

## Product datasheet

# Mouse VEGFA Matched Antibody Pair Kit ab220118

Recombinant

## 2 Images

### Overview

<b>Product name</b>	Mouse VEGFA Matched Antibody Pair Kit
<b>Detection method</b>	Colorimetric
<b>Assay type</b>	ELISA set
<b>Sensitivity</b>	1.42 pg/ml
<b>Range</b>	7.81 pg/ml - 500 pg/ml
<b>Species reactivity</b>	<b>Reacts with:</b> Mouse
<b>Product overview</b>	Mouse VEGFA Matched Antibody Pair Kits include a capture and a biotinylated detector antibody pair, along with a calibrated protein standard, suitable for sandwich ELISA. The Matched Antibody Pair Kit can be used to quantify native and recombinant mouse VEGFA.

Optimization of the kit reagents to sample type, immunoassay format or instrumentation may be required. Guidelines for use of this kit in a standard 96-well microplate sandwich ELISA using HRP/TMB system of colorimetric detection is described in this assay procedure for the purposes of quantification.

Protocol information and tips on the use of the Matched Antibody Pair kits for sandwich ELISA can be found on our [website](#). An accessory pack can be purchased which includes buffer reagents required to perform 10 x 96-well plate sandwich ELISAs ([ab210905](#)).

For additional information on the performance of the antibody pair used in this kit, please see our equivalent SimpleStep ELISA kit [ab209882](#). Please note that while the antibody pair is the same provided in the corresponding SimpleStep ELISA Kit, due to differences in their formulation, this antibody pair cannot be used with the consumables provided with our SimpleStep ELISA Kits.

### Tested applications

**Suitable for:** ELISA

### Platform

Reagents

### Properties

#### Storage instructions

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

Components	
Mouse VEGFA Capture Antibody	2 x 50µg
Mouse VEGFA Detector Antibody	2 x 12.5µg
Mouse VEGFA Lyophilized Protein	2 x 1 vial
Mouse VEGFA Capture Antibody	1 x 50µg
Mouse VEGFA Detector Antibody	1 x 12.5µg
Mouse VEGFA Lyophilized Protein	1 x 1 vial

<b>Function</b>	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.
<b>Tissue specificity</b>	Isoform VEGF189, isoform VEGF165 and isoform VEGF121 are widely expressed. Isoform VEGF206 and isoform VEGF145 are not widely expressed.
<b>Involvement in disease</b>	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.
<b>Sequence similarities</b>	Belongs to the PDGF/VEGF growth factor family.
<b>Cellular localization</b>	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.

## Applications

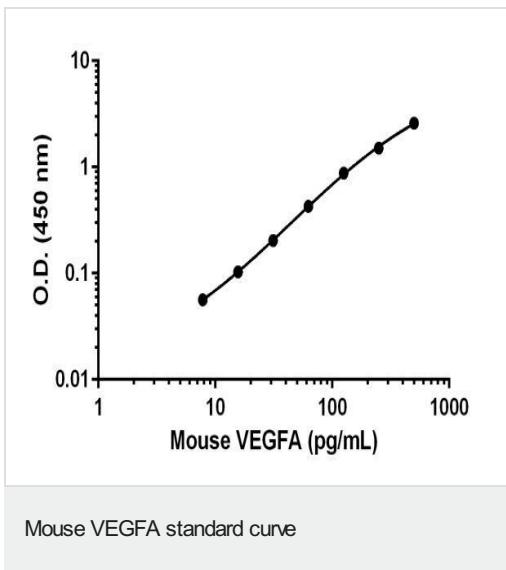
### The Abpromise guarantee

Our **Abpromise guarantee** covers the use of ab220118 in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
ELISA		Use at an assay dependent concentration.

## Images



Standard calibration curve. Background subtracted values are graphed.



To learn more about the advantages of recombinant antibodies see [here](#).

**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.com/abpromise> or contact our technical team.

## **Terms and conditions**

---

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