

Human PDGF BB ELISA Kit ab100624

9 References 2 Images

Overview

Product name	Human PDGF BB ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 1 pg/ml
Range	0.549 pg/ml - 400 pg/ml
Recovery	93 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	95.43	84% - 104%
Serum	93.57	82% - 103%
Plasma	91.99	83% - 104%

Assay duration Multiple steps standard assay

Species reactivity Reacts with: Human

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

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

Get results in 90 minutes with Human PDGF BB ELISA Kit ([ab184860](#)) from our SimpleStep

ELISA® range.

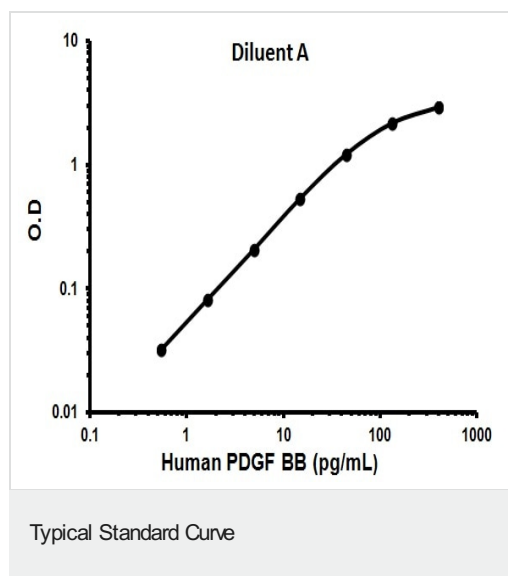
<b>Notes</b>	<b>Note:</b> Optimization may be required with urine samples, and very low level are detectable in serum
<b>Platform</b>	Microplate

## Properties

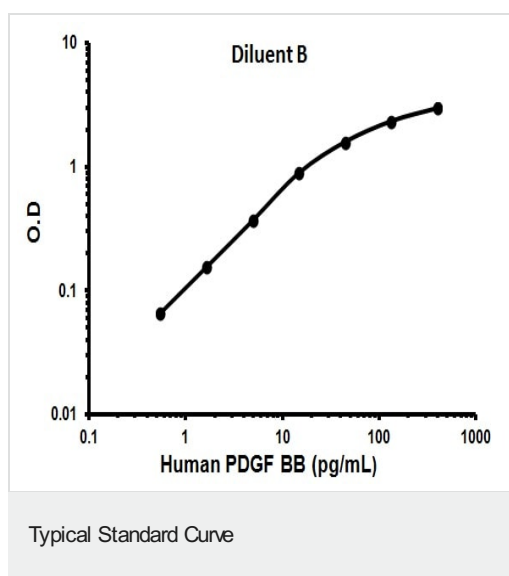
**Storage instructions** Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
800X HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent A	1 x 30ml
Biotinylated anti-Human PDGF BB	2 vials
PDGF BB Microplate (12 x 8 wells)	1 unit
Recombinant Human PDGF BB Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

## Images



Representative Standard Curve using ab100624.



Representative Standard Curve using ab100624.

**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