

Human VEGF ELISA Kit ab100662

30 References 画像数 4

製品の概要

|         |   |
|---------|---|
| 製品名     | Human VEGF ELISA Kit                    |
| 検出方法    | Colorimetric                            |
| サンプルの種類 | Cell culture supernatant, Serum, Plasma |
| アッセイタイプ | Sandwich (quantitative)                 |
| 検出感度    | < 10 pg/ml                              |
| 検出範囲    | 8.23 pg/ml - 6000 pg/ml                 |
| 添加回収試験  | > 100 %                                 |

特定サンプルでの回収試験

| サンプルの種類                  | 平均 %  | 測定範囲       |
|--------------------------|-------|------------|
| Cell culture supernatant | 103.5 | 92% - 113% |
| Serum                    | 104.4 | 92% - 115% |
| Plasma                   | 105.7 | 93% - 114% |

|            |  |
|------------|--|
| ステップ       | Multiple steps standard assay  |
| 種交差性       | 交差種: Human   |
| 製品の概要      | <p>Abcam's VEGF Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Human VEGF in serum, plasma and cell culture supernatants.</p> <p>This assay employs an antibody specific for Human VEGF coated on a 96-well plate. Standards and samples are pipetted into the wells and VEGF present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human VEGF 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 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.</p> |
| 特記事項       | Optimization may be required with urine samples.   |
| 試験プラットフォーム | Microplate   |

## 製品の特性

### 保存方法

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

| 内容  | 1 x 96 tests |
|---|--------------|
| 20X Wash Buffer                               | 1 x 25ml     |
| 300X HRP-Streptavidin Concentrate             | 1 x 200µl    |
| 5X Assay Diluent B                            | 1 x 15ml     |
| Assay Diluent A                               | 1 x 30ml     |
| Biotinylated anti-Human VEGF                  | 2 vials      |
| Recombinant Human VEGF Standard (lyophilized) | 2 vials      |
| Stop Solution                                 | 1 x 8ml      |
| TMB One-Step Substrate Reagent                | 1 x 12ml     |
| VEGF Microplate (12 x 8 wells)                | 1 unit       |

### 機能

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.

### 組織特異性

Isoform VEGF189, isoform VEGF165 and isoform VEGF121 are widely expressed. Isoform VEGF206 and isoform VEGF145 are not widely expressed.

### 関連疾患

Defects in VEGFA are a cause of susceptibility to microvascular complications of diabetes type 1 (MVCD1) [MIM:603933]. These are pathological conditions that develop in numerous tissues and organs as a consequence of diabetes mellitus. They include diabetic retinopathy, diabetic nephropathy leading to end-stage renal disease, and diabetic neuropathy. Diabetic retinopathy remains the major cause of new-onset blindness among diabetic adults. It is characterized by vascular permeability and increased tissue ischemia and angiogenesis.

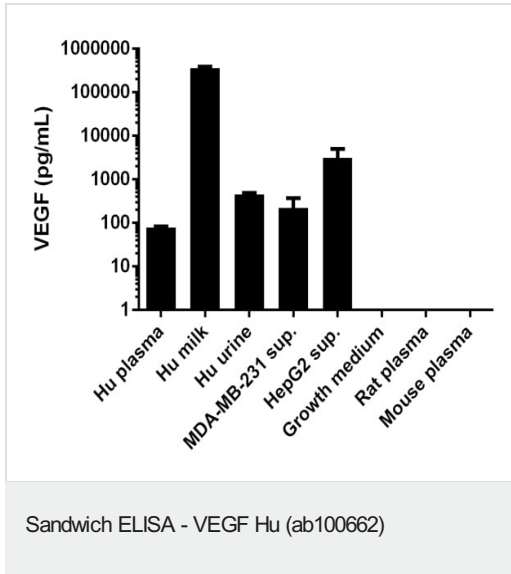
### 配列類似性

Belongs to the PDGF/VEGF growth factor family.

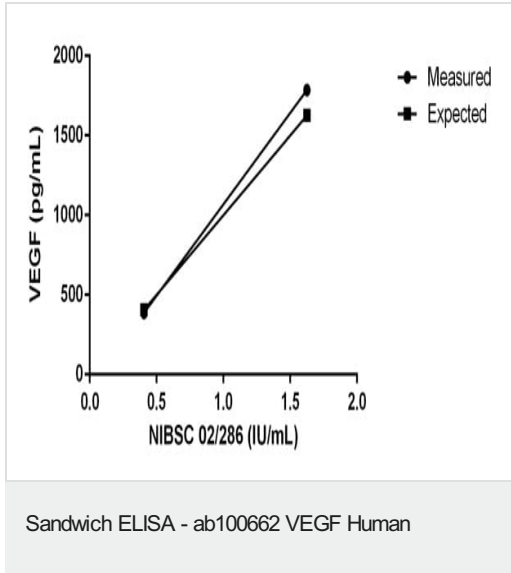
### 細胞内局在

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 plasmin.

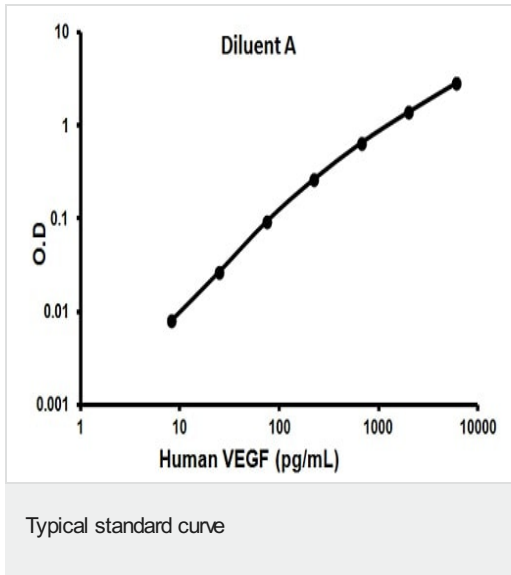
## 画像



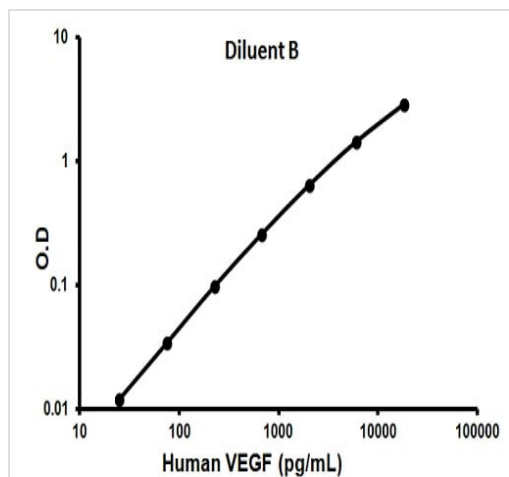
VEGF measured in biological fluids showing quantity (pg) per milliliter of sample tested



Curves showing concentrations huVEGF165 in NIBSC 02/286 using ab100662 VEGF human



Representative standard curve using ab100662



Representative standard curve using ab100662

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.co.jp/abpromise> or contact our technical team.

### Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors