

Human PDGF AA ELISA Kit ab100622

3 References 1 Image

Overview

| | |
|------------------|---|
| Product name | Human PDGF AA ELISA Kit |
| Detection method | Colorimetric |
| Sample type | Cell culture supernatant, Serum, Plasma |
| Assay type | Sandwich (quantitative) |
| Sensitivity | < 40 pg/ml |
| Range | 40 pg/ml - 30000 pg/ml |
| Recovery | |

Sample specific recovery

| Sample type | Average % | Range |
|--------------------------|-----------|------------|
| Cell culture supernatant | 92.54 | 82% - 102% |
| Serum | 76.71 | 70% - 83% |
| Plasma | 78.06 | 72% - 85% |

Assay duration Multiple steps standard assay

Species reactivity Reacts with: Human

Product overview Abcam's PDGF AA Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human PDGF AA in serum, plasma, and cell culture supernatants.

This assay employs an antibody specific for Human PDGF AA coated on a 96- well plate. Standards and samples are pipetted into the wells and PDGF AA present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human PDGF AA 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 PDGF AA bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

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 PDGF AA | 2 vials |
| PDGF AA Microplate (12 x 8 wells) | 1 unit |
| Recombinant Human PDGF AA Standard (lyophilized) | 2 vials |
| Stop Solution | 1 x 8ml |
| TMB One-Step Substrate Reagent | 1 x 12ml |

Function

Platelet-derived growth factor is a potent mitogen for cells of mesenchymal origin. Binding of this growth factor to its receptor elicits a variety of cellular responses. It is released by platelets upon wounding and plays an important role in stimulating adjacent cells to grow and thereby heals the wound.

Sequence similarities

Belongs to the PDGF/VEGF growth factor family.

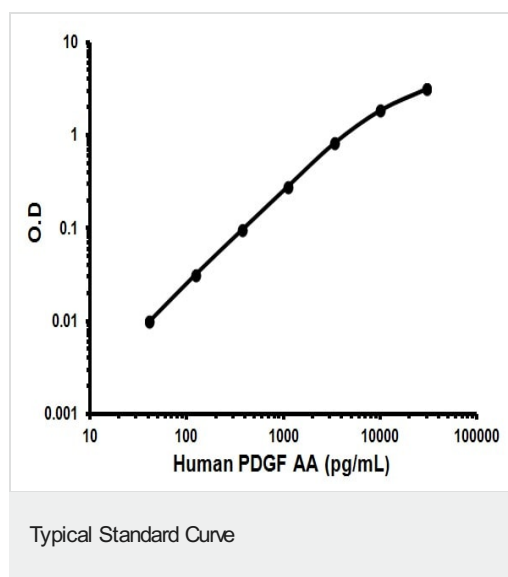
Domain

The long form contains a basic insert which acts as a cell retention signal.

Cellular localization

Secreted.

Images



Representative Standard Curve using ab100622.

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