

# 15-deoxy- $\Delta^{12,14}$ -PGJ<sub>2</sub> ELISA kit

**First-to-market 15-d-PGJ<sub>2</sub> ELISA  
commercially available for eicosanoid  
research.**

The 15-deoxy- $\Delta^{12,14}$ -PGJ<sub>2</sub> ELISA kit is a colorimetric competitive enzyme immunoassay kit with results in 5 hours. Absorbance is read at 405 nm. This kit saves time and is much less time consuming than GC/MS analysis. Ready-to-use liquid color-coded reagents reduce laboratory error.

15-deoxy- $\Delta^{12,14}$ -Prostaglandin J<sub>2</sub> (15-d-PGJ<sub>2</sub>) is one of the ultimate dehydration products of PGD<sub>2</sub>, formed via the intermediate  $\Delta^{12}$ -PGJ<sub>2</sub>. 15-d-PGJ<sub>2</sub> is physiologically present in body fluids at concentration of 10<sup>-12</sup> to 10<sup>-9</sup> M, but increases dramatically under stress such as infection and inflammation. 15-d-PGJ<sub>2</sub> exerts both pro- and anti-inflammatory effects in a cell-type and concentration-dependent manner, with many of these effects mediated via binding to the  $\gamma$  isoform of the peroxisome proliferators-activated receptor (PPAR- $\gamma$ ). 15-d-PGJ<sub>2</sub> appears to induce HO-1 expression by a PPAR- $\gamma$ -independent mechanism mediated by the generation of reactive oxygen species (ROS).

Citations: 32

[View Online »](#)

## Ordering Information

[Order Online »](#)

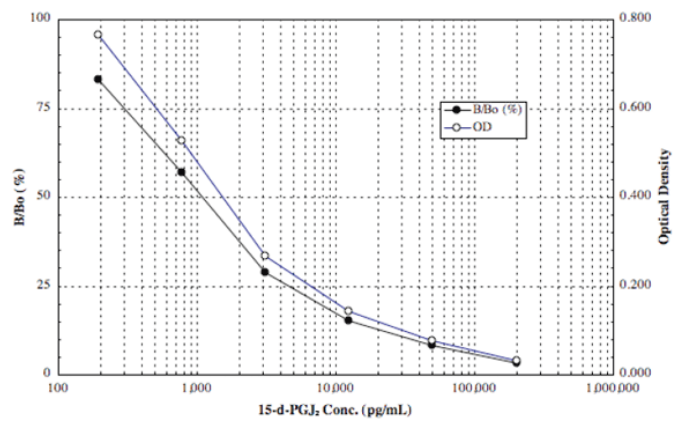
ADI-900-023

96 wells

## Manuals, SDS & CofA

[View Online »](#)

- High sensitivity, measuring as low as 36.8 pg/ml
- Fast results, much less time consuming than GC/MS analysis
- Simple and ready-to-use liquid color-coded reagents reduce error
- 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** 15-deoxy- $\Delta^{12,14}$ -Prostaglandin J2, 15-d-PGJ2

**Application** Colorimetric detection, ELISA

**Application Notes** For the quantitative determination of 15-deoxy- $\Delta^{12,14}$ -PGJ<sub>2</sub> in culture supernatants, plasma, saliva, and urine from any species.

**Assay Time** 5 hours

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

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

**Sensitivity** 36.8 pg/ml (range 195 – 200,000 pg/ml)

**Species Reactivity** Species independent

**Wavelength** 405 nm

**Last modified: August 23, 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-be@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-fr@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)