

## AB315039 – Human Brevican SimpleStep ELISA Kit

For the quantitative measurement of Brevican in Human serum, plasma (EDTA), plasma (heparin), cell culture supernatant, and CSF.

For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit: [www.abcam.com/AB315039](http://www.abcam.com/AB315039)

**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	Storage Condition
Human Brevican Capture Antibody 10X	600 µL	+4°C
Human Brevican Detector Antibody 10X	600 µL	+4°C
Human Brevican 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
TMB Development Solution	12 mL	+4°C
Stop Solution	12 mL	+4°C
SimpleStep Pre-Coated 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:

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.

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

**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.

**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.

### 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 Brevican standard sample by adding the volume indicated on the protein vial label. For **serum, plasma (EDTA), and plasma (heparin) samples measurements**, use Sample Diluent NS + 1X Enhancer. For cell culture supernatant **samples measurements**, use Sample Diluent NS. For **CSF samples measurements**, use 1X Wash Buffer PT. Hold at room temperature for 10 minutes. Mix thoroughly and gently. This is the 16,000 pg/mL **Stock Standard** Solution.
1. Label eight tubes, Standards 1–8.
2. Use the same Sample Diluent as used to resuspend the Stock Standard to prepare the standard curve. Add 375 µL of Sample Diluent into tube number 1 and 150 µL of Sample Diluent into numbers 2-8.
3. 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	<b>Stock Standard</b>	25	375	16,000	1,000
2	Standard#1	150	150	1,000	500
3	Standard#2	150	150	500	250
4	Standard#3	150	150	250	125
5	Standard#4	150	150	125	62.5
6	Standard#5	150	150	62.5	31.25
7	Standard#6	150	150	31.25	15.63
8	Blank Control	0	150	N/A	N/A

## Sample Preparation

Typical Sample Dynamic Range	
Sample Type	Range
Serum	≤25%
Plasma – EDTA	≤50%
Plasma – Heparin	≤25%
SH-SY5Y Cell Culture Supernatant	≤100%
Cerebrospinal Fluid	≤2.25%

**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 + 1X Enhancer and assay. Store un-diluted serum at -20°C or below. Avoid repeated freeze-thaw cycles.

**Plasma** Collect plasma using EDTA or Heparin. Centrifuge samples at 2,000 x g for 10 minutes. Dilute samples at least 1:2 for plasma (EDTA) or 1:4 for plasma (heparin) into Sample Diluent NS + 1X Enhancer 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 (citrate) samples.

**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.

**Cerebrospinal Fluid (CSF)** Dilute cerebrospinal fluid at least 1:44.4 into 1X Wash Buffer PT and assay. Store un-diluted 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, 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 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](http://www.abcam.com/protocols/the-complete-elisa-guide)

**For technical support contact information, visit:** [www.abcam.com/contactus](http://www.abcam.com/contactus)

# AB315039 – Human Brevican SimpleStep ELISA® Kit

## Additional information

### ASSAY SPECIFICITY

This kit is designed for the quantification of Human Brevican.

The standard protein in this kit is an N-terminal fragment representing ~38% of mature Human Brevican.

Native signal was detected in serum, plasma (EDTA), plasma (heparin), cell culture supernatant, and CSF.

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

This kit is incompatible with plasma (citrate).

### SPECIES REACTIVITY

Other species reactivity was determined by measuring 100% 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.

Reactivity < 3% was determined for the following species: Cow

No signal was observed in undiluted serum samples from the following species: Monkey, Mouse, and Rat

Other species reactivity not determined.

### 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	0.041	0.042	0.041
15.63	0.095	0.092	0.093
31.25	0.141	0.138	0.139
62.5	0.239	0.222	0.230
125	0.437	0.430	0.433
250	0.769	0.757	0.763
500	1.544	1.527	1.535
1,000	2.961	2.810	2.886

Table 1. Example of Human Brevican standard curve in Sample Diluent NS + 1X Enhancer. The Brevican standard curve was prepared as described in the Standard Preparation section. The table shows raw data values.

Standard Curve Measurements			
Concentration (pg/mL)	O.D 450 nm		Mean O.D
	1	2	
0	0.041	0.041	0.041
15.63	0.093	0.091	0.092
31.25	0.147	0.142	0.145
62.5	0.261	0.263	0.262
125	0.465	0.446	0.456
250	0.856	0.824	0.840
500	1.671	1.648	1.659
1,000	3.158	3.206	3.182

Table 2. Example of Human Brevican standard curve in Sample Diluent NS. The Brevican standard curve was prepared as described in the Standard Preparation section. The table shows raw data values.

Standard Curve Measurements			
Concentration (pg/mL)	O.D 450 nm		Mean O.D
	1	2	
0	0.043	0.043	0.043
15.63	0.087	0.089	0.088
31.25	0.131	0.136	0.133
62.5	0.199	0.210	0.205
125	0.381	0.373	0.377
250	0.720	0.766	0.743
500	1.361	1.469	1.415
1,000	2.905	2.905	2.905

Table 3. Example of Human Brevican standard curve in 1X PT Wash Buffer. The Brevican standard curve was prepared as described in the Standard Preparation section. The table shows raw data values.

### TYPICAL SAMPLE VALUES

#### Sensitivity:

The minimal detectable dose (MDD) was determined by calculating the mean of zero standard replicates and adding 2 standard deviations then extrapolating the corresponding concentration.

Sample Diluent Buffer	N=	Minimal Detectable Dose
Sample Diluent NS + 1X Enhancer	24	2.63 pg/mL
Sample Diluent NS	32	2.28 pg/mL
1X Wash Buffer PT	32	2.32 pg/mL

#### Recovery

Three concentrations of Brevican 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% Human Serum	87	86 - 88
25% Human Plasma (EDTA)	92	90 - 95
25% Human Plasma (Heparin)	83	78 - 89
50% Cell Culture Supernatant	106	87 - 119
0.38% CSF	100	94 - 107

RPMI 1640 containing 10% fetal calf 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 Brevican was measured in the following biological samples in a 2-fold dilution series. Sample dilutions are made in Sample Diluent NS + 1X Enhancer for serum and plasma; Sample Diluent NS for cell culture supernatant; and 1X Wash Buffer PT for CSF.

Dilution Factor	Interpolated value	25% Human Serum	50% Human Plasma (EDTA)	25% Human Plasma (Heparin)	100% SH-SY5Y Supernatant	2.25% CSF
Undiluted	pg/mL	146.1	176.8	129.4	245.5	1056.9
	% Expected value	100	100	100	100	100
2	pg/mL	80.5	97.3	72.3	127.5	535.5
	% Expected value	110	110	112	104	101
4	pg/mL	46.4	50.8	38.3	70.4	260.9
	% Expected value	127	115	118	115	99
8	pg/mL	20	25.6	20.1	35.8	131
	% Expected value	120	116	124	117	99
16	pg/mL	ND	ND	ND	ND	64
	% Expected value	ND	ND	ND	ND	97

ND – Not Detected, below product dynamic range

#### Precision

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

	Intra-assay	Inter-assay
N=	8	3
CV (%)	2.17	0.76

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

[www.abcam.com/protocols/the-complete-elisa-guide](http://www.abcam.com/protocols/the-complete-elisa-guide)

For technical support contact information, visit: [www.abcam.com/contactus](http://www.abcam.com/contactus)