

# **ab133025 – 8-iso-PGF<sub>2</sub> alpha ELISA Kit**

## Instructions for Use

For quantitative detection of 8-iso-PGF<sub>2</sub> alpha in tissue culture media and urine.

This product is for research use only and is not intended for diagnostic use.

# Table of Contents

---

## INTRODUCTION

1. BACKGROUND	2
2. ASSAY SUMMARY	3

## GENERAL INFORMATION

3. PRECAUTIONS	4
4. STORAGE AND STABILITY	5
5. MATERIALS SUPPLIED	5
6. MATERIALS REQUIRED, NOT SUPPLIED	6
7. LIMITATIONS	6
8. TECHNICAL HINTS	7

## ASSAY PREPARATION

9. REAGENT PREPARATION	8
10. STANDARD PREPARATIONS	9
11. SAMPLE COLLECTION AND STORAGE	11
12. PLATE PREPARATION	13

## ASSAY PROCEDURE

13. ASSAY PROCEDURE	14
---------------------	----

## DATA ANALYSIS

14. CALCULATIONS	16
15. TYPICAL DATA	17
16. TYPICAL SAMPLE VALUES	19
17. ASSAY SPECIFICITY	21

## RESOURCES

18. TROUBLESHOOTING	22
19. NOTES	23

## 1. BACKGROUND

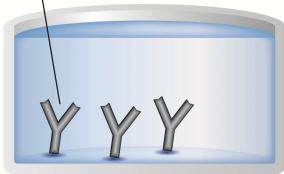
Abcam's 8-iso-PGF<sub>2</sub> alpha *in vitro* competitive ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of 8-iso-PGF<sub>2</sub> alpha in tissue culture media and urine.

A goat anti-rabbit IgG antibody has been precoated onto 96-well plates. Standards or test samples are added to the wells, along with an alkaline phosphatase (AP) conjugated-8-iso-PGF<sub>2</sub> alpha antigen and a polyclonal rabbit antibody specific to 8-iso-PGF<sub>2</sub> alpha. After incubation the excess reagents are washed away. pNpp substrate is added and after a short incubation the enzyme reaction is stopped and the yellow color generated is read at 405 nm. The intensity of the yellow coloration is inversely proportional to the amount of 8-iso-PGF<sub>2</sub> alpha captured in the plate.

The 8-epimer of Prostaglandin F<sub>2</sub> alpha (8-iso-PGF<sub>2</sub> alpha) is produced *in vivo* by both non-cyclooxygenase and cyclooxygenase dependant mechanisms from arachidonic acid. 8-iso-PGF<sub>2</sub> alpha has been shown to be a potent vasoconstrictor, a potential mediator of hepatorenal syndrome and atherosclerosis and a mutagen in 3T3 cells and in vascular smooth muscle cells. It has also been postulated to participate as a pathophysiological mediator and is able to modify the fluidity and integrity of membranes. 8-iso-PGF<sub>2</sub> alpha has been shown to circulate in plasma and is excreted in urine.

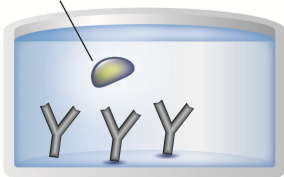
## 2. ASSAY SUMMARY

**Capture Antibody**



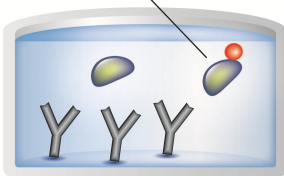
Prepare all reagents and samples as instructed.

**Sample**



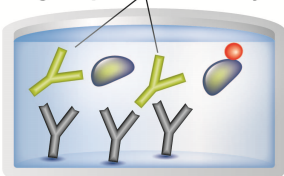
Add standards and samples to appropriate wells.

**Labeled AP-Conjugate**



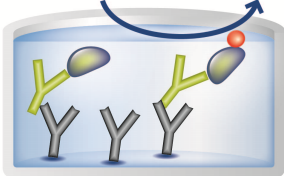
Add prepared labeled AP-conjugate to appropriate wells.

**Target Specific Antibody**



Add 8-iso-PGF<sub>2</sub> alpha antibody to appropriate wells. Incubate at room temperature.

**Substrate**      **Colored Product**



Add pNpp substrate to each well. Incubate at room temperature. Add Stop Solution to each well. Read immediately.

### 3. PRECAUTIONS

**Please read these instructions carefully prior to beginning the assay.**

- 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
- Stop Solution is a solution of trisodium phosphate. This solution is caustic; care should be taken in use
- The activity of the alkaline phosphatase conjugate is dependent on the presence of  $Mg^{2+}$  and  $Zn^{2+}$  ions. The activity of the conjugate is affected by concentrations of chelators ( $>10$  mM) such as EDTA and EGTA
- 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
- The 8-iso-PGF<sub>2</sub> alpha Standard provided is supplied in ethanolic buffer at a pH optimized to maintain 8-iso-PGF<sub>2</sub> alpha integrity. Care should be taken handling this material because of the known and unknown effects of prostaglandin.

#### 4. STORAGE AND STABILITY

**Store kit at +4°C immediately upon receipt. Avoid multiple freeze-thaw cycles.**

Refer to list of materials supplied for storage conditions of individual components.

#### 5. MATERIALS SUPPLIED

Item	Amount	Storage Condition (Before Preparation)
Goat anti-rabbit IgG Microplate (12 x 8 wells)	96 Wells	+4°C
8-iso-PGF <sub>2</sub> alpha Alkaline Phosphatase Conjugate	5 mL	+4°C
8-iso-PGF <sub>2</sub> alpha EIA Antibody	5 mL	+4°C
8-iso-PGF <sub>2</sub> alpha Standard	500 µL	+4°C
10X Assay Buffer	27 mL	+4°C
20X Wash Buffer Concentrate	27 mL	+4°C
pNpp Substrate	20 mL	+4°C
Stop Solution	5 mL	+4°C

### **6. MATERIALS REQUIRED, NOT SUPPLIED**

These materials are not included in the kit, but will be required to successfully utilize this assay:

- Standard microplate reader - capable of reading at 405 nm, preferably with correction between 570 and 590 nm.
- Automated plate washer (optional)
- Adjustable pipettes and pipette tips. Multichannel pipettes are recommended when large sample sets are being analyzed
- Eppendorf tubes
- Microplate Shaker
- Absorbent paper for blotting
- 200 mg C18 Reverse Phase Extraction Columns (only required for extraction of samples containing low levels of 8-iso-PGF<sub>2</sub> alpha)
- 2M hydrochloric acid (only required for extraction of samples containing low levels of 8-iso-PGF<sub>2</sub> alpha)
- Deionized water
- Ethanol
- Hexane
- Ethyl acetate

### **7. LIMITATIONS**

- Assay kit intended for research use only. Not for use in diagnostic procedures
- Do not mix or substitute reagents or materials from other kit lots or vendors. Kits are QC tested as a set of components and performance cannot be guaranteed if utilized separately or substituted

### 8. TECHNICAL HINTS

- Standards can be made up in either glass or plastic tubes.
- Pre-rinse the pipette tip with the reagent, use fresh pipette tips for each sample, standard and reagent.
- Pipette standards and samples to the bottom of the wells.
- Add the reagents to the side of the well to avoid contamination.
- 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.
- 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.
- Prior to addition of substrate, ensure that there is no residual wash buffer in the wells. Any remaining wash buffer may cause variation in assay results.
- **This kit is sold based on number of tests. A ‘test’ simply refers to a single assay well. The number of wells that contain sample, control or standard will vary by product. Review the protocol completely to confirm this kit meets your requirements. Please contact our Technical Support staff with any questions**



## 9. REAGENT PREPARATION

Equilibrate all reagents and samples to room temperature (18 - 25°C) prior to use.

### 9.1 **1X Assay Buffer**

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

### 9.2 **Conjugate 1:10 Dilution for Total Activity Measurement**

Prepare the conjugate 1:10 Dilution by diluting 50µL of the supplied conjugate with 450 µL of the 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.

### 9.3 **1X Wash Buffer**

Prepare the 1X Wash Buffer by diluting 5 mL of the 20X Wash Buffer Concentrate in 95 mL of deionized water. Mix thoroughly and gently.

## 10. STANDARD PREPARATIONS

Prepare serially diluted standards immediately prior to use. Always prepare a fresh set of standards for every use. Preparation of the 8-iso-PGF<sub>2α</sub> standard should be prepared no more than 1 hour prior to the experiment.

### 10.1 For:

10.1.1 **Urine** samples dilute the 8-iso-PGF<sub>2</sub> alpha standard with Assay Buffer.

10.1.2 **Cell culture supernatant** samples dilute the 8-iso-PGF<sub>2</sub> alpha standard with tissue culture media.

10.2 Allow the 1,000,000 pg/mL 8-iso-PGF<sub>2α</sub> Stock Standard solution to warm to room temperature. The standard solution should be stored at -20°C. Avoid repeated freeze-thaw cycles.

10.3 Label eight tubes with #1 –# 8

10.4 Add 900 µL of appropriate diluent (Assay Buffer or Tissue Culture Media) to tube #1.

10.5 Add 750 µL appropriate diluent to tubes #2 through #8.

10.6 Prepare a 100,000 pg/mL **Standard 1** by adding 100 µL of the 1,000,000 pg/mL Stock Standard to tube 1. Vortex thoroughly.

10.7 Prepare **Standard 2** by transferring 250 µL from Standard 1 to tube 2. Vortex thoroughly.

10.8 Prepare **Standard 3** by transferring 250 µL from Standard 2 to tube 3. Vortex thoroughly.

10.9 Using the table below as a guide, repeat for tubes 4 through 8.

## ASSAY PREPARATION

Standard	Sample to Dilute	Volume to Dilute (μL)	Volume of Diluent (μL)	Starting Conc. (pg/mL)	Final Conc. (pg/mL)
1	Standard	100	900	1,000,000	100,000
2	Standard 1	250	750	100,000	25,000
3	Standard 2	250	750	25,000	6,250
4	Standard 3	250	750	6,250	1562.5
5	Standard 4	250	750	1562.5	390.6
6	Standard 5	250	750	390.6	97.7
7	Standard 6	250	750	97.7	24.4
8	Standard 7	250	750	24.4	6.1



## 11. SAMPLE COLLECTION AND STORAGE

- The 8-iso-PGF<sub>2</sub> alpha kit is compatible with 8-iso-PGF<sub>2</sub> alpha samples in a wide range of matrices after dilution in Assay Buffer. However, the end user must verify that the recommended dilutions are appropriate for their samples. Samples containing rabbit IgG may interfere with the assay.
- Samples in the majority of tissue culture media, can be read in the assay, provided the standards have been diluted into the tissue culture media instead of Assay Buffer. There will be a small change in binding associated with running the standards and samples in media.
- For tissue homogenates or urine, prostaglandin synthetase inhibitors such as indomethacin or meclofenamic acid at concentrations up to 10 µg/mL should be added to either the tissue homogenate or urine samples. Some samples normally have very low levels of 8-iso-PGF<sub>2</sub> alpha present and extraction may be necessary for accurate measurement. A suitable extraction procedure is outlined below:
  - 11.1. Acidify the sample by addition of 2M HCl to pH of 3.5. Allow to sit at 4°C for 15 minutes. Centrifuge samples in a microcentrifuge for 2 minutes to remove any precipitate.
  - 11.2. Prepare the C18 reverse phase column (see Section 6) by washing with 10 mL of ethanol followed by 10 mL of deionized water.
  - 11.3. 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.
  - 11.4. 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 sample. Mix well and hold at room temperature for 5 minutes. Repeat twice more. If analysis is to be delayed, store samples as the eluted ethyl acetate

solutions at  $-80^{\circ}\text{C}$  until the immunoassay is to be run. Evaporate the organic solvent under a stream of nitrogen prior to running assay and reconstitute as above.

- Store samples to be assayed within 24 hours at  $2 - 8^{\circ}\text{C}$ . For long-term storage, aliquot and freeze samples at  $-20^{\circ}\text{C}$ . Avoid repeated freeze-thaw cycles.

## 12. PLATE PREPARATION

- The 96 well plate strips included with this kit are supplied ready to use. It is not necessary to rinse the plate prior to adding reagents
- Unused well strips should be returned to the plate packet and stored at 4°C
- For statistical reasons, we recommend each sample should be assayed with a minimum of two replicates (duplicates)
- Well effects have not been observed with this assay. Contents of each well can be recorded on the template sheet included in the Resources section

	1	2	3	4
A	B <sub>s</sub>	Std 1	Std 5	Sample 1
B	B <sub>s</sub>	Std 1	Std 5	Sample 1
C	TA	Std 2	Std 6	Sample 2
D	TA	Std 2	Std 6	Sample 2
E	NSB	Std 3	Std 7	etc
F	NSB	Std 3	Std 7	etc
G	B <sub>0</sub>	Std 4	Std 8	
H	B <sub>0</sub>	Std 4	Std 8	

Plate layout shows controls, blanks and standards required for each assay. Use additional strips of wells to assay all your samples.

### Key:

**B<sub>s</sub>** = Blank; contains substrate only.

**TA** = Total Activity; contains conjugate (5 µL) and substrate.

**NSB** = Non-specific binding; contains standard diluent, assay buffer, conjugate and substrate.

**B<sub>0</sub>** = 0 pg/mL standard; contains standard diluent, conjugate, antibody and substrate

## 13. ASSAY PROCEDURE

- Equilibrate all materials and prepared reagents to room temperature prior to use
- It is recommended to assay all standards, controls and samples in duplicate
- Refer to the recommended plate layout in Section 12 before proceeding with the assay

13.1 Add 150  $\mu$ L appropriate diluent\* into the NSB (non-specific binding) wells. (\*Use the same diluent used to prepare standards in section 10, either Assay Buffer or Tissue Culture Media).

13.2 Add 100  $\mu$ L appropriate diluent (Assay Buffer or tissue culture media) into the B<sub>0</sub> (0 pg/mL standard) wells

13.3 Add 100  $\mu$ L of prepared standards and diluted samples to appropriate wells.

13.4 Add 50  $\mu$ L of 8-iso-PGF<sub>2</sub> alpha Alkaline Phosphatase Conjugate (blue) into NSB, B<sub>0</sub>, standard and sample wells, i.e. not the Total Activity (TA) and B<sub>s</sub> wells.

13.5 Add 50  $\mu$ L of 8-iso-PGF<sub>2</sub> alpha antibody (yellow) into B<sub>0</sub>, standard and sample wells, i.e. not B<sub>s</sub>, TA and NSB wells.

**NOTE:** Every well used should be green in color except the NSB wells which should be blue. The B<sub>s</sub> and TA wells are empty at this point and have no color.

13.6 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.

13.7 Empty the contents of the wells and wash by adding 400  $\mu$ L of 1X Wash Buffer to every well. Repeat the wash 2 more times for a total of 3 washes. After the final wash, empty or aspirate the wells, and firmly tap the plate on a lint free paper towel to remove any remaining wash buffer.

13.8 Add 5  $\mu$ L of the 8-iso-PGF<sub>2</sub> alpha Alkaline Phosphatase Conjugate to the TA wells.

## ASSAY PROCEDURE

- 13.9 Add 200  $\mu$ L of the pNpp Substrate solution to every well. Incubate at room temperature for 45 minutes without shaking.
- 13.10 Add 50  $\mu$ L Stop Solution into each well. The plate should be read immediately.
- 13.11 Read the O.D. absorbance at 405 nm, preferably with correction between 570 and 590 nm.



## 14. CALCULATIONS

- 14.1 Calculate the average net absorbance measurement (Average Net OD) for each standard and sample by subtracting the average NSB absorbance measurement from the average absorbance measurement (Average OD) for each standard and sample.

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

- 14.2 Calculate the binding of each pair of standard wells as a percentage of the maximum binding wells ( $B_0$ ), using the following formula

$$\text{Percent Bound} = \frac{\text{Average Net OD}}{\text{Average Net } B_0 \text{ OD}} \times 100$$

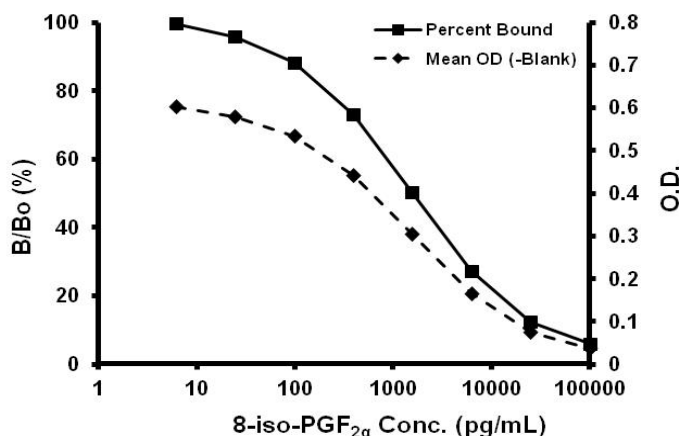
- 14.3 Plot the Percent Bound ( $B/B_0$ ) and the net OD versus concentration of 8-iso-PGF<sub>2</sub> alpha for the standards. The concentration of 8-iso-PGF<sub>2</sub> alpha in the unknowns can be determined by interpolation of net OD values.

A four parameter algorithm (4PL) provides the best fit, though other equations can be examined to see which provides the most accurate (e.g. linear, semi-log, log/log, 4 parameter logistic). Interpolate protein concentrations for unknown samples from the standard curve plotted.

Samples producing signals greater than that of the highest standard should be further diluted and reanalyzed, then multiplying the concentration found by the appropriate dilution factor

## 15. TYPICAL DATA

**TYPICAL STANDARD CURVE** – Data provided for **demonstration purposes only**. A new standard curve must be generated for each assay performed.



Sample	Mean OD (- B <sub>s</sub> )	Percent Bound	8-iso-PGF <sub>2</sub> alpha (pg/mL)
B <sub>s</sub>	(0.089)		
TA	0.401		
NSB	0.001	0	
(B <sub>0</sub> )	0.605	100	0
S1	0.037	6	100,000
S2	0.076	12.4	25,000
S3	0.166	27.3	6,250
S4	0.305	50.3	1,562.50
S5	0.442	73	390.6
S6	0.534	88.2	97.7
S7	0.579	95.7	24.4
S8	0.602	99.5	6.1
Unknown 1	0.171	28.2	5,944
Unknown 2	0.355	58.6	965

### TYPICAL QUALITY CONTROL PARAMETERS –

Total Activity Added	= $0.401 \times 10 \times 10 = 40.1$
%B <sub>0</sub> /TA	= 15.06
Quality of Fit	= 1.000 (Calculated from 4 parameter logistic curve fit)
20% Intercept	= 11,286 pg/mL
50% Intercept	= 1,563 pg/mL
80% Intercept	= 229 pg/mL

## 16. TYPICAL SAMPLE VALUES

### SENSITIVITY –

The sensitivity, minimum detectable dose of 8-iso-PGF<sub>2</sub> alpha using this Abcam ELISA kit was found to be 16.3 pg/mL. This was determined by the average optical density of the 0 pg/mL Standard and comparing to the average optical density for Standard 7. The detection limit was determined as the concentration of 8-iso-PGF<sub>2</sub> alpha measured at two standard deviations from the zero along the standard curve.

### SAMPLE RECOVERY –

Recovery was determined by 8-iso-PGF<sub>2</sub> alpha into tissue culture media, and Human urine. Mean recoveries are as follows:

Sample Type	Average % Recovery	Recommended Dilution
Tissue Culture Media	129.9	1:10
Human Urine	101.58	Neat - 1:60

### LINEARITY OF DILUTION –

A sample containing 10,000 pg/mL 8-iso-PGF<sub>2</sub> alpha was diluted 8 times 1:2 in the kit Assay Buffer and measured in the assay. The data was plotted graphically as actual 8-iso-PGF<sub>2</sub> alpha concentration versus measured 8-iso-PGF<sub>2</sub> alpha concentration.

The line obtained had a slope of 1.01 and a correlation coefficient of 0.991.

## PRECISION –

### Intra-Assay

	8-iso-PGF <sub>2</sub> alpha (pg/mL)	%CV
Low	232	11.0
Medium	1,128	4.4
High	8,263	5.8

### Inter-Assay

	8-iso-PGF <sub>2</sub> alpha (pg/mL)	%CV
Low	183	11.0
Medium	855	8.8
High	6,239	5.0

## 17. ASSAY SPECIFICITY

### CROSS REACTIVITY –

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 100,000 to 6 pg/ml. These samples were then measured in the 8-iso-PGF<sub>2</sub>α assay and the measured 8-iso-PGF<sub>2</sub>α concentration at 50% B/Bo was calculated. The % cross reactivity was calculated by comparison with the actual concentration of cross reactant in the sample and expressed as a percentage.

8-iso-PGF <sub>2</sub> α	100 %
PGF <sub>1</sub> α	4.6 %
PGF <sub>2</sub> α	1.85 %
PGE <sub>1</sub>	0.19 %
TXB <sub>2</sub>	0.023 %
PGB <sub>1</sub>	0.02 %
PGE <sub>3</sub>	0.012 %
6-keto-PGF <sub>1</sub> α	0.008 %
13,14-dihydro-15-keto-PGF <sub>2</sub> α	0.008 %
6,15-keto-13,14-dihydro-PGF <sub>1</sub> α	0.005 %
8-iso-PGE <sub>1</sub>	<0.001 %
PGA <sub>2</sub>	<0.001 %
PGJ <sub>2</sub>	<0.001 %
2-Arachidonoylglycerol	<0.001 %
Anandamide	<0.001 %

## 18. TROUBLESHOOTING

Problem	Cause	Solution
Poor standard curve	Inaccurate pipetting	Check pipettes
	Improper standards dilution	Prior to opening, briefly spin the stock standard tube and dissolve the powder thoroughly by gentle mixing
Low Signal	Incubation times too brief	Ensure sufficient incubation times; change to overnight standard/sample incubation
	Inadequate reagent volumes or improper dilution	Check pipettes and ensure correct preparation
Samples give higher value than the highest standard	Starting sample concentration is too high.	Dilute the specimens and repeat the assay
Large CV	Plate is insufficiently washed	Review manual for proper wash technique. If using a plate washer, check all ports for obstructions
	Contaminated wash buffer	Prepare fresh wash buffer
Low sensitivity	Improper storage of the kit	Store the all components as directed.

### 19. NOTES









## **Technical Support**

Copyright © 2022 Abcam, All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

**For all technical or commercial enquiries please go to:**

[www.abcam.com/contactus](http://www.abcam.com/contactus)

[www.abcam.cn/contactus](http://www.abcam.cn/contactus) (China)

[www.abcam.co.jp/contactus](http://www.abcam.co.jp/contactus) (Japan)