

PGF2 alpha High Sensitivity ELISA Kit ab133056

2 References 1 Image

Overview

Product name PGF2 alpha High Sensitivity ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
107.5pmol/ml	24			6.53%
33.79pmol/ml	24			4.94%
11.5pmol/ml	24			7.16%

Inter-assay

Sample	n	Mean	SD	CV%
117.8pmol/ml				10.97%
35.81pmol/ml				8.88%
10.1pmol/ml				10.14%

Sample type Cell culture supernatant, Saliva, Milk, Urine, Serum, Plasma

Assay type Competitive

Sensitivity 0.98 pg/ml

Range 1.95 pg/ml - 2000 pg/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Saliva	103.5	% - %
Urine	107	% - %
Serum	100.3	% - %

Sample type	Average %	Range
Plasma	96.9	% - %


Assay time

2h 00m

Assay duration

Multiple steps standard assay

Species reactivity

Predicted to work with: a wide range of other species 

Product overview

Abcam's PGF_{2α} High Sensitivity *in vitro* competitive ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of PGF_{2α} in milk, culture supernatants, saliva, urine, serum and plasma.

A donkey anti-sheep 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-PGF_{2α} antigen and a polyclonal sheep antibody specific to PGF_{2α}. 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 PGF_{2α} captured in the plate.

Notes

Prostaglandin F_{2α} (PGF_{2α}) is formed in a variety of cells from PGH₂, which itself is synthesized from arachidonic acid by the enzyme prostaglandin synthetase. PGF_{2α} is often viewed as an antagonist to PGE₂ due to their opposing effects on various tissues. PGF_{2α} is a potent bronchoconstrictor and has been implicated in asthma attacks. PGF_{2α} is also involved in reproductive functions including corpus luteum regulation, uterine contractions, and sperm motility. This has led to its use in terminating pregnancies and inducing labor at term. High levels of PGF_{2α} have also been associated with preeclampsia.

Compound	Cross Reactivity
PGF _{2α}	100%
PGF _{1α}	11.82%
PGD ₂	3.62%
6-keto-PGF _{1α}	1.38%
PGI ₂	1.25%
PGE ₂	0.77%
Thromboxane B ₂	0.77%
8-iso PGF _{2α}	0.73%

PGE ₁	0.39%
PGA ₂	<0.1%
6, 15-keto-13, 14-dihydro-PGF _{1α}	<0.01%
2-Arachidonoylglycerol	<0.01%
Anandamide	<0.01%

Platform

Microplate

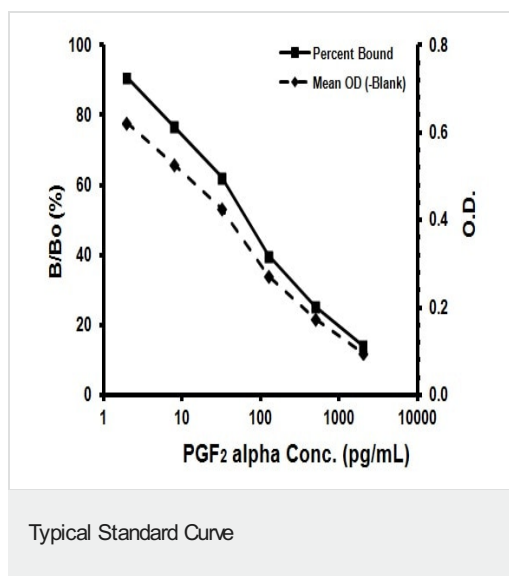
Properties

Storage instructions

Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
20X Wash Buffer Concentrate	1 x 27ml
Assay Buffer	1 x 27ml
Donkey anti-sheep IgG Microplate (12 x 8 wells)	1 unit
PGF _{2α} Alkaline Phosphatase Conjugate	1 x 2.5ml
PGF _{2α} Antibody	1 x 2.5ml
PGF _{2α} Standard	1 x 0.5ml
Plate Sealer	1 unit
pNpp Substrate	1 x 20ml
Stop Solution	1 x 5ml

Images



Representative Standard Curve using ab133056.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors