatants, tissue, and urine from any species. Cited sample types include plasma.

Assay Time <3 hours

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

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

Sensitivity 16.3 pg/ml (range 6.1 – 100,000 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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com