



# KNG1 (Human) ELISA Kit

Catalog Number KA1040

96 assays

Version: 14

Intended for research use only

[www.abnova.com](http://www.abnova.com)

## Table of Contents

<b>Introduction .....</b>	<b>3</b>
Background .....	3
Principle of the Assay .....	3
<b>General Information .....</b>	<b>4</b>
Materials Supplied .....	4
Storage Instruction .....	4
Materials Required but Not Supplied .....	4
Precautions for Use .....	5
<b>Assay Protocol .....</b>	<b>6</b>
Reagent Preparation .....	6
Sample Preparation .....	6
Assay Procedure .....	7
<b>Data Analysis.....</b>	<b>9</b>
Calculation of Results.....	9
Performance Characteristics .....	10
<b>Resources.....</b>	<b>12</b>
Troubleshooting.....	12
References .....	13
Plate Layout .....	14

## **Introduction**

### **Background**

High molecular weight kininogen (HK) is a plasma protein coagulation cofactor serving for the activation of zymogens (prekallikrein, factor XI, and factor XII) and is a substrate of each of their proteolytic forms. It circulates as a complex with these zymogens and links the plasma coagulation, fibrinolysis, complement activation, and blood pressure control. HK is produced by the liver and weighs 120 kDa with 626 amino acids (1-5). HK exhibits anticoagulant properties and is a strong inhibitor of cysteine proteases. Upon cleavage by kallikrein, the released active peptide, bradykinin, mediates NO release, vasodilation, hypotension, and pain. The remaining cleaved HK (HKa) exhibits antiadhesive and antiangiogenic activity, enhancing cell-associated fibrinolysis and releasing cytokines and chemokines to enhance inflammation. Patients with HK deficiency exhibit abnormal surface-mediated activation of fibrinolysis (6-7).

### **Principle of the Assay**

The KNG1 (Human) ELISA Kit is designed for detection of kininogen in human plasma and serum samples. This assay employs a quantitative competitive enzyme immunoassay technique that measures human kininogen in approximately 3 hours. A polyclonal antibody specific for human kininogen has been pre-coated onto a 96-well microplate with removable strips. Kininogen in standards and samples is competed with a biotinylated human kininogen protein sandwiched by the immobilized antibody and streptavidin-peroxidase (SP) conjugate. All unbound material is washed away and a peroxidase enzyme substrate is added. The color development is stopped and the intensity of the color is measured.

## General Information

### Materials Supplied

List of component

Component	Amount
Human Kininogen (HMW) Microplate: A 96-well polystyrene microplate (12 strips of 8 wells) coated with a polyclonal antibody against human kininogen.	96 (8x12) wells
Sealing Tapes: Each kit contains 3 precut, pressure sensitive sealing tapes that can be cut to fit the format of the individual assay.	3 slices
Human Kininogen (HMW) Standard: Human kininogen in a buffered protein base, lyophilized.	1.8 µg/vial, 2 vials
Biotinylated Human Kininogen (HMW) Protein (1x): Lyophilized.	2 vials
MIX Diluent Concentrate (10x): A 10-fold concentrated buffered protein base.	30 mL
Wash Buffer Concentrate (20x): A 20-fold concentrated buffered surfactant.	30 mL
SP Conjugate (100x): A 100-fold concentrate.	80 µL
Chromogen Substrate (1x): A stabilized peroxidase chromogen substrate tetramethylbenzidine.	7 mL
Stop Solution (1x): A 0.5 N hydrochloric acid solution to stop the chromogen substrate reaction.	11 mL

### Storage Instruction

- ✓ Upon arrival, immediately store components of the kit at recommended temperatures up to the expiration date.
- ✓ Store SP Conjugate at -20°C.
- ✓ Store Microplate, Diluent Concentrate (10x), Wash Buffer, Stop Solution, and Chromogen Substrate at 2-8°C.
- ✓ Unused microplate wells may be returned to the foil pouch with the desiccant packs and resealed. May be stored for up to 30 days in a vacuum desiccator.
- ✓ Store Standard and Biotinylated Protein at 2-8°C before reconstituting with Diluent.

### Materials Required but Not Supplied

- ✓ 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

**Precautions for Use**

- ✓ This product is for Research Use Only and is not intended for use in diagnostic procedures.
- ✓ Prepare all reagents (diluent buffer, wash buffer, standard, biotinylated protein, and SP conjugate) as instructed, prior to running the assay.
- ✓ Prepare all samples prior to running the assay. The dilution factors for the samples are suggested in this insert. However, the user should determine the optimal dilution factor.
- ✓ Spin down the SP conjugate vial before opening and using contents.
- ✓ The Stop Solution is an acidic solution.
- ✓ The kit should not be used beyond the expiration date.

## Assay Protocol

### Reagent Preparation

- ✓ Freshly dilute all reagents and bring all reagents to room temperature before use.
- ✓ MIX Diluent Concentrate (10x): Dilute the MIX Diluent Concentrate 10- fold with reagent grade water to produce a 1x solution. When diluting the concentrate, make sure to rinse the bottle thoroughly to extract any precipitates left in the bottle. Mix the 1x solution gently until the crystals have completely dissolved. Store for up to 30 days at 2-8°C.
- ✓ Human Kininogen (HMW) Standard: Reconstitute the Human Kininogen (HMW) Standard (1.8 µg) with 0.45 mL of MIX Diluent to generate a 4 µg/mL standard stock solution. Allow the vial to sit for 10 minutes with gentle agitation prior to making dilutions. Prepare duplicate or triplicate standard points by serially diluting from the standard stock solution (4 µg/mL) 2-fold with equal volume of MIX Diluent to produce 2, 1, 0.5, 0.25, 0.125 and 0.063 µg/mL solutions. MIX Diluent serves as the zero standard (0 µg/mL). Reconstitute a new vial for each assay.

Standard Point	Dilution	[Kininogen] (µg/mL)
P1	1 part Standard (4 µg/mL)	4.0
P2	1 part P1 + 1 part MIX Diluent	2.0
P3	1 part P2 + 1 part MIX Diluent	1.0
P4	1 part P3 + 1 part MIX Diluent	0.5
P5	1 part P4 + 1 part MIX Diluent	0.25
P6	1 part P5 + 1 part MIX Diluent	0.125
P7	1 part P6 + 1 part MIX Diluent	0.063
P8	MIX Diluent	0.000

- ✓ Biotinylated Human Kininogen (HMW) Protein (1x): Reconstitute the Biotinylated Human Kininogen (HMW) Protein with 3 mL of MIX Diluent to generate a stock solution. Allow the vial to sit for 10 minutes with gentle agitation prior to use. Reconstitute a new vial for each assay.
- ✓ Wash Buffer Concentrate (20x): Dilute the Wash Buffer Concentrate 20-fold with reagent grade water to produce a 1x solution. When diluting the concentrate, make sure to rinse the bottle thoroughly to extract any precipitates left in the bottle. Mix the 1x solution gently until the crystals have completely dissolved.
- ✓ SP Conjugate (100x): Spin down the SP Conjugate briefly and dilute the desired amount of the conjugate 100-fold with MIX Diluent to produce a 1x solution. The undiluted conjugate should be stored at -20°C.

### 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 200-fold sample dilution is suggested into MIX Diluent; 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 200-fold sample dilution is suggested into MIX Diluent; 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.

*Applicable samples may also include biofluids, cell culture, and tissue homogenates. If necessary, user should determine optimal dilution factor depending on application needs.*

- ✓ Refer to Dilution Guidelines for further instruction.

Guidelines for Dilutions of 100-fold or Greater (for reference only; please follow the insert for specific dilution suggested)	
100x	10000x
A. 4 µL sample: 396 µL buffer (100x) = 100-fold dilution  Assuming the needed volume is less than or equal to 400 µL.	A. 4 µL sample: 396 µL buffer (100x) B. 4 µL of A : 396 µL buffer (100x) = 10000-fold dilution  Assuming the needed volume is less than or equal to 400 µL.
1000x	100000x
A. 4 µL sample: 396 µL buffer (100x) B. 24 µL of A : 216 µL buffer (10x) = 1000-fold dilution  Assuming the needed volume is less than or equal to 240 µL.	A. 4 µL sample: 396 µL buffer (100x) B. 4 µL of A : 396 µL buffer (100x) C. 24 µL of B : 216 µL buffer (10x) = 100000-fold dilution  Assuming the needed volume is less than or equal to 240 µL.

### **Assay Procedure**

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 desiccant inside. Reseal the pouch securely to minimize exposure to water vapor and store in a vacuum desiccator.
3. Add 25 µL of Human Kininogen (HMW) Standard or sample to each well, and immediately add 25 µL of Biotinylated Human Kininogen (HMW) Protein to each well (on top of the standard or sample). Gently tap plate to ensure thorough mixing. 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 the microplate manually or automatically using a microplate washer. Invert the plate and decant the contents; hit 4-5 times on absorbent material to completely remove the liquid. If washing manually, wash

five times with 200  $\mu$ L of Wash Buffer per well. Invert the plate each time and decant the contents; hit 4-5 times on absorbent material to completely remove the liquid. If using a microplate washer, wash six times with 300  $\mu$ L of Wash Buffer per well; invert the plate and hit 4-5 times on absorbent material to completely remove the liquid.

5. Add 50  $\mu$ L 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.
6. Wash a microplate as described above.
7. Add 50  $\mu$ L of Chromogen Substrate to each well. Gently tap plate to thoroughly coat the wells. Break any bubbles that may have formed. Incubate in ambient light for 30 minutes or until the optimal blue color density develops.
8. Add 50  $\mu$ L of Stop Solution to each well. The color will change from blue to yellow. Gently tap plate to ensure thorough mixing. Break any bubbles that may have formed.
9. 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 low concentration points after stopping the reaction for about 10 minutes, which will reduce the readings.

✓ Assay Summary

1. Add 25  $\mu$ L of Standard or Sample and 25  $\mu$ L of Biotinylated Protein per well. Incubate 2 hours.
2. Wash, then add 50  $\mu$ L of SP Conjugate per well. Incubate 30 minutes.
3. Wash, then add 50  $\mu$ L of Chromogen Substrate per well. Incubate 30 minutes.
4. Add 50  $\mu$ L of Stop Solution per well. Read at 450 nm immediately.



## Data Analysis

### Calculation of Results

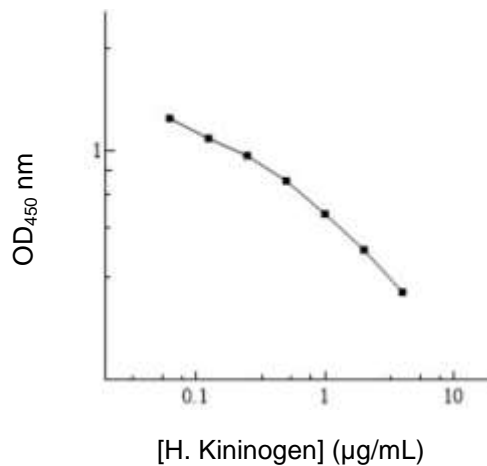
- ✓ Calculate the mean value of the duplicate or triplicate readings for each standard and sample.
- ✓ To generate a Standard Curve, plot the graph using the standard concentrations on the x-axis and the corresponding mean 450 nm absorbance (OD) on the y-axis. The best fit line can be determined by regression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

The typical data is provided for reference only. Individual laboratory means may vary from the values listed. Variations between laboratories may be caused by technique differences.

Standard Point	µg/mL	OD	Average OD
P1	4.0	0.242 0.241	0.242
P2	2.0	0.370 0.370	0.370
P3	1.0	0.536 0.526	0.531
P4	0.5	0.747 0.733	0.740
P5	0.25	0.961 0.943	0.952
P6	0.125	1.135 1.125	1.130
P7	0.063	1.388 1.379	1.384
P8	0.0	1.788 1.747	1.768
Sample: Pooled Normal Sodium Citrate Plasma (200x)		0.816 0.800	0.808

- ✓ Standard Curve

The curve is provided for illustration only. A standard curve should be generated each time the assay is performed.



✓ **Reference Value**

Plasma and serum samples from healthy adults were tested (n=40). On average, human kininogen (HMW) level was 102 µg/mL.

Sample	n	Average Value (µg/mL)
Pooled Normal Plasma	10	107.4
Normal Plasma	20	99.6
Pooled Normal Serum	10	97.7

**Performance Characteristics**

- ✓ The minimum detectable dose of human kininogen as calculated by 2SD from the mean of a zero standard was established to be 50 ng/mL.
- ✓ Intra-assay precision was determined by testing three plasma samples twenty times in one assay.
- ✓ Inter-assay precision was determined by testing three plasma samples in twenty assays.

	Intra-Assay Precision			Inter-Assay Precision		
Sample	1	2	3	1	2	3
n	20	20	20	20	20	20
CV (%)	4.7%	4.9%	5.3%	10.0%	9.8%	9.9%
Average CV (%)	5.0%			9.9%		

✓ **Recovery**

Standard Added Value	0.25 – 2 µg/mL
Recovery %	91-113%
Average Recovery %	97%

✓ Linearity

Plasma and serum samples were serially diluted to test for linearity.

Average Percentage of Expected Value (%)		
Sample Dilution	Plasma	Serum
100x	107%	106%
200x	101%	98%
400x	92%	96%

✓ Cross-Reactivity

Species	Cross Reactivity (%)
Canine	None
Bovine	None
Monkey	None
Mouse	None
Rat	None
Swine	None
Rabbit	None
Protein	Cross Reactivity (%)
Kininogen (LMW)	10%

## Resources

### Troubleshooting

Issue	Causes	Course of Action
Low Precision	Use of improper components	Check the expiration date listed before use. Do not interchange components from different lots.
	Improper wash step	Check that the correct wash buffer is being used. Check that all wells are empty after aspiration. Check that the microplate washer is dispensing properly. If washing by pipette, check for proper pipetting technique.
	Splashing of reagents while loading wells	Pipette properly in a controlled and careful manner.
	Inconsistent volumes loaded into wells	Pipette properly in a controlled and careful manner. Check pipette calibration. Check pipette for proper performance.
	Insufficient mixing of reagent dilutions	Thoroughly agitate the lyophilized components after reconstitution. Thoroughly mix dilutions.
	Improperly sealed microplate	Check the microplate pouch for proper sealing. Check that the microplate pouch has no punctures. Check that three desiccants are inside the microplate pouch prior to sealing.
Unexpectedly Low or High Signal Intensity	Microplate was left unattended between steps	Each step of the procedure should be performed uninterrupted.
	Omission of step	Consult the provided procedure for complete list of steps.
	Step performed in incorrect order	Consult the provided procedure for the correct order.
	Insufficient amount of reagents added to wells	Check pipette calibration. Check pipette for proper performance.
	Wash step was skipped	Consult the provided procedure for all wash steps.
	Improper wash buffer	Check that the correct wash buffer is being used.
	Improper reagent preparation	Consult reagent preparation section for the correct dilutions of all reagents.
	Insufficient or prolonged incubation periods	Consult the provided procedure for correct incubation time.

Deficient Standard Curve Fit	Non-optimal sample dilution	<p>Sandwich ELISA: If samples generate OD values higher than the highest standard point (P1), dilute samples further and repeat the assay.</p> <p>Competitive ELISA: If samples generate OD values lower than the highest standard point (P1), dilute samples further and repeat the assay.</p> <p>User should determine the optimal dilution factor for samples.</p>
	Contamination of reagents	A new tip must be used for each addition of different samples or reagents during the assay procedure.
	Contents of wells evaporate	Verify that the sealing film is firmly in place before placing the assay in the incubator or at room temperature.
	Improper pipetting	<p>Pipette properly in a controlled and careful manner.</p> <p>Check pipette calibration.</p> <p>Check pipette for proper performance.</p>
	Insufficient mixing of reagent dilutions	<p>Thoroughly agitate the lyophilized components after reconstitution.</p> <p>Thoroughly mix dilutions.</p>

## **References**

1. Reddigari SR et al. (1993) Blood 81:1306-1311
2. Maier M et al. (1983) Proc. Natl. Acad. Sci. USA 80:3928-3932
3. Kellermann J et al. (1986) Eur. J. Biochem. 154:471-478
4. Kitamura N et al. (1985) J. Biol. Chem. 260:8610-8617
5. Kerbiriou DM and Griffin JH (1979) J. Biol. Chem. 254:12020-12027
6. Coffman LG et al. (2009) Proc. Natl. Acad. Sci. USA 106:570-575
7. Hassan S et al. (2007) Am. J. Physiol Heart Circ Physiol. 292:H2959-H2965

**Plate Layout**

12								
11								
10								
9								
8								
7								
6								
5								
4								
3								
2								
1								
	A	B	C	D	E	F	G	H