

Human VEGF Receptor 2 ELISA Kit ab100665

2 References 2 图像

概述

产品名称	人VEGF Receptor 2 ELISA试剂盒
检测方法	Colorimetric
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 70 pg/ml
范围	34.3 pg/ml - 25000 pg/ml
回收率	95 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	96.23	86% - 104%
Serum	96.46	85% - 105%
Plasma	94.78	84% - 103%

实验步骤	Multiple steps standard assay
种属反应性	与反应: Human
产品概述	Abcam's VEGF Receptor 2 Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Human VEGF Receptor 2 in serum, plasma and cell culture supernatants. (Genecards: KDR)

This assay employs an antibody specific for Human VEGF Receptor 2 coated on a 96-well plate. Standards and samples are pipetted into the wells and VEGF Receptor 2 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human VEGF Receptor 2 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 Receptor 2 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

说明	Optimization may be required with urine samples.
平台	Microplate

性能

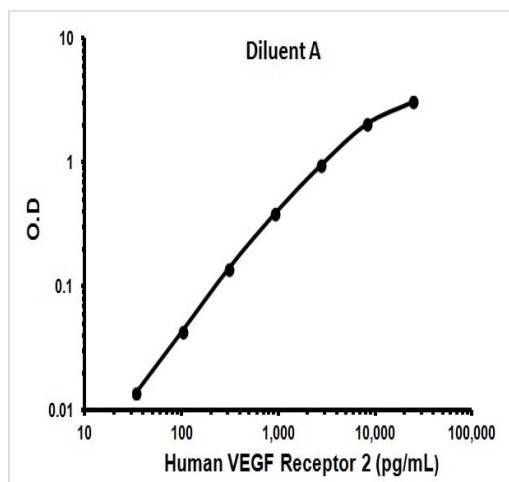
存放说明 Store at -20°C. Please refer to protocols.

组件	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
600X HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent A	1 x 30ml
Biotinylated anti-Human VEGF Receptor 2	2 vials
Recombinant Human VEGF Receptor 2 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml
VEGF Receptor 2 Microplate (12 x 8 wells)	1 unit

相关性 VEGF receptor 2 is a member of a receptor tyrosine kinase family whose activation plays an essential role in a large number of biological processes such as embryonic development, wound healing, cell proliferation, migration and differentiation. Like other growth factor receptors, upon ligand binding VEGF receptor 2 dimerises and is autophosphorylated on multiple tyrosine residues. These sites can be involved in the regulation of kinase activity or serve as binding sites for SH2 and phosphotyrosine binding containing signalling proteins. Phosphorylation of Tyrosines 1054 and 1059 in the activation loop is required for activation of VEGF receptor 2 and its intrinsic tyrosine kinase activity. In case of HIV-1 infection, the interaction with extracellular viral Tat protein seems to enhance angiogenesis in Kaposi's sarcoma lesions.

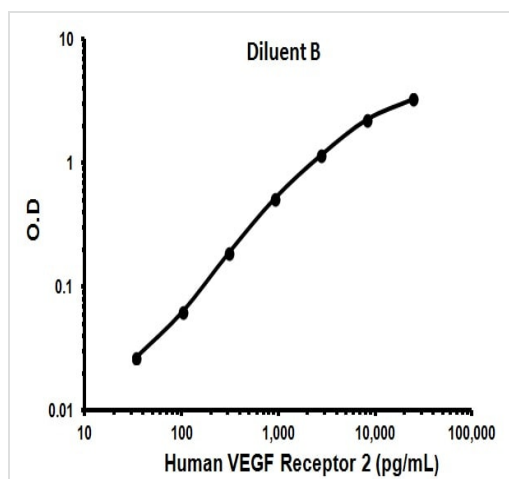
细胞定位 Cell Membrane, Cytoplasmic, Nuclear and Secreted

图片



Representative standard curve using ab100665

Typical standard curve



Representative standard curve using ab100665

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.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