

15-deoxy- $\Delta^{12,14}$ -PGJ₂ ELISA kit

**First-to-market 15-d-PGJ₂ ELISA
commercially available for eicosanoid
research.**

The 15-deoxy- $\Delta^{12,14}$ -PGJ₂ EIA 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 J₂ (15-d-PGJ₂) is one of the ultimate dehydration products of PGD₂, formed via the intermediate Δ^{12} -PGJ₂. 15-d-PGJ₂ is physiologically present in body fluids at concentration of 10⁻¹² to 10⁻⁹ M, but increases dramatically under stress such as infection and inflammation. 15-d-PGJ₂ 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 γ isoform of the peroxisome proliferators-activated receptor (PPAR- γ). 15-d-PGJ₂ appears to induce HO-1 expression by a PPAR- γ -independent mechanism mediated by the generation of reactive oxygen species (ROS).

Citations: 32

[View Online »](#)

Ordering Information

[Order Online »](#)

ADI-901-023

5x96 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
- Save with bulk kit package!

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- $\Delta^{12,14}$ -PGJ₂ 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 Microtiter plate, Conjugate, Antibody, Assay buffer concentrate, 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 17, 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)