

ab324923 – Human Complement C4 Binding Protein Beta SimpleStep ELISA® Kit (C4BPB)

For the quantitative measurement of Complement C4 Binding Protein Beta in human serum, plasma (citrate), plasma (EDTA), plasma (heparin), and cell culture media.
For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit: www.abcam.com/ab324923

This kit is available in a 384-well plate format. This plate utilises smaller volumes of standards and samples per well. Directions for using this format can be found on pages 5-6.

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.

Materials Supplied

Item	Quantity 1 x 96 tests	Quantity 10 x 96 tests	Storage Condition
Human Complement C4 Binding Protein Beta Capture Antibody 10X	600 µL	10 x 600 µL	+4°C
Human Complement C4 Binding Protein Beta Detector Antibody 10X	600 µL	10 x 600 µL	+4°C
Human Complement C4 Binding Protein Beta Lyophilized Recombinant Protein	2 Vials	10 x 2 Vials	+4°C
Antibody Diluent CPR2	6 mL	10 x 6 mL	+4°C
Cell Extraction Buffer SSW	10 mL	2 x 50 mL	+4°C
Sample Diluent NS	12 mL	N/A	+4°C
Wash Buffer PT 10X	20 mL	200 mL	+4°C
TMB Development Solution	12 mL	120 mL	+4°C
Stop Solution	12 mL	120 mL	+4°C
SimpleStep Pre-Coated 96-Well Microplate	96 wells	10 x 96 wells	+4°C
Plate Seal	1	10	+4°C

Sample Diluent NS is provided but not necessary for this product.

Materials Required, Not Supplied

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

- Microplate reader capable of measuring absorbance at 450 or 600 nm.
- Deionized water.
- Multi- and single-channel pipettes.
- Tubes for standard dilution.
- Plate shaker for all incubation steps.
- Heat block
- Optional: Phenylmethylsulphonyl 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 100 mL 1X Wash Buffer PT combine 10 mL Wash Buffer PT 10X with 90 mL deionized water. Mix thoroughly and gently.

1X Sample Diluent SSW: Prepare 1X Sample Diluent SSW by diluting Cell Extraction Buffer SSW Stock to 1X with 1X Wash Buffer PT. To make 10 mL 1X Sample Diluent SSW, combine 9.6 mL 1X Wash Buffer PT and 0.4 mL Cell Extraction Buffer SSW Stock. Mix thoroughly and gently.

Antibody Cocktail: Prepare Antibody Cocktail by diluting the capture and detector antibodies in Antibody Diluent CPR2. 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 CPR2. Mix thoroughly and gently.

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 Complement C4 Binding Protein Beta standard sample by adding the volume of 1X Sample Diluent SSW indicated on the protein vial label. Hold at room temperature for 10 minutes. Mix thoroughly and gently. This is the 32 ng/mL **Stock Standard** Solution.
2. Label eight tubes, Standards 1–8.
3. Add 294 µL of 1X Sample Diluent SSW into tube number 1 and 150 µL of 1X Sample Diluent SSW into numbers 2-8.
4. Use the **Stock Standard** to prepare the following dilution series. Standard #8 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	294	32,000	7,500
2	Standard#1	150	150	7,500	3,750
3	Standard#2	150	150	3,750	1,875
4	Standard#3	150	150	1,875	937.5
5	Standard#4	150	150	937.5	468.8
6	Standard#5	150	150	468.8	234.4
7	Standard#6	150	150	234.4	117.2
8	Blank Control	0	150	0	0

Sample Preparation

Typical Sample Dynamic Range	
Sample Type	Range
Serum	0.31 - 5%
Plasma – Citrate	0.31 - 5%
Plasma – EDTA	0.31 - 2.5%
Plasma – Heparin	0.31 - 5%
Cell Culture Media* without Fetal Bovine Serum	≤ 50%
Cell Culture Media* containing 10% Fetal Bovine Serum	≤ 50%

*Media is RPMI 1640.

Note: Heat treatment of biological samples is required before use in assay.

- 1. Dilute to appropriate concentration.
- 2. Heat sample for 30 minutes at 50°C in a heat block.
- 3. Cool to room temperature, mix gently, spin down and assay.

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 1:20 into 1X Sample Diluent SSW. Heat samples as described above 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. Dilute plasma (citrate or heparin) samples at least 1:20 and plasma (EDTA) samples at least 1:40 into 1X Sample Diluent SSW. Heat samples as described above 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. Dilute samples at least 1:2 into 1X Sample Diluent SSW. Heat samples as described above and assay. Store un-diluted samples at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.

Note: Cell culture media containing 10% fetal bovine serum (FBS) was tested and ~0.15 O.D. was observed. It is recommended to use either serum-free media or to normalize data using an appropriate media-only control.

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 1 hour at room temperature on a plate shaker set to 400 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 10 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 100 µL of TMB Development Solution to each well and incubate for 10 minutes in the dark on a plate shaker set to 400 rpm.

Given variability in laboratory environmental conditions, optimal incubation time may vary between 5 and 20 minutes.

Note: The addition of Stop Solution will change the color from blue to yellow and enhance the signal intensity about 3X. To avoid signal saturation, proceed to the next step before the high concentration of the standard reaches a blue color of O.D.600 equal to 1.0.

- 8. Add 100 µL of Stop Solution to each well. Shake plate on a plate shaker for 1 minute to mix. Record the OD at 450 nm. This is an endpoint reading.
- 9. Alternative to 7 – 8: Instead of the endpoint reading at 450 nm, record the development of TMB Substrate kinetically. Immediately after addition of TMB Development Solution begin recording the blue color development with elapsed time in the microplate reader prepared with the following settings:

Mode	Kinetic
Wavelength:	600 nm
Time:	up to 20 min
Interval:	20 sec - 1 min
Shaking:	Shake between readings

Note that an endpoint reading can also be recorded at the completion of the kinetic read by adding 100 µL Stop Solution to each well and recording the OD at 450 nm.

Download our ELISA guide for technical hints, results, calculation, and troubleshooting tips:
www.abcam.com/protocols/the-complete-elisa-guide

For technical support contact information, visit: www.abcam.com/contactus

ab324923 – Human Complement C4 Binding Protein Beta SimpleStep ELISA® Kit (C4BPB)

Additional information

ASSAY SPECIFICITY

This kit is designed for the quantification of human Complement C4 Binding Protein Beta.

The standard protein in this kit is an N-terminal fragment representing approximately 75% of the mass of the mature beta chain of human Complement C4 Binding Protein Beta.

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

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

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

INTERFERENCE

7.5 ng/mL of recombinant human Complement Factor 4 alpha, human Complement Factor 4 Binding Protein alpha, and human Protein S were tested for interference with 7.5 ng/mL of recombinant human Complement Factor 4 Binding Protein beta. No interference was observed.

SPECIES REACTIVITY

Other species reactivity was determined by measuring 1:20 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.

All values were below the detectable range of the assay for the following species: Monkey, Mouse, Rat, Cow

Other species reactivity not determined.

Note: Cell culture media containing 10% fetal bovine serum (FBS) was tested and ~0.15 O.D. was observed. It is recommended to use either serum-free media or to normalize data using an appropriate media-only control.

CALCULATION

- Calculate the average absorbance value for the blank control (zero) standards. Subtract the average blank control standard absorbance value from all other absorbance values.
- Create a standard curve by plotting the average blank control subtracted absorbance 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.

Δ Note: Most microplate 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 absorbance 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 absorbance values greater than that of the highest standard should be further diluted and reanalysed. Similarly, samples which measure at absorbance 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)	O.D. 450 nm		Mean O.D.
	1	2	
0	0.099	0.099	0.099
117.2	0.160	0.157	0.158
234.4	0.214	0.214	0.214
468.8	0.320	0.310	0.315
937.5	0.551	0.547	0.549
1,875	0.964	0.922	0.943
3,750	1.808	1.682	1.745
7,500	3.193	3.260	3.226

Table 1. Example of human Complement C4 Binding Protein Beta standard curve in 1X Sample Diluent SSW. The Complement C4 Binding Protein Beta 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 54.5 pg/mL. The MDD was determined by calculating the mean of zero standard replicates (n=24) and adding 2 standard deviations then extrapolating the corresponding concentration.

Recovery

Three concentrations of Complement C4 Binding Protein Beta were spiked into the indicated biological matrix to evaluate signal recovery in the working range of the assay.

Sample Type	Average % Recovery	Range (%)
0.63% Serum	91	85 - 97
0.63% Plasma – Citrate	94	91 - 97
0.31% Plasma – EDTA	90	87 - 94
0.63% Plasma – Heparin	84	81 - 88
50% Cell Culture Media*	95	90 - 100
50% Cell Culture Media* + 10% FBS	98	96 - 102

*Media is RPMI 1640.

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 Complement C4 Binding Protein Beta was measured in the following biological samples in a 2-fold dilution series. Sample dilutions are made in 1X Sample Diluent SSW.

Dilution Factor	Interpolated value	5% Human Serum	5% Human Plasma (Citrate)	2.5% Human Plasma (EDTA)	5% Human Plasma (Heparin)
Undiluted	pg/mL	2,833.9	2,417.6	5,020	1,388.7
	% Expected value	100	100	100	100
2	pg/mL	1,434.4	1,104.2	2,451	674.6
	% Expected value	101	91	98	97
4	pg/mL	705.5	509.6	1,227.7	316
	% Expected value	100	84	98	91
8	pg/mL	327.7	261.1	498.4	169.2
	% Expected value	93	86	79	97
16	pg/mL	162.3	128.8	NL	86.9
	% Expected value	92	85	NL	100

NL – Non-Linear

Recombinant Complement C4 Binding Protein Beta was spiked into the following biological samples and then diluted in a 2-fold dilution series. Sample dilutions are made in 1X Sample Diluent SSW.

Dilution Factor	Interpolated value	50% Cell Culture Media*	50% Cell Culture Media* with 10% FBS
Undiluted	pg/mL	3,682.3	4,487.6
	% Expected value	100	100
2	pg/mL	1,842.8	2,183.5
	% Expected value	100	97
4	pg/mL	1,009.9	1,086.0
	% Expected value	110	97
8	pg/mL	517.2	572.2
	% Expected value	112	102
16	pg/mL	246.1	285.8
	% Expected value	107	102

*Media is RPMI 1640.

Precision

Mean coefficient of variations of interpolated values of Complement C4 Binding Protein Beta from a single concentration of human serum within the working range of the assay.

	Intra-assay	Inter-assay
N=	8	3
CV (%)	6.3	9.3

Download our ELISA guide for technical hints, results, calculation, and troubleshooting tips: www.abcam.com/protocols/the-complete-elisa-guide

For technical support contact information, visit: www.abcam.com/contactus

Copyright © 2025 Abcam, All Rights Reserved. All information / detail is correct at time of going to print.
Version 1 | 2025-11-05

DIRECTIONS FOR 384-WELL PLATE FORMAT:

Materials Supplied for 384-well Format

Item	Quantity	Storage Condition
Human Complement C4 Binding Protein Beta Capture Antibody 10X	600 µL	+4°C
Human Complement C4 Binding Protein Beta Detector Antibody 10X	600 µL	+4°C
Human Complement C4 Binding Protein Beta Lyophilized Recombinant Protein	2 Vials	+4°C
Antibody Diluent CPR2	6 mL	+4°C
Cell Extraction Buffer SSW	50 mL	+4°C
Wash Buffer PT 10X	2 x 20 mL	+4°C
TMB Development Solution	2 x 12 mL	+4°C
Stop Solution	2 x 12 mL	+4°C
SimpleStep Pre-Coated 384-Well Microplate	384 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:

- Microplate reader capable of measuring absorbance at 450 or 600 nm in a 384-well plate.
- Deionized water.
- Multi- and single-channel pipettes.
- Tubes for standard dilution.
- Heat block
- Plate shaker for all incubation steps.
- Optional: Phenylmethylsulphonyl Fluoride (PMSF) (or other protease inhibitors).
- Optional: Automated liquid handler.

Reagent Preparation

Equilibrate all reagents to room temperature (18-25°C) prior to use. The kit contains enough reagents for one full plate. The sample volumes below are sufficient for running all 384 wells; adjust volumes as needed for the number of samples and dilution scheme for 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 200 mL 1X Wash Buffer PT combine 20 mL Wash Buffer PT 10X with 180 mL deionized water. Mix thoroughly and gently.

1X Sample Diluent SSW: Prepare 1X Sample Diluent SSW by diluting Cell Extraction Buffer SSW Stock to 1X with 1X Wash Buffer PT. To make 100 mL 1X Sample Diluent SSW combine 96 mL 1X Wash Buffer PT and 4 mL Cell Extraction Buffer SSW Stock. Mix thoroughly and gently. If required protease inhibitors can be added.

Antibody Cocktail: Prepare Antibody Cocktail by diluting the capture and detector antibodies in Antibody Diluent CPR2. To make 6 mL of the Antibody Cocktail combine 600 µL 10X Capture Antibody and 600 µL 10X Detector Antibody with 4.8 mL Antibody Diluent CPR2. Mix thoroughly and gently.

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 Complement C4 Binding Protein Beta standard sample by adding the volume of 1X Sample Diluent SSW indicated on the protein vial label. Hold at room temperature for 10 minutes. Mix thoroughly and gently. This is the 32 ng/mL **Stock Standard** Solution.
2. Label eight tubes, Standards 1– 8.
3. Add 147 µL of 1X Sample Diluent SSW into tube number 1 and 75 µL of 1X Sample Diluent SSW into numbers 2-8.
4. Use the **Stock Standard** to prepare the following dilution series. Standard #8 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	45	147	32,000	7,500
2	Standard#1	75	75	7,500	3,750
3	Standard#2	75	75	3,750	1,875
4	Standard#3	75	75	1,875	937.5
5	Standard#4	75	75	937.5	468.8
6	Standard#5	75	75	468.8	234.4
7	Standard#6	75	75	234.4	117.2
8	Blank Control	0	75	0	0

Plate Preparation

The 384-well plate included with this kit are supplied ready to use. It is not necessary to rinse the plate prior to adding reagents.

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 for 384-well Plate Format

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. Add 12.5 µL of all sample or standard to appropriate wells.
3. Add 12.5 µL of the Antibody Cocktail to each well.
4. Seal the plate and incubate for 1 hour at room temperature on a plate shaker set to 700 rpm.
5. Wash each well with 3 x 100 µL 1X Wash Buffer PT. Wash by aspirating or decanting from wells then dispensing 100 µL 1X Wash Buffer PT into each well. Wash Buffer PT should remain in wells for at least 10 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.
6. Add 25 µL of TMB Development Solution to each well and incubate for 10 minutes in the dark on a plate shaker set to 700 rpm.
Given variability in laboratory environmental conditions, optimal incubation time may vary between 5 and 20 minutes.
Note: The addition of Stop Solution will change the color from blue to yellow and enhance the signal intensity about 3X. To avoid signal saturation, proceed to the next step before the high concentration of the standard reaches a blue color of O.D.600 equal to 1.0.
7. Add 25 µL of Stop Solution to each well. Shake plate on a plate shaker for 1 minute to mix. Record the OD at 450 nm. This is an endpoint reading. Proper mixing of the Stop Solution is required for proper measurement.
8. Alternative to 6 – 7: Instead of the endpoint reading at 450 nm, record the development of TMB Substrate kinetically. Immediately after addition of TMB Development Solution begin recording the blue color development with elapsed time in the microplate reader prepared with the following settings:

Mode	Kinetic
Wavelength:	600 nm
Time:	up to 20 min
Interval:	20 sec – 1 min
Shaking:	Shake between readings

Note that an endpoint reading can also be recorded at the completion of the kinetic read by adding 25 µL Stop Solution to each well and recording the OD at 450 nm.

Download our ELISA guide for technical hints, results, calculation, and troubleshooting tips:
www.abcam.com/protocols/the-complete-elisa-guide

For technical support contact information, visit: www.abcam.com/contactus