

ab284009 – Human KAP1 ELISA Kit

For the quantitative measurement of KAP1 in human plasma, serum, saliva, cell culture, cell lysate, and tissue samples.

For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit:

<http://www.abcam.com/ab284009>

Storage and Stability: Store kit at +4°C immediately upon receipt, apart from the Standard, Biotinylated Antibody and SP Conjugate which should be stored at -20°C.

Materials Supplied

Item	Quantity	Storage Condition
100X Streptavidin-Peroxidase Conjugate	80 µL	-20°C
10X Diluent N Concentrate	30 mL	4°C
1X Standard Diluent	2 mL	4°C
20X Wash Buffer Concentrate	2 x 30 mL	4°C
KAP1 Standard	1 vial	-20°C
Chromogen Substrate	7 mL	4°C
Human KAP1 Microplate	96 wells	4°C
Sealing Tapes	3	4°C
Stop Solution	11 mL	4°C
50X Biotinylated Human KAP1 Antibody	1 vial	-20°C

Materials Required, Not Supplied

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

- Microplate reader capable of measuring absorbance at 450 nm.
- Pipettes (1-20 µL, 20-200 µL, 200-1000 µL, and multiple channel).
- Deionized or distilled reagent grade water.

Reagents Preparation

Equilibrate all reagents to room temperature (18-25°C) prior to use. The kit contains enough reagents for 96 wells.

Prepare only as much reagent as is needed on the day of the experiment.

If crystals have formed in the concentrate, mix gently until the crystals have completely dissolved.

Δ Note: Concentration of the kit components are lot-specific, and the end user should always refer to the vial label.

10X Diluent N Concentrate: If crystals have formed in the concentrate, mix gently until the crystals have completely dissolved. Dilute the Diluent N Concentrate 1:10 with reagent grade water to produce a 1x solution. Store for up to 30 days at 2-8°C.

50X Biotinylated Human KAP1 Antibody: Spin down the antibody briefly and dilute the desired amount of the antibody 1:50 with 1X Diluent N to produce a 1x solution. The undiluted antibody should be stored at -20°C.

20X Wash Buffer Concentrate: If crystals have formed in the concentrate, mix gently until the crystals have completely dissolved. Dilute the Wash Buffer Concentrate 1:20 with reagent grade water to produce a 1x solution.

100X Streptavidin-Peroxidase Conjugate: Spin down the SP Conjugate briefly and dilute the desired amount of the conjugate 1:100 with Diluent N to produce a 1x solution. The undiluted conjugate should be stored at -20°C.

Standard Preparation

Prepare serially diluted standards immediately prior to use. Always prepare a fresh set of standards for every use.

Any remaining standard should be stored at -20°C after reconstitution and used within 48 hours. This procedure prepares sufficient standard dilutions for duplicate wells.

- 1) Reconstitution of the KAP1 Standard vial with 1X Standard Diluent to prepare the 100 ng/mL Factor XI **Stock Standard**.

1.1) First consult the KAP1 Standard vial to determine the mass of protein in the vial.

1.2) Calculate the appropriate volume of 1X Standard Diluent to add when resuspending the KAP1 Standard vial to produce a 100 ng/mL KAP1 **Stock Standard** by using the following equation:

C_s = Starting mass of KAP1 Standard (see vial label) (ng)

C_f = 100 ng/mL KAP1 **Stock Standard** final required concentration

V_d = Required volume of Standard Diluent for reconstitution (µL)

Calculate total required volume Standard Diluent for resuspension:

$$(C_s / C_f) \times 1,000 = V_d$$

Example:

NOTE: This example is for demonstration purposes only. Please remember to check your standard vial for the actual amount of standard provided.

C_s = 50 ng of KAP1 Standard in vial

C_f = 100 ng/mL KAP1 **Stock Standard** final concentration

V_d = Required volume of 1X Standard Diluent for reconstitution

$$(50 \text{ ng} / 100 \text{ ng/mL}) \times 1,000 = 500 \text{ µL}$$

1.3) First briefly spin the KAP1 Standard Vial to collect the contents on the bottom of the tube.

1.4) Reconstitute the KAP1 Standard vial by adding the appropriate calculated amount V_d of 1X Standard Diluent to the vial to generate the 100 ng/mL Factor XI **Stock Standard**. Mix gently and thoroughly.

- 2) Allow the reconstituted 100 ng/mL KAP1 **Stock Standard** to sit for 10 minutes with gentle agitation prior to making subsequent dilutions
- 3) Label tubes #1 – 8.
- 4) Add 360 µL of **1X Diluent N** to tube #1.
- 5) Prepare the 25 ng/mL **Standard #1** by adding 120 µL of 100 ng/mL **Stock Standard** into tube #1. Mix thoroughly and gently.
- 6) Add 120 µL of **1X Diluent N** to tubes #2 – 8.
- 7) To prepare Standard #2, add 120 µL of the Standard #1 into tube #2 and mix gently.
- 8) To prepare Standard #3, add 120 µL of the Standard #2 into tube #3 and mix gently.
- 9) Using the table below as a guide, prepare subsequent serial dilutions. 1X Diluent N serves as the zero standard (0 ng/mL)

Standard #	Volume to dilute (µL)	Volume 1X Diluent N (µL)	KAP1 (ng/ml)
1	120 (Stock Standard)	360	25.00
2	120 (Standard #1)	120	12.50
3	120 (Standard #2)	120	6.250
4	120 (Standard #3)	120	3.125
5	120 (Standard #4)	120	1.563
6	120 (Standard #5)	120	0.781
7	120 (Standard #6)	120	0.391
8	-	120	0.0

Sample Preparation

Plasma: Collect plasma using one-tenth volume of 0.1 M sodium citrate as an anticoagulant. Centrifuge samples at 3000 x g for 10 minutes and collect plasma. A 4-fold sample dilution is suggested into Diluent N or within the range of 1x – 10x; however, user should determine optimal dilution factor depending on application needs. The undiluted samples can be stored at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles (EDTA or Heparin can also be used as an anticoagulant).

Serum: Samples should be collected into a serum separator tube. After clot formation, centrifuge samples at 3000 x g for 10 minutes and remove serum. A 4-fold sample dilution is suggested into Diluent N or within the range of 1x – 10x; however, user should determine optimal dilution factor depending on application needs. The undiluted samples can be stored at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.

Saliva: Collect saliva using sample tube. Centrifuge samples at 800 x g for 10 minutes. The sample is suggested for use at 1x; however, user should determine optimal dilution factor depending on application needs. The undiluted samples can be stored at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.

Cell Culture Supernatants: Centrifuge cell culture media at 1500 rpm for 10 minutes at 4°C to remove debris and collect supernatants. Samples can be stored at -80°C. Avoid repeated freeze-thaw cycles.

Cell Lysate: Rinse cell with cold PBS and then scrape the cell into a tube with 5 ml of cold PBS and 0.5 M EDTA. Centrifuge suspension at 1500 rpm for 10 minutes at 4°C and aspirate supernatant. Re-suspend pellet in ice-cold Lysis Buffer (10 mM Tris, pH8.0, 130 mM NaCl, 1% Triton X-100, protease inhibitor cocktail). For every 1 x 10⁶ cells, add approximately 100 µL of ice-cold Lysis Buffer. Incubate on ice for 60 minutes. Centrifuge at 13000 rpm for 30 minutes at 4°C and collect supernatant.

Tissue: Extract tissue samples with 0.1 M phosphate-buffered saline (pH 7.4) containing 1% Triton X-100 and centrifuge at 14000 x g for 20 minutes. Collect the supernatant and measure the protein concentration. If necessary, dilute samples into Diluent N; user should determine optimal dilution factor depending on application needs. Store remaining extract at -80°C. Avoid repeated freeze-thaw cycles.

Plate Preparation

The 96 well plate strips included with this kit are supplied ready to use. It is not necessary to rinse the plate prior to adding reagents. Unused plate strips should be immediately returned to the foil pouch containing the desiccant pack, resealed and stored at 4°C.

For each assay performed, a minimum of two wells must be used as the zero control. For statistical reasons, we recommend each sample should be assayed with a minimum of two replicates (duplicates).

Assay Procedure

Equilibrate all materials and prepared reagents to room temperature prior to use. We recommend that you assay all standards, controls and samples in duplicate.

1. Prepare all reagents, standard solutions, and samples as instructed. Bring all reagents to room temperature before use. The assay is performed at room temperature (20-25°C).
2. Remove excess microplate strips from the plate frame and return them immediately to the foil pouch with desiccants inside. Reseal the pouch securely to minimize exposure to water vapor and store in a vacuum desiccator.
3. Add 50 µl of KAP1 Standard or sample to each well. Gently tap plate to thoroughly coat the wells. Break any bubbles that may have formed. Cover wells with a sealing tape and incubate for 2 hours. Start the timer after the last addition.
4. Wash five times with 200 µl of Wash Buffer manually. Invert the plate each time and decant the contents; hit 4-5 times on absorbent material to completely remove the liquid. If using a machine, wash six times with 300 µl of Wash Buffer and then invert the plate, decanting the contents; hit 4-5 times on absorbent material to completely remove the liquid.
5. Add 50 µl of Biotinylated Human KAP1 Antibody to each well. Gently tap plate to thoroughly coat the wells. Break any bubbles that may have formed. Cover wells with a sealing tape and incubate for 2 hours.
6. Wash the microplate as described above.
7. Add 50 µl of SP Conjugate to each well. Gently tap plate to thoroughly coat the wells. Break any bubbles that may have formed. Cover wells with a sealing tape and incubate for 30 minutes. Turn on the microplate reader and set up the program in advance.
8. Wash the microplate as described above.
9. Add 50 µl of Chromogen Substrate to each well. Gently tap plate to thoroughly coat the wells. Break any bubbles that may have formed. Incubate for 20 minutes or until the optimal blue colour density develops.
10. Add 50 µl of Stop Solution to each well. The colour will change from blue to yellow. Gently tap plate to ensure thorough mixing. Break any bubbles that may have formed.
11. Read the absorbance on a microplate reader at a wavelength of 450 nm immediately. If wavelength correction is available, subtract readings at 570 nm from those at 450 nm to correct optical imperfections. Otherwise, read the plate at 450 nm only. Please note that some unstable black particles may be generated at high concentration points after stopping the reaction for about 10 minutes, which will reduce the readings.

Calculation

1. Calculate the mean value of the duplicate or triplicate readings for each standard and sample.
2. To generate a standard curve, plot the graph using the standard concentrations on the x-axis and the corresponding mean 450 nm absorbance (OD) on the y-axis. The best fit line can be determined by regression analysis using log-log or four-parameter logistic curve fit.
3. Determine the unknown sample concentration from the Standard Curve and multiply the value by the dilution factor.

PERFORMANCE CHARACTERISTICS

- 1. This assay recognizes both natural and recombinant human KAP-1.
- 2. The minimum detectable dose of human KAP1 as calculated by 2SD from the mean of a zero standard was established to be 0.25 ng/ml.
- 3. Intra-assay precision was determined by testing three samples twenty times in one assay.
- 4. Inter-assay precision was determined by testing three samples in twenty assays.

	Intra-Assay Precision	Inter-Assay Precision
CV (%)	4.9	10.2

RECOVERY

Standard Added Value	0.781 - 12.5 ng/ml
Recovery %	90 - 108
Average Recovery %	97

CROSS REACTIVITY

Species	Cross-Reactivity (%)
Canine	30
Bovine	None
Monkey	80
Mouse	10
Rat	50
Swine	50
Rabbit	None
Protein	Cross-Reactive (%)
Livin	None
PPP1R14A	None
PPP1R3B	None
Ro52	None

Download our ELISA guide for technical hints, results, calculation, and troubleshooting tips:
www.abcam.com/protocols/the-complete-elisa-guide

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)