

ab326079 – Human alpha 1 Acid Glycoprotein SimpleStep ELISA® Kit, Chemiluminescent

For the quantitative measurement of alpha 1 Acid Glycoprotein in human serum, plasma (heparin), plasma (EDTA), plasma (citrate), cell culture supernatant, saliva, urine, and milk.
For research use only - not intended for diagnostic use.
Patent pending.

For overview, typical data and additional information please visit: www.abcam.com/ab326079

Storage and Stability: Store kit at 2-8°C immediately upon receipt. Refer to list of materials supplied for storage conditions of individual components. Observe the storage conditions for individual prepared components in the Standard Preparation and Reagent preparation sections.

Limitations: All data, except typical standard curve and sensitivity were collected using the colorimetric version of this kit (ab243675).

Materials Supplied

Item	Quantity 1 x 96 tests	Storage Condition
Human alpha 1 Acid Glycoprotein Capture Antibody 10X	600 µL	+4°C
Human alpha 1 Acid Glycoprotein Detector Antibody 10X	600 µL	+4°C
Human alpha 1 Acid Glycoprotein Lyophilized Recombinant Protein	2 Vials	+4°C
Antibody Diluent 5BI	6 mL	+4°C
Sample Diluent NS	2 x 50 mL	+4°C
Wash Buffer PT 10X	20 mL	+4°C
ChemiHRP Reagent A	3 mL	+4°C
ChemiHRP Reagent B	3 mL	+4°C
SimpleStep Pre-Coated Black 96-Well Microplate	96 Wells	+4°C
Plate Seal	1	+4°C

Materials Required, Not Supplied

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

Luminometer with the following settings: 0.5-1 second/well read time; summation mode (all wavelengths).

Deionized water.

Multi- and single-channel pipettes.

Tubes for standard dilution.

Orbital microplate shaker for all incubation steps: capable of 750 rpm shaking speed.

Optional: Phenylmethylsulfonyl Fluoride (PMSF) (or other protease inhibitors).

Reagent Preparation

Equilibrate all reagents to room temperature (18-25°C) prior to use. The kit contains enough reagents for 96 wells. The sample volumes below are sufficient for 48 wells (6 x 8-well strips); adjust volumes as needed for the number of strips in your experiment.

Prepare only as much reagent as is needed on the day of the experiment. Capture and Detector Antibodies have only been tested for stability in the provided 10X formulations.

1X Wash Buffer PT: Prepare 1X Wash Buffer PT by diluting Wash Buffer PT 10X with deionized water. To make 50 mL 1X Wash Buffer PT combine 5 mL Wash Buffer PT 10X with 45 mL deionized water. Mix thoroughly and gently.

Antibody Cocktail: Prepare Antibody Cocktail by diluting the capture and detector antibodies in Antibody Diluent 5BI. To make 3 mL of the Antibody Cocktail combine 300 µL 10X Capture Antibody and 300 µL 10X Detector Antibody with 2.4 mL Antibody Diluent 5BI. Mix thoroughly and gently.

Lumi HRP Development Solution: Just prior to use, prepare Lumi HRP Development Solution by mixing equal volume of the ChemiHRP Reagent A and the ChemiHRP Reagent B. To make 3 mL of the Lumi HRP Development Solution combine 1.5 mL of ChemiHRP Reagent A and 1.5 mL of ChemiHRP Reagent B. Mix thoroughly and gently by inversion or slow pipetting (Avoid shaking or vortexing). Protect the prepared solution from light until use.

Standard Preparation

Always prepare a fresh set of standards for every use. Discard working standard dilutions after use as they do not store well. The following section describes the preparation of a standard curve for duplicate measurements (recommended).

1. Reconstitute the alpha 1 Acid Glycoprotein standard sample by adding the volume of Sample Diluent NS indicated on the protein vial label. Hold at room temperature for 10 minutes. Mix thoroughly and gently. This is the 192 ng/mL **Stock Standard #1** Solution.
2. Label nine tubes, Standards 1–9.
3. Add 150 µL of Sample Diluent NS into numbers 2-9.
4. Use the **Stock Standard #1** to prepare the following dilution series. Standard #9 contains no protein and is the Blank control:

Standard #	Dilution Sample	Volume to Dilute (µL)	Volume of Diluent (µL)	Starting Conc. (ng/mL)	Final Conc. (ng/mL)
1	Stock Standard #1	150	0	192	192
2	Stock Standard #1	75	150	192	64
3	Standard#2	75	150	64	21.33
4	Standard#3	75	150	21.33	7.11
5	Standard#4	75	150	7.11	2.37
6	Standard#5	75	150	2.37	0.79
7	Standard#6	75	150	0.79	0.26
8	Standard#7	75	150	0.26	0.09
9	Blank Control	0	150	0	0

Sample Preparation

Typical Sample Dynamic Range	
Sample Type	Range
Serum	1:80,000 - 1:1,280,000
Plasma – Heparin	1:80,000 - 1:1,280,000
Plasma – EDTA	1:80,000 - 1:1,280,000
Plasma – Citrate	1:40,000 - 1:640,000
HepG2 Cell Culture Supernatant	1:28 – 1:200
Urine	1:800 – 1:12,800
Saliva	1:800 – 1:12,800
Milk	1:500 – 1:8,000

Note: Due to the high dilutions required for biological samples, we recommend to pre-dilute the samples sequentially. As an example, the tables below show the steps suggested to pre-dilute the samples. Final dilutions should be carried in Sample Diluent NS.

Serum and plasma (citrate, EDTA, and heparin) samples:

Tube #	Sample to Dilute	Volume to Dilute (µL)	Volume of Diluent (µL)	Diluent	Starting Dilution	Ending Dilution
1	Neat	5	245	Wash Buffer PT	Neat	1:50
2	Tube #1	5	195	Wash Buffer PT	1:50	1:2,000

Urine, saliva, and milk samples:

Tube #	Sample to Dilute	Volume to Dilute (µL)	Volume of Diluent (µL)	Diluent	Starting Dilution	Ending Dilution
1	Neat	5	95	Wash Buffer PT	Neat	1:20

Serum Samples should be collected into a serum separator tube. After clot formation, centrifuge samples at 2,000 x g for 10 minutes and collect serum. Pre-dilute serum samples 1: 2,000 into 1X Wash Buffet PT. Further dilute samples at least to 1:80,000 into Sample Diluent NS and assay. Store un-diluted serum at -20°C or below. Avoid repeated freeze-thaw cycles.

Plasma Collect plasma using citrate, EDTA or heparin. Centrifuge samples at 2,000 x g for 10 minutes. Pre-dilute plasma samples 1: 2,000 into 1X Wash Buffet PT. Further dilute plasma (citrate) samples at least to 1: 40,000 into Sample Diluent NS and assay.

Further dilute plasma (EDTA and heparin) samples at least to 1: 80,000 into Sample Diluent NS and assay. Store un-diluted plasma samples at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.

Cell Culture Supernatants Centrifuge cell culture media at 2,000 x g for 10 minutes to remove debris. Collect supernatants and assay. Or dilute samples into Sample Diluent NS and assay. Store un-diluted samples at -20°C or below. Avoid repeated freeze-thaw cycles.

Urine Centrifuge urine at 2,000 x g for 10 minutes to remove debris. Pre-dilute urine samples 1: 20 into 1X Wash Buffet PT. Further dilute samples at least to 1:800 into Sample Diluent NS and assay. Store un-diluted urine samples at -20°C or below. Avoid repeated freeze-thaw cycles.

Saliva Centrifuge saliva at 800 x g for 10 minutes to remove debris. Collect supernatants. Pre-dilute urine samples 1: 20 into 1X Wash Buffet PT. Further dilute samples at least to 1:800 into Sample Diluent NS and assay. Store un-diluted samples at -20°C or below. Avoid repeated freeze-thaw cycles.

Milk De-fat milk samples as follows. Centrifuge milk samples at 500 x g for 15 minutes at 4°C and collect the aqueous fraction using syringe attached to needle. Centrifuge the aqueous fraction at 3,000 x g for 15 minutes at 4°C and collect the final aqueous fraction (de-fatted milk) using syringe attached to needle. Pre-dilute milk samples 1: 20 into 1X Wash Buffet PT. Further dilute samples at least to 1:500 into Sample Diluent NS and assay. Store un-diluted de-fatted milk at -20°C or below. 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).

Differences in well absorbance or "edge effects" have not been observed with this assay.

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, working standards, and samples as directed in the previous sections.
2. Remove excess microplate strips from the plate frame, return them to the foil pouch containing the desiccant pack, reseal and return to 4°C storage.
3. Add 50 µL of all sample or standard to appropriate wells.
4. Add 50 µL of the Antibody Cocktail to each well.
5. Seal the plate and incubate for 30 minutes at room temperature on a plate shaker set to 750 rpm.
6. Wash each well with 3 x 350 µL 1X Wash Buffer PT. Wash by aspirating or decanting from wells then dispensing 350 µL 1X Wash Buffer PT into each well. Wash Buffer PT should remain in wells for at least 30 seconds. Complete removal of liquid at each step is essential for good performance. After the last wash invert the plate and tap gently against clean paper towels to remove excess liquid.
7. Add 50 µL of prepared Lumi HRP Development Solution to each well and incubate for 1 minute in the dark on a plate shaker set to 750 rpm. Further optimization of incubation time vs signal strength can be performed if needed. Avoid introducing bubbles into the wells.
8. Measure the produced light of each well using a microplate luminometer with the following settings: 1 second/well read time in summation mode (all wavelengths). Relative light unit (RLU) readings may vary between luminometer models. It is recommended to configure instrument settings according to the manufacturer's specifications. Note: Relative light unit (RLU) values may change over the course of the 15-minute reading window.
9. Analyze the data as described below.

Mode:	Luminescence
Instrument settings:	Endpoint
Detection Mode:	All wavelengths
Read Time:	0.5-1 sec
Read:	Top

Note For microplate readers with Pre-Read Optimization option, the Read Height as well as Microplate Optimization is recommended before the first read.

Download our ELISA guide for technical hints, results, calculation, and troubleshooting tips:

<https://www.abcam.com/en-us/technical-resources/guides/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.

Version 1a | 2026-03-11

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)

ab326079 – Human alpha 1 Acid Glycoprotein SimpleStep ELISA® Kit, Chemiluminescent

Additional information

ASSAY SPECIFICITY

This kit is designed for the quantification of human alpha 1 Acid Glycoprotein.

Native signal was detected in serum, plasma (citrate), plasma (EDTA), plasma (heparin), cell culture supernatant, urine, saliva, and milk sample types.

CSF, cell extract, and tissue extract samples have not been tested with this kit.

CROSS REACTIVITY

2 ng/mL of recombinant human alpha 1 Acid Glycoprotein 2 was tested for cross reactivity. No cross reactivity was observed.

INTERFERENCE

50 ng/mL of recombinant human alpha 1 Acid Glycoprotein 2 and 2 ng/mL of recombinant human alpha 1 Acid Glycoprotein 2 were tested for interference. No interference was observed.

SPECIES REACTIVITY

Other species reactivity was determined by measuring 1:160,000 serum samples of various species, interpolating the protein concentrations from the human standard curve, and expressing the interpolated concentrations as a percentage of the protein concentration in human serum assayed at the same dilution.

Interpolated values were below the detectable range of the assay for the following species: Monkey (Rhesus macaque), Mouse, Rat, Cow.

Other species reactivity not determined.

CALIBRATION

This immunoassay is calibrated against a highly purified human alpha 1 Acid Glycoprotein. European Reference Material ERM®-DA470k/IFCC was evaluated in this kit. The dose response curve of the standard alpha 1 Acid Glycoprotein parallels the SimpleStep standard curve.

To convert sample values obtained with the SimpleStep Human alpha 1 Acid Glycoprotein kit to approximate ERM®-DA470k/IFCC units, use the equation below.

$ERM®-DA470k/IFCC \text{ approximate value (g/L)} = 1.14 \times \text{SimpleStep Human alpha 1 Acid Glycoprotein value (g/L)}$.

CALCULATION

- Preconfigured protocols are available when using SoftMax Pro software from Molecular Devices.
- Calculate the average chemiluminescence value for the blank control (zero) standards. Subtract the average blank control standard chemiluminescence value from all other chemiluminescence values.

- Create a standard curve by plotting the average blank control subtracted chemiluminescence value for each standard concentration (y-axis) against the target protein concentration (x-axis) of the standard. Use graphing software to draw the best smooth curve through these points to construct the standard curve.
- Note: Most chemiluminescence reader software or graphing software will plot these values and fit a curve to the data. A four-parameter curve fit (4PL) is often the best choice; however, other algorithms (e.g. linear, semi-log, log/log, 4-parameter logistic) can also be tested to determine if it provides a better curve fit to the standard values.
- Determine the concentration of the target protein in the sample by interpolating the blank control subtracted chemiluminescence values against the standard curve. Multiply the resulting value by the appropriate sample dilution factor, if used, to obtain the concentration of target protein in the sample.
- Samples generating chemiluminescence values greater than that of the highest standard should be further diluted and reanalyzed. Similarly, samples which measure at chemiluminescence values less than that of the lowest standard should be retested in a less dilute form.

TYPICAL DATA

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

Standard Curve Measurements			
Concentration (ng/mL)	RLU		Mean RLU
	1	2	
0	2,276	2,276	2,276
0.09	3,438	5,379	4,409
0.26	19,729	8,829	14,279
0.79	29,090	18,184	23,637
2.37	74,143	79,792	76,968
7.11	249,797	221,663	235,730
21.33	791,175	760,107	775,641
64	2,232,691	2,232,500	2,232,596
192	5,324,559	5,398,165	5,361,362

Table 1. Example of human alpha 1 Acid Glycoprotein standard curve in Sample Diluent NS. The alpha 1 Acid Glycoprotein standard curve was prepared as described in the Standard Preparation section. The table shows raw data values.

TYPICAL SAMPLE VALUES

Sensitivity:

The calculated minimal detectable dose (MDD) is 52.54 pg/mL. The MDD was determined by calculating the mean of zero standard replicates (n=16) and adding 2 standard deviations then extrapolating the corresponding concentration.

Recovery

Three concentrations of alpha 1 Acid Glycoprotein recombinant protein were spiked in duplicate to the indicated biological matrix to evaluate signal recovery in the working range of the assay.

Sample Type	Average % Recovery	Range (%)
1:333,333 Serum	90	85 - 97
1:333,333 Plasma – Heparin	106	100 - 112
1:333,333 Plasma – EDTA	107	105 - 108
1:333,333 Plasma – Citrate	95	88 - 105
1% HepG2 Cell Culture Supernatant	99	93 - 106
1:1,000 Urine	103	93 - 112
1:2,000 Saliva	112	110 - 117
1:2,000 Milk	94	79 - 104
100% Cell Culture Media*	84	78 - 89

*Media is RPMI 1640 containing 10% fetal bovine serum.

Linearity of Dilution

Linearity of dilution is determined based on interpolated values from the standard curve. Linearity of dilution defines a sample concentration interval in which interpolated target concentrations are directly proportional to sample dilution.

Native alpha 1 Acid Glycoprotein was measured in the following biological samples in a 2-fold dilution series. Sample dilutions are made in Sample Diluent NS.

Dilution Factor	Interpolated value	1:80,000 Human Serum	1:40,000 Human Plasma (Citrate)	1:80,000 Human Plasma (EDTA)	1:80,000 Human Plasma (Heparin)
Undiluted	ng/mL	10.476	11.628	11.425	10.004
	% Expected	100	100	100	100
2	ng/mL	4.905	5.984	5.412	4.693
	% Expected	96	103	95	94
4	ng/mL	2.559	2.879	2.721	2.404
	% Expected	98	99	95	96
8	ng/mL	1.321	1.501	1.412	1.191
	% Expected	101	103	99	95
16	ng/mL	0.625	0.801	0.667	0.591
	% Expected	95	110	93	94

Native alpha 1 Acid Glycoprotein was measured in the following biological samples in a 2-fold dilution series. Sample dilutions are made in Sample Diluent NS.

Dilution Factor	Interpolated value	1:800 Human Urine	1:500 Human Milk	1:800 Human Saliva	3.5% Human HepG2 Cell Culture Supernatant
Undiluted	ng/mL	9.099	11.052	11.614	8.666
	% Expected	100	100	100	100
2	ng/mL	4.268	5.621	5.329	4.075
	% Expected	94	102	92	94
4	ng/mL	2.105	2.616	2.549	1.972
	% Expected	93	95	88	91
8	ng/mL	1.047	1.307	1.253	0.949
	% Expected	92	95	86	88
16	ng/mL	0.526	0.636	0.632	0.491
	% Expected	92	92	87	91

Precision

Mean coefficient of variations of interpolated values of alpha 1 Acid Glycoprotein from three concentrations of human serum within the working range of the assay.

	Intra-assay	Inter-assay
N=	8	3
CV (%)	5.9	5.4

Download our ELISA guide for technical hints, results, calculation, and troubleshooting tips:

<https://www.abcam.com/en-us/technical-resources/guides/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.

Version 1a | 2026-03-11

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)