

Human PDGF BB ELISA Kit ab184860

SimpleStep ELISA

★★★★★ [1 Abreviews](#) [4 References](#) [7 图像](#)

概述

产品名称 人PDGF BB ELISA试剂盒

检测方法 Colorimetric

精确度 批次内

样品	n	Mean	SD	CV%
Human Serum	8			3.3%

批次间

样品	n	Mean	SD	CV%
Human Serum	3			5.3%

样品类型 Serum, Plasma, Cell culture media

检测类型 Sandwich (quantitative)

灵敏度 2.9 pg/ml

范围 11.7 pg/ml - 750 pg/ml

回收率 特定样本回收率

样品类型	平均%	范围
Serum	95	85% - 100%
Cell culture media	104	101% - 106%
Hep Plasma	104	92% - 114%
EDTA Plasma	99	91% - 107%
Cit plasma	109	100% - 114%

检测时间 1h 30m

实验步骤 One step assay

种属反应性**与反应:** Human**产品概述**

Human PDGF BB ELISA kit (ab184860) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of PDGF BB protein in human serum, plasma, and cell culture supernatant. It uses our proprietary SimpleStep ELISA® technology. Quantitate human PDGF BB with 2.9 pg/mL sensitivity.

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate ([ab203359](#)) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

说明

Platelet derived growth factor B chain plays an essential role in the regulation of embryonic development, cell proliferation, cell migration, survival and chemotaxis. As a homodimer it is a potent mitogen for cells of mesenchymal origin and is required for normal proliferation and recruitment of pericytes and vascular smooth muscle cells in the central nervous system, skin, lung, heart and placenta. PDGF-BB is essential for normal blood vessel development, for normal development of kidney glomeruli and in wound healing. Signaling is modulated by the formation of heterodimers with PDGF A chain to yield PDGF-AB.

平台

Microplate

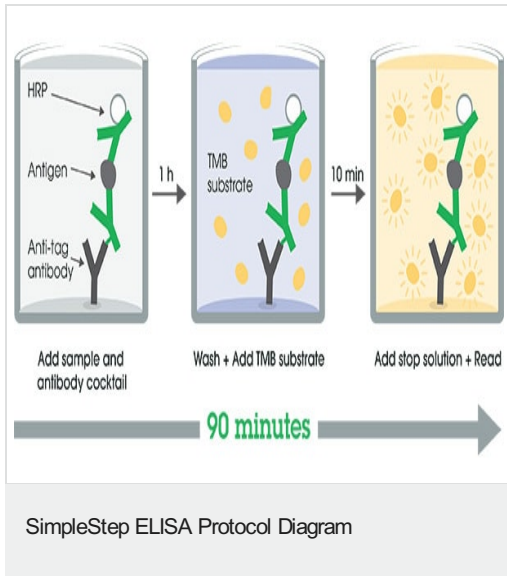
性能**存放说明**

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

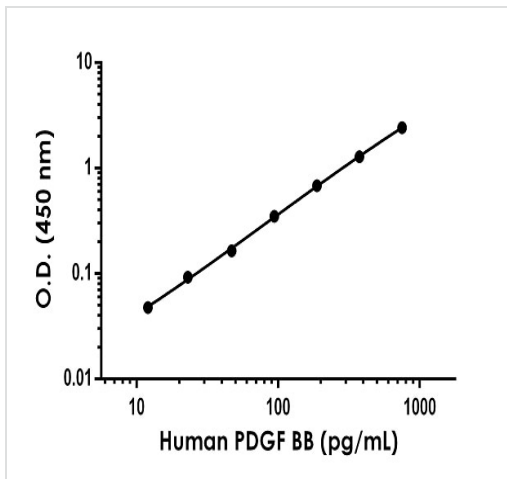
组件	1 x 96 tests	10 x 96 tests
10X Human PDGF BB Capture Antibody	1 x 600µl	10 x 600µl
10X Human PDGF BB Detector Antibody	1 x 600µl	10 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml	1 x 200ml
Antibody Diluent 4BI	1 x 6ml	10 x 6ml
Human PDGF BB Lyophilized Recombinant Protein	2 vials	20 vials
Plate Seals	1 unit	10 units
Sample Diluent NS (ab193972)	1 x 50ml	2 x 250ml

组件	1 x 96 tests	10 x 96 tests
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit	10 units
Stop Solution	1 x 12ml	1 x 120ml
TMB Development Solution	1 x 12ml	1 x 120ml

图片

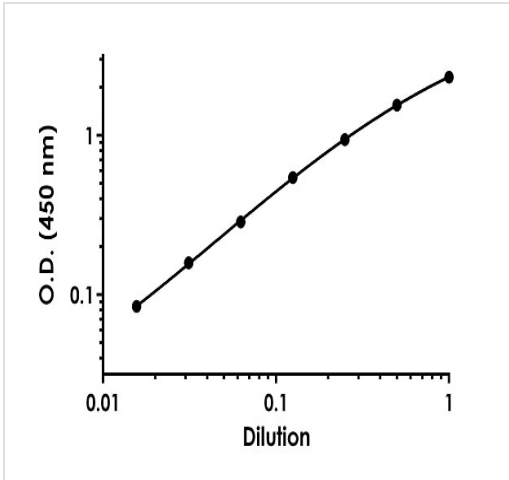


SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.



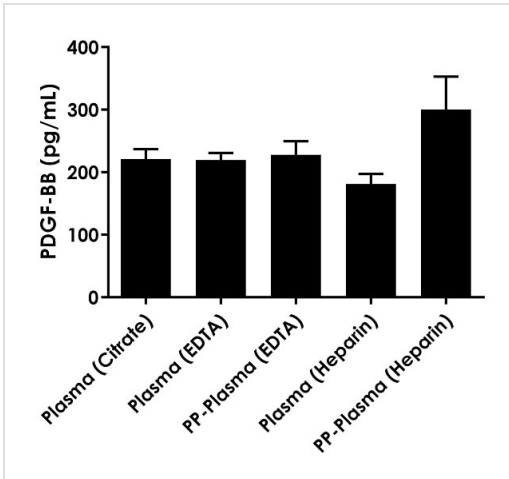
The PDGF BB standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.

Example of human PDGF BB standard curve in Sample Diluent NS.

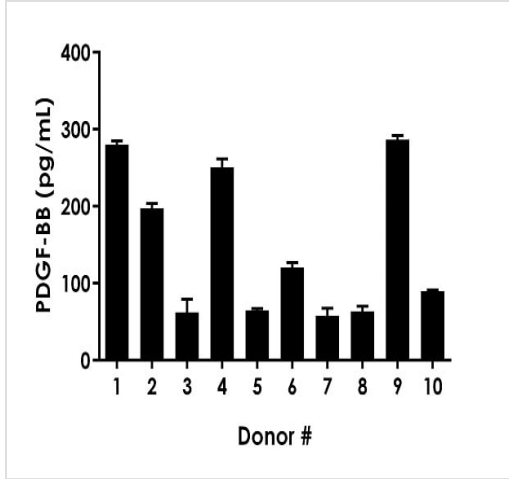


Background subtracted data from duplicate measurements are plotted.

Titration of human serum within the working range of the assay.

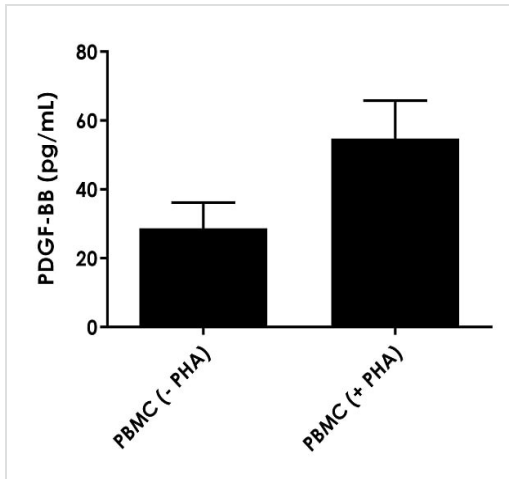


Observed PDGF-BB levels in pooled donor normal human plasma and normal human platelet-poor plasma (PP-Plasma).



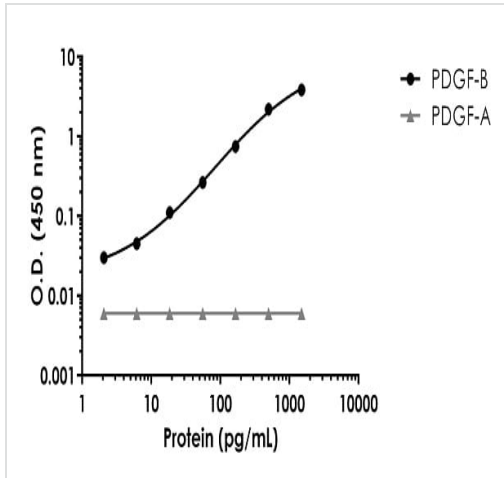
Mean human PDGF-BB is 145 pg/mL and values fall within expected normal reference ranges (WHO).

Observed PDGF-BB levels in individual donor normal human serum (n=10).



Cells were cultured unstimulated or stimulated with 10 µg/mL PHA. Conditioned media was harvested after 48 hours aliquoted and assayed for endogenous PDGF-BB levels.

Human peripheral blood cells (1x10⁶ cells/mL) were cultured in RPMI media supplemented with 10% fetal calf serum, 100 U/mL penicillin, and 100 µg/mL streptomycin sulfate.



The PDGF-BB SimpleStep ELISA was validated for assay specificity against PDGF-A protein.

Standard curve was prepared as described in Section 10.

Background-subtracted data values (mean +/- SD) are graphed. In addition, PDGF-A was prepared at 15 ng/mL and assayed for cross reactivity. No cross reactivity or interference was observed.

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.cn/abpromise> or contact our technical team.

Terms and conditions

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