



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

Catalog No. ADI-901-023

480 Well (5 by 96 Well) Kit

Table of Contents

Description	Page	2
Introduction		2
Precautions		2
Materials Supplied		3
Storage		3
Materials Needed but Not Supplied		3
Sample Handling		4
Procedural Notes		5
Reagent Preparation		5
Assay Procedure		6
Calculation of Results		7
Typical Results		7
Typical Standard Curve		8
Typical Quality Control Parameters		8
Performance Characteristics		9
Sample Dilution Recommendations		11
References		11
Limited Warranty		12

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.

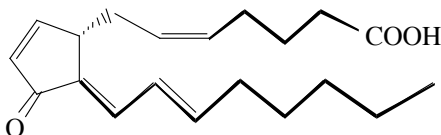
Description

The 15-deoxy- $\Delta^{12,14}$ -PGJ₂ ELISA kit is a competitive immunoassay for the quantitative determination of 15-deoxy- $\Delta^{12,14}$ -Prostaglandin J₂ (15-d-PGJ₂) in biological fluids. Please read the complete kit insert before performing this assay. The kit uses a polyclonal antibody to 15-d-PGJ₂ to bind, in a competitive manner, the 15-d-PGJ₂ in the sample or an alkaline phosphatase molecule which has 15-d-PGJ₂ covalently attached to it. After a simultaneous incubation the excess reagents are washed away and substrate is added. After a short incubation time the enzyme reaction is stopped and the yellow color generated read on a microplate reader at 405nm. The intensity of the bound yellow color is inversely proportional to the concentration of 15-d-PGJ₂ in either standards or samples. The measured optical density is used to calculate the concentration of 15-d-PGJ₂. For further explanation of the principles and practice of immunoassays please see the excellent books by Chard¹³ or Tijssen¹⁴

Introduction

15-deoxy- $\Delta^{12,14}$ -Prostaglandin J₂ is one of the ultimate dehydration products of PGD₂. PGD₂ is formed from PGH₂ which itself is synthesized from arachidonic acid by the enzyme prostaglandin synthase¹⁻⁴. In the aqueous solution PGD₂ forms PGJ₂⁵. Albumin or other serum proteins present in the sample will convert PGD₂ to the isomeric compound, Δ^{12} -PGJ₂⁶⁻⁸. In males about 152 ng of Δ^{12} -PGJ₂ is excreted in urine per day, and in women about half this amount is excreted⁹. Excellent recent reviews^{10,11} of the biological actions of these compounds have been published. 15-d-PGJ₂ has been shown as an inducer of adipogenesis and an activator of the γ isoform of the peroxisome proliferated receptor (PPAR γ)¹².

15-deoxy- $\Delta^{12,14}$ -Prostaglandin J₂



Precautions

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.

1. Some kit components contain azide, which may react with lead or copper plumbing. When disposing of reagents always flush with large volumes of water to prevent azide build-up.
2. Stop solution is a solution of trisodium phosphate. This solution is caustic; care should be taken in use.
3. The activity of the alkaline phosphatase conjugate is dependent on the presence of Mg²⁺ and Zn²⁺ ions. The activity of the conjugate is affected by concentrations of chelators (>10 mM) such as EDTA and EGTA.
4. We test this kit's performance with a variety of samples, however it is possible that high levels of interfering substances may cause variation in assay results.
5. The 15-deoxy- $\Delta^{12,14}$ -PGJ₂ standard provided, Catalog No. 80-0244, is supplied in ethanolic buffer at a pH optimized to maintain 15-d-PGJ₂ integrity. Care should be taken handling this material because of the known and unknown effects of prostaglandins.

Materials Supplied

- 1. Goat anti-Rabbit IgG Microtiter Plate, Five Plates of 96 Wells, Catalog No. 80-0060**
Plates using break-apart strips coated with goat antibody specific to rabbit IgG.
- 2. 15-d-PGJ₂ ELISA Conjugate, 25 mL, Catalog No. 80-0245**
A blue solution of alkaline phosphatase conjugated with 15-d-PGJ₂.
- 3. 15-d-PGJ₂ ELISA Antibody, 25 mL, Catalog No. 80-0246**
A yellow solution of a polyclonal antibody to 15-d-PGJ₂.
- 4. Assay Buffer Concentrate, 27 mL, Catalog No. 80-1545**
Tris buffered saline, containing proteins and sodium azide as preservative.
- 5. Wash Buffer Concentrate, 100 mL, Catalog No. 80-1287**
Tris buffered saline containing detergents.
- 6. 15-deoxy- $\Delta^{12,14}$ -Prostaglandin J₂ Standard, 3 x 0.5 mL, Catalog No. 80-0244**
A solution of 1,000,000 pg/mL 15-d-PGJ₂.
- 7. pNpp Substrate, 100 mL, Catalog No. 80-0076**
A solution of p-nitrophenyl phosphate in buffer. Ready to use.
- 8. Stop Solution, 27 mL, Catalog No. 80-0248**
A solution of trisodium phosphate in water. Keep tightly capped. Caution: **Caustic**.
- 9. 15-d-PGJ₂ Assay Layout Sheet, 1 each, Catalog No. 30-0062**
- 10. Plate Sealer, 10 each, Catalog No. 30-0012**

Storage

All components of this kit, **except the conjugate and standard**, are stable at 4°C until the kit's expiration date. The conjugate and standard **must** be stored frozen at -20°C.

Materials Needed but Not Supplied

1. Deionized or distilled water.
2. Precision pipets for volumes between 5 μ L and 1,000 μ L.
3. Repeater pipets for dispensing 50 μ L and 200 μ L.
4. Disposable beaker for diluting buffer concentrates.
5. Graduated cylinders.
6. A Microplate shaker.
7. Adsorbent paper for blotting.
8. A 37°C Incubator.
9. Microplate reader capable of reading at 405 nm, preferably with correction between 570 and 590 nm.

Sample Handling

The 15-d-PGJ₂ ELISA kit is compatible with 15-d-PGJ₂ samples in a wide range of matrices. Samples diluted sufficiently into Assay Buffer can be read directly from the standard curve. Please refer to the Sample Recovery recommendations on page 11 for details of suggested dilutions. However, the end user **must verify** that the recommended dilutions are appropriate for their samples. **Samples containing rabbit IgG may interfere with the assay.** Some samples normally have very low levels of 15-d-PGJ₂ present and extraction may be necessary for accurate measurement. A suitable extraction procedure is outlined below:

Materials Needed

1. 15-d-PGJ₂ Standard to allow extraction efficiency to be accurately determined.
2. 2M hydrochloric acid, deionized water, ethanol, hexane and ethyl acetate.
3. 200 mg C₁₈ Reverse Phase Extraction Columns.

Procedure

1. Acidify the plasma, urine or tissue homogenate by addition of 2M HCl to pH of 3.5. Approximately 50 µL of HCl will be needed per mL of plasma. Allow to sit at 4°C for 15 minutes. Centrifuge samples in a microcentrifuge for 2 minutes to remove any precipitate.
2. Prepare the C₁₈ reverse phase column by washing with 10 mL of ethanol followed by 10 mL of deionized water.
3. Apply the sample under a slight positive pressure to obtain a flow rate of about 0.5 mL/minute. Wash the column with 10 mL of water, followed by 10 mL of 15% ethanol, and finally 10 mL hexane. Elute the sample from the column by addition of 10 mL ethyl acetate.
4. If analysis is to be carried out immediately, evaporate samples under a stream of nitrogen. Add at least 250 µL of Assay Buffer to the dried samples. Vortex well then allow to sit 5 minutes at room temperature. Repeat twice more. If analysis is to be delayed, store samples as the eluted ethyl acetate solutions at -80°C until the immunoassay is to be run. Evaporate the organic solvent under a stream of nitrogen prior to running assay and reconstitute as above.

Please refer to references 15-18 for details of extraction protocols.

Procedural Notes

1. Do not mix components from different kit lots or use reagents beyond the kit expiration date.
2. Allow all reagents to warm to room temperature for at least 30 minutes before opening.
3. Standards can be made up in either glass or plastic tubes.
4. Pre-rinse the pipet tip with reagent, use fresh pipet tips for each sample, standard and reagent.
5. Pipet standards and samples to the bottom of the wells.
6. Add the reagents to the side of the well to avoid contamination.
7. This kit uses break-apart microtiter strips, which allow the user to measure as many samples as desired. Unused wells must be kept desiccated at 4°C in the sealed bag provided. The wells should be used in the frame provided.

8. **Care must be taken to minimize contamination by endogenous alkaline phosphatase.** Contaminating alkaline phosphatase activity, especially in the substrate solution, may lead to high blanks. Care should be taken not to touch pipet tips and other items that are used in the assay with bare hands.
9. **Prior to the addition of substrate, ensure that there is no residual wash buffer in the wells. Any remaining wash buffer may cause variation in the assay results.**

Reagent Preparation

1. Assay Buffer

Just before use, prepare the Assay Buffer by diluting 10 mL of the supplied concentrate with 90 mL of deionized water. Discard unused assay buffer or add up to 0.09% sodium azide (w/v) for storage.

2. 15-d-PGJ₂ Standard

Allow the 1,000,000 pg/mL 15-d-PGJ₂ standard solution to warm to room temperature. Label six 12 x 75 mm glass tubes #1 through #6. Pipet 800 µL of standard diluent (Assay Buffer or Tissue Culture Media) into tube #1. Pipet 750 µL of standard diluent into tubes #2 through #6. Add 200 µL of the 1,000,000 pg/mL standard to tube #1. Vortex thoroughly. Add 250 µL of tube #1 to tube #2 and vortex thoroughly. Continue this for tubes #2 through #6.

The concentration of 15-d-PGJ₂ in tubes #1 through #6 will be 200,000, 50,000, 12,500, 3,125, 781 and 195 pg/mL respectively. See 15-d-PGJ₂ Assay Layout Sheet for dilution details.

Diluted Standards should be used within 60 minutes of preparation.

3. 15-d-PGJ₂ Conjugate

Allow the conjugate to warm to room temperature. Any unused conjugate should be aliquoted and re-frozen at or below -20°C. Avoid repeated freeze/thaws of the aliquots.

4. Conjugate 1:10 Dilution for Total Activity Measurement

Prepare the Conjugate 1:10 Dilution by diluting 50 µL of the supplied Conjugate with 450 µL of Assay Buffer. The dilution should be used within 3 hours of preparation. **This 1:10 dilution is intended for use in the Total Activity wells ONLY.**

5. Wash Buffer

Prepare the Wash Buffer by diluting 5 mL of the supplied concentrate with 95 mL of deionized water. This can be stored at room temperature until the kit expiration date, or for 3 months, whichever is earlier.

Assay Procedure

Bring all reagents to room temperature for at least 30 minutes prior to opening.

All standards and samples should be run in duplicate.

1. Refer to the Assay Layout Sheet to determine the number of wells to be used and put any remaining wells with the desiccant back into the pouch and seal the ziploc. Store unused wells at 4°C.
2. Pipet 100 µL of standard diluent (Assay Buffer or Tissue Culture Media) into the NSB and the Bo (0 pg/mL Standard) wells.
3. Pipet 100 µL of Standards #1 through #6 into the appropriate wells.
4. Pipet 100 µL of the Samples into the appropriate wells.
5. Pipet 50 µL of Assay Buffer into the NSB wells.
6. Pipet 50 µL of blue Conjugate into each well, except the Total Activity (TA) and Blank wells.
7. Pipet 50 µL of yellow Antibody into each well, except the Blank, TA and NSB wells.

NOTE: Every well used should be **Green** in color except the NSB wells which should be **Blue**. The Blank and TA wells are empty at this point and have no color.

8. Incubate the plate at room temperature on a plate shaker for 2 hours at ~500 rpm. The plate may be covered with the plate sealer provided, if so desired.
9. Empty the contents of the wells and wash by adding 400 µL of wash solution to every well. Repeat the wash 2 more times for a total of **3 Washes**.
10. After the final wash empty or aspirate the wells and firmly tap the plate dry on a lint free paper towel to remove any remaining wash buffer.
11. Add 5 µL of the light blue Conjugate 1:10 Dilution (**see Step 4, Reagent Preparation, on Page 5**) to the TA wells.
12. Add 200 µL of the pNpp Substrate solution to every well. Cover plate with the plate sealer provided and incubate at 37°C for 3 hours.
13. Add 50 µL of Stop Solution to every well. This stops the reaction and the plate should be read immediately.
14. Blank the plate reader against the Blank wells, read the optical density at 405 nm, preferably with correction between 570 and 590 nm. If the plate reader is not able to be blanked against the Blank wells, manually subtract the mean optical density of the Blank wells from all readings.

Calculation of Results

Several options are available for the calculation of the concentration of 15-d-PGJ₂ in the samples. We recommend that the data be handled by an immunoassay software package utilizing a 4 parameter logistic curve fitting program. If this sort of data reduction software is not readily available, the concentration of 15-d-PGJ₂ can be calculated as follows:

1. Calculate the average net Optical Density (OD) bound for each standard and sample by subtracting the average NSB OD from the average OD bound:

$$\text{Average Net OD} = \text{Average Bound OD} - \text{Average NSB OD}$$

2. Calculate the binding of each pair of standard wells as a percentage of the maximum binding wells (Bo), using the following formula:

$$\text{Percent Bound} = \frac{\text{Net OD}}{\text{Net Bo OD}} \times 100$$

3. Plot Percent Bound versus Concentration of 15-d-PGJ₂ for the standards. Approximate a straight line through the points. The concentration of 15-d-PGJ₂ in the unknowns can be determined by interpolation.

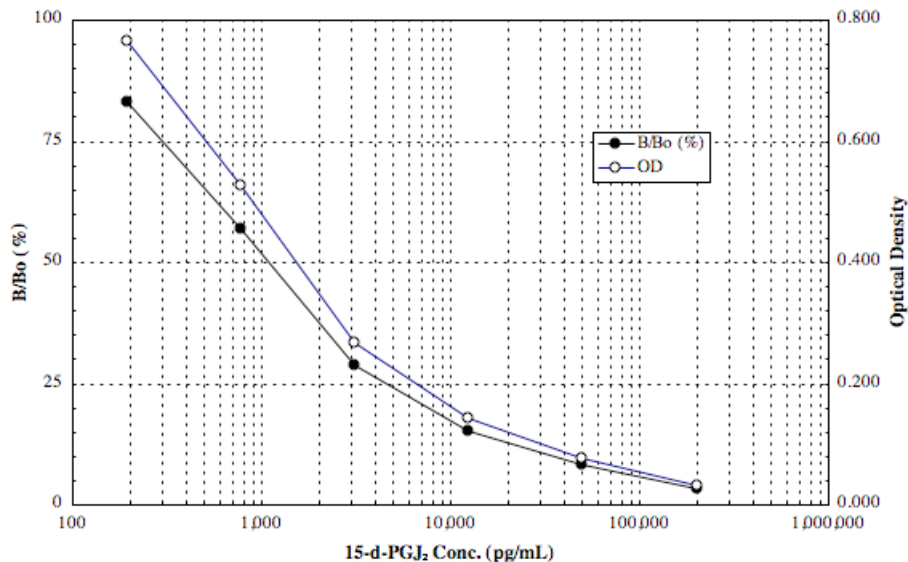
Typical Results

The results shown below are for illustration only and **should not** be used to calculate results from another assay.

<u>Sample</u>	<u>Mean OD (-Blank)</u>	<u>Average Net OD</u>	<u>Percent Bound</u>	<u>15-d-PGJ₂ - (pg/mL)</u>
Blank OD	(0.134)			
TA	2.696			
NSB	0.009	0.000	0.00%	
Bo	0.932	0.923	100%	0
S1	0.034	0.031	3.4%	200,000
S2	0.086	0.077	8.3%	50,000
S3	0.151	0.142	15.4%	12,500
S4	0.276	0.267	28.9%	3,125
S5	0.535	0.526	57.0%	781
S6	0.775	0.766	83.0%	195
Unknown 1	0.624	0.615	66.6%	486
Unknown 2	0.102	0.093	10.1%	23,638

Typical Standard Curve

A typical standard curve is shown below. This curve **must not** be used to calculate 15-d-PGJ₂ concentrations; each user must run a standard curve for each assay.



Typical Quality Control Parameters

Total Activity Added	=	2.696 x 10 x 10 = 269.6
%NSB	=	0.0%
%Bo/TA	=	0.34%
Quality of Fit	=	0.9997 (Calculated from 4 parameter logistic curve fit)

20% Intercept	=	7,636 pg/mL
50% Intercept	=	1,104 pg/mL
80% Intercept	=	247 pg/mL

Performance Characteristics

The following parameters for this kit were determined using the guidelines listed in the National Committee for Clinical Laboratory Standards (NCCLS) Evaluation Protocols¹⁹.

Sensitivity

Sensitivity was calculated by determining the average optical density bound for sixteen (16) wells run as Bo, and comparing to the average optical density for sixteen (16) wells run with Standard #6. The detection limit was determined as the concentration of 15-d-PGJ₂ measured at two (2) standard deviations from the zero along the standard curve.

Average Optical Density for the Bo = 1.300 ± 0.030 (2.3%)

Average Optical Density for Standard #6 = 0.983 ± 0.050 (5.1%)

Delta Optical Density (0-195 pg/mL) = 0.318

2 SD's of the Zero Standard = 2 x 0.030 = 0.060

Sensitivity = $\frac{0.060}{0.318} \times 195 \text{ pg/mL} = \mathbf{36.8 \text{ pg/mL}}$

Linearity

A sample containing 13,011 pg/mL 15-d-PGJ₂ was diluted 4 times 1:2 in the kit Assay Buffer and measured in the assay. The data was plotted graphically as actual 15-d-PGJ₂ concentration versus measured 15-d-PGJ₂ concentration.

The line obtained had a slope of 0.974 and a correlation coefficient of 0.998.

Precision

Intra-assay precision was determined by taking samples containing low, medium and high concentrations of 15-d-PGJ₂ and running these samples multiple times (n=16) in the same assay. Inter-assay precision was determined by measuring three samples with low, medium and high concentrations of 15-d-PGJ₂ in multiple assays (n=8).

The precision numbers listed below represent the percent coefficient of variation for the concentrations of 15-d-PGJ₂ determined in these assays as calculated by a 4 parameter logistic curve fitting program.

	15-d-PGJ ₂ (pg/mL)	Intra-assay %CV	Inter-assay %CV
Low	1,136	5.6	
Medium	3,062	5.7	
High	4,394	7.4	
Low	1,195		15.7
Medium	4,406		13.0
High	7,219		14.5

Cross Reactivities

The cross reactivities for a number of related eicosanoid compounds was determined by dissolving the cross reactant (purity checked by N.M.R. and other analytical methods) in Assay Buffer at concentrations from 500,000 to 5 pg/mL. These samples were then measured in the 15-d-PGJ₂ assay, and the measured 15-d-PGJ₂ concentration at 50% B/Bo calculated. The % cross reactivity was calculated by comparison with the actual concentration of cross reactant in the sample and expressed as a percentage.

<u>Compound</u>	<u>Cross Reactivity</u>
15-deoxy - $\Delta^{12,14}$ -PGJ ₂	100%
PGJ ₂	49.2%
Δ^{12} -PGJ ₂	5.99%
PGD ₂	4.92%
Arachidonic Acid	0.03%
PGE ₂	<0.01%
PGF _{2α}	<0.01%
9 α ,11 β -PGF _{2α}	<0.01%
Thromboxane B ₂	<0.01%
2-Arachidonoylglycerol	<0.01%
Anandamide	<0.01%

Sample Recoveries

Please refer to pages 4 and 5 for Sample Handling recommendations and Standard preparation.

15-d-PGJ₂ concentrations were measured in a variety of different samples including Tissue Culture Media, human saliva, human urine and porcine plasma. For samples in Tissue Culture Media, ensure that the standards have been diluted into the same media. For saliva, urine and plasma samples 15-d-PGJ₂ was spiked into samples of these media which were diluted with the kit Assay Buffer and assayed in the kit. The following results were obtained:

<u>Sample</u>	<u>% Recovery*</u>	<u>Recommended Dilution*</u>
Tissue Culture Media	97.4	1:2
Human Saliva	94.4	None
Human Urine	113.9	1:16
Porcine Plasma	105.0	≥1:4

* See Sample Handling instructions on page 4 for details.

Reference

1. T. Chard, "An Introduction to Radioimmunoassay & Related Techniques, 4th Ed.", (1990), Amsterdam: Elsevier.
2. P. Tijssen, "Practice & Theory of Enzyme Immunoassays", (1985) Amsterdam: Elsevier.
3. P.W. Ramwell, Biol. Reprod., (1977) 16: 70.
4. R.J. Flower & G.J. Blackwell, Biochem. Pharm., (1976) 25: 285.
5. S. Moncada & J.R. Vane, Pharm. Rev., (1979) 30: 293.
6. B. Samuelsson, et al., Ann. Rev. Biochem., (1978) 47: 997.
7. M. Fukushima, et al., Biochem. Biophys. Res. Comm., (1982) 109: 626-633.
8. F.A. Fitzpatrick & M.A. Wynalda, J. Biol. Chem., (1983) 258: 11713-11718.
9. Y. Kikawa, et al., Proc. Natl. Acad. Sci., USA, (1984) 81: 1317-1321.
10. S. Narumiya, Biochem. Biophys. Res. Comm., (1985) 127: 739-745.
11. Y. Hirata, et al., J. Biol. Chem., (1988) 263: 16619-16625.
12. M. Fukushima, Prostaglandins, Leukotrienes & Ess. Fatty Acids, (1992) 47: 1-12.
13. M. Negishi, et al., J. Lipid Med. Cell Signal., (1995) 12: 443-448.
14. B.F. Forman, et al., Cell, (1995) 83, 803-812.
15. K. Green, et al., Anal. Biochem., (1973) 54: 434.
16. J. Frolich, et al., J. Clin. Invest., (1975) 55: 763.
17. J.E. Shaw & P.W. Ramwell, Meth. Biochem. Anal., (1969) 17: 325.
18. K. Green, et al., Adv. Prostaglandin & Thromboxane Res., (1978) 5: 15.
19. National Committee for Clinical Laboratory Standards Evaluation Protocols, SC1, (1989) Villanova, PA: NCCLS.



USE FOR RESEARCH PURPOSES ONLY

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products

www.enzolifesciences.com
Enabling Discovery in Life Science®

LIMITED WARRANTY; DISCLAIMER OF WARRANTIES

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NONINFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

TRADEMARKS AND PATENTS

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending.

Global Headquarters

Enzo Life Sciences Inc.

10 Executive Blvd
Farmingdale, NY 11735
(p) 1-800-942-0430
(f) 1-631-694-7501
(e) info-usa@enzolifesciences.com

Enzo Life Sciences (ELS) AG

Industriestrasse 17, Postfach
CH-4415 Lause / Switzerland
(p) +41/0 61 926 89 89
(f) +41/0 61 926 89 79
(e) info-ch@enzolifesciences.com

Please visit our website at www.enzolifesciences.com for additional contact information.

