

Porcine PDGF-BB ELISA ab218297

1 图像

概述

产品名称	Porcine PDGF-BB ELISA				
检测方法	Colorimetric				
精确度	批次内				
	样品	n	Mean	SD	CV%
	1				< 10%
	批次间				
	样品	n	Mean	SD	CV%
	1				< 12%
样品类型	Cell culture supernatant, Serum, Plasma				
检测类型	Sandwich (quantitative)				
灵敏度	10 pg/ml				
范围	10 pg/ml - 2500 pg/ml				
回收率	特定样本回收率				
	样品类型	平均%		范围	
	Cell culture supernatant	120.5		108% - 130%	
	Serum	108.6		87% - 123%	
	Plasma	112.3		98% - 115%	
检测时间	4h 45m				
实验步骤	Multiple steps standard assay				
种属反应性	与反应: Human, Pig				
产品概述	The Porcine PDGF-BB ELISA kit (ab218297) is an <i>in vitro</i> enzyme-linked immunosorbent assay (ELISA) for the quantitative measurement of porcine PDGF-BB in serum, plasma and cell culture supernatants.				

This assay employs an antibody specific for porcine 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-porcine 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.

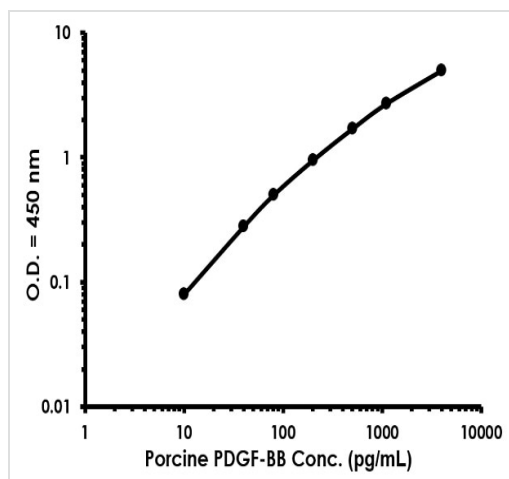
平台 Pre-coated microplate (12 x 8 well strips)

性能

存放说明 Store at -20°C. Please refer to protocols.

组件	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent	1 x 15ml
800X Streptavidin-Conjugated HRP	1 x 200μl
Anti-Human PDGF-BB coated Microplate (12 x 8 wells)	1 x 96 tests
Biotinylated (PDGF-BB) Detection Antibody	2 vials
Lyophilized (PDGF-BB) Standard	2 vials
Stop Solution	1 x 8ml
TMB Substrate	1 x 12ml

图片



Porcine PDGF-BB ELISA kit (ab218297) Standard curve.

Porcine PDGF-BB ELISA kit (ab218297) 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.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