

Version 1c Last updated 21 April 2026

ab264501

Ramucirumab ELISA Kit

For the measurement of Ramucirumab in human serum and plasma.

This product is for research use only and is not intended for diagnostic use.

Table of Contents

1. Overview	1
2. Protocol Summary	2
3. Precautions	3
4. Storage and Stability	3
5. Limitations	3
6. Materials Supplied	4
7. Materials Required, Not Supplied	4
8. Technical Hints	5
9. Reagent Preparation	6
10. Standard and Control Preparation	6
11. Sample Preparation	7
12. Assay Procedure	8
13. Calculations	9
14. Typical Data	10
15. Typical Sample Values	10
16. Notes	11

1. Overview

Ramucirumab ELISA Kit (ab264501) is a highly specific and sensitive kit designed for the *in vitro* determination of Ramucirumab in biological matrices such as human serum and plasma. The density of color is proportional to the amount of Ramucirumab captured from the samples and can be quantified when compared with standard curve.

Ramucirumab is a human monoclonal antibody (IgG1) against vascular endothelial growth factor receptor 2 (VEGFR2), a type II trans-membrane tyrosine kinase receptor expressed on endothelial cells. By binding to VEGFR2, Ramucirumab prevents binding of its ligands (VEGF-A, VEGF-C, and VEGF-D), thereby preventing VEGF-stimulated receptor phosphorylation and downstream ligand-induced proliferation, permeability, and migration of human endothelial cells.

2. Protocol Summary

Prepare all reagents, samples, and standards as instructed



Add 50 μ L Assay Buffer to each well.



Add 50 μ L standard or sample to appropriate wells. Cover and incubate for 30 minutes at room temperature



Discard incubation solution and wash plate 3 times with 300 μ L diluted Wash Buffer



Add 100 μ L HRP-conjugate to each well. Cover and incubate for 30 minutes at room temperature



Discard the solution and wash plate 3 times with 300 μ L diluted Wash Buffer



Add 100 μ L 1X TMB Substrate and incubate the plate in the dark at room temperature for 10 minutes.



Add 100 μ L Stop Solution and read OD at 450 nm within 30 minutes.

3. Precautions

Please read these instructions carefully prior to beginning the assay.

- Observe good laboratory practices. Gloves, lab coat, and protective eyewear should always be worn. Never pipet by mouth. Do not eat, drink or smoke in the laboratory areas.
- For general guidelines, precautions, limitations on the use of our assay kits and general assay troubleshooting tips, particularly for first time users, please consult our guide: <https://www.abcam.com/en-us/products/biochemical-assays>
- If applicable, please refer to the current Safety Data Sheet (SDS) provided with this product for safety, handling, and disposal information. The most up to date and current versions are available on our website <https://www.abcam.com/en-us>.

4. Storage and Stability

Store kit at +4°C immediately upon receipt.

Refer to list of materials supplied for storage conditions of individual components.

5. Limitations

- Assay kit intended for research use only. Not for use in diagnostic procedures.
- Do not mix or substitute reagents or materials from other kit lots or vendors.

6. Materials Supplied

Item	Quantity	Storage condition
Micro ELISA Plate	1 unit	+4°C
Ramucirumab Standard S1	0.3 mL	+4°C
Ramucirumab Standard S2	0.3 mL	+4°C
Ramucirumab Standard S3	0.3 mL	+4°C
Ramucirumab Standard S4	0.3 mL	+4°C
Ramucirumab Standard S5	0.3 mL	+4°C
Ramucirumab Standard S6	0.3 mL	+4°C
Ramucirumab Standard S7	0.3 mL	+4°C
Assay Buffer	50 mL	+4°C
HRP-conjugate Probe	12 mL	+4°C
TMB Substrate	12 mL	+4°C
Stop Solution	12 mL	+4°C
Wash Buffer (20X)	50 mL	+4°C
Plate sealers	2 units	+4°C

7. Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully perform this assay:

- Microplate reader capable of measuring absorbance at OD 450 nm
- Deionized water.
- Multi- and single-channel pipettes.
- Tubes for sample dilution.
- Plate shaker for all incubation steps.
- Absorbent paper

8. Technical Hints

- Samples generating values higher than the highest standard should be further diluted.
- Avoid foaming or bubbles when mixing or reconstituting components.
- Avoid cross contamination of samples or reagents by changing tips between sample, standard and reagent additions.
- Ensure plates are properly sealed or covered during incubation steps.
- Complete removal of all solutions and buffers during wash steps is necessary to minimize background.
- All samples should be mixed thoroughly and gently.
- Avoid multiple freeze/thaw of samples.
- Incubate ELISA plates on a plate shaker during all incubation steps.
- When generating positive control samples, it is advisable to change pipette tips after each step.

9. Reagent Preparation

- Equilibrate all reagents to room temperature (18-25°C) prior to use. Before using the kit, spin tubes and bring down all components to the bottom of tubes.
- Prepare only as much reagent as is needed on the day of the experiment.

9.1 20X Wash Buffer:

Dilute the 20X Wash Buffer to 1X solution in ddH₂O (10 mL of Wash Buffer stock to 190 mL of ddH₂O). Mix the 1X solution thoroughly by vortex manually. The working stock can be stable for 2 weeks after preparation at 4°C.

10. Standard and Control Preparation

Standard and controls, S1 – S7, are ready to use, please see table below for concentrations:

Name	S1	S2	S3	S4	S5	S6	S7
Conc. ng/mL	600	200	60	20	0	High control	Low control

Concentration for high and low controls are indicated on vials.

11. Sample Preparation

General sample information:

- We recommend performing several dilutions of your sample to ensure the readings are within the standard value range.
- We recommend that you use fresh samples for the most reproducible assay.

11.1 Serum/plasma:

1. First dilute samples at 1:10 (10 μ L serum/plasma + 90 μ L ddH₂O).
2. Then dilute samples 1:100 (5 μ L diluted sample + 495 μ L ddH₂O) (Dilution factor of 1000).
3. Diluted samples should further be diluted if the concentration of Ramucirumab is higher than the measuring range.
4. Samples are stable at 4°C for 2 days and -20°C for 6 months. Avoid freeze-and-thaw cycle.

▲ Note: The usual precautions for venipuncture should be observed.

12. Assay Procedure

- Prepare reagents within 30 minutes before the experiment.
- Equilibrate all materials and prepared reagents to room temperature 15 minutes prior to use.
- We recommend that you assay all standards, controls and samples in duplicate.
- A standard curve must be run with each assay.

- 12.1** Add 50 μL of Assay Buffer to each well. Then add 50 μL standards and diluted-samples into appropriate wells. Cover wells and incubate for 30 minutes at room temperature (RT).
- 12.2** Discard incubation solution. Wash plate 3 times each with 300 μL of diluted Wash Buffer. Remove excess solution by tapping the inverted plate on a paper towel.
- 12.3** Add 100 μL of HRP-conjugate into each well. Cover wells with adhesive plate sealer and incubate at RT for 30 minutes.
- 12.4** Discard the solution and wash the wells as step 12.2.
- 12.5** Add 100 μL of 1X TMB substrate solution and incubate the plate in dark at RT for 10 minutes.
- 12.6** Add 100 μL of Stop solution to stop the reaction.
- 12.7** Read the absorbance in micro plate reader set to 450 nm within 30 minutes. (reference wavelength is 650 nm).

13. Calculations

- 13.1 Calculate the average absorbance value for the blank control (0 ng/mL) standards. Subtract the average blank control standard absorbance value from all other absorbance values.
- 13.2 **Create a standard curve** by plotting the average blank control subtracted absorbance value for each standard concentration (y axis) against the target protein concentration (x axis) of the standard.
- 13.3 Construct a standard curve of difference data using software capable of generating four-parameter logistic (4PL) or point-to-point calculation curve fit.
- 13.4 To obtain the exact values of the samples, the concentration determined from the standard curve should be multiplied by the dilution factor.
- 13.5 Determine the concentration of the target protein in the sample by interpolating the blank control subtracted **absorbance values against the standard curve**. Multiply the resulting value by the appropriate dilution factor to obtain the concentration of target protein in the sample.
- 13.6 Samples generating absorbance values greater than that of the highest standard should be further diluted and reanalyzed. Similarly, samples which measure at an absorbance values less than that of the lowest standard should be retested in a less dilute form.

14. Typical Data

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

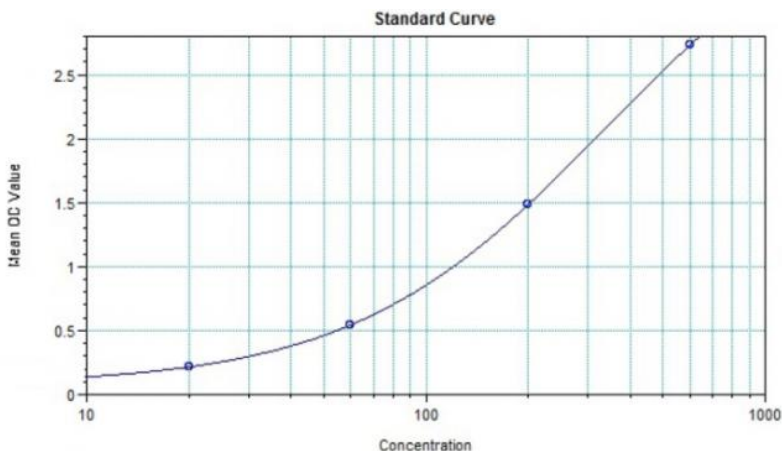


Figure 1. Typical Standard Curve: This standard curve is for demonstration only. A standard curve must be run with each assay.

15. Typical Sample Values

Detection Range: 20 - 600 ng/mL.

Sensitivity: 20 ng/mL.

Assay Precision: Intra-Assay: CV < 15%; Inter-Assay: CV < 15%

(CV (%) = SD/mean X 100)

Cross Reactivity: No significant cross-reactivity or interference with other proteins present in native human serum or other therapeutic immunoglobulins.

Recovery rate: 85 – 115% with normal human serum samples with known concentrations.

16. Notes

Technical Support

Copyright © 2026 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

<https://www.abcam.com/en-us/contact-us>

<https://www.abcam.cn/contact-us> (China)

<https://www.abcam.co.jp/contact-us> (Japan)