

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

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).

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

ADI-915-043-100	100 $\mu$ l
-----------------	-------------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** 15-deoxy- $\Delta$ 12,14-Prostaglandin J2

**Application** ELISA

**Formulation** Liquid. In PBS containing 1.0mg/ml Rabbit IgG and 0.1% sodium azide.

**Host** Rabbit

**Immunogen** 15-deoxy- $\Delta$ 12,14-PGJ2 conjugated to a protein carrier molecule.

**Purity Detail** Sodium sulfate precipitation purified.

**Source** Purified from rabbit serum.

**Species Reactivity** Species independent

**Worry-free Guarantee** This antibody is covered by our [Worry-Free Guarantee](#).



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-  
fr@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)