15-deoxy-Δ12,14-PGJ2 ELISA kit

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

The 15-deoxy- $\Delta^{12,14}$ -PGJ $_2$ 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 J_2 (15-d-PG J_2) is one of the ultimate dehydration products of PGD $_2$, formed via the intermediate Δ^{12} -PG J_2 . 15-d-PG J_2 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-PG J_2 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-PG J_2 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-Δ12,14-Prostaglandin J2, 15-d-PGJ2

Application Colorimetric detection, ELISA

For the quantitative determination of 15-deoxy-Δ^{12,14}-PGJ **Application Notes**

 $_{
m 2}$ 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

info-



eu@enzolifesciences.com