

Porcine VEGF-A ELISA Kit ab218298

★★★★★ [1 Abreviews](#) [2 References](#) [1 图像](#)

概述

产品名称 Porcine VEGF-A ELISA试剂盒

检测方法 Colorimetric

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 40 pg/ml

范围 40 pg/ml - 10000 pg/ml

特定样本回收率		
样品类型	平均%	范围
Cell culture supernatant	104.8	97% - 113%
Serum	99.39	86% - 112%
Plasma	80.15	67% - 104%

检测时间 4h 45m

实验步骤 Multiple steps standard assay

种属反应性 与反应: Pig

产品概述 The Porcine VEGF-A ELISA kit (ab218298) is an *in vitro* enzyme-linked immunosorbent assay (ELISA) for the quantitative measurement of Porcine VEGF-A in serum (Porcine VEGF-A concentration is low in normal serum and plasma samples and may not be detectable in this assay), plasma and cell culture supernatants.

This assay employs an antibody specific for VEGF-A coated on a 96-well plate. Standards and samples are pipetted into the wells and VEGF-A present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Porcine VEGF-A 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 VEGF-A 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
300X Streptavidin-Conjugated HRP	1 x 200µl
5X Assay Diluent B	1 x 15ml
Porcine VEGF-A coated Microplate (12 x 8 wells)	1 x 96 tests
Assay Diluent C	1 x 30ml
Biotinylated Anti-Pig VEGF-A Detection Antibody	2 vials
Lyophilized Pig VEGF-A Standard	2 vials
Stop Solution	1 x 8ml
TMB Substrate	1 x 12ml

**功能** Growth factor active in angiogenesis, vasculogenesis and endothelial cell growth. Induces endothelial cell proliferation, promotes cell migration, inhibits apoptosis and induces permeabilization of blood vessels. Binds to the FLT1/VEGFR1 and KDR/VEGFR2 receptors, heparan sulfate and heparin. NRP1/Neuropilin-1 binds isoforms VEGF-165 and VEGF-145. Isoform VEGF165B binds to KDR but does not activate downstream signaling pathways, does not activate angiogenesis and inhibits tumor growth.

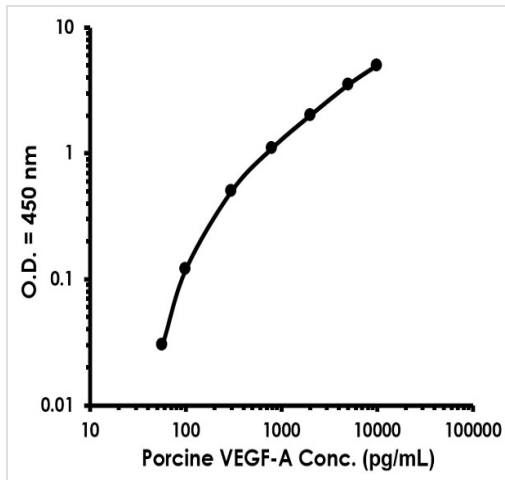
**组织特异性** Isoform VEGF189, isoform VEGF165 and isoform VEGF121 are widely expressed. Isoform VEGF206 and isoform VEGF145 are not widely expressed. A higher level expression seen in pituitary tumors as compared to the pituitary gland.

**疾病相关** Microvascular complications of diabetes 1

**序列相似性** Belongs to the PDGF/VEGF growth factor family.

**细胞定位** Secreted. VEGF121 is acidic and freely secreted. VEGF165 is more basic, has heparin-binding properties and, although a significant proportion remains cell-associated, most is freely secreted. VEGF189 is very basic, it is cell-associated after secretion and is bound avidly by heparin and the extracellular matrix, although it may be released as a soluble form by heparin, heparinase or

## 图片



Porcine VEGF-A ELISA kit (ab218298) Standard curve.

Porcine VEGF-A ELISA kit (ab218298) 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