

Rat PDGF-AA ELISA Kit ab155464

[2 References](#) [2 Images](#)

Overview

Product name Rat PDGF-AA ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
overall				< 10%

Inter-assay

Sample	n	Mean	SD	CV%
overall				< 12%

Sample type Serum, Plasma, Cell culture media

Assay type Sandwich (quantitative)

Sensitivity < 90 pg/ml

Range 93.75 pg/ml - 6000 pg/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Serum	100	91% - 108%
Plasma	101.6	84% - 114%
Cell culture media	130.4	122% - 138%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Rat

Product overview

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

This assay employs an antibody specific for rat 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-rat 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.

Platform Microplate

Properties

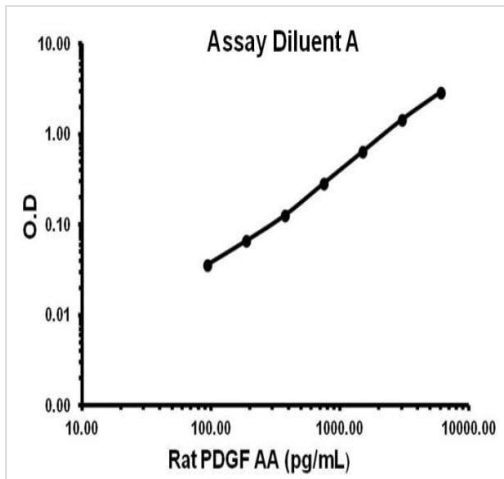
Storage instructions Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
280X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-rat PDGF-AA	2 vials
PDGF-AA Microplate (12 x 8 wells)	1 unit
Recombinant Rat PDGF-AA Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Relevance

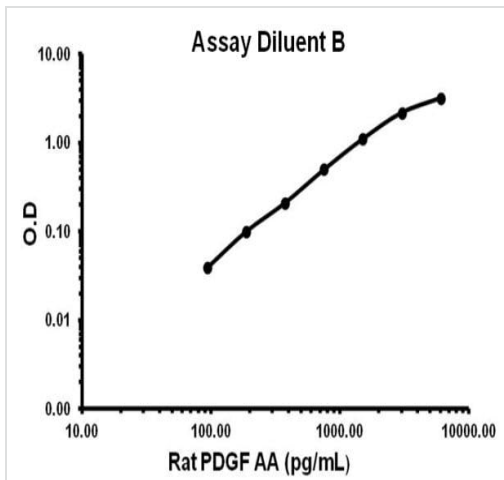
Function: Growth factor that plays an essential role in the regulation of embryonic development, cell proliferation, cell migration, survival and chemotaxis. Potent mitogen for cells of mesenchymal origin. Required for normal lung alveolar septum formation during embryogenesis, normal development of the gastrointestinal tract, normal development of Leydig cells and spermatogenesis. Required for normal oligodendrocyte development and normal myelination in the spinal cord and cerebellum. Plays an important role in wound healing. Signaling is modulated by the formation of heterodimers with PDGFB. Subunit structure: Homodimer (AA); antiparallel disulfide-linked dimer. Heterodimer with PDGFB (AB); antiparallel disulfide-linked dimer. The PDGFA homodimer interacts with PDGFRA homodimers, and with heterodimers formed by PDGFRA and PDGFRB. The heterodimer composed of PDGFA and PDGFB interacts with PDGFRA homodimers, and with heterodimers formed by PDGFRA and PDGFRB. Interacts with CSPG4.

Images



Representative standard curve using ab155464 - Assay Diluent A.

Typical Standard Curve



Representative standard curve using ab155464 - Assay Diluent B.

Typical 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