

Human PLGF ELISA Kit ab100629

7 References 1 Image

Overview

Product name	Human PLGF ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 2 pg/ml
Range	1.372 pg/ml - 1000 pg/ml
Recovery	89 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	103.4	75% - 115%
Serum	88.09	79% - 96%
Plasma	75.69	68% - 93%

Assay duration	Multiple steps standard assay
Species reactivity	<b>Reacts with:</b> Human
Product overview	<p>Abcam's PLGF Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Human PLGF in serum, plasma, and cell culture supernatants.</p> <p>This assay employs an antibody specific for Human PLGF coated on a 96-well plate. Standards and samples are pipetted into the wells and PLGF present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human PLGF antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of PGF bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.</p>
Notes	Optimization may be required with urine samples.
Platform	Microplate

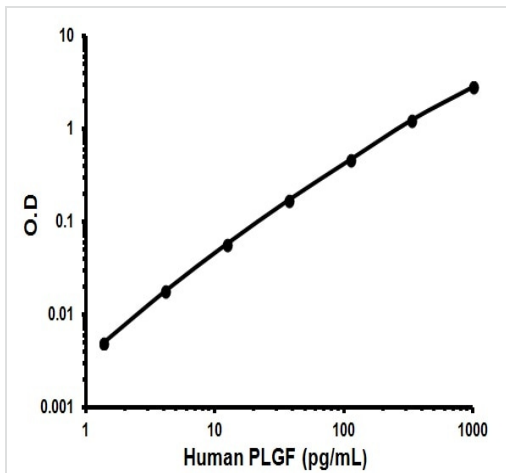
## Properties

**Storage instructions** Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent	1 x 15ml
Biotinylated anti-Human PLGF	2 vials
PLGF Microplate (12 x 8 wells)	1 unit
Recombinant Human PLGF Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

<b>Function</b>	Growth factor active in angiogenesis and endothelial cell growth, stimulating their proliferation and migration. It binds to the receptor FLT1/VEGFR-1. Isoform PIGF-2 binds NRP1/neuropilin-1 and NRP2/neuropilin-2 in a heparin-dependent manner.
<b>Tissue specificity</b>	While the three isoforms are present in most placental tissues, PIGF-2 is specific to early (8 week) placenta and only PIGF-1 is found in the colon and mammary carcinomas.
<b>Sequence similarities</b>	Belongs to the PDGF/VEGF growth factor family.
<b>Domain</b>	Isoform PIGF-2 contains a basic insert which acts as a cell retention signal.
<b>Post-translational modifications</b>	N-glycosylated.
<b>Cellular localization</b>	Secreted. The three isoforms are secreted but PIGF-2 appears to remain cell attached unless released by heparin.

## Images



Representative Standard Curve using ab100629.

Typical Standard Curve

**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