

Product datasheet

Human PDGF AB ELISA Kit ab100623

2 Images

Overview

Product name	Human PDGF AB ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 10 pg/ml
Range	10.24 pg/ml - 2500 pg/ml
Recovery	74 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	74.29	68% - 78%
Serum	73.96	67% - 78%
Plasma	74.85	67% - 79%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

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

This assay employs an antibody specific for Human PDGF AB coated on a 96- well plate. Standards and samples are pipetted into the wells and PDGF AB present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human PDGF AB 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 AB 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

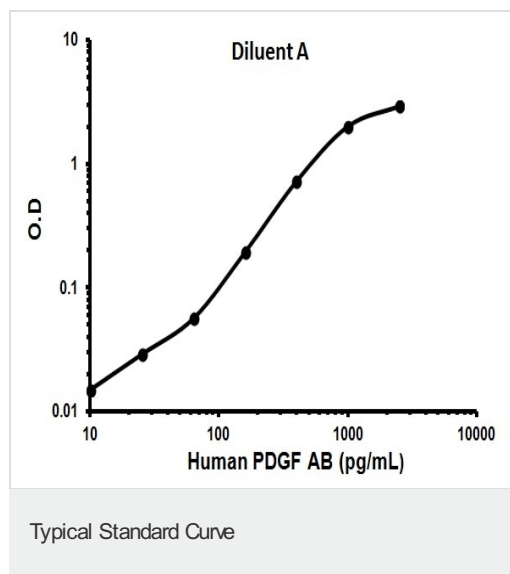
Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
500X HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent A	1 x 30ml
Biotinylated anti-Human PDGF AB	2 vials
PDGF AB Microplate (12 x 8 wells)	1 unit
Recombinant Human PDGF AB Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

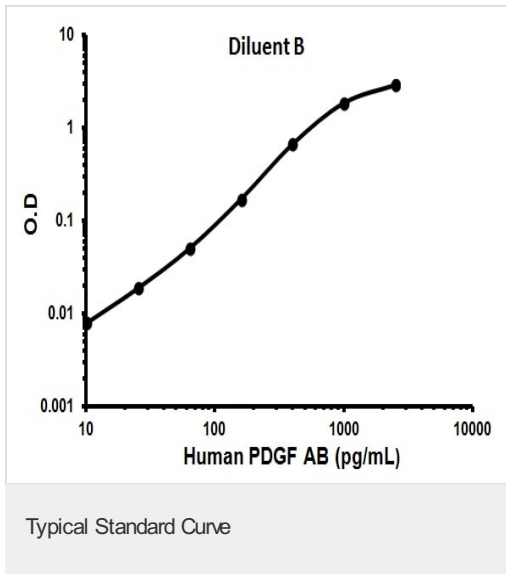
Relevance

Platelet Derived Growth Factor (PDGF) is a potent mitogen for a wide range of cell types including fibroblasts, smooth muscle and connective tissue. PDGF, which is composed of a dimer of two chains joined by disulphide bonds, termed the A chain and B chain, can be present as AA or BB homodimers or as an AB heterodimer. Human PDGF-AB is a 25.5 kDa heterodimer protein consisting of a 13.3 kDa A chain and a 12.2 kDa B chain.

Images



Representative Standard Curve using ab100623.



Representative Standard Curve using ab100623.

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