

Human VEGF-165 ELISA Kit ab273164

画像数 1

医薬用外毒物

製品の概要

製品名	Human VEGF-165 ELISA Kit				
検出方法	Colorimetric				
再現性	Intra-Assay (同時再現性)				
	サンプル	N	平均値	SD	CV%
	Various				3.8%
	Inter-Assay (日差再現性)				
	サンプル	N	平均値	SD	CV%
	Various				7.6%
サンプルの種類	Cell culture supernatant, Serum, Plasma, Other biological fluids, Hep Plasma, EDTA Plasma, Cit plasma				
アッセイタイプ	Sandwich (quantitative)				
検出感度	15.9 pg/ml				
検出範囲	31.25 pg/ml - 1000 pg/ml				
ステップ	Multiple steps standard assay				
種交差性	交差種: Human				
製品の概要	Human VEGF <sub>165</sub> ELISA kit (ab273164) is for the in-vitro qualitative and quantitative determination of VEGF <sub>165</sub> in supernatants, buffered solutions or serum and plasma samples. This assay will recognise both natural and recombinant human VEGF <sub>165</sub> .				

A capture antibody highly specific for VEGF-A has been coated to the wells of the microtiter strip plate provided during manufacture. Binding of VEGF165 in samples and known standards to the capture antibodies is completed and then any excess unbound analyte is removed.

During the next incubation period the binding of the biotinylated anti-VEGF165 secondary antibody to the analyte occurs. Any excess unbound secondary antibody is then removed.

The HRP conjugate solution is then added to every well including the zero wells, following incubation excess conjugate is removed by careful washing.

A chromogen substrate is added to the wells resulting in the progressive development of a blue coloured complex with the conjugate. The colour development is then stopped by the addition of acid turning the resultant final product yellow. The intensity of the produced coloured complex is directly proportional to the concentration of VEGF165 present in the samples and standards.

The absorbance of the colour complex is measured and the generated OD values for each standard are plotted against expected concentration forming a standard curve. This standard curve can then be used to accurately determine the concentration of VEGF165 in any sample tested.

#### 試験プラットフォーム

Microplate (12 x 8 well strips)

#### 製品の特性

##### 保存方法

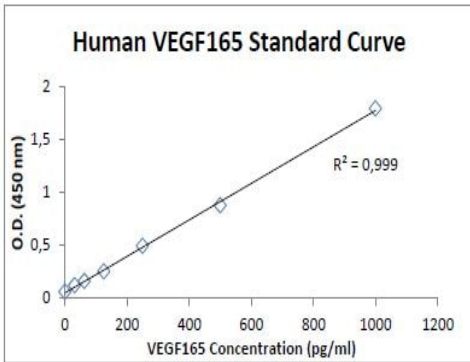
Store at +4°C. Please refer to protocols.

内容	1 x 96 tests
Biotinylated Antibody diluent	1 x 7ml
Biotinylated Antibody VEGF-165	1 x 400µl
HRP Diluent	1 x 12ml
Plastic Plate Cover	2 units
Standard diluent Buffer	1 x 15ml
Stop Reagent	1 x 11ml
Streptavidin-HRP	2 x 5µl
TMB Substrate	1 x 11ml
VEGF-165 Coated Microwell strips	1 unit
VEGF-165 Standard	2 vials
Wash Buffer	1 x 10ml

#### 細胞内局在

VEGF 165A: 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



Example data

**Typical standard curve** – data provided **for demonstration purposes only**. A new standard curve must be generated for each assay performed.

Example of HumanVEGF<sub>165</sub> ELISA kit 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