

15-deoxy- $\Delta^{12,14}$ -PGJ₂ polyclonal antibody

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).

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

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ADI-915-043-100	100 μ 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- Δ 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- Δ 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](#)

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