

Human VEGFC ELISA Kit ab100664

1 [References](#) 2 [图像](#)

概述

产品名称	人VEGFC 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	< 15 pg/ml
范围	13.72 pg/ml - 10000 pg/ml
回收率	98 %

特定样本回收率

样品类型	平均 %	范围
Cell culture supernatant	90.46	83% - 99%
Serum	109.4	100% - 118%
Plasma	96.02	86% - 104%

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

产品概述 Abcam's VEGFC Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human VEGFC in serum, plasma, cell culture supernatants.

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

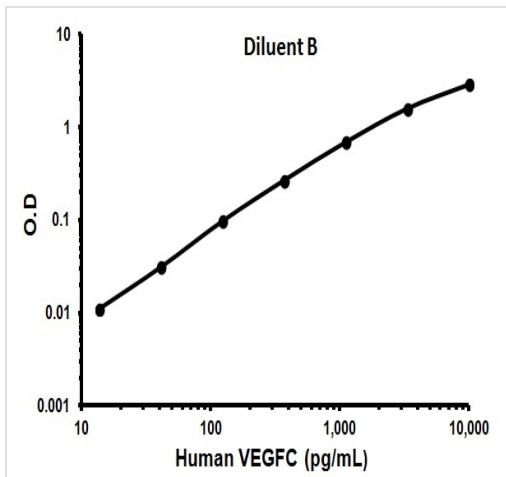
存放说明 Store at -20°C. Please refer to protocols.

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

相关性

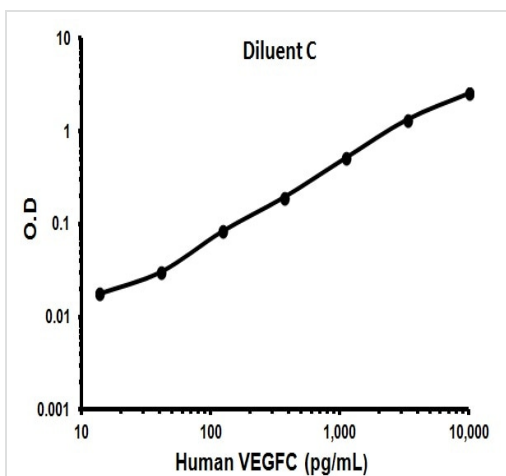
Vascular endothelial growth factors (VEGFs), also known as vasculotropins, are a family of closely related growth factors having a conserved pattern of eight cysteine residues and sharing common VEGF receptors. VEGFs stimulate the proliferation of endothelial cells, induce angiogenesis, promote cell migration, increase vascular permeability, and inhibit apoptosis. The mitogenic activity of VEGFs appears to be mediated by specific VEGF receptors. The target cell specificity of VEGF is restricted to vascular endothelial cells. Vascular Endothelial Growth Factor C (VEGFC) is a member of the VEGF subfamily of PDGF-related growth factors. It is the ligand for Flt4 (VEGFR3) and KDR (VEGFR2). VEGFC binds Flt4 and induces tyrosine autophosphorylation of VEGFR3 and VEGFR2. VEGFC also stimulates the migration of bovine capillary endothelial cells in collagen gel. It is a specific growth factor for the lymphatic vascular system and mediates lymphangiogenesis. VEGFC is abundantly expressed in heart and skeletal muscle. Other tissues such as lung and kidney also express VEGFC.

图片



Representative standard curve using ab100664

Typical standard curve



Representative standard curve using ab100664

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