

Product datasheet

Mouse VEGF R3 Matched Antibody Pair Kit ab221433

Recombinant

2 Images

Overview

Product name	Mouse VEGF R3 Matched Antibody Pair Kit
Detection method	Colorimetric
Assay type	ELISA set
Sensitivity	31.47 pg/ml
Range	125 pg/ml - 8000 pg/ml
Species reactivity	Reacts with: Mouse
Product overview	<p>Mouse VEGF R3 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 VEGF R3.</p> <p>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.</p> <p>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).</p> <p>For additional information on the performance of the antibody pair used in this kit, please see our equivalent SimpleStep ELISA kit ab203367. 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.</p>
Tested applications	Suitable for: ELISA
Platform	Reagents

Properties

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

Components	10 x 96 tests	5 x 96 tests
Mouse VEGF R3 Capture Antibody	2 x 50µg	1 x 50µg
Mouse VEGF R3 Detector Antibody	2 x 12.5µg	1 x 12.5µg
Mouse VEGF R3 Lyophilized Protein	2 vials	1 vial

Function	Receptor for VEGFC. Has a tyrosine-protein kinase activity.
Tissue specificity	Placenta, lung, heart, and kidney, does not seem to be expressed in pancreas and brain.
Involvement in disease	Defects in FLT4 are the cause of lymphedema hereditary type 1A (LMPH1A) [MIM:153100]; also known as Nonne-Milroy lymphedema or Milroy disease. Hereditary lymphedema is a chronic disabling condition which results in swelling of the extremities due to altered lymphatic flow. Patients with lymphedema suffer from recurrent local infections and physical impairment. Note=Defects in FLT4 are found in juvenile hemangioma. Juvenile hemangiomas are the most common tumors of infancy, occurring as many as 10% of all births. These benign vascular lesions enlarge rapidly during the first year of life by hyperplasia of endothelial cells and attendant pericytes, and then spontaneously involute over a period of years, leaving loose fibrofatty tissue.
Sequence similarities	Belongs to the protein kinase superfamily. Tyr protein kinase family. CSF-1/PDGF receptor subfamily. Contains 7 Ig-like C2-type (immunoglobulin-like) domains. Contains 1 protein kinase domain.
Cellular localization	Membrane.

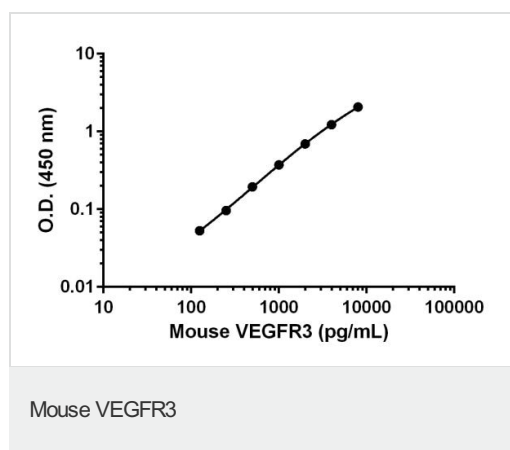
Applications

Our [Abpromise guarantee](#) covers the use of **ab221433** 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.

Powered by
recombinant antibodies



Research with confidence
Consistent and reproducible results



Long-term and scalable supply
Recombinant technology



Success from the first experiment
Confirmed specificity



Ethical standards compliant
Animal-free production

Sandwich ELISA - Mouse VEGF R3 Matched
Antibody Pair Kit (ab221433)

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