

ab327316 – Monkey CRP SimpleStep ELISA® Kit (C-Reactive Protein), Chemiluminescent

For the quantitative measurement of CRP in monkey serum, plasma (heparin), plasma (EDTA), plasma (citrate), cell culture supernatant, and CSF.
For research use only - not intended for diagnostic use.
Patent pending.

For overview, typical data and additional information please visit: www.abcam.com/ab327316

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 (ab260062).

Materials Supplied

Item	Quantity 1 x 96 tests	Quantity 10 x 96 tests	Storage Condition
Monkey CRP Capture Antibody 10X	600 µL	10 x 600 µL	+4°C
Monkey CRP Detector Antibody 10X	600 µL	10 x 600 µL	+4°C
Monkey CRP Lyophilized Recombinant Protein	2 Vials	10 x 2 Vials	+4°C
Antibody Diluent 4BI	6 mL	10 x 6 mL	+4°C
Sample Diluent NS	50 mL	2 x 250 mL	+4°C
Wash Buffer PT 10X	20 mL	200 mL	+4°C
ChemiHRP Reagent A	3 mL	50 mL	+4°C
ChemiHRP Reagent B	3 mL	50 mL	+4°C
SimpleStep Pre-Coated Black 96-Well Microplate	96 Wells	10 x 96 Wells	+4°C
Plate Seal	1	10	+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.
Hydrochloric acid (HCl)
Sodium hydroxide (NaOH)
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 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).

- Reconstitute the CRP standard sample by adding the volume of NS indicated on the protein vial label. Hold at room temperature for 10 minutes. Mix thoroughly and gently. This is the 800 ng/mL **Stock Standard** Solution.
Acid Treatment Protocol: For dissociation of CRP in standard.
 - Dilute CRP Stock Standard to 40 ng/mL: Add 10 µL of CRP Stock Standard to 190 µL of Sample Diluent NS.
 - Add 100 µL of 0.25N HCl to the diluted CRP. Incubate for 15 minutes at room temperature while rotating.
 - Add 100 µL of 0.25N NaOH to acid treated sample. Incubate for 3 minutes at room temperature while rotating.
 - CRP **Treated Stock Standard #1** after treatment is 20,000 pg/mL. Dilute CRP **Treated Stock Standard #1** in Sample Diluent NS to prepare dilution series.
- Label nine tubes, Standards 1–9.
- Add 150 µL of NS into numbers 2-9.
- Use the **Treated 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. (pg/mL)	Final Conc. (pg/mL)
------------	-----------------	-----------------------	------------------------	------------------------	---------------------

1	Treated Stock Standard #1	150	0	20,000	20,000
2	Treated Stock Standard #1	75	150	20,000	6,666.6
3	Standard#2	75	150	6,666.6	2,222.2
4	Standard#3	75	150	2,222.2	740.74
5	Standard#4	75	150	740.74	246.91
6	Standard#5	75	150	246.91	82.30
7	Standard#6	75	150	82.30	27.43
8	Standard#7	75	150	27.43	9.14
9	Blank Control	0	150	0	0

Sample Preparation

Typical Sample Dynamic Range	
Sample Type	Range
Serum	1:8,000 - 1:500
Plasma – Citrate	1:8,000 - 1:500
Plasma – EDTA	1:16,000 - 1:2,000
Plasma – Heparin	1:8,000 - 1:500
Cell Culture Media*	≤ 25%
Cerebrospinal fluid	0.39 – 3.13%

*Based on spiked sample

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. Perform acid treatment protocol before assaying.

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. Perform acid treatment protocol before assaying.

Acid Treatment Protocol: For dissociation of CRP in serum/plasma samples.

- 1) Dilute serum/plasma to 10%: Add 10 µL of sample to 90 µL of Sample Diluent NS.
- 2) Add 50 µL of 0.25N HCl to 10% serum/plasma. Incubate for 15 minutes at room temperature while rotating.
- 3) Add 50 µL of 0.25N NaOH to acid treated sample. Incubate for 3 minutes at room temperature while rotating.
- 4) Serum/plasma concentration after treatment is 5%, dilute treated sample in Sample Diluent NS to in assay concentration (see typical sample dynamic range).

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.

Acid Treatment Protocol: For dissociation of CRP in cell culture supernatant.

- 1) Add 50 µL of 0.25N HCl to 100 µL neat cell culture supernatant. Incubate for 15 minutes at room temperature while rotating.
- 2) Add 50 µL of 0.25N NaOH to acid treated sample. Incubate for 3 minutes at room temperature while rotating.
- 3) Cell culture supernatant concentration after treatment is 50%, dilute treated sample in Sample Diluent NS to in assay concentration (see typical sample dynamic range).

Cerebrospinal Fluid (CSF) Treat cerebrospinal fluid according to the below protocol and assay. Store un-diluted samples at -20°C or below. Avoid repeated freeze-thaw cycles.

Acid Treatment Protocol: For dissociation of CRP in cerebrospinal fluid samples.

- 1) Add 50 µL of 0.25N HCl to 100 µL neat cerebrospinal fluid samples. Incubate for 15 minutes at room temperature while rotating.
- 2) Add 50 µL of 0.25N NaOH to acid treated sample. Incubate for 3 minutes at room temperature while rotating.
- 3) Cerebrospinal fluid concentration after treatment is 50%, dilute treated sample in Sample Diluent NS to in assay concentration (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 1 | 2026-04-29

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)

ab327316 – Monkey CRP SimpleStep ELISA® Kit (C-Reactive Protein), Chemiluminescent

Additional information

ASSAY SPECIFICITY

This kit is designed for the quantification of rhesus macaque CRP.

The standard protein in this kit is full length rhesus macaque CRP.

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

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

Saliva, urine, milk, cell extract, and tissue extract samples have not been tested with this kit.

CROSS REACTIVITY

Recombinant mouse CRP and rat CRP were prepared at 50 ng/mL and assayed for cross reactivity. No cross-reactivity was observed.

INTERFERENCE

Recombinant mouse CRP and rat CRP were prepared at 50 ng/mL and assayed for interference. No interference was observed.

SPECIES REACTIVITY

Other species reactivity was determined by measuring 1:500 serum samples of various species, interpolating the protein concentrations from the monkey standard curve, and expressing the interpolated concentrations as a percentage of the protein concentration in monkey serum assayed at the same dilution.

Reactivity was determined for the following species: Mouse (< 3%), Rat (< 3%), Cow (< 3%).

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	74	74	74
9.14	409	355	382
27.43	1,514	1,384	1,449
82.30	6,679	5,956	6,318
246.91	22,116	22,499	22,308
740.74	72,730	68,951	70,841
2,222.2	224,330	220,159	222,245
6,666.6	683,101	658,382	670,742
20,000	1,922,988	1,886,111	1,904,550

Table 1. Example of monkey CRP standard curve in NS. The CRP 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 1.65 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 CRP 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: 2,000 Serum	115	110 - 119

1: 2,000 Plasma - Citrate	115	113 - 117
1: 10,000 Plasma - EDTA	115	111 - 117
1: 2,000 Plasma - Heparin	113	112 - 115
25% Cell culture media*	100	97 - 103
1.56% Cerebrospinal fluid	104	99 - 110

*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 CRP was measured in the following biological samples in a 2-fold dilution series. Sample dilutions are made in NS.

Dilution Factor	Interpolated value	1:500 Monkey Serum	1:500 Monkey Plasma (Citrate)	1:2,000 Monkey Plasma (EDTA)	1:500 Monkey Plasma (Heparin)	3.13% Monkey CSF
Undiluted	pg/mL	1,780	1,597	1,471	1,714	353
	% Expected value	100	100	100	100	100
2	pg/mL	915	805	693	884	165
	% Expected value	103	101	94	103	93
4	pg/mL	429	377	332	433	79
	% Expected value	96	94	90	101	89
8	pg/mL	184	177	149	193	40
	% Expected value	83	89	81	90	90
16	pg/mL	91	81	NL	94	ND
	% Expected value	81	81	NL	88	ND

NL – Non-Linear

ND – Not Detected – below product dynamic range

Recombinant CRP was spiked in in the following biological samples in a 2-fold dilution series. Sample dilutions are made in Sample Diluent NS.

Dilution Factor	Interpolated value	25% Cell culture media
Undiluted	pg/mL	958
	% Expected value	100

2	pg/mL	454
	% Expected value	95
4	pg/mL	233
	% Expected value	97
8	pg/mL	123
	% Expected value	102
16	pg/mL	58
	% Expected value	98

Precision

Mean coefficient of variations of interpolated values of CRP from three concentrations of macaque serum within the working range of the assay.

	Intra-assay	Inter-assay
n=	8	3
CV (%)	2.4	0.9

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 1 | 2026-04-29

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)