

ab326752 – Human AMH SimpleStep ELISA® Kit, Chemiluminescent

For the quantitative measurement of AMH in human serum, plasma (heparin), plasma (EDTA), plasma (citrate), and cell culture supernatant.

For research use only - not intended for diagnostic use.

Patent pending.

For overview, typical data and additional information please visit: www.abcam.com/ab326752

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 (ab318930).

Materials Supplied

Item	Quantity 1 x 96 tests	Storage Condition
Human AMH Capture Antibody 10X	600 µL	+4°C
Human AMH Detector Antibody 10X	600 µL	+4°C
Human AMH Lyophilized Recombinant Protein	2 Vials	+4°C
Antibody Diluent 4BI	6 mL	+4°C
Cell Extraction Enhancer Solution 50X	1 mL	+4°C
Sample Diluent NS	12 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 capable of reading strip plates 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.

The provided Cell Extraction Enhancer Solution 50X may precipitate when stored at + 4°C. To dissolve, warm briefly at + 37°C and mix gently. The Cell Extraction Enhancer Solution 50X can be stored at room temperature to avoid precipitation.

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.

Sample Diluent NS + 1X Enhancer: Prepare **Sample Diluent NS + 1X Enhancer** by combining Sample Diluent NS and 50X Cell Extraction Enhancer Solution. To make 5 mL **Sample Diluent NS + 1X Enhancer**, combine 4.9 mL Sample Diluent NS and 100 µL Cell Extraction Enhancer Solution 50X. Mix thoroughly and gently.

Antibody Cocktail: Prepare Antibody Cocktail by diluting the capture and detector antibodies in Antibody Diluent 4BI. 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 4BI. 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 AMH standard sample by adding the volume of **Sample Diluent NS + 1X Enhancer** indicated on the protein vial label. Hold at room temperature for 10 minutes. Mix thoroughly and gently. This is the 28.8 ng/mL **Stock Standard** Solution.
- 2. Label nine tubes, Standards 1– 9.
- 3. Add 203 µL of **Sample Diluent NS + 1X Enhancer** into tube number 1 and 150 µL of **Sample Diluent NS + 1X Enhancer** into numbers 2-9.
- 4. Use the **Stock Standard** 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. (pg/mL)	Final Conc. (pg/mL)
1	Stock Standard	85	203	28,800	8,500
2	Standard#1	75	150	8,500	2,833
3	Standard#2	75	150	2,833	944.4
4	Standard#3	75	150	944.4	314.8
5	Standard#4	75	150	314.8	104.9
6	Standard#5	75	150	104.9	34.98
7	Standard#6	75	150	34.98	11.66
8	Standard#7	75	150	11.66	3.89
9	Blank Control	0	150	0	0

Sample Preparation

Typical Sample Dynamic Range	
Sample Type	Range
Serum	2.8 – 45%
Plasma – Citrate	5.6 – 45%
Plasma – EDTA	2.8 – 45%
Plasma – Heparin	2.8 – 45%
Cell Culture Supernatant	≤ 50%

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. 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. 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. Store un-diluted samples at -20°C or below. Avoid repeated freeze-thaw cycles.

Biological Sample Treatment Protocol To activate latent AMH to the immunoreactive form, biologicals must be treated with 50X Enhancer prior to sample dilution. Do not treat the kit standard as it already contains the immunoreactive form of human AMH.

- 1. Add 90 µL of sample to a tube.
- 2. Add 10µL of 50X Enhancer. Mix well.
- 3. Incubate for 10 minutes at room temperature.
- 4. Biological sample concentration after treatment is 90%. Dilute biological to desired concentration in **Sample Diluent NS + 1X Enhancer** (see typical sample dynamic range).

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: 0.5-1 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 | 2026-05-08

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)

ab326752 – Human AMH SimpleStep ELISA® Kit, Chemiluminescent

Additional information

ASSAY SPECIFICITY

This kit is designed for the quantification of human AMH.

The standard protein in this kit is an N-terminal fragment of the human AMH propeptide.

Native signal was detected in serum, plasma (citrate), plasma (EDTA), and plasma (heparin) sample types.

Spiked protein experiments were used to validate cell culture supernatant sample types.

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

CROSS REACTIVITY

50 ng/mL of recombinant human AMHR2 (MIS RII) was tested for cross reactivity. No cross reactivity was observed.

INTERFERENCE

50 ng/mL of recombinant human AMHR2 (MISR II) was tested for interference with 700 pg/mL of recombinant Human AMH. No interference was observed.

SPECIES REACTIVITY

Other species reactivity was determined by measuring serum samples of various species. Sample dilutions were 45% cow serum, 22.5% rat serum, and 2.25% monkey serum.

No signal was observed for the following species: Cow

Reactivity was determined for the following species: Monkey, Rat

This kit is not recommended for the quantification of AMH in mouse samples.

Other species reactivity not determined.

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 (pg/mL)	RLU		Mean RLU
	1	2	
0	828	828	828
3.89	2,596	2,590	2,593
11.66	7,602	7,702	7,652
34.98	18,089	18,994	18,542
104.9	52,107	55,179	53,643
314.8	206,660	223,410	215,035
944.4	457,570	558,080	507,825
2,833	1,169,300	1,386,300	1,277,800
8,500	4,545,400	3,759,700	4,152,550

Table 1. Example of human AMH standard curve in **Sample Diluent NS + 1X Enhancer**. The AMH 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 1.12 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 AMH 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 (%)
11% Serum	100	87 - 108
22.5% Plasma (citrate)	107	100 – 117
11% Plasma (EDTA)	99	94 – 105
22.5% Plasma – Heparin	94	86 – 105
50% Cell Culture Media	98	82 - 106

*Media is DMEM 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 AMH was measured in the following biological samples in a 2-fold dilution series. Sample dilutions are made in **Sample Diluent NS + 1X Enhancer**.

Dilution Factor	Interpolated value	45% Human Serum	45% Human Plasma (Citrate)	45% Human Plasma (EDTA)	45% Human Plasma (Heparin)
Undiluted	pg/mL	935.7	460.7	980.3	618.2
	% Expected value	100	100	100	100
2	pg/mL	514.1	253.9	521.4	329.2
	% Expected value	110	110	106	107
4	pg/mL	263.5	142.7	292.2	177.0
	% Expected value	113	124	119	115
8	pg/mL	112.9	58.4	134.1	92.1
	% Expected value	97	101	109	119
16	pg/mL	52.4	ND	64.5	41.7
	% Expected value	90	ND	105	108

ND – Not Detected – below product dynamic range

Recombinant AMH was spiked into the following biological samples and then diluted in a 2-fold dilution series. Sample dilutions are made in **Sample Diluent NS + 1X Enhancer**.

Dilution Factor	Interpolated value	50% Cell Culture Media
Undiluted	pg/mL	1,306.8
	% Expected value	100
2	pg/mL	668.7
	% Expected value	102
4	pg/mL	315.7
	% Expected value	97
8	pg/mL	155.3
	% Expected value	95
16	pg/mL	77.4
	% Expected value	95

Precision

Mean coefficient of variations of interpolated values of AMH from a single concentration of human serum within the working range of the assay.

	Intra-assay	Inter-assay
N=	8	3
CV (%)	6.1	10.6

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 | 2026-05-08

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)