

ab326005 – Human IL-11 SimpleStep ELISA® Kit, Chemiluminescent

For the quantitative measurement of IL-11 in human serum, 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/ab326005

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 (ab324512).

Materials Supplied

Item	Quantity 1 x 96 tests	Storage Condition
Human IL-11 Capture Antibody 10X	600 µL	+4°C
Human IL-11 Detector Antibody 10X	600 µL	+4°C
Human IL-11 Lyophilized Recombinant Protein	2 Vials	+4°C
Antibody Diluent CPI2	6 mL	+4°C
Sample Diluent NS	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 CPI2. 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 CPI2. 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 IL-11 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 40,000 pg/mL **Stock Standard** Solution.
2. Label nine, Standards 1– 9.
3. Add 150 µL of Sample Diluent NS into tube number 1 and 150 µL of Sample Diluent NS 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	90	150	40,000	15,000
2	Standard#1	75	150	15,000	5,000
3	Standard#2	75	150	5,000	1,666.7
4	Standard#3	75	150	1,666.7	555.6
5	Standard#4	75	150	555.6	185.2
6	Standard#5	75	150	185.2	61.7
7	Standard#6	75	150	61.7	20.6
8	Standard#7	75	150	20.6	6.9
9	Blank Control	0	150	0	0

Sample Preparation

Typical Sample Dynamic Range	
Sample Type	Range
Serum	≤ 25%
Plasma – Citrate	≤ 25%
Plasma – EDTA	≤ 25%
A549 Cell Culture Supernatant*	12.5 – 100%
Untreated MRC-5 Cell Culture Supernatant**	3.75 – 60%
PMA + IL-1 alpha Treated MRC-5 Cell Culture Supernatant**	0.94 – 15%

*Media is FK12 containing 10% fetal bovine serum

**Media is DMEM containing 10% fetal bovine serum

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. Dilute samples at least 1:4 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 or EDTA. Centrifuge samples at 2,000 x g for 10 minutes. Dilute samples at least 1:4 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. **Note:** This kit is incompatible with plasma (heparin) samples.

Cell Culture Supernatants Centrifuge cell culture media at 2,000 x g for 10 minutes to remove debris. Collect supernatants. Assay, or dilute samples into Sample Diluent NS and assay. Store undiluted samples 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, resealed 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 | 2026-03-13

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)

ab326005 – Human IL-11 SimpleStep ELISA® Kit, Chemiluminescent

Additional information

ASSAY SPECIFICITY

This kit is designed for the quantification of human IL-11.

The standard protein in this kit is mature full-length human IL-11.

Native signal was detected in cell culture supernatant sample types.

Spiked protein experiments were used to validate serum, plasma (citrate), and plasma (EDTA) sample types.

25% pooled serum and plasma (citrate and EDTA) samples from healthy donors was measured in duplicate. All values were below the detectable range of the assay.

25% serum from ten individual healthy human female donors was measured in duplicate. All values were below the detectable range of the assay.

25% pooled rheumatoid arthritis serum was measured in duplicate. Values were below the detectable range of the assay.

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

This kit is incompatible with plasma (heparin) samples.

CROSS REACTIVITY

44,800 pg/mL of recombinant human IL-11 RA was tested for cross reactivity. No cross reactivity was observed.

INTERFERENCE

44,800 pg/mL of recombinant human IL-11 RA was tested for interference with 2,800 pg/mL of recombinant human IL-11. No interference was observed.

SPECIES REACTIVITY

5,600 pg/mL of recombinant mouse IL-11 and 5,600 pg/mL of recombinant monkey IL-11 were tested for reactivity. No reactivity was observed with recombinant mouse IL-11. 14% reactivity was observed with recombinant monkey IL-11.

Other species reactivity not determined.

CALIBRATION

This immunoassay is calibrated against a highly purified human IL-11. The NIBSC/WHO unclassified purified human IL-11 preparation 92/788 was evaluated in this kit.

The dose response curve of the unclassified standard IL-11 parallels the SimpleStep standard curve. To convert sample values obtained with the SimpleStep Human IL-11 kit to approximate NIBSC 92/788 units, use the equation below.

NIBSC (92/788) approximate value (units/mL) = 0.02029 x SimpleStep Human IL-11 value (pg/mL).

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	1,218	1,218	1,218
6.9	2,240	2,101	2,171
20.6	3,811	3,789	3,800
61.7	12,441	12,832	12,637
185.2	31,274	30,531	30,903
555.6	97,924	99,026	98,475
1,666.7	377,340	372,760	375,050
5,000	931,330	881,170	906,250
15,000	3,424,200	3,237,200	3,330,700

Table 1. Example of human IL-11 standard curve in Sample Diluent NS. The IL-11 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 4.5 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 IL-11 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 (%)
25% Serum	96	93 - 100
25% Plasma – Citrate	98	88 - 110
25% Plasma – EDTA	84	83 - 85
50% A549 Cell Culture Supernatant*	96	93 - 99
30% Untreated MRC-5 Cell Culture Supernatant**	93	86 - 95
7.5% PMA + IL-1 alpha Treated MRC-5 Cell Culture Supernatant**	100	98 - 102

*Media is FK12 containing 10% fetal bovine serum

**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 IL-11 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	100% A549 Cell Supernatant	60% Untreated MRC-5 Cell Supernatant	15% Treated MRC-5 Cell Supernatant
Undiluted	pg/mL	434.4	906.99	1,128.8
	% Expected value	100%	100%	100%
2	pg/mL	238.1	470.15	607.2
	% Expected value	110%	104%	108%
4	pg/mL	114.9	268.30	303.6
	% Expected value	106%	118%	108%
8	pg/mL	60.3	130.95	155.9
	% Expected value	111%	116%	111%
16	pg/mL	ND	68.91	74.2
	% Expected value	ND	122%	105%

ND – Not Detected – below product dynamic range

Recombinant IL-11 was spiked into the following biological samples and then diluted in a 2-fold dilution series. Sample dilutions are made in Sample Diluent NS.

Dilution Factor	Interpolated value	25% Serum	25% Plasma (Citrate)	25% Plasma (EDTA)
Undiluted	pg/mL	609.0	763.0	688.1
	% Expected value	100%	100%	100%
2	pg/mL	341.2	312.2	309.3
	% Expected value	112%	82%	90%
4	pg/mL	177.7	158.5	156.0
	% Expected value	117%	83%	91%
8	pg/mL	75.8	70.5	67.1
	% Expected value	100%	74%	78%
16	pg/mL	45.3	42.0	50.0
	% Expected value	119%	88%	116%

Precision

Mean coefficient of variations of interpolated values of IL-11 from two concentrations of PMA + IL-1 alpha treated MRC-5 cell culture supernatant within the working range of the assay.

	Intra-assay	Inter-assay
N=	8	3
CV (%)	6.1	1.0

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 | 2026-03-13

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)