

Rat PDGF-AB ELISA Kit ab213906

1 Image

Overview

Product nameRat PDGF-AB ELISA Kit

Detection methodColorimetric

PrecisionIntra-assay

Sample	n	Mean	SD	CV%
Sample 1	16	246pg/ml	14.27	5.8%
Sample 2	16	632pg/ml	34.76	5.5%
Sample 3	16	1068pg/ml	50.2	4.7%

Inter-assay

Sample	n	Mean	SD	CV%
Sample 1	24	275pg/ml	20.35	7.4%
Sample 2	24	723pg/ml	45.55	6.3%
Sample 3	24	1249pg/ml	73.7	5.9%

Sample typeCell culture supernatant, Serum, Cell Lysate, Hep Plasma, EDTA Plasma, Tissue Homogenate

Assay typeSandwich (quantitative)

Sensitivity< 3 pg/ml

Range31.2 pg/ml - 2000 pg/ml

Assay time3h 30m

Assay durationMultiple steps standard assay

Species reactivityReacts with: Rat

Product overviewThe Rat PDGF-AB Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213906) is designed for the quantitative measurement of Rat PDGF-AB in cell culture supernatants, cell lysates, tissue homogenates, serum and plasma (heparin, EDTA).

The ELISA kit is based on standard sandwich enzyme-linked immune-sorbent assay technology.

A monoclonal antibody from mouse specific for PDGF-AB has been pre-coated onto 96-well plates. Standards and test samples are added to the wells; a biotinylated detection polyclonal antibody from goat specific for PDGF-AB is added subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex is added and unbound conjugates are washed away with PBS or TBS buffer. HRP substrate TMB is used to visualize HRP enzymatic reaction. TMB is catalyzed by HRP to produce a blue color product that changed into yellow after adding acidic stop solution. The density of yellow is proportional to the rat PDGF-AB amount of sample captured in plate.

Notes

The platelet-derived growth factor (PDGF) is a mitogen derived from human platelets consisting of two related polypeptides termed A and B chains. The genes for PDGF A chain, B chain/c-sis, and the PDGF receptor are expressed in human malignant glioma cell lines. Normal human endothelial cells in culture express the B chain of PDGF, and that endothelial-derived PDGF B chain is synthesized as a predicted precursor polypeptide of Mr 27,281. The entire B chain of PDGF is highly (96%) homologous to a portion of p28sis, the transforming protein of simian sarcoma virus (SSV). It has been suggested that p28sis exerts its transforming potential by mimicking the growth promoting activity of PDGF and stimulating the cell in an autocrine manner. PDGF A-chain precursor polypeptide is assigned to the proximal long arm of chromosome 7, band q11.23. The human homolog (PDGF Bchain/c-sis) of the transforming gene of simian sarcoma virus is assigned to chromosome 22.

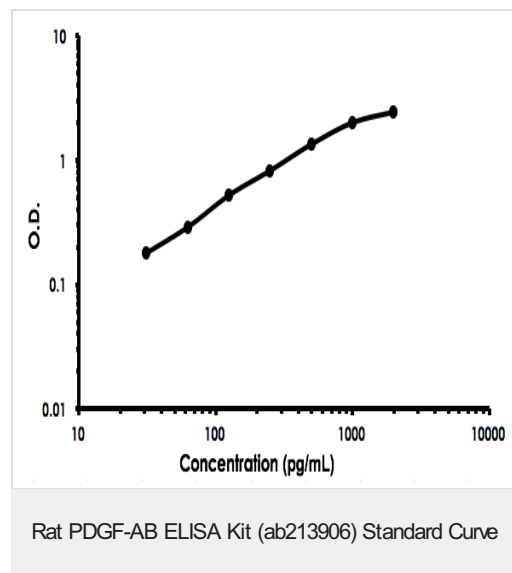
Platform

Pre-coated microplate (12 x 8 well strips)

Properties

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

Components	Identifier	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml
Adhesive Plate Seal		4 units
Antibody Diluent Buffer	Green Cap	1 x 12ml
Anti-rat PDGF-AB coated Microplate (12 x 8 wells)		1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl
Biotinylated anti- Rat PDGF-AB antibody		1 x 100µl
Lyophilized recombinant Rat PDGF-AB standard		2 vials
Sample Diluent Buffer	Green Cap	1 x 30ml
TMB Color Developing Agent	Black Cap	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml
Wash Buffer (25X)		1 x 20ml



Rat PDGF-ABELISA Kit (ab213906) 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