

Human PLGF ELISA Kit ab100629

7 References 画像数 1

製品の概要

製品名	Human PLGF ELISA Kit
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Serum, Plasma
アッセイタイプ	Sandwich (quantitative)
検出感度	< 2 pg/ml
検出範囲	1.372 pg/ml - 1000 pg/ml
添加回収試験	89 %

特定サンプルでの回収試験

サンプルの種類	平均 %	測定範囲
Cell culture supernatant	103.4	75% - 115%
Serum	88.09	79% - 96%
Plasma	75.69	68% - 93%

ステップ	Multiple steps standard assay
種交差性	交差種: Human
製品の概要	<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>
特記事項	Optimization may be required with urine samples.
試験プラットフォーム	Microplate

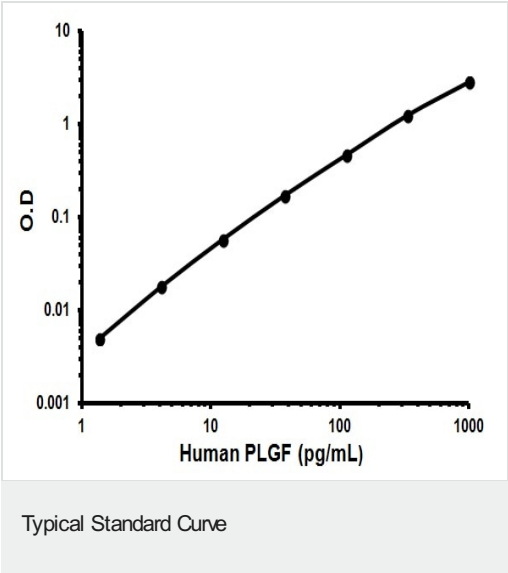
製品の特性

保存方法Store at -20°C. Please refer to protocols.

内容	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

機能	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.
組織特異性	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.
配列類似性	Belongs to the PDGF/VEGF growth factor family.
ドメイン	Isoform PIGF-2 contains a basic insert which acts as a cell retention signal.
翻訳後修飾	N-glycosylated.
細胞内局在	Secreted. The three isoforms are secreted but PIGF-2 appears to remain cell attached unless released by heparin.

画像



Representative Standard Curve using ab100629.

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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